



---

---

# SUPELCO

## Air Methods Guide

---

---

**2nd edition**

T411091D

 **SUPELCO** ISO 9001  
REGISTERED

©1999 Sigma-Aldrich Co.

# Supelco Air Methods Guide

The Supelco Air Methods Guide is a resource of sampling and analytical methods for more than 1000 compounds in air. Air monitoring of many matrices – workplace, indoor, ambient, and source – are integrated into this one easy-to-read reference.

## The guide lists:

- compound (alphabetically)
- method reference
- sample volume and flow rate
- sampling media description and Supelco catalog number
- analytical technique

Note that a compound of interest may be listed by more than one chemical name and/or category classification (e.g., VOCs).

## The guide also features:

- NIOSH, OSHA, EPA, and ASTM validated methods
- cross reference to SKC part numbers
- latest updates through May 1999

### **Please note:**

This guide is for general information only. It is not intended to replace careful review of applicable agency methods, equipment operating instructions, or national or regional legal regulations.

## Key to Abbreviations

### Sampling Devices

Abbrev.	Description
1-2PP	1-(2-pyridyl)piperazine
2-HMP c/w	2-(hydroxymethyl)piperadine coated with
DMBA	3,4-dimethoxybenzylamine
DNPH	2,4-dinitrophenylhydrazine
DOP	di-n-octyl phthalate
MCE	mixed cellulose ester
NBDC	7-chloro-4-nitrobenzo-2-oxa-1,3-diazole
NITC	1-naphthylisothiocyanate
NPMPA	N-[(4-nitrophenyl)methyl]propylamine
OVS	OSHA versatile sampler
PVC t/w	polyvinyl chloride treated with
TBC	4-tert-butylcatechol
TEA	triethanolamine

### Other

Abbrev.	Description
NA	not available

### Analytical Techniques

Abbrev.	Description
AACV	atomic adsorption cold vapor
AAFAG	atomic adsorption flame arsine generation
AAGF	atomic adsorption graphite furnace
AAS	atomic adsorption spectrophotometry
AAS-HGA	flame atomic adsorption spectrophotometry - heated graphite furnace atomizer
ASV	anodic stripping voltametry
DPP	differential pulse polarography
EGA-TOA	evolved gas analysis - thermal optical sensor
ELISA	enzyme linked immunosorbent assay
FAAS	flame atomic adsorption spectrophotometry
GC-AED	gas chromatography - atomic emission detector
GC-ECD	gas chromatography - electron capture detector
GC-ECN	gas chromatography - electrolytic conductivity detector
GC-ELCD	gas chromatography - electrochemical detector
GC-FID	gas chromatography - flame ionization detector
GC-FPD	gas chromatography - flame photometric detector
GC-MS	gas chromatography - mass spectrometry
GC-NPD	gas chromatography - nitrogen/phosphorus detector
GC-PID	gas chromatography - photoionization detector
GC-SCLD	gas chromatography - sulfur chemiluminescence detector
GC-TCD	gas chromatography - thermal conductivity detector
GC-TEA	gas chromatography - thermal energy analyzer
HPLC-AAGF	high performance liquid chromatography - atomic adsorption graphite furnace
HPLC-FL	high performance liquid chromatography - fluorescence
HPLC-TEA/UV	high performance liquid chromatography - thermal energy analyzer/ultraviolet
HPLC-UV	high performance liquid chromatography - ultraviolet
HPLC-UV/ELCD	high performance liquid chromatography - ultraviolet/electrochemical detector
HPLC-UV/FL	high performance liquid chromatography - ultraviolet/fluorescence
HPLC-VIS	high performance liquid chromatography - visible absorption spectrophotometry
IC	ion chromatography
IC-AAS	ion chromatography - atomic absorption spectrophotometry
IC-CD	ion chromatography - conductivity detector
IC-ELCD	ion chromatography - electrochemical detector
IC-PCR	ion chromatography - post-column reactor
IC-PED	ion chromatography - pulsed electrochemical detector
IC-UV	ion chromatography - ultraviolet
ICP-AES	inductively coupled plasma/atomic emission spectroscopy
IR	infrared spectrophotometry
ISE	ion-specific electrode
NDIR	nondispersive infrared spectrophotometry
PCD	post-column derivatization
PCM	phase contrast microscopy
PES	plasma emission spectrometry
PLM	polarized light microscopy
TD-GC/MS	thermal desorption - gas chromatography - mass spectrometry
TD-IR	thermally desorbed infrared spectrophotometry
TEM	transmission electron microscopy
UV	ultraviolet
VIS	visible absorption spectrophotometry
XPD	x-ray powder diffraction
XRD	x-ray diffraction
XRF	x-ray fluorescence
XRF/XRD	x-ray fluorescence/x-ray diffraction

### Trademarks

Anasorb – SKC, Inc.  
Cellosolve – Union Carbide Corp.  
Chromosorb – Celite Corp.  
Florisil – U.S. Silica Co.  
Freon – E.I. duPont de Nemours & Co., Inc.  
Gas Chrom – Applied Science Laboratories, Inc.  
Hopcalite – Mine Safety Appliances Company.  
Porapak – Waters Associates, Inc.  
Tedlar – E.I. du Pont de Nemours & Co., Inc.  
Teflon – E.I. du Pont de Nemours & Co., Inc.  
Tenax – Enka Research Institute Arnhem  
XAD – Rohm and Haas Co.  
Carbosieve, Carbotrap, Carboxen, ORBO,  
Supelpak – Sigma-Aldrich Co.

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equip	Analytical Technique
Acenaphthene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Acenaphthene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	GC-FID
Acenaphthylene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Acenaphthylene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	GC-FID
Acetaldehyde	NIOSH 2538	1-12	0.01-0.05	XAD-2 c/w 2-HMP	ORBO-25	20357	226-27	GC-FID
Acetaldehyde	NIOSH 2539	5	0.01-0.05	XAD-2 c/w 2 HMP	ORBO-23	20257-U	226-118	GC-FID
Acetaldehyde	NIOSH 3507	6-60	0.1-0.5	fritted bubbler	-----	64835-U	225-36-2	HPLC-UV
Acetaldehyde	OSHA 68	3	0.05	XAD-2 c/w 2-HMP	ORBO-25	20357	226-27	GC-NPD
Acetic Acid	NIOSH 1603	20-300	0.01-0.1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Acetic Acid	US Army	10-25		Chromosorb P c/w Na2CO3	ORBO-70	20256-U	NA	GC-FID
Acetic Anhydride	NIOSH 3506	25-100	0.2-1	fritted bubbler	-----	64835-U	225-36-2	VIS
Acetic Anhydride	OSHA 102	7.5	0.05-0.5	glass fiber filter c/w DMBA & DOP	894	custom	225-9010	GC-NPD
Acetic Anhydride	OSHA 82	0.75	0.05	glass fiber filter c/w 1-2PP	821	custom	225-9009	GC-NPD
Acetone	NIOSH 1300	0.50-3	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Acetone	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Acetone	OSHA 69	3	0.05	Carbosieve S-III (Anasorb CMS)	ORBO-91	20360	226-121	GC-FID
Acetone Cyanohydrin	NIOSH 2506	0.30-12	0.2	Porapak QS	ORBO-1101	20061	226-59-09	GC-NPD
Acetonitrile	NIOSH 1606	1-25	0.01-0.2	coconut charcoal	ORBO-32L	20228	226-09	GC-FID
Acetylene Dichloride	NIOSH 1003	0.2-5	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Acetylene Tetrabromide	NIOSH 2003	50-100	0.2-1	silica gel	ORBO-52S	20229	226-10	GC-FID
Acetylene Tetrachloride	NIOSH 1019	3-30	0.01-0.2	petroleum charcoal	ORBO-303	20040-U	226-38	GC-FID
Acid Mist	OSHA ID 165SG	96	0.20	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Acids, Inorganic	NIOSH 7903	3-100	0.2-0.5	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Acridine	OSHA 58	960	2	glass fiber filter	-----	23378	225-7	HPLC-UV
Acrolein	NIOSH 2501	1.50-48	0.01-0.1	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-NPD
Acrolein	NIOSH 2539	5	0.01-0.05	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-FID
Acrolein	OSHA 52	48	0.1	XAD-2 c/w 2-HMP	ORBO-24	20231	226-117	GC-NPD
Acrylamide	OSHA 21	120	1	silica gel glass fiber filter	ORBO-52S ----	20229 23376	226-10 225-16	GC-NPD
Acrylic Acid	OSHA 28	24	0.1	XAD-8	ORBO-601	20048	226-30-08	HPLC-UV
Acrylic acid	OSHA 312	24	0.1	(2) XAD-8	ORBO-601	20048	226-30-08	HPLC-UV
Acrylonitrile	NIOSH 1604	3.50-20	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Acrylonitrile	OSHA 37	20	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-NPD
Aerobic Bacteria by GC-FAME	NIOSH 0801	50-300	28.30	Andersen Impactor				GC-FID
Air Exchange Rate	EPA IP-4A			perfluorocarbon tracer				GC-ECD
Air Exchange Rate	EPA IP-4B			tracer gas				GC-ECD
Alachlor	NIOSH 5603	70-1750	1	SPE disk				ELISA
Aldehydes	EPA 0011			DNPH absorbing solution				HPLC-UV
Aldehydes (screening)	NIOSH 2539	5	0.01-0.05	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-FID
Aldehydes/Ketones	EPA 0100		1	silica gel c/w DNPH	LpDNPH S10	21026-U	226-119	HPLC-UV
Aldehydes/Ketones	EPA IP-6A		0.5-1.2	silica gel c/w DNPH	LpDNPH S10	21026-U	226-119	HPLC-UV
Aldehydes/Ketones	EPA IP-6B			continuous colorimetric analyzer				colorimetric
Aldehydes/Ketones	EPA IP-6C			diffusive sampler	-----	NA		HPLC-UV
Aldehydes/Ketones	EPA TO-11A	280	0.1-2	silica gel c/w DNPH ozone scrubber	LpDNPH S10 -----	21026-U 505285	226-119 NA	HPLC-UV
Aldehydes/Ketones	EPA TO-5	<80	0.1-1	DNPH absorbing solution				HPLC-UV
Aldicarb (Temik)	OSHA 74	480	1	XAD-2/glass fiber filter (OVS)	ORBO-49P	20350	226-30-16	GC-NPD
Aldrin	NIOSH 5502	18-240	0.2-1	glass fiber filter fritted bubbler	----- -----	23378 64835-U	225-7 225-36-2	GC-ECD
Alkaline Dusts	NIOSH 7401	70-1000	1-4	PTFE filter	-----	23383	225-1710	titration
Allyl Alcohol	NIOSH 1402	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Allyl Chloride	NIOSH 1000	16-100	0.01-1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Allyl Chloride	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Allyl Glycidyl Ether	NIOSH 2545	1.5-8	0.01-0.2	Tenax	ORBO-402	20832-U	226-35-03	GC-FID
Allyl Trichloride	NIOSH 1003	2-60	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Alumina	NIOSH 0500	7-133	1-2	PVC filter	-----	23387	225-8-01	gravimetric
Aluminum	NIOSH 7300	5-100	1-4	MCE filter	-----	23381	225-5	ICP-AES
Aluminum, Al compounds (as Al)	NIOSH 7013	10-400	1-3	MCE filter	-----	23381	225-5	FAAS
Amines, Aliphatic	NIOSH 2010	3-30	0.01-1	silica gel	ORBO-52S	20229	226-10	GC-FID
Amines, Aliphatic	Supelco			Carbotrap	ORBO-100	20255-U	NA	GC-NPD
Amines, Aromatic	NIOSH 2002	30-150	0.2-1	silica gel	ORBO-52S	20229	226-10	GC-FID
Aminobenzene	NIOSH 2002	30-150	0.2-1	silica gel	ORBO-52S	20229	226-10	GC-FID
4-Aminodiphenyl (4-Aminobiphenyl)	NIOSH 4/269	0.2		silica gel glass fiber filter	ORBO-502 -----	20030-U 23376	226-51 225-16	GC-FID
4-Aminodiphenyl (4-Aminobiphenyl)	OSHA 93	100	1.0	glass fiber filter c/w H2SO4	822	20065	225-9004	GC-ECD
Aminoethanol Compounds II	NIOSH 3509	5-300	0.5-1	impinger	-----	20270-U	225-36-1	IC
2-Aminoethanol (Ethanolamine)	NIOSH 2007	4-24	0.01-0.2	silica gel	ORBO-506	20032	226-10-04	GC-FID
2-Aminoethanol (Ethanolamine)	NIOSH 3509	5-300	0.5-1	impinger	-----	20270-U	225-36-1	IC
p-Aminophenylarsonic Acid	NIOSH 5022	50-1000	1-3	PTFE filter	-----	23383	225-1710	IC-AAS
2-Aminopyridine	NIOSH 4/S158	12	0.01-0.2	Tenax	ORBO-405	custom	226-35	GC-FID
2-Aminotoluene	NIOSH 2002	10-150	0.02-1	silica gel	ORBO-52S	20229	226-10	GC-FID
Ammonia	NIOSH 6015	0.1-96	0.1-0.2	silica gel c/w H2SO4	ORBO-555	custom	226-10-06	VIS
Ammonia	NIOSH 6016	0.1-96	0.1-0.5	silica gel c/w H2SO4	ORBO-555	custom	226-10-06	IC-CD
Ammonia	OSHA ID 188	7.50-24	0.1-0.5	Purosieve c/w H2SO4	ORBO-77	20036	226-29	IC
Ammonium Sulfamate	NIOSH 5/S348	90	1.5-2	MCE filter	-----	23381	225-5	IC-CD
Amyl Acetate	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
n-Amyl Acetate	NIOSH 1450	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Amyl Acetate	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
sec-Amyl Acetate	NIOSH 1450	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
sec-Amyl Acetate	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Aniline	NIOSH 2002	5-30	0.02-0.2	silica gel	ORBO-52S	20229	226-10	GC-FID
Aniline	NIOSH 2017	5-50	0.2	silica gel + glass fiber filter c/w H2SO4	ORBO-507 822	20870-U 20065	226-15 225-9004	GC-FID
Anisidine	NIOSH 2514	24-320	0.5-1	XAD-2	ORBO-608	20050-U	226-30-05	HPLC-UV
Anthracene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Anthracene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	GC-FID
Anthracene	OSHA 58	960	2	glass fiber filter	-----	23378	225-7	HPLC-UV
Antimony	NIOSH 1/173	180	1.5	MCE filter	-----	23381	225-5	AAS
Antimony	NIOSH 1/189		1.5-2	MCE filter	-----	23381	225-5	ASV
Antimony	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Antimony, Sb Compounds (as Sb)	NIOSH 2/S2		1.5	MCE filter	-----	23381	225-5	AAS
Antimony, Sb Compounds (as Sb)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Antimony, Sb Compounds (as Sb)	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Antimony Particulates	NIOSH 4/261	45	1.5	MCE filter	-----	23381	225-5	AAGF
ANTU (Alphanaphthyl thiourea)	NIOSH 5/S276	480	1.5-2	PTFE filter	-----	23383	225-1710	HPLC-UV
Aqua Fortis	NIOSH 7903	3-100	0.2-0.5	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
p-Arsanilic acid	NIOSH 5022	50-1000	1-3	PTFE filter	-----	23383	225-1710	IC-AAS
Arsenic	NIOSH 7300	5-2000	1-4	MCE filter	-----	23381	225-5	ICP-AES
Arsenic, As Compounds (as As)	NIOSH 7900	30-1000	1-3	MCE filter	-----	23381	225-5	AAFAG
Arsenic, Organo	NIOSH 5022	50-1000	1-3	PTFE filter	-----	23383	225-1710	IC-AAS
Arsenic, Particulate & Volatiles	OSHA ID 105	480-960	2	MCE filter + support pad c/w Na2CO3	845	custom	225-9005 225-9001	AAGF

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equip	Analytical Technique
Arsenic Trioxide (as As)	NIOSH 7901	30-1000	1-3	MCE filter w/ support pad c/w Na2CO3	824	20067	NA	AAGF
Arsine	NIOSH 6001	0.1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	AAGF
Arsine	OSHA ID 105	120-240	0.5	MCE filter + support pad c/w Na2CO3	845	custom	225-9005 225-9001	AAGF
Asbestos	NIOSH 9000			bulk		-----		XRD
Asbestos	NIOSH 9002			bulk		-----		PLM
Asbestos	OSHA ID 191			bulk		-----		PLM
Asbestos Fibers	NIOSH 7400	400-	0.5-16	MCE filter	-----	23380-U	225-19	PCM
Asbestos Fibers	NIOSH 7402	400-	0.5-16	MCE filter	-----	23380-U	225-19	TEM
Asbestos in Air	OSHA ID 160	25-1200	0.5-2.5	MCE filter	-----	23380-U	225-19	PCM
Aspartame	NIOSH 5031	70-1200	1-3	PTFE filter	-----	23383	225-1710	HPLC-UV
Asphalt Fumes (Petroleum)	OSHA 58	960	2	glass fiber filter	-----	23378	225-7	HPLC-UV
Azelaic Acid	NIOSH 5019	200-1000	1-3	PVC filter	-----	23387	225-8-01	GC-FID
Azinphos, Methyl	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Barium	NIOSH 1/173	180	1.5	MCE filter	-----	23381	225-5	AAS
Barium (soluble compounds)	NIOSH 7056	50-2000	1-4	MCE filter	-----	23381	225-5	AAS
Barium (soluble compounds)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Benz(a)anthracene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----- 23390-U	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Benz(a)anthracene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----- 23390-U	20258 23390-U	226-30-04 225-17-07	GC-FID
Benzaldehyde	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Benzene	NIOSH 1500	2-30	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Benzene	NIOSH 1501	5-30	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Benzene	NIOSH 2549	1-6	0.01-0.5	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Benzene	NIOSH 3700	<80% vol	0.02-5	5L sampling bag	-----	24655	232-05	GC-PID
Benzene	OSHA 12	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Benzene-Soluble Fraction & Total Particulate (Asphalt Fume)	NIOSH 5042	28-400	1-4	PTFE filter	-----	23390-U	225-17-07	gravimetric
Benzene-Soluble Particulate Matter	ASTM D4600	960	2	glass fiber filter	-----	23378	225-7	gravimetric
Benzidine	NIOSH 5509	20-100	0.2	glass fiber filter	-----	23376	225-16	HPLC-UV
Benzidine	OSHA 65	100	1	glass fiber filter c/w H2SO4	822	20065	225-9004	GC-ECD
Benzidine Dyes	NIOSH 5013	150-500	1-3	PTFE filter	-----	23389	225-17	HPLC-UV
Benzin (petroleum ether)	NIOSH 1550	1.3-20	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Benzo(a)pyrene (BAP)	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----- 23390-U	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Benzo(a)pyrene (BAP)	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----- 23390-U	20258 23390-U	226-30-04 225-17-07	GC-FID
Benzo(a)pyrene (BAP)	OSHA 58	960	2	glass fiber filter	-----	23378	225-7	HPLC-UV
Benzo(b)fluoranthene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----- 23390-U	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Benzo(b)fluoranthene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----- 23390-U	20258 23390-U	226-30-04 225-17-07	GC-FID
Benzo(e)pyrene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----- 23390-U	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Benzo(e)pyrene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----- 23390-U	20258 23390-U	226-30-04 225-17-07	GC-FID
Benzo(ghi)perylene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----- 23390-U	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Benzo(ghi)perylene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----- 23390-U	20258 23390-U	226-30-04 225-17-07	GC-FID
Benzo(k)fluoranthene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----- 23390-U	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Benzo(k)fluoranthene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----- 23390-U	20258 23390-U	226-30-04 225-17-07	GC-FID
Benzothiazole in Asphalt Fume	NIOSH 2550	480-960	1-2	XAD-2 PTFE filter	ORBO-42L -----	20264-U 23390-U	NA 225-17-07	GC-SCLD
Benzoyl Peroxide	NIOSH 5009	40-400	1-3	MCE filter	-----	23381	225-5	HPLC-UV

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equip	Analytical Technique
Benzyl Chloride	NIOSH 1003	6-50	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Benzyl Chloride	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Beryllium	NIOSH 7300	1250-2000	1-4	MCE filter	----	23381	225-5	ICP-AES
Beryllium	OSHA ID 125G	480	2	MCE filter	----	23381	225-5	ICP-AES
Beryllium, Be Compounds (as Be)	NIOSH 7102	25-1000	1-4	MCE filter	----	23381	225-5	AAGF
BHT (Butylated Hydroxytoluene)	NIOSH 1/226	10	0.2	silica gel	ORBO-504	20031	NA	GC-FID
Bioaerosol Sampling	NIOSH 0800			Andersen cascade impactor Andersen N-6 sampler				
Biphenyl (Diphenyl)	NIOSH 2530	15-30	0.01-0.5	Tenax	ORBO-404	custom	226-35-01	GC-FID
bis-Chloromethyl Ether	NIOSH 1/213	5	0.08	Chromosorb 101	ORBO-703	custom	226-102	GC-MS
bis-Chloromethyl Ether	OSHA 10	50	0.5	fritted bubbler	----	64835-U	225-36-2	GC-ECD
Bismuth	NIOSH 1/173	180	1.5	MCE filter	----	23381	225-5	AAS
Bisphenol A (BPA)	NIOSH 6/333	288	1.6	glass fiber filter	----	23378	225-7	HPLC-UV
Boroethane (Diborane)	NIOSH 6006	60-260	0.5-1	charcoal c/w oxidizer PTFE filter	ORBO-352 ----	20043 special order	226-151	PES
Boron Carbide	NIOSH 7506	100-1000	1.7 or 2.2	PVC filter	----	23387	225-8-01	XPD
Boron Oxide	NIOSH 0500	25-133	1.5-2	PVC filter	----	23387	225-8-01	gravimetric
BPA (Bisphenol A)	NIOSH 6/333	288	1.6	glass fiber filter	----	23378	225-7	HPLC-UV
Bromine	NIOSH 6011	8-360	0.3-1	PTFE filter + silver membrane	885	custom	225-9006	IC-CD
Bromine	OSHA ID 108	7.5-30	0.5	fritted bubbler	----	64835-U	225-36-2	IC
Bromoethane (Ethyl Bromide)	NIOSH 1011	0.5-4	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Bromoform	NIOSH 1003	4-70	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Bromoform	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Bromotrifluoromethane	NIOSH 1017	0.1-1	0.01-0.05	(2) coconut charcoal	ORBO-32S + ORBO-32L	20267-U 20228	226-01 226-09	GC-FID
Bromoxynil	NIOSH 5010	2-400	1-3	PTFE filter	----	23390-U	225-17-07	HPLC-UV
Bromoxynil Octanoate	NIOSH 5010	90-400	1-3	PTFE filter	----	23390-U	225-17-07	HPLC-UV
1,3-Butadiene	NIOSH 1024	5-25	0.01-0.5	coconut charcoal	ORBO-32L	20228	226-09	GC-FID
1,3-Butadiene	OSHA 56	3	0.05	charcoal c/w TBC	ORBO-351	20042	226-73	GC-FID
1-Butanethiol (n-Butyl Mercaptan)	NIOSH 2525	1-4	0.01-0.05	Chromosorb 104	----	NA	226-109	GC-FPD
2-Butanone (MEK)	NIOSH 2500	.25-12	0.01-0.2	Carboxen 564	ORBO-90	20358	226-81 226-121	GC-FID
2-Butanone (MEK)	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
2-Butanone (MEK)	OSHA 84	3	0.05	Carbosieve S-III	ORBO-91	20360	226-121	GC-FID
2-Butoxyethanol (Butyl Cellosolve)	NIOSH 1403	2-10	0.01-0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2-Butoxyethanol	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
2-Butoxyethanol	OSHA 83	48	0.1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2-Butoxyethyl Acetate (Butyl Cellosolve Acetate)	OSHA 83	48	0.1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Butyl Acetate	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Butyl Acetate	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Butyl Acetate, sec- & tert-	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Butyl Acetate	NIOSH 1450	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
sec-Butyl Acetate	NIOSH 1450	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
tert-Butyl Acetate	NIOSH 1450	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Butyl Alcohol	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Butyl Alcohol	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Butyl Alcohol, n- & sec-	NIOSH 1401	2-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
sec-Butyl Alcohol	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
tert-Butyl Alcohol	NIOSH 1400	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
tert-Butyl Alcohol	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Butylamine	NIOSH 2012	2-100	0.01-1	silica gel c/w H2SO4	ORBO-554	20033	226-53	GC-FID

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Butyl Cellosolve - see 2-Butoxyethanol								
Butyl Cellosolve Acetate - see 2-Butoxyethyl Acetate								
n-Butyl Glycidyl Ether	NIOSH 1616	15-30	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Butyl Glycidyl Ether	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Butyl Mercaptan	NIOSH 2525	1-4	0.01-0.05	Chromosorb 104	-----	NA	226-109	GC-FPD
n-Butyl Mercaptan	NIOSH 2542	10-150	0.01-0.2	glass fiber filter c/w mercuric acetate	826	20068	225-9007	GC-FPD
p-tert-Butyltoluene	NIOSH 1501	1-29	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
p-tert-Butyltoluene	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Butylated Hydroxytoluene (BHT)	NIOSH 1/226	10	0.2	silica gel	ORBO-504	20031	NA	GC-FID
1,3-Butylene Glycol	NIOSH 5523	5-60	0.5-2	XAD-7/glass fiber filter (OVS)	ORBO-47P	custom	226-57	GC-FID
Butyraldehyde	NIOSH 2539	5	0.01-0.05	XAD-2 c/w 2 HMP	ORBO-23	20257-U	226-118	GC-FID
Cadmium	NIOSH 7300	13-2000	1-4	MCE filter	-----	23381	225-5	ICP-AES
Cadmium	OSHA ID 189	960	2.0	MCE filter	-----	23381	225-5	AAS or AAS-HGA
Cadmium, Cd Compounds (as Cd)	NIOSH 7048	25-1500	1-3	MCE filter	-----	23381	225-5	AAS
Cadmium Dust (as Cd)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Cadmium Dust (as Cd)	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Cadmium Fume (as Cd)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Cadmium Fume (as Cd)	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Calcium	NIOSH 7300	5-200	1-4	MCE filter	-----	23381	225-5	ICP-AES
Calcium & compounds	NIOSH 7020	20-400	1-3	MCE filter	-----	23381	225-5	AAS
Camphor	NIOSH 1301	1-25	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Camphor	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Capsaicin	NIOSH 5041	5-1000	1-3	glass fiber filter	-----	23376	225-16	HPLC-FL
Carbaryl (Sevin)	NIOSH 5006	20-400	1-3	glass fiber filter	-----	23378	225-7	VIS
Carbaryl (Sevin)	OSHA 63	60	1	XAD-2/glass fiber filter (OVS)	ORBO-49P	20350	226-30-16	HPLC-UV
Carbitol	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Carbon, Elemental (Diesel Particulate)	NIOSH 5040	142-19,000	2-4	quartz fiber filter (37mm)	-----	special order	225-1809	EGA-TOA
Carbon Black	NIOSH 5000	30-570	1-2	PVC filter	-----	23387	225-8-01	gravimetric
Carbon Black	OSHA ID 196	480-960	2	PVC filter	-----	23387	225-8-01	gravimetric
Carbon Dioxide	EPA IP-3A			nondispersive infrared				NDIR
Carbon Dioxide	EPA IP-3B			continuous gas filter correlation				CFC Spectrometry
Carbon Dioxide	EPA IP-3C			electrochemical CO monitor				Electrochemical Oxidation
Carbon Dioxide	NIOSH 6603	80% vol	0.02-1	5L sampling bag	-----	24655	232-05	GC-TCD
Carbon Dioxide	OSHA ID 172	2-5	0.01-0.05	5L sampling bag	-----	24655	232-05	GC-TCD
Carbon Disulfide	NIOSH 1600	2-25	0.01-0.2	coconut charcoal in-line trap	ORBO-32S	20267-U	226-01	GC-FPD
Carbon Monoxide	EPA IP-3A			nondispersive infrared				NDIR
Carbon Monoxide	EPA IP-3B			continuous gas filter correlation				CFC Spectrometry
Carbon Monoxide	EPA IP-3C			electrochemical CO monitor				Electrochemical Oxidation sensor
Carbon Monoxide	NIOSH 6604			portable instrument				
Carbon Monoxide	OSHA ID 209			direct reading instrument				
Carbon Monoxide	OSHA ID 210	2-5	0.01-0.05	5L sampling bag	-----	24655	232-05	GC-TCD
Carbon Tetrachloride	NIOSH 1003	3-150	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Carbon Tetrachloride	OSHA 7	15	0.1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Carbonyls - see Aldehydes/Ketones								

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
delta-3-Carene	NIOSH 1552	2-30	0.01-0.02	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Chlordane	ASTM D4947	250	1-5	polyurethane foam (PUF)	ORBO-1000	20557	226-92	GC-ECD
Chlordane	NIOSH 5510	10-200	0.5-1	XAD-2 MCE filter stainless steel screen	ORBO-706 ----- -----	20057 23381 custom	226-107 225-5	GC-ECD
Chlordane	OSHA 67	480	1	XAD-2/glass fiber filter (OVS)	ORBO-49P	20350	226-30-16	GC-ECD
Chlorides	EPA 0050			filter + absorbing solutions				IC
Chlorides	EPA 0051	120	2	filter + absorbing solutions				IC
Chlorinated Camphene (Toxaphene)	NIOSH 5039	2-30	0.2-1	MCE filter	-----	23381	225-5	GC-ECD
Chlorinated Diphenyl Ether	NIOSH 5025	8-200	0.5-1.5	MCE filter	-----	23381	225-5	GC-ECD
Chlorinated Terphenyl (60% chlorine)	NIOSH 5014	100-1500	1-3	glass fiber filter	-----	23378	225-7	GC-ECD
Chlorine	NIOSH 6011	2-90	0.3-1	PTFE filter + silver membrane	885	custom	225-9006	IC-CD
Chlorine	OSHA ID 101	15-240	1	fritted bubbler	-----	64835-U	225-36-2	ISE
Chlorine Dioxide	OSHA ID 202	7.50-120	0.5	fritted bubbler	-----	64835-U	225-36-2	IC
Chloroacetaldehyde	NIOSH 2015	3-16	0.05-0.2	silica gel	ORBO-507	20870-U	226-15	GC-ECD
Chloroacetaldehyde	OSHA 76	7.5	0.5	silica gel	ORBO-507	20870-U	226-15	GC-ECD
Chloroacetic Acid	NIOSH 2008	1-100	0.05-0.2	silica gel	ORBO-505	custom	226-47-01	IC
α-Chloroacetophenone	NIOSH 5/291	12	0.01-0.2	Tenax	ORBO-405	custom	226-35	GC-FID
Chlorobenzene	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Chlorobenzene (Monochlorobenzene)	NIOSH 1003	1.50-40	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
o-Chlorobenzylidene Malonitrile	NIOSH 5/304	90	1.5	Tenax PTFE stainless steel screen	ORBO-408 ----- -----	custom 23383 custom	226-35-07 225-1710	HPLC-UV
Chlorobromomethane	NIOSH 1003	0.5-8	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Chlorodifluoromethane	NIOSH 1018	1-4	0.01-0.05	(2) coconut charcoal	ORBO-32S + ORBO-32L	20267-U 20228	226-01 226-09	GC-FID
Chlorodiphenyl (42% & 54% Cl)	NIOSH 5503	1-50	0.05-0.2	glass fiber filter florisil	----- ORBO-60	23376 20351	225-16 226-39	GC-ECD
Chloroethane (Ethyl Chloride)	NIOSH 2519	0.3-3	0.02-0.05	(2) coconut charcoal	ORBO-32L	20228	226-09	GC-FID
2-Chloroethanol	NIOSH 2513	2-35	0.01-0.2	petroleum charcoal	ORBO-303	20040-U	226-38	GC-FID
Chloroform (Trichloromethane)	NIOSH 1003	1-50	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Chloroform (Trichloromethane)	OSHA 5	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Chloromethyl Methyl Ether	NIOSH 1/220		1.5	fritted bubbler	-----	64835-U	225-36-2	GC-ECD
Chloromethyl Methyl Ether	OSHA 10	50	0.5	fritted bubbler	-----	64835-U	225-36-2	GC-ECD
4-Chloronitrobenzene	NIOSH 2005	1-150	0.01-1	silica gel	ORBO-52S	20229	226-10	GC-FID
1-Chloro-1-nitropropane	NIOSH 5/S211	12	0.04-0.2	Chromosorb 108	ORBO-710	20058	226-113	GC-FID
p-Chlorophenol	NIOSH 2014	1.5-40	0.05-0.2	silica gel	ORBO-52S	20229	226-10	HPLC-UV
Chloroprene	NIOSH 1002	1.5-8	0.01-0.1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Chloroprene	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
beta-Chloroprene	OSHA 112	6	0.05	Chromosorb 106	ORBO-711	20059	226-111	GC-ECD
Chlorpyrifos (Dursban)	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Chlorpyrifos (Dursban)	OSHA 62	480	1	XAD-2/glass fiber filter (OVS)	ORBO-49P	20350	226-30-16	GC-FPD
Chromic Acid	NIOSH 7600	8-400	1-4	PVC filter	-----	23387	225-8-01	VIS
Chromic Acid	NIOSH 7604	100-1000	1-4	PVC filter	-----	23387	225-8-01	IC-CD
Chromic Acid, Chromates (as CrO3)	OSHA ID 103	30-960	2	PVC filter	-----	23387	225-8-01	DPP
Chromium	NIOSH 7300	5-1000	1-4	MCE filter	-----	23381	225-5	ICP-AES
Chromium, Cr Compounds (as Cr)	NIOSH 7024	10-1000	1-3	MCE filter	-----	23381	225-5	AAS

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Chromium (metal & insoluble compounds)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Chromium (metal & insoluble compounds)	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Chromium Acetate (listed as II & III compounds)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Chromium Carbonate (listed as II & III compounds)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Chromium, Hexavalent	EPA 0061			KOH absorbing solution				IC/PCR
Chromium, Hexavalent	NIOSH 7600	8-400	1-4	PVC filter	-----	23387	225-8-01	VIS
Chromium, Hexavalent	NIOSH 7604	100-1000	1-4	PVC filter	-----	23387	225-8-01	IC-CD
Chromium, Hexavalent	NIOSH 9101			chromate test kit				
Chromium, Hexavalent	OSHA ID 215	960	2	PVC filter	-----	23387	225-8-01	IC-UV, PCD
Chromium Phosphate (listed as II & III compounds)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Chromium, Soluble Salts (IC/OUS) (listed as II & III compounds)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Chrysene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----- 23390-U	20258 226-30-04 225-17-07	226-30-04 226-30-04	HPLC-UV/FI
Chrysene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----- 23390-U	20258 226-30-04 225-17-07	226-30-04 225-17-07	GC-FID
Chrysene	OSHA 58	960	2	glass fiber filter	-----	23378	225-7	HPLC-UV
Chrysotile Fibers	NIOSH 7400	400- varies	0.5-16	MCE filter	-----	23380-U	225-19	PCM
Coal Tar Naphtha	NIOSH 1550	1.3-20	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Coal Tar Pitch Volatiles	OSHA 58	960	2	glass fiber filter	-----	23378	225-7	HPLC-UV
Cobalt	NIOSH 7300	25-2000	1-4	MCE filter	-----	23381	225-5	ICP-AES
Cobalt	OSHA ID 213	30-480	2	MCE filter	-----	23381	225-5	ICP-AES
Cobalt, Co Compounds (as Co)	NIOSH 7027	30-1500	1-3	MCE filter	-----	23381	225-5	AAS
Cobalt, Metal, Dust & Fume	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Cobalt, Metal, Dust & Fume	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Coke Oven Emissions	OSHA 58	960	2	glass fiber filter	-----	23378	225-7	HPLC-UV
Copper	NIOSH 7300	5-1000	1-4	MCE filter	-----	23381	225-5	ICP-AES
Copper, Dust & Fume	NIOSH 7029	50-1500	1-3	MCE filter	-----	23381	225-5	AAS
Copper, Dusts & Mists	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Copper, Dusts & Mists	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Copper Fume	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Copper Fume	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Crag Herbicide (total dust)	NIOSH 5/S356	90	1-1.5	MCE filter	-----	23381	225-5	VIS
Cresol	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Cresol	OSHA 32	24	0.1	XAD-7 (Supelpak 70)	ORBO-47	20349	226-95	HPLC-UV
Cresol, all isomers	NIOSH 2546	1-24	0.01-0.1	XAD-7 (Supelpak 70)	ORBO-47	20349	226-95	GC-FID
Cresols (Methylphenols)	EPA TO-8	<80	0.1-1	NaOH absorbing solution				HPLC UV/ELCD/FI
Cristobalite	OSHA ID 142	816	1.7	PVC filter 10mm nylon cyclone	----- ----- NA	23387 NA	225-8-01	XRD
Crocidolite Fibers	NIOSH 7400	400-	0.5-16	MCE filter	-----	23380-U	225-19	PCM
Crotonaldehyde	NIOSH 2539	5	0.01-0.05	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-FID
Crotonaldehyde	NIOSH 3516	1-49	0.1-0.2	fritted bubbler	-----	64835-U	225-36-2	DPP
Crotonaldehyde	OSHA 81	6	0.1	glass fiber filter c/w DNPH	827	20069	225-9003	HPLC-UV
Cryofluorane	NIOSH 1018	1-4	0.01-0.05	(2) coconut charcoal	ORBO-32S + ORBO-32L	20267-U 20228	226-01 226-09	GC-FID
Cryolite	NIOSH 7902	12-800	1-2	MCE filter + support pad c/w Na <sub>2</sub> CO <sub>3</sub>	845	custom	225-9005 225-9001	ISE
Cumene	NIOSH 1501	1-30	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Cumene	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Cyanide (as CN)	OSHA ID 120	90-120	1	MCE filter impinger	-----	23381 20270-U	225-5 225-36-1	ISE
Cyanides	NIOSH 6010	2-90	0.05-0.2	soda lime	ORBO-1006	custom	226-28	VIS
Cyanides, Aerosol & Gas	NIOSH 7904	10-180	0.5-1	MCE filter fritted bubbler	-----	23381 64835-U	225-5 225-36-2	ISE
Cyanuric Acid	NIOSH 5030	10-1000	1-3	PVC filter	-----	23387	225-8-01	HPLC-UV
Cyclohexane	NIOSH 1500	2.50-5	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Cyclohexane	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Cyclohexanol	NIOSH 1402	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Cyclohexanol	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Cyclohexanone	NIOSH 1300	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Cyclohexanone	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Cyclohexanone	OSHA 1	10	0.05-0.2	Chromosorb 106	ORBO-704	custom	226-110	GC-FID
Cyclohexene	NIOSH 1500	5-7	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Cyclohexene	OSHA 7	5	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,3-Cyclopentadiene	NIOSH 2523	1-5	0.01-0.05	Chromosorb 104 c/w maleic anhydride	-----	NA	NA	GC-FID
2,4-D	NIOSH 5001	15-200	1-3	glass fiber filter	-----	23378	225-7	HPLC-UV
DDT	NIOSH 3/S274	90	1.5	glass fiber filter	-----	23378	225-7	GC-ELCD
n-Decane	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Demeton	NIOSH 5514	30-500	0.2-1	XAD-2 MCE filter stainless steel screen	ORBO-608 ----- -----	20050-U 23381 custom	226-30-05 225-5	GC-FPD
Desflurane	OSHA 106	3	0.05	Carbosieve S-III (Anasorb 747)	ORBO-91	20360	226-81A	GC-FID
Di-(2-ethylhexyl) phthalate	NIOSH 5020	10-200	1-3	MCE filter	-----	23381	225-5	GC-FID
Di-(2-ethylhexyl) phthalate	OSHA 104	240	1	Tenax/glass fiber filter (OVS)	ORBO-410P	custom	226-56	GC-FID
Diacetone Alcohol	NIOSH 1402	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Diacetone Alcohol	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
o-Dianisidine	NIOSH 5013	150-500	1-3	PTFE filter	-----	23389	225-17	HPLC-UV
o-Dianisidine	OSHA 71	100	1	glass fiber filter c/w H2SO4	822	20065	225-9004	GC-ECD
Diatomaceous Earth	NIOSH 7501	50-400	1-3	PVC filter	-----	23387	225-8-01	XRD
Diazinon	NIOSH 5600	12-240	0.2-1.0	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Diazinon	OSHA 62	480	1	XAD-2/glass fiber filter (OVS)	ORBO-49P	20350	226-30-16	GC-FPD
Diazomethane	NIOSH 2515	6-30	0.2	XAD-2 c/w octanoic acid	ORBO-651	custom	226-23	GC-FID
Dibenz(a,h)anthracene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Dibenz(a,h)anthracene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	GC-FID
Diborane	NIOSH 6006	60-260	0.5-1	charcoal c/w oxidizer PTFE filter	ORBO-352 -----	20043 special order	226-151	PES
Dibromodifluoromethane	NIOSH 1012	2.5-10	0.01-0.2	(2) coconut charcoal	ORBO-32S or ORBO-301	20267-U or 20039	226-01 NA	GC-FID
1,2-Dibromoethane (Ethylene Dibromide)	NIOSH 1008	0.1-25	0.02-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-ECD
1,2-Dibromoethane	OSHA 2	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-ECD
2-Dibutylaminoethanol	NIOSH 2007	4-24	0.01-0.2	silica gel	ORBO-506	20032	226-10-04	GC-FID
Dibutyl Phosphate	NIOSH 5017	50-250	1-3	PTFE filter	-----	23383	225-1710	GC-FPD
Dibutyl Phthalate	NIOSH 5020	6-200	1-3	MCE filter	-----	23381	225-5	GC-FID
Dibutyl Phthalate	OSHA 104	240	1	Tenax/glass fiber filter (OVS)	ORBO-410P	custom	226-56	GC-FID
Dibutyltin bis (Isooctyl Mercaptoacetate)	NIOSH 5504	50-500	1-1.5	XAD-2 glass fiber filter	ORBO-614 -----	custom 23378	226-30 225-7	HPLC-AAGF
o-Dichlorobenzene	NIOSH 1003	1-60	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
o-Dichlorobenzene	OSHA 7	3	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
o,p-Dichlorobenzene	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
p-Dichlorobenzene	NIOSH 1003	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
p-Dichlorobenzene	OSHA 7	3	0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
3,3'-Dichlorobenzidine	NIOSH 5509	20-100	0.2	glass fiber filter	-----	23376	225-16	HPLC-UV
3,3'-Dichlorobenzidine	OSHA 65	100	1	glass fiber filter c/w H2SO4	822	20065	225-9004	GC-ECD
Dichlorodifluoromethane (F-12)	NIOSH 1018	1-4	0.01-0.05	(2) coconut charcoal	ORBO-32S + ORBO-32L	20267-U 20228	226-01 226-09	GC-FID
1,1-Dichloroethane	NIOSH 1003	0.5-15	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,1-Dichloroethane	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,2-Dichloroethane – see Ethylene dichloride								
1,1-Dichloroethene – see Vinylidene Chloride								
1,2-Dichloroethylene	NIOSH 1003	0.2-5	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,2-Dichloroethylene	OSHA 7	3	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Dichloroethyl Ether	NIOSH 1004	2-15	0.01-1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Dichloroethyl Ether	OSHA 7	15	1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,1-Dichloro-1-fluoroethane (Freon 141b)	OSHA 113	1	0.05	(2) Carbosieve S-III (Anasorb CMS)	ORBO-91	20360	226-121	GC-FID
Dichlorofluoromethane	NIOSH 2516	0.25-3	0.01-0.05	(2) coconut charcoal	ORBO-32L	20228	226-09	GC-FID
Dichloromethane (Methylene Chloride)	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Dichloromethane	OSHA 80	3	0.05	Carbosieve S-III	ORBO-91	20360	226-121	GC-FID
Dichloromethane	NIOSH 1005	0.5-2.5	0.01-0.2	(2) coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,1-Dichloro-1-nitroethane	NIOSH 1601	1.5-15	0.01-1	petroleum charcoal	ORBO-303	20040-U	226-38	GC-FID
1,1-Dichloro-1-nitroethane	OSHA 7	15	1	petroleum charcoal	ORBO-303	20040-U	226-38	GC-FID
1,2-Dichloropropane	NIOSH 1013	0.10-3.50	0.01-0.2	petroleum charcoal	ORBO-303	20040-U	226-38	GC-ECN
1,2-Dichlorotetrafluoroethane	NIOSH 1018	1-4	0.01-0.05	(2) coconut charcoal	ORBO-32S + ORBO-32L	20267-U 20228	226-01 226-09	GC-FID
Dichlorvos	NIOSH 5/295	120	0.5-1	XAD-2	ORBO-603	custom	226-30-03	GC-FPD
Dichlorvos	OSHA 62	480	1	XAD-2/glass fiber filter (OVS)	ORBO-49P	20350	226-30-16	GC-FPD
Dicrotophos	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Dieldrin	NIOSH 3/S283	180	1.5	glass fiber filter quartz fiber filter (37mm)	-----	23378	225-7	GC-ELCD
Diesel Exhaust	NIOSH 5040	106-4300	1-4	impinger	-----	special order	225-1809	EGA-TOA
Diethanolamine	NIOSH 3509	5-300	0.5-1	silica gel	ORBO-52S	20229	226-10	GC-FID
Diethylamine	NIOSH 2010	3-30	0.01-1	XAD-7 c/w 10% NBDC	ORBO-653	custom	226-96	HPLC-VIS
Diethylamine	OSHA 41	10	0.2	silica gel	ORBO-506	20032	226-10-04	GC-FID
2-Diethylaminoethanol	NIOSH 2007	4-24	0.01-0.2	Porapak P	ORBO-1102	20062	226-114	GC-FID
Diethyl Carbamoyl Chloride	NIOSH 6/317	12	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Diethyl Ether	NIOSH 1610	0.25-3	0.01-0.2	Tenax/glass fiber filter (OVS)	ORBO-410P	custom	226-56	GC-FID
Diethyl Phthalate	OSHA 104	240	1					
Diethylene Dioxide - see Dioxane								
Diethylene Glycol	NIOSH 5523	5-60	0.5-2	XAD-7/glass fiber filter (OVS)	ORBO-47P	custom	226-57	GC-FID
Diethylene Glycol Ether	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Diethylenetriamine (DETA)	NIOSH 2540	1-20	0.01-0.1	XAD-2 c/w 10% NITC	ORBO-654	custom	226-30-18	HPLC-UV
Diethylenetriamine (DETA)	OSHA 60	10	0.1	XAD-2 c/w 10% NITC	ORBO-654	custom	226-30-18	HPLC-UV
Difluorodibromomethane	NIOSH 1012	2.5-10	0.01-0.2	(2) coconut charcoal	ORBO-32S or ORBO-301	20267-U or 20039	226-01 NA	GC-FID
Difluorodibromomethane	OSHA 7	10	0.2	(2) coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Difluorodichloromethane	NIOSH 1018	1-4	0.01-0.05	(2) coconut charcoal	ORBO-32S + ORBO-32L	20267-U 20228	226-01 226-09	GC-FID
Diglycidyl Ether of Bisphenyl A (DGEBA)	NIOSH 6/333	288	1.6	glass fiber filter	-----	23378	225-7	HPLC-UV
Dihydrocapsaicin	NIOSH 5041	7-1000	1-3	glass fiber filter	-----	23376	225-16	HPLC-FL

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Diisobutyl Ketone	NIOSH 1300	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Diisobutyl Ketone	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Dimethoxymethane (Methylal)	NIOSH 1611	1-3	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Dimethylacetamide	NIOSH 2004	15-80	0.01-1	silica gel	ORBO-52S	20229	226-10	GC-FID
Dimethylamine	NIOSH 2010	3-30	0.01-1	silica gel	ORBO-52S	20229	226-10	GC-FID
Dimethylamine	OSHA 34	10	0.2	XAD-7 c/w 10% NBDC	ORBO-653	custom	226-96	HPLC-VIS
N,N-Dimethylaniline	NIOSH 2002	3-30	0.02-1	silica gel	ORBO-52S	20229	226-10	GC-FID
N,N-Dimethylaniline	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Dimethylarsinic Acid	NIOSH 5022	50-1000	1-3	PTFE filter	-----	23383	225-1710	IC-AAS
Dimethylformamide	NIOSH 2004	15-80	0.01-1	silica gel	ORBO-52S	20229	226-10	GC-FID
N,N-Dimethylformamide	OSHA 66	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-NPD
1,1-Dimethylhydrazine	NIOSH 3515	2-100	0.2-1	fritted bubbler	-----	64835-U	225-36-2	VIS
Dimethyl Phthalate	OSHA 104	240	1	Tenax/glass fiber filter (OVS)	ORBO-410P	custom	226-56	GC-FID
Dimethyl Sulfate	NIOSH 2524	0.25-12	0.01-0.2	Porapak P	ORBO-1102	20062	226-114	GC-ECD
Dimethyl Sulfate	NIOSH 5/301	12	0.05	Porapak P	ORBO-1102	20062	226-114	GC-ECN
N,N-Dimethyl-p-toluidine	NIOSH 2002	20-100	0.02-1	silica gel	ORBO-52S	20229	226-10	GC-FID
Dinitrobenzene (all isomers)	NIOSH 4/S214	90	1.5	MCE filter fritted bubbler	----- -----	23381 64835-U	225-5 225-36-2	HPLC-UV
Dinitro-o-cresol	NIOSH 5/S166	180	1.5	MCE filter stainless steel screen fritted bubbler	----- -----	23381 custom 64835-U	225-5 NA 225-36-2	HPLC-UV
2,4-Dinitrotoluene	OSHA 44	60	1	Tenax/glass fiber filter	ORBO-79	custom	NA	GC-TEA
Di-n-octyl Phthalate	OSHA 104	240	1	Tenax/glass fiber filter (OVS)	ORBO-410P	custom	226-56	GC-FID
Dioxane (Diethylene Dioxide)	NIOSH 1602	0.5-15	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Dioxane (Diethylene Dioxide)	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Dioxins	EPA TO-9A	325,000	225	quartz fiber filter polyurethane foam (PUF)	----- ORBO-2000	custom 20037	226-131	GC-MS
Dioxins	EPA 0023A			glass fiber filter XAD-2	----- Supelpak 2	custom 20279	NA	GC-MS
Diphenyl (Biphenyl)	NIOSH 2530	15-30	0.01-0.5	Tenax	ORBO-404	custom	226-35-01	GC-FID
Diphenylamine	OSHA 78	100	1	(2) glass fiber filter c/w H2SO4	822	20065	225-9004	HPLC-UV
Dipropylene Glycol Methyl Ether	NIOSH 2/S69	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Dipropylene Glycol Methyl Ether	OSHA 101	10	0.1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Dipropylene Glycol Methyl Ether	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Disul	NIOSH 5/S356	90	1-1.5	MCE filter	-----	23381	225-5	VIS
Disulfoton	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
2,6-Di-t-butyl-p-cresol (DBPC)	NIOSH 1/226	10	0.2	silica gel	ORBO-504	20031	NA	GC-FID
Divinylbenzene	OSHA 89	12	0.05	charcoal c/w TBC	ORBO-356	20047	NA	GC-FID
Dursban	OSHA 62	480	1	XAD-2/glass fiber filter (OVS)	ORBO-49P	20350	226-30-16	GC-FPD
Dust, Respirable Nuisance	NIOSH 0600	75-1000	1.7	PVC filter	-----	23387	225-8-01	gravimetric
Dust, Total Nuisance	NIOSH 0500	7-133	1-2	PVC filter	-----	23387	225-8-01	gravimetric
Dyes, Benzidine	NIOSH 5013	150-500	1-3	PTFE filter	-----	23389	225-17	HPLC-UV
Dyes, o-Dianisidine	NIOSH 5013	150-500	1-3	PTFE filter	-----	23389	225-17	HPLC-UV
Dyes, o-Toluidine	NIOSH 5013	150-500	1-3	PTFE filter	-----	23389	225-17	HPLC-UV
EDA – see Ethylenediamine								
Elements	NIOSH 7300	varies	1-4	MCE filter	-----	23381	225-5	ICP-AES
Endrin	NIOSH 5519	12-400	0.5-1	Chromosorb 102 MCE filter	ORBO-706 -----	20057 23381	226-107 225-5	GC-ECD

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Enflurane (Ethrane)	OSHA 103	12	0.05	Carbosieve S-III (Anasorb 747)	ORBO-91	20360	226-81A	GC-FID
Epichlorohydrin	NIOSH 1010	2-30	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Epichlorohydrin	OSHA 7	20	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
EPN – see o-Ethyl-o-(p-nitro-phenyl) phenylphosphorothioate								
Ethanol	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Ethanol (Ethyl Alcohol)	NIOSH 1400	0.1-1	0.01-0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethanol (Ethyl Alcohol)	OSHA 100	12	0.05	Carbosieve S-III (Anasorb 747)	ORBO-91	20360	226-82	GC-FID
Ethanol (Ethyl Alcohol)	OSHA 7	1	0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethanolamine - see 2-Aminoethanol								
3-Ethenylpyridine	ASTM D5075	90	<1.5	XAD-4	ORBO-613	20052	226-93	GC-NPD
Ethion	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Ethoprop	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
2-Ethoxyethanol (Cellosolve)	NIOSH 1403	1-6	0.01-0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2-Ethoxyethanol (Cellosolve)	OSHA 79	15-48	0.1-1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2-Ethoxyethyl Acetate (Cellosolve Acetate)	NIOSH 1450	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2-Ethoxyethyl Acetate	OSHA 79	15-48	0.1-1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2-Ethoxyethyl Acetate	OSHA 7	10	0.1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethrane (Enflurane)	OSHA 103	12	0.05	Carbosieve S-III (Anasorb 747)	ORBO-91	20360	226-81A	GC-FID
Ethyl Acetate	NIOSH 1457	0.1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethyl Acetate	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Ethyl Acetate	OSHA 7	5	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethyl Acrylate	NIOSH 1450	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethyl Acrylate	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethyl Acrylate	OSHA 92	12	0.05	charcoal c/w TBC	ORBO-356	20047	NA	GC-FID
Ethylamine	NIOSH 3/S144	30	0.2	silica gel	ORBO-52S	20229	226-10	GC-FID
Ethylamine	OSHA 36	10	0.2	XAD-7 c/w 10% NBDC	ORBO-653	custom	226-96	HPLC-VIS
Ethyl Amyl Ketone (5-Methyl-3-heptanone)	NIOSH 1301	1-25	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethylbenzene	NIOSH 1501	1-24	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethylbenzene	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethyl Bromide	NIOSH 1011	0.5-4	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethyl Bromide	OSHA 7	5	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethyl Butyl Ketone (3-Heptanone)	NIOSH 1301	1-25	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethyl Butyl Ketone (3-Heptanone)	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethyl Chloride	NIOSH 2519	0.3-3	0.02-0.05	(2) coconut charcoal	ORBO-32L	20228	226-09	GC-FID
Ethyl-2-cyanoacrylate	OSHA 55	12	0.1	XAD-7 t/w H3PO4	ORBO-655	20054	226-98	HPLC-UV
Ethyl Ether	NIOSH 1610	0.25-3	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethyl Ether	OSHA 7	3	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethyl Formate	NIOSH 1452	0.3-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethyl Formate	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethyl Mercaptan	NIOSH 2542	10-150	0.1-0.2	glass fiber filter c/w mercuric acetate	826	20068	225-9007	GC-FPD
Ethyl Silicate	NIOSH 3/S264	9	0.05	XAD-2	ORBO-605	20049	NA	GC-FID
Ethylene Chlorohydrin	NIOSH 2513	2-35	0.01-0.2	petroleum charcoal	ORBO-303	20040-U	226-38	GC-FID
Ethylene Chlorohydrin	OSHA 7	20	0.2	petroleum charcoal	ORBO-303	20040-U	226-38	GC-FID
Ethylenediamine (EDA)	NIOSH 2540	1-20	0.01-0.1	XAD-2 c/w 10% NITC	ORBO-654	custom	226-30-18	HPLC-UV
Ethylenediamine (EDA)	OSHA 60	10	0.1	XAD-2 c/w 10% NITC	ORBO-654	custom	226-30-18	HPLC-UV

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equip	Analytical Technique
Ethylene Dibromide – see 1,2-Dibromoethane								
Ethylene Dichloride (1,2-Dichlorethane)	NIOSH 1003	1-50	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethylene Dichloride (1,2-Dichlorethane)	OSHA 3	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-ECD
Ethylene Glycol	NIOSH 5523	5-60	0.5-2	XAD-7/glass fiber filter (OVS)	ORBO-47P	custom	226-57	GC-FID
Ethylene Glycol Dinitrate	NIOSH 2507	3-100	0.2-1	Tenax	ORBO-402	20832-U	226-35-03	GC-ECD
Ethylene Glycol Dinitrate	OSHA 43	15	1	Tenax	ORBO-402	20832-U	226-35-03	HPLC-TEALUV
Ethylene Oxide	ASTM D4413	3-20	0.01-0.2	coconut charcoal	ORBO-32J	custom	226-16	GC-FID
Ethylene Oxide	ASTM D5578	1-10	0.02-0.1	Carboxen-564 c/w HBr	ORBO-78	20355	NA	GC-ECD
Ethylene Oxide	NIOSH 1614	1-24	0.05-0.15	petroleum charcoal c/w HBr	ORBO-353	20044	226-38-03	GC-ECD
Ethylene Oxide	NIOSH 3702	<80% vol	>0.02	5L sampling bag	----	24655	232-05	GC-PID
Ethylene Oxide	OSHA 49		0.0493	3M #3551 diffusive sampler	----	20253-U		GC-ECD
Ethylene Oxide	OSHA 50	24	0.1	petroleum charcoal c/w HBr	ORBO-353	20044	226-38-03	GC-ECD
Ethylene Oxide (Qazi-Ketcham)	AIHAJ 11/77	4.8-9.6	0.2	petroleum charcoal	ORBO-33	20259	226-36	GC-FID
Ethylene Thiourea	NIOSH 5011	200-800	1-3	MCE filter	----	23381	225-5	VIS
Ethylene Thiourea	OSHA 95	480	2	(2) glass fiber filter	----	23376	225-16	HPLC-UV
Ethylenimine	NIOSH 3514	1-48	0.2	fritted bubbler	----	64835-U	225-36-2	HPLC-UV
N-Ethylmorpholine	NIOSH 3/S146	10	0.2	silica gel	ORBO-52S	20229	226-10	GC-FID
o-Ethyl-o-(p-nitrophenyl)-phenylphosphorothioate (EPN)	NIOSH 5012	15-700	1-2	glass fiber filter	----	23378	225-7	GC-FPD
Ethyl sec-Amyl Ketone (5-Methyl-3-heptanone)	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ethylvinylbenzene	OSHA 89	12	0.05	charcoal c/w TBC	ORBO-356	20047	NA	GC-FID
Fenamiphos	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Ferrovandium Dust	NIOSH 1/173	180	1.5	MCE filter	----	23381	225-5	AAS
Ferrovandium Dust	OSHA ID 121	480-960	2	MCE filter	----	23381	225-5	AAS
Ferrovandium Dust	OSHA ID 125G	480	2	MCE filter	----	23381	225-5	ICP-AES
Fluoranthene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Fluoranthene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----	20258 23390-U	226-30-04 225-17-07	GC-FID
Fluorene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Fluorene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----	20258 23390-U	226-30-04 225-17-07	GC-FID
Fluoride (F- & HF)	OSHA ID 110	22.5-90	1.5	MCE filter + support pad c/w Na2CO3	845	custom	225-9005 225-9001	ISE
Fluoride Particles	OSHA ID 110	22.5-90	1.5	MCE filter	----	23381	225-5	ISE
Fluorides	ASTM D4765	varies	2	MCE filter + support pad c/w Na2CO3	845	custom	225-9005 225-9001	ISE
Fluorides, Aerosol & Gas	NIOSH 7906	1-120	1-2	MCE filter + support pad c/w Na2CO3	845	custom	225-9005 225-9001	IC-CD
Fluorides, Aerosol & Gas	NIOSH 7902	12-800	1-2	MCE filter + support pad c/w Na2CO3	845	custom	225-9005 225-9001	ISE
Fluorotrichloromethane	NIOSH 1006	0.3-7	0.01-0.05	coconut charcoal	ORBO-32L	20228	226-09	GC-FID
Fonofos	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Formaldehyde	ASTM D5197	see method	0.5-1.5	silica gel c/w DNPH	LpDNPH S10	21026-U	226-119	HPLC-UV
Formaldehyde	EPA 0100		1	silica gel c/w DNPH	LpDNPH S10	21026-U	226-119	HPLC-UV
Formaldehyde	NIOSH 2016	1-15	0.1-1.5	silica gel c/w DNPH	LpDNPH S10	21026-U	226-119	HPLC-UV
Formaldehyde	NIOSH 2539	5	0.01-0.05	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-FID
Formaldehyde	NIOSH 2541	1-36	0.01-0.1	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-FID
Formaldehyde	NIOSH 3500	1-100	0.2-1	PTFE filter (2) impingers	----	23383 20270-U	225-1710 225-36-1	VIS
Formaldehyde	OSHA 52	3-24	0.1-0.2	XAD-2 c/w 2-HMP	ORBO-24	20231	226-117	GC-NPD

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equip	Analytical Technique
Formaldehyde	OSHA ID 102	30-160	0.5-1	MCE filter fritted bubbler	----- -----	23381 64835-U	225-5 225-36-2	DPP
Formaldehyde	OSHA ID 205		0.0614	3M #3721 diffusive sampler	-----	NA		colorimetric
Formaldehyde (on dust)	NIOSH 5700	240-1050	2	PVC filter inhalable sampler	----- -----	special order		HPLC-UV
Formic Acid	NIOSH 2011	1-24	0.05-0.2	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Formic Acid	US Army	10-25		Chromosorb P c/w Na <sub>2</sub> CO <sub>3</sub>	ORBO-70	20256-U	NA	GC-FID
Furfural	NIOSH 2529	1-12	0.01-0.05	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-FID
Furfural	NIOSH 2539	5	0.05	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-FID
Furfural	OSHA 72	180	1	petroleum charcoal	ORBO-303	20040-U	226-38	GC-FID
Furfuryl Alcohol	NIOSH 2505	3-25	0.01-0.05	Porapak Q	ORBO-1103	20063	226-115	GC-FID
Gallium	NIOSH 1/177	100	1.5-2	MCE filter	-----	23381	225-5	ASV
Glass, Fibrous	NIOSH 7400	400-	0.5-16	MCE filter	-----	23380-U	225-19	PCM
Glutaraldehyde	NIOSH 2532	1-30	0.05-0.5	silica gel c/w DNPH	LpDNPH H10	505315	226-119	HPLC-UV
Glutaraldehyde	OSHA 64	15	1	glass fiber filter c/w DNPH	827	20069	225-9003	HPLC-UV
Glycerin Mist	NIOSH 0500	7-133	1-2	PVC filter	-----	23387	225-8-01	gravimetric
Glycidol (2,3-Epoxy-1-propanol)	NIOSH 1608	5-100	0.01-1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Glycidol (2,3-Epoxy-1-propanol)	OSHA 7	50	1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Glycols	NIOSH 5523	5-60	0.5-2	XAD-7/glass fiber filter (OVS)	ORBO-47P	custom	226-57	GC-FID
Hafnium	NIOSH 5/S194	90	1.5-2	MCE filter	-----	23381	225-5	PES
Halothane	OSHA 103	12	0.05	Carbosieve S-III (Anasorb 747)	ORBO-91	20360	226-81A	GC-FID
Heptachlor	ASTM D4947	250	1-5	polyurethane foam (PUF)	ORBO-1000	20557	226-92	GC-ECD
Heptachlor	NIOSH 5/S287	60	0.01-1	Chromosorb 102	ORBO-706	20057	226-107	GC-ECD
Heptanal	NIOSH 2539	5	0.01-0.05	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-FID
n-Heptane	NIOSH 1500	4	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Heptane	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
n-Heptane	OSHA 7	5	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
3-Heptanone	NIOSH 1301	1-25	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
3-Heptanone	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Herbicides, Chlorinated and Organonitrogen (Hand Wash)	NIOSH 9200			polyethylene bag				GC-ECD
Herbicides, Chlorinated and Organonitrogen (Patch)	NIOSH 9201			PUF pads				GC-ECD
Herbicides, Chlorinated, Organonitrogen, and Acid	NIOSH 5602	12-480	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-ECD
Hexachlorobutadiene	NIOSH 2543	1-100	0.05-0.2	XAD-2	ORBO-608	20050-U	226-30-05	GC-ECD
Hexachloro-1,3- cyclopentadiene	NIOSH 2518	0.25-90	0.01-0.2	(2) Porapak T	ORBO-1105	custom	226-116	GC-ECD
Hexachloroethane	NIOSH 1003	3-70	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Hexachloroethane	OSHA 7	30	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Hexachloronaphthalene	NIOSH 2/S100	30	1	MCE filter	-----	23381	225-5	GC-ECD
1,6-Hexamethylene Diisocyanate (HDI)	OSHA 42	15	1	glass fiber filter c/w 1-2PP	ORBO-80	20811	225-9002	HPLC-UV
1,6-Hexamethylene- diisocyanate (HDI)	NIOSH 5521	5-500	1	impinger	-----	20270-U	225-36-1	HPLC-UV ELCD
Hexamethylene Diisocyanate	NIOSH 5522	15-360	1-2	impinger	-----	20270-U	225-36-1	HPLC-FL
Hexamethylenetetramine (HMTA)	NIOSH 4/263	15	1.5	MCE filter fritted bubbler	----- -----	23381 64835-U	225-5 225-36-2	VIS
Hexanal	NIOSH 2539	5	0.01-0.05	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-FID
Hexanal	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
n-Hexane	NIOSH 1500	4	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Hexane	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
n-Hexane	OSHA 7	5	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2-Hexanone (MBK)	NIOSH 1300	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2-Hexanone (MBK)	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Hexavalent Chromium - see Chromium								

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equip	Analytical Technique
Hexone (MIBK)	NIOSH 1300	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Hexone (MIBK)	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Hexone (MIBK)	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
sec-Hexyl Acetate	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Hydrazine	NIOSH 3503	7-100	0.2-1	fritted bubbler	-----	64835-U	225-36-2	VIS
Hydrazine	OSHA 108	240	1	(2) glass fiber filters c/w H2SO4	840	custom	NA	HPLC-UV
Hydrazoic Acid	OSHA ID 211	>5	1.0	silica gel c/w NaOH PVC filter	ORBO-560 -----	custom 23387	226-55 225-8-01	IC-UV
Hydrazoic Acid	US Army	10-25		Chromosorb P c/w Na2CO3	ORBO-70	20256-U	NA	VIS
Hydrobromic Acid	NIOSH 7903	3-100	0.2-0.5	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Hydrocarbons, General	Supelco			Carbotrap	ORBO-100	20255-U	NA	GC-FID
Hydrocyanic Acid	NIOSH 6010	2-90	0.05-0.2	soda lime	ORBO-1006	custom	226-28	VIS
Hydrocyanic Acid	NIOSH 7904	7-100	0.5-1	MCE filter fritted bubbler	----- -----	23381 64835-U	225-5 225-36-2	ISE
Hydrocyanic Acid	OSHA ID 120	90-120	1	MCE filter impinger	----- -----	23381 20270-U	225-5 225-36-1	ISE
Hydrofluoric Acid	NIOSH 7902	12-800	1-2	MCE filter + support pad c/w Na2CO3	845	custom	225-9005 225-9001	ISE
Hydrofluoric Acid	NIOSH 7903	3-100	0.2-0.5	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Hydrogen Bromide	NIOSH 7903	3-100	0.2-0.5	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Hydrogen Bromide	OSHA ID 165SG	96	0.20	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Hydrogen Chloride	NIOSH 7903	3-100	0.2-0.5	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Hydrogen Chloride	OSHA ID 174	7.5	0.5	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Hydrogen Cyanide	NIOSH 6010	2-90	0.05-0.2	soda lime	ORBO-1006	custom	226-28	VIS
Hydrogen Cyanide	NIOSH 7904	7-100	0.5-1	MCE filter fritted bubbler	----- -----	23381 64835-U	225-5 225-36-2	ISE
Hydrogen Cyanide	OSHA ID 120	90-120	1	MCE filter impinger	----- -----	23381 20270-U	225-5 225-36-1	ISE
Hydrogen Fluoride	NIOSH 7906	1-800	1-2	MCE filter + support pad c/w Na2CO3	845	custom	225-9005 225-9001	IC
Hydrogen Fluoride – see Hydrofluoric Acid								
Hydrogen Sulfide	NIOSH 6013	1.2-40	0.1-1.5	coconut charcoal	ORBO-34	20211	NA	IC-CD
Hydrogen Sulfide	OSHA ID 141	2-6	0.1-0.2	MCE filter c/w AgNO3	828	20070-U	225-9008	DPP
Hydroquinone	NIOSH 5004	30-180	1-4	MCE filter	-----	23381	225-5	HPLC-UV
Indeno(1,2,3-cd)pyrene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Indeno(1,2,3-cd)pyrene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	GC-FID
Indium	NIOSH 1/173	180	1.5	MCE filter	-----	23381	225-5	AAS
Indium	NIOSH 1/190	100	1.5-2	MCE filter	-----	23381	225-5	ASV
Iodine	NIOSH 6005	15-225	0.5-1	charcoal c/w KOH	ORBO-354	20045	226-27	IC
Iodine	OSHA ID 212	2.5-7.5	0.5	beaded carbon c/w KOH	ORBO-	custom	226-80	IC-PED
Iron	NIOSH 7300	5-100	1-4	MCE filter	-----	23381	225-5	ICP-AES
Iron	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Iron Oxide Fume	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Iron Oxide Fume	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Isoamyl Acetate	NIOSH 1450	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isoamyl Acetate	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isoamyl Alcohol	NIOSH 1402	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isobutyl Acetate	NIOSH 1450	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isobutyl Acetate	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isobutyl Alcohol	NIOSH 1401	2-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isobutyl Alcohol	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isobutyraldehyde	NIOSH 2539	5	0.01-0.05	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-FID
Isocyanates	NIOSH 5521	5-500	1	impinger	-----	20270-U	225-36-1	HPLC-UV ELCD
Isocyanates	NIOSH 5522	15-360	1-2	impinger	-----	20270-U	225-36-1	HPLC-FL
Isocyanates – see HDI, MDI, TDI								
Isoflurane	OSHA 103	12	0.05	Carbosieve S-III (Anasorb 747)	ORBO-91	20360	226-81A	GC-FID

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Isophorone	NIOSH 2508	2-25	0.01-1	petroleum charcoal	ORBO-303	20040-U	226-38	GC-FID
Isophorone	OSHA 7	10	0.2	petroleum charcoal	ORBO-303	20040-U	226-38	GC-FID
Isopropyl Acetate	NIOSH 1454	0.1-9	0.02-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isopropyl Acetate	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isopropyl Alcohol	NIOSH 1400	0.2-3	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isopropyl Alcohol	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Isopropyl Alcohol	OSHA 7	3	0.1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isopropylamine	NIOSH 3/S147	100	1	fritted bubbler	-----	64835-U	225-36-2	GC-FID
N-Isopropylaniline	OSHA 78	100	1	(2) glass fiber filter c/w H2SO4	822	20065	225-9004	HPLC-UV
Isopropyl Ether	NIOSH 1618	0.1-3	0.01-0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isopropyl Ether	OSHA 7	3	0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isopropyl Glycidyl Ether (IGE)	NIOSH 1620	1-30	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isopropyl Glycidyl Ether (IGE)	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Isovaleraldehyde	NIOSH 2539	5	0.01-0.05	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-FID
Kepone	NIOSH 5508	50-600	0.5-1	MCE filter impinger	----- -----	23381 20270-U	225-5 225-36-1	GC-ECD
Kerosene	NIOSH 1550	1.3-20	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Ketene	NIOSH 2/S92	50	1	fritted bubbler	-----	64835-U	225-36-2	colorimetric
Ketones	EPA 0011			DNPH absorbing solution				HPLC-UV
Ketones – see Aldehydes								
Lead	NIOSH 7082	200-1500	1-4	MCE filter	-----	23381	225-5	AAS
Lead	NIOSH 7105	1-1500	1-4	MCE filter	-----	23381	225-5	AAGF
Lead	NIOSH 7300	1250-	1-4	MCE filter	-----	23381	225-5	ICP-AES
Lead	NIOSH 7700	10-240	2	MCE filter	-----	23381	225-5	Spot
Lead	NIOSH 7701	20-1500	1-4	MCE filter	-----	23381	225-5	ASV
Lead	NIOSH 9100			wipes	-----			AAS, ICP
Lead by Field Portable XRF	NIOSH 7702	570-1900	1-4	MCE filter	-----	23381	225-5	XRF
Lead, Inorganic Fumes & Dusts (as Pb)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Lead, Inorganic Fumes & Dusts (as Pb)	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Lead Sulfide	NIOSH 7505	600-1000	1.7 or 2.2	PVC filter	-----	23387	225-8-01	XPD
Limonene	NIOSH 1552	2-30	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Limonene	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Lindane (γ-BHC)	NIOSH 5502	18-240	0.2-1	glass fiber filter impinger	----- -----	23378 64835-U	225-7 225-36-2	GC-ECD
Lithium	NIOSH 7300	100-2000	1-4	MCE filter	-----	23381	225-5	ICP-AES
Lithium Hydroxide	NIOSH 7401	70-1000	1-4	PTFE filter	-----	23383	225-1710	titration
Magnesium	NIOSH 7300	5-67	1-4	MCE filter	-----	23381	225-5	ICP-AES
Magnesium	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Magnesium Oxide Fume	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Malathion	NIOSH 5012	15-130	1-2	glass fiber filter	-----	23378	225-7	GC-FPD
Malathion	NIOSH 5600	12-60	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Malathion	OSHA 62	60	1	XAD-2/glass fiber filter (OVS)	ORBO-49P	20350	226-30-16	GC-FPD
Maleic Anhydride	NIOSH 3512	40-500	0.2-1.5	fritted bubbler	-----	64835-U	225-36-2	HPLC-UV
Maleic Anhydride	OSHA 86	60	0.5	(2) glass fiber filter c/w DMBA	834	20072	NA	HPLC-UV
Maneb	OSHA 107	500	2	MCE filter	-----	23380-U	225-19	HPLC-UV
Manganese	NIOSH 7300	5-200	1-4	MCE filter	-----	23381	225-5	ICP-AES
Manganese, Mn Compounds (as Mn)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Manganese, Mn Compounds (as Mn)	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Manganese Oxide	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Manganese Oxide	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
MBK – see 2-Hexanone								

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
MDI – see Methylene Bisphenyl Isocyanate and Isocyanates								
Mercaptans	NIOSH 2542	10-150	0.1-0.2	glass fiber filter c/w mercuric acetate	826	20068	225-9007	GC-FPD
Mercury	NIOSH 6009	2-100	0.15-0.25	Hopcalite	ORBO-1002	20863	226-17-1A	AAS
Mercury, Organic	NIOSH 6/S342	3-12	0.01-0.2	Carbosieve B	ORBO-910	custom	NA	AAS
Mercury, Particulate	OSHA ID 145	10	2	MCE filter	-----	23381	225-5	AACV
Mercury, Vapor	OSHA ID 140	3-100	0.2	Hopcalite	ORBO-1002	20863	226-17-1A	AACV
Mesityl Oxide	NIOSH 1301	1-25	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Mesityl Oxide	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Metal Particulates	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Metal Particulates	OSHA ID 125G	30-480	2	MCE filter	-----	23381	225-5	ICP-AES
Metal/Metalloid Particulates	OSHA ID 206	480	2	MCE filter	-----	23381	225-5	ICP-AES
Metals	ASTM D4185	varies	2	MCE filter	-----	23381	225-5	AAS
Metals	EPA 0060			filter + absorbing solutions				ICP-AES
Metals	NIOSH 7300	varies-2000	1-4	MCE filter	-----	23381	225-5	ICP-AES
Methacrylic Acid	OSHA 312	24	0.1	(2) XAD-8	ORBO-601	20048	226-30-08	HPLC-UV
Methamidophos	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Methanol	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Methanol (Methyl Alcohol)	NIOSH 2000	1-5	0.02-0.2	silica gel	ORBO-502	20030-U	226-51	GC-FID
Methanol (Methyl Alcohol)	OSHA 91	3-5	0.05	Carbosieve S-III (Anasorb 747)	ORBO-91	20360	226-82	GC-FID
Methoxychlor	NIOSH 4/S371	100	1.5	glass fiber filter	-----	23378	225-7	GC-ELCD
2-Methoxyethanol (Methyl Cellosolve)	NIOSH 1403	6-50	0.01-0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2-Methoxyethanol	OSHA 79	15-48	0.1-1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2-Methoxyethyl Acetate (Methyl Cellosolve Acetate)	NIOSH 1451	0.2-20	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2-Methoxyethyl Acetate	OSHA 79	15-48	0.1-1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1-Methoxy-2-propanol (1M2P)	OSHA 99	10	0.1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2-Methoxy-1-propanol (2M1P)	OSHA 99	10	0.1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1-Methoxy-2-propyl Acetate (1M2PA)	OSHA 99	10	0.1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2-Methoxy-1-propyl Acetate (2M1PA)	OSHA 99	10	0.1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methyl Acetate	NIOSH 1458	0.2-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methyl Acetate	OSHA 7	5	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methyl Acetylene	NIOSH 5/S84	3-4	0.01-0.05	5L sampling bag	-----	24655	232-05	GC-FID
Methyl Acetylene Propadiene Mix	OSHA 7	2	0.05	(2) coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methyl Acetylene Propadiene Mix (MAPP)	NIOSH 6/S85	3-4	0.01-0.05	5L sampling bag	-----	24655	232-05	GC-FPD
Methyl Acrylate	NIOSH 1459	1-5	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methyl Acrylate	OSHA 7	5	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methyl Acrylate	OSHA 92	12	0.05	charcoal c/w TBC	ORBO-356	20047	NA	GC-FID
Methyl Amine	NIOSH 4/277		0.2	silica gel	ORBO-508	custom	226-14	GC-FID
Methyl Amine	NIOSH 6/S148	24	0.02-0.2	silica gel	ORBO-508	custom	226-14	IC-ELCD
Methyl Amine	OSHA 40	10	0.2	XAD-7 c/w 10% NBDC	ORBO-653	custom	226-96	HPLC-VIS
4-Methylamino-azobenzene	NIOSH 4/284	72	0.2	Gas Chrom P glass fiber filter	ORBO-1001	custom	226-52	GC-FID
Methyl(n-amy)ketone	NIOSH 1301	1-25	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Methyl Bromide	NIOSH 2520	1-5	0.01-0.1	Na2SO4 drying tube (2) petroleum charcoal	ORBO-971 ORBO-310	custom custom	226-44-02 226-38-02	GC-AED
Methyl-n-butyl Ketone – see 2-Hexanone								
Methyl Chloride	NIOSH 1001	0.40-3	0.01-0.1	(2) coconut charcoal	ORBO-32S ORBO-32L	20267-U 20228	226-01 226-09	GC-FID
Methyl Chloroform	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Methyl Chloroform – see 1,1,1-Trichloroethane								
Methyl-2-cyanoacrylate	OSHA 55	12	0.1	XAD-7 t/w H3PO4	ORBO-655	20054	226-98	HPLC-UV
Methyl Cyclohexane	NIOSH 1500	4	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methyl Cyclohexane	OSHA 7	5	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methyl Ethyl Ketone	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Methyl Ethyl Ketone (MEK) – see 2-Butanone								
Methyl Ethyl Ketone Peroxide	NIOSH 3508	52-520	0.5-2	impinger	-----	20270-U	225-36-1	VIS
Methyl Ethyl Ketone Peroxide	OSHA 77	15	1	XAD-4	ORBO-613	20052	226-93	HPLC-UV PCD
Methyl Formate	NIOSH 5/S291	3	0.01-0.05	Carbosieve G	ORBO-901	custom	NA	GC-FID
5-Methyl-3-heptanone	NIOSH 1301	1-25	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
5-Methyl-3-heptanone	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methyl Iodide	NIOSH 1014	15-50	0.01-1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methyl Isoamyl Acetate	NIOSH 1450	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methyl Isobutyl Carbinol	NIOSH 1402	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methyl Isobutyl Carbinol	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methyl Isobutyl Ketone (MIBK) – see Hexone								
Methyl Isocyanate	OSHA 54	15	0.05	XAD-7 c/w H2SO4 & 1-2PP	ORBO-657	20055	NA	HPLC-UV
Methyl Mercaptan	NIOSH 2542	10-150	0.1-0.2	glass fiber filter c/w mercuric acetate	826	20068	225-9007	GC-FPD
Methyl Mercaptan	OSHA 26	20	0.2	glass fiber filter c/w mercuric acetate	826	20068	225-9007	GC-FPD
Methyl Methacrylate	NIOSH 2537	1-8	0.01-0.05	XAD-2	ORBO-609	20051	226-30-06	GC-FID
Methyl Methacrylate	OSHA 94	3	0.05	charcoal c/w TBC	ORBO-356	20047	NA	GC-FID
Methyl Parathion	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Methyl Phenol	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
N-Methyl-2-pyrrolidinone	NIOSH 1302	0.5-125	0.05-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-NPD/FID
α-Methyl Styrene	NIOSH 1501	1-30	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
α-Methyl Styrene	OSHA 7	3	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methyl tert-Butyl Ether (MTBE)	NIOSH 1615	2-96	0.1-0.2	(2) coconut charcoal	ORBO-32L	20228	226-09	GC-FID
N-Methyl-N-2,4,6- tetranitrobenzamine	NIOSH 3/S225	1	1.5	MCE filter	-----	23381	225-5	VIS
Methylal	NIOSH 1611	1-3	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methylal	OSHA 7	2	0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methylarsonic Acid	NIOSH 5022	50-1000	1-3	PTFE filter	-----	23383	225-1710	IC-AAS
Methylcyclohexanol	NIOSH 1404	1-15	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Methylcyclohexanone	NIOSH 2521	1-6	0.01-0.05	Porapak Q	ORBO-1103	20063	226-115	GC-FID
4,4'-Methylene Bis(2- chloroaniline) (MOCA)	NIOSH 1/236	3-480	0.2-1	silica gel glass fiber filter	ORBO-501 -----	custom 23376	226-63 225-16	HPLC-UV
4,4'-Methylene Bis(2- chloroaniline)	OSHA 71	100	1	glass fiber filter c/w H2SO4	822	20065	225-9004	GC-ECD
Methylene Bisphenyl Isocyanate (MDI)	NIOSH 5521	5-500	1	impinger	-----	20270-U	225-36-1	HPLC-UV ELCD
Methylene Bisphenyl Isocyanate (MDI)	OSHA 47	15	1	glass fiber filter c/w 1-2PP	ORBO-80	20811	225-9002	HPLC-UV
Methylene Chloride	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Methylene Chloride	OSHA 59	10	0.05	coconut charcoal (3-section, 350mg)	ORBO-32TB	custom	226-09-02	GC-FID
Methylene Chloride – see Dichloromethane								
4,4'-Methylene Dianiline (MDA)	NIOSH 5029	10-1000	1-2	filter gasket glass fiber filter c/w H2SO4	----- 822	NA 20065	225-9004	HPLC-UV ELCD
4,4'-Methylene Dianiline (MDA)	OSHA 57	100	1	glass fiber filter c/w H2SO4	822	20065	225-9004	GC-ECD
Mevinphos (Phosdrin)	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Mineral Spirits	NIOSH 1550	1.3-20	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Mineral Spirits	OSHA 48	3	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
MOCA - see 4,4'-Methylene Bis(2-chloroaniline)								
Molybdenum	NIOSH 7300	5-67	1-4	MCE filter	-----	23381	225-5	ICP-AES
Molybdenum Insolubles (as Mo)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Molybdenum Insolubles (as Mo)	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Molybdenum Solubles (as Mo)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Monochloroacetic Acid	NIOSH 2008	1-100	0.05-0.2	silica gel	ORBO-505	custom	226-47-01	IC
Monochlorobenzene	NIOSH 1003	1.5-40	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Monochlorobenzene	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Monocrotophos	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Monomethylaniline	NIOSH 3511	11-100	0.2-1	fritted bubbler	-----	64835-U	225-36-2	GC-FID
Monomethylhydrazine	NIOSH 3510	3-20	0.5-1.5	fritted bubbler	-----	64835-U	225-36-2	VIS
Morpholine	NIOSH 3/S150	20	0.2	silica gel	ORBO-52S	20229	226-10	GC-FID
Mycobacterium Tuberculosis Airborne	NIOSH 0900		>4	PTFE filter	-----	23391	225-1705	Polymerase Chain
Naphtha	OSHA 48	3	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Naphtha (Coal Tar)	NIOSH 1550	1.3-20	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Naphthalene	NIOSH 1501	100-200	0.01-1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Naphthalene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Naphthalene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	GC-FID
Naphthalene	OSHA 35	10	0.2	Chromosorb 106	ORBO-704	custom	226-110	GC-FID
α-Naphthalthiourea	NIOSH 5/S276	480	1.5-2	PTFE filter	-----	23383	225-1710	HPLC-UV
α-Naphthylamine (1-Naphthylamine)	NIOSH 5518	30-100	0.2-0.8	silica gel glass fiber filter	ORBO-502 -----	20030-U 23376	226-51 225-16	GC-FID
α-Naphthylamine (1-Naphthylamine)	OSHA 93	100	1	glass fiber filter c/w H2SO4	822	20065	225-9004	GC-ECD
β-Naphthylamine (2-Naphthylamine)	NIOSH 5518	30-100	0.2-0.8	silica gel glass fiber filter	ORBO-502 -----	20030-U 23376	226-51 225-16	GC-FID
β-Naphthylamine (2-Naphthylamine)	OSHA 93	100	1	glass fiber filter c/w H2SO4	822	20065	225-9004	GC-ECD
Nickel	NIOSH 7300	25-1000	1-4	MCE filter	-----	23381	225-5	ICP-AES
Nickel Carbonyl	NIOSH 6007	7-80	0.05-0.2	low Ni charcoal MCE filter	ORBO-304 -----	20041 23381	NA 225-5	AAGF
Nickel, Insoluble Compounds (as Ni)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Nickel, Insoluble Compounds (as Ni)	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Nickel, Metal & Soluble Compounds (as Ni)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Nickel, Metal & Soluble Compounds (as Ni)	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equip	Analytical Technique
Nicotine	ASTM D5075	90	<1.5	XAD-4	ORBO-613	20052	226-93	GC-NPD
Nicotine	EPA IP-2A		1	XAD-4	ORBO-613	20052	226-93	GC-NPD
Nicotine	EPA IP-2B		1.7-3	2 Teflon coated glass fiber filters, 1 t/w sodium	----	custom		GC-NPD
Nicotine	NIOSH 2544	60-400	1	XAD-2	ORBO-43	20258	226-30-04	GC-NPD
Nicotine	NIOSH 2551	0.5-600	0.1-1	XAD-4	ORBO-613	20052	226-93	GC-NPD
Nicotine	NIOSH 3/S293	100	1	XAD-2	ORBO-605	20049	NA	GC-FID
Nitric Acid	NIOSH 7903	3-100	0.2-0.5	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Nitric Acid	OSHA ID 127	120	1	fritted bubbler	----	64835-U	225-36-2	IC
Nitric Acid	OSHA ID 165SG	96	0.20	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Nitric Oxide	NIOSH 6014	1.5-6	0.025	molecular sieve c/w TEA + oxidizer	ORBO-904	custom	226-40	VIS
Nitric Oxide	OSHA ID 190	6	0.025	(2) molecular sieve c/w TEA + (1) oxidizer tube	ORBO-905 ORBO-905	custom custom	226-40-02 NA	IC-CD
p-Nitroaniline	NIOSH 5033	16-350	1-3	MCE filter	----	23381	225-5	HPLC-UV
Nitrobenzene	NIOSH 2005	10-150	0.01-1	silica gel	ORBO-52S	20229	226-10	GC-FID
Nitrobenzene	NIOSH 2017	5-50	0.2	silica gel + glass fiber filter c/w H2SO4	ORBO-507 822	20870-U 20065	226-15 225-9004	GC-FID
4-Nitrobiphenyl	NIOSH 4/273	48	0.2	silica gel glass fiber filter	ORBO-502	20030-U 23376	226-51 225-16	GC-FID
p-Nitrochlorobenzene	NIOSH 2005	1-150	0.01-1	silica gel	ORBO-52S	20229	226-10	GC-FID
Nitroethane	NIOSH 2526	1.5-3	0.01-0.05	XAD-2	ORBO-610	custom	226-30-02	GC-FID
Nitrogen Dioxide	EPA IP-5A			continuous luminol monitor				
Nitrogen Dioxide	EPA IP-5B			Palmer diffusion tube	----	NA		VIS
Nitrogen Dioxide	EPA IP-5C			diffusive sampler	----	NA		IC
Nitrogen Dioxide	NIOSH 6014	1.5-6	0.025-0.2	molecular sieve c/w TEA + oxidizer	ORBO-904	custom	226-40	VIS
Nitrogen Dioxide	NIOSH 6700	0.25-8hr		Palmer diffusion tube	----	NA		VIS
Nitrogen Dioxide	OSHA ID 182	3	0.2	(2) molecular sieve c/w TEA + (1) oxidizer tube	ORBO-906 ORBO-905	custom custom	226-40-02 NA	IC
Nitrogen Dioxide	OSHA ID 190	6	0.025	(2) molecular sieve c/w TEA + (1) oxidizer tube	ORBO-906 ORBO-905	custom custom	226-40-02 NA	IC-CD
Nitroglycerin	NIOSH 2507	3-100	0.2-1	Tenax	ORBO-402	20832-U	226-35-03	GC-ECD
Nitroglycerin	OSHA 43	15	1	Tenax	ORBO-402	20832-U	226-35-03	HPLC-
Nitromethane	NIOSH 2527	1.2-3	0.01-0.05	Chromosorb 106	ORBO-711	20059	226-111	GC-FPD
1-Nitropropane	OSHA 46	4	0.1	XAD-4	ORBO-613	20052	226-93	GC-FID
2-Nitropropane	NIOSH 2528	0.1-2	0.01-0.05	Chromosorb 106	ORBO-704	custom	226-110	GC-FID
2-Nitropropane	OSHA 46	4	0.1	XAD-4	ORBO-613	20052	226-93	GC-FID
Nitrosamines	NIOSH 2522	15-1000	0.2-2	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosodiamylamine (NDAMA)	OSHA 38	75	1	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosodibutylamine	NIOSH 2522	15-1000	0.2-2	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosodi-n-butylamine	OSHA 27	75	1	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosodiethanolamine	OSHA 31	480	2	glass fiber filter	----	23378	225-7	GC-TEA
N-Nitrosodiethylamine	NIOSH 2522	15-1000	0.2-2	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosodiethylamine	OSHA 27	75	1	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosodi-isopropylamine (NDIPA)	OSHA 38	75	1	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosodimethylamine	EPA TO-7	<300	0.1-2	Thermosorb/N	----	NA	NA	GC-MS
N-Nitrosodimethylamine	NIOSH 2522	15-1000	0.2-2	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosodimethylamine	OSHA 27	75	1	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosodiphenylamine	OSHA 23	250	1	fritted bubbler	----	64835-U	225-36-2	HPLC-UV
N-Nitrosodi-n-propylamine	OSHA 27	75	1	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosodipropylamine	NIOSH 2522	15-1000	0.2-2	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosoethyl-n-butylamine (NEBA)	OSHA 38	75	1	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosomethyl-n-butylamine (NMBA)	OSHA 38	75	1	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosomethyl-ethylamine (NMEA)	OSHA 38	75	1	Thermosorb/N	----	NA	NA	GC-TEA

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equip	Analytical Technique
N-Nitrosomorpholine	NIOSH 2522	15-1000	0.2-2	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosomorpholine	OSHA 27	75	1	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosopiperidine	NIOSH 2522	15-1000	0.2-2	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosopiperidine	OSHA 27	75	1	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitroso-n-propyl-n-butylamine (NPBA)	OSHA 38	75	1	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosopyrrolidine	NIOSH 2522	15-1000	0.2-2	Thermosorb/N	----	NA	NA	GC-TEA
N-Nitrosopyrrolidine	OSHA 27	75	1	Thermosorb/N	----	NA	NA	GC-TEA
Nitrotoluene	NIOSH 2005	1-30	0.01-0.02	silica gel	ORBO-52S	20229	226-10	GC-FID
Nitrous Oxide	NIOSH 6600	80% vol		5L sampling bag	----	24655	232-05	IR
Nitrous Oxide	OSHA ID 166			Landauer diffusive sampler				TD-IR
Non-Methane Organic Compounds	EPA TO-12			glass bead trap canister		custom NA	NA	FID
Octachloronaphthalene	NIOSH 2/S97	30	1.30	MCE filter	----	23381	225-5	GC-ECD
Octamethylcyclotetrasiloxane	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Octane	OSHA 7	5	0.1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Octane	NIOSH 1500	4	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Octane	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
1-Octanethiol	NIOSH 2510	1-15	0.01-0.2	Tenax	ORBO-403	20034	NA	GC-FPD
Oil Mist, Mineral	NIOSH 5026	20-500	1-3	MCE filter or	----	23381	225-5	IR
				glass fiber filter or	----	23378	225-7	
				PVC filter or	----	23387	225-8-01	
				PTFE filter	----	23390-U	225-17-07	
Oil of Vitriol (Sulfuric Acid)	NIOSH 7903	3-100	0.2-0.5	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Organic Vapors	ASTM D3686	see method	0.001-0.2	coconut charcoal	ORBO-32S ORBO-32L	20267-U 20228	226-01 226-09	GC-FID
Organics	EPA 0040	20	0.5	Tedlar bag	----	24656	231-15	GC-MS
Organotin Compounds (as Sn)	NIOSH 5504	50-500	1-1.5	XAD-2 glass fiber filter	ORBO-614	custom 23378	226-30 225-7	HPLC-AAGF
Oxygen	NIOSH 6601	1		portable O2 monitor				sensor
Ozone	NIOSH 1/153	variable	1-2	impinger	----	20270-U	225-36-1	colorimetric
Ozone	NIOSH 1/154	45	1	impinger	----	20270-U	225-36-1	colorimetric
Ozone	NIOSH 2/S8		1	impinger	----	20270-U	225-36-1	colorimetric
Ozone	OSHA ID 214	22.5-90	0.25-1.5	(2) glass fiber filter c/w NaNO <sub>2</sub> , K <sub>2</sub> CO <sub>3</sub> , glycerol	850	custom	NA	IC-UV
PAHs – see Polynuclear Aromatic Hydrocarbons								
Palladium	NIOSH 1/173	180	1.5	MCE filter	----	23381	225-5	AAS
Paraquat	NIOSH 5003	40-1000	1-4	PTFE filter	----	23383	225-1710	HPLC-UV
Parathion	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Parathion	OSHA 62	480	1	XAD-2/glass fiber filter (OVS)	ORBO-49P	20350	226-30-16	GC-FPD
Particulate Matter, Respirable	EPA IP-10A			size specific impaction				gravimetric
Particulate Matter, Respirable	EPA IP-10B			continuous particulate monitor				gravimetric
Particulates	NIOSH 0600	20-400	1.7 or 2.2	PVC filter cyclone	----	23387 NA	225-8-01	gravimetric
PCBs – Aroclor 1242 (42% Cl)	NIOSH 5503	1-50	0.05-0.2	glass fiber filter Florisil	ORBO-60	23376 20351	225-16 226-39	GC-ECD
PCBs – Aroclor 1254 (54% Cl)	NIOSH 5503	1-50	0.05-0.2	glass fiber filter Florisil	ORBO-60	23376 20351	225-16 226-39	GC-ECD
Pentachlorobenzene	NIOSH 5517	3-12	0.01-0.2	XAD-2 PTFE filter	ORBO-44	20260-U special order	226-30-04	GC-ECD
Pentachloroethane	NIOSH 2517	1-10	0.01-0.2	Porapak R	ORBO-1104	20064	226-59-04	GC-ECD
Pentachloronaphthalene	NIOSH 2/S96	250	1.3	glass fiber filter	----	23378	225-7	GC-ECD
				fritted bubbler	----	64835-U	225-36-2	
Pentachlorophenol	NIOSH 5512	48-480	0.5-1	MCE filter	----	23381	225-5	HPLC-UV
				stainless steel screen fritted bubbler	----	custom 64835-U	225-36-2	

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Pentachlorophenol	OSHA 39	48	0.2	XAD-7 XAD-7/glass fiber filter (set of 3)	ORBO-602	custom	226-97	HPLC-UV
Pentamidine Isethionate	NIOSH 5032	50-1500	1-2	PVC filter	-----	23387	225-8-01	HPLC-FL
Pentane	OSHA 7	2	0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Pentane	NIOSH 1500	2	0.01-0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Pentane	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
2-Pentanone	NIOSH 1300	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2-Pentanone	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Perchloroethylene	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Perchloroethylene in exhaled breath and air	NIOSH 3704		0.02-5	1L sampling bag	-----	24633	232-01	portable GC-PID
Perchloroethylene – see also Tetrachloroethylene								
Pesticides	ASTM D4861	240-7200	1-5	polyurethane foam (PUF)	ORBO-1000	20557	226-92	GC-ECD
Pesticides	US Air Force		0.6-1	XAD-2	ORBO-42S	20262	NA	GC-ECD or HPLC-UV
Pesticides, Organochlorine	EPA IP-8		1-5	polyurethane foam (PUF) glass fiber filter cartridge	ORBO-1000 -----	20557 21031	226-126	GC-ECD
Pesticides, Organonitrogen	NIOSH 5601	<480	0.1-1	XAD-2/glass fiber filter XAD-2/quartz filter (OVS)	ORBO-49P or ORBO-49Q	20350 custom	226-30-16 226-58	HPLC-UV
Pesticides, Organophosphorus	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Pesticides/PCBs	EPA TO-10A		1-5	polyurethane foam (PUF) quartz filter cartridge	ORBO-1000 -----	20557 21031	226-126	GC/ECD/NPD/FPD/MS
Pesticides/PCBs	EPA TO-4A	320,000	225	polyurethane foam (PUF)	ORBO-2000	20037	226-131	GC/ECD/NPD/FPD/MS
Petroleum Distillates (Naphtha)	NIOSH 1550	1.3-20	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Petroleum Distillates (Naphtha)	OSHA 48	3	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Petroleum Ether (Benzin)	NIOSH 1550	1.3-20	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Phenanthrene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Phenanthrene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	GC-FID
Phenanthrene	OSHA 58	960	2	glass fiber filter	-----	23378	225-7	HPLC-UV
Phenol	EPA TO-8	<80	0.1-1	NaOH absorbing solution				HPLC UV/ELCD/FI
Phenol	NIOSH 2546	1-24	0.01-0.1	XAD-7 (Supelpak 70)	ORBO-47	20349	226-95	GC-FID
Phenol	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Phenol	OSHA 32	24	0.1	XAD-7 (Supelpak 70)	ORBO-47	20349	226-95	HPLC-UV
Phenyl Ether	NIOSH 1617	1-50	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Phenyl Ether	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Phenyl Ether – Biphenyl Mixture	NIOSH 2013	1-40	0.01-0.2	silica gel	ORBO-52S	20229	226-10	GC-FID
Phenyl Glycidyl Ether	NIOSH 1619	80-150	0.01-1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Phenyl Glycidyl Ether	OSHA 7	50	0.1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Phenylhydrazine	NIOSH 3518	25-120	0.2-1	fritted bubbler	-----	64835-U	225-36-2	VIS
N-Phenyl-1-naphthylamine	OSHA 96	240	1	glass fiber filter c/w ascorbic acid	893	custom	NA	HPLC-FI
N-Phenyl-2-naphthylamine	OSHA 96	240	1	glass fiber filter c/w ascorbic acid	893	custom	NA	HPLC-FI
o-, m-, p-Phenylenediamine	OSHA 87	100	1	glass fiber filter c/w H2SO4	822	20065	225-9004	HPLC-UV
Phorate	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Phosdrin (Mevinphos)	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Phosgene	EPA TO-6	<50	<1	aniline in toluene absorbing solution				HPLC-UV
Phosgene	NIOSH 1/219	25-50	1	impinger	----	20270-U	225-36-1	colorimetric
Phosgene	OSHA 61	240	1	XAD-2 c/w 2-HMP	ORBO-24	20231	226-117	GC-NPD
Phosphine	NIOSH 6002	1-16	0.01-0.2	silica gel c/w mercuric cyanide	ORBO-553	NA	NA	VIS
Phosphine	OSHA ID 180	36	0.05-0.15	carbon c/w KOH	ORBO-1004	20035	226-31	IC
Phosphoric Acid	OSHA ID 111	960	2	MCE filter	----	23381	225-5	IC
Phosphoric acid	OSHA ID 165SG	96	0.20	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
m-Phosphoric Acid	NIOSH 7903	3-100	0.2-0.5	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
o-Phosphoric Acid	NIOSH 7903	3-100	0.2-0.5	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Phosphorus	NIOSH 7300	50-2000	1-4	MCE filter	----	23381	225-5	ICP-AES
Phosphorus	NIOSH 7905	5-100	0.01-0.2	Tenax	ORBO-402	20832-U	226-35-03	GC-FPD
Phosphorus Pentachloride	NIOSH 5/S257	48	0.2	PVC filter fritted bubbler	---- ----	custom 64835-U	225-36-2	colorimetric
Phosphorus Trichloride	NIOSH 6402	11-100	0.05-0.2	fritted bubbler	----	64835-U	225-36-2	VIS
Phthalic Anhydride	NIOSH 3/S179	100	1.5	MCE filter	----	23381	225-5	HPLC-UV
Phthalic Anhydride	OSHA 90	75	1	(2) glass fiber filter c/w DMBA	835	custom	NA	HPLC-UV
Picric Acid	NIOSH 4/S228	180	1.5	MCE filter	----	23381	225-5	HPLC-UV
Pinene	NIOSH 1552	2-30	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Pinene	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Platinum	NIOSH 7300	13-2000	1-4	MCE filter	----	23381	225-5	ICP-AES
Polychlorinated Dibenzofurans	EPA 23			XAD-2	Supelpak 2	20279	NA	GC-MS
Polychlorinated Dibenzop-dioxins – see Dioxins	EPA 23			XAD-2	Supelpak 2	20279	NA	GC-MS
Polychlorobenzenes	NIOSH 5517	3-12	0.01-0.2	XAD-2 PTFE filter	ORBO-44 ----	20260-U special order	226-30-04	GC-ECD
Polychlorobiphenyls (also see PCBs)	ASTM D4861	240-7200	1-5	polyurethane foam (PUF)	ORBO-1000	20557	226-92	GC-ECD
Polycyclic Aromatic Compounds, Total (PACs)	NIOSH 5800	5-1000	1-2	XAD-2 PTFE filter	ORBO-43 ----	20258 23390-U	226-30-04 225-17-07	Flow injection, FI
Polymethylsiloxane Mist	NIOSH 1/227	408	1.7	MCE filter	----	23381	225-5	AAS
Polynuclear Aromatic Hydrocarbons (PAHs)	EPA IP-7	30,000	20	polyurethane foam (PUF)	ORBO-2000	20037	226-131	GC-FID HPLC-UV
Polynuclear Aromatic Hydrocarbons	EPA TO-13A	<320,000	225	polyurethane foam (PUF) or XAD-2	ORBO-2000 Supelpak 2	20037 or 20279	226-131	GC-MS
Polynuclear Aromatic Hydrocarbons	ASTM D6209	<350,000	225	polyurethane foam (PUF) or XAD-2	ORBO-2000 Supelpak 2	20037 or 20279	226-131	GC-MS
Polynuclear Aromatic Cmpds by HPLC	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 or ORBO-42L ----	20258 20264-U 23390-U	226-30-04 NA 225-17-07	HPLC-UV/FI
Polynuclear Aromatic Hydrocarbons	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 ----	20258 23390-U	226-30-04 225-17-07	GC-FID
Polynuclear Aromatic Hydrocarbons, Selected	OSHA 58	960	2	glass fiber filter	----	23378	225-7	HPLC-UV
Portland Cement (Total Dust)	OSHA ID 207	240	1	PVC filter	----	23387	225-8-01	gravimetric, XRF
Potassium	NIOSH 1/173	180	1.5	MCE filter	----	23381	225-5	AAS
Potassium, K	OSHA ID 121	480-960	2	MCE filter	----	23381	225-5	AAS
Potassium Hydroxide	NIOSH 7401	70-1000	1-4	PTFE filter	----	23383	225-1710	titration
Propargyl Alcohol	OSHA 97	6	0.05	petroleum charcoal c/w HBr	ORBO-353	20044	226-38-03	GC-ECD
Propionaldehyde	NIOSH 2539	5	0.01-0.05	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-FID
n-Propyl Acetate	NIOSH 1450	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Propyl Acetate	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Propyl Alcohol	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Propyl Alcohol	NIOSH 1401	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Propyl Nitrate	NIOSH 3/S227	70	1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
n-Propyl Nitrate	OSHA 7	50	0.1	petroleum charcoal	ORBO-303	20040-U	226-38	GC-FID

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Propylene Dichloride	NIOSH 1013	0.1-3.5	0.01-0.2	petroleum charcoal	ORBO-303	20040-U	226-38	GC-ECN
Propylene Dichloride	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Propylene Glycol	NIOSH 5523	5-60	0.5-2	XAD-7/glass fiber filter (OVS)	ORBO-47P	custom	226-57	GC-FID
Propylene Glycol Methomethyl Ethers/Acetates - see Methoxypropanols, Methoxypropyl Acetates								
Propylene Oxide	NIOSH 1612	0.5-5	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Propylene Oxide	OSHA 88	5	0.1	Carbosieve S-III (Anasorb 747)	ORBO-91	20360	226-81A	GC-FID
Prussic Acid	NIOSH 7904	7-100	0.5-1	MCE filter fritted bubbler	----- -----	23381 64835-U	225-5 225-36-2	ISE
Prussic Acid	OSHA ID 120	90-1280	1	MCE filter impinger	----- -----	23381 20270-U	225-5 225-36-1	ISE
Prussic Acid (Hydrogen Cyanide)	NIOSH 6010	2-90	0.05-0.2	soda lime	ORBO-1006	custom	226-28	VIS
Pyrene	NIOSH 5506	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	HPLC-UV/FI
Pyrene	NIOSH 5515	200-1000	2	XAD-2 PTFE filter	ORBO-43 -----	20258 23390-U	226-30-04 225-17-07	GC-FID
Pyrene	OSHA 58	960	2	glass fiber filter	-----	23378	225-7	HPLC-UV
Pyrethrum	NIOSH 5008	20-400	1-4	glass fiber filter	-----	23378	225-7	HPLC-UV
Pyrethrum	OSHA 70	60	1	XAD-2/glass fiber filter (OVS)	ORBO-49P	20350	226-30-16	GC-ECD
Pyridine	NIOSH 1613	18-150	0.01-1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Pyridine	OSHA 7	50	1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Qualitative X-ray Fluorescence Analysis of Workplace Substances	OSHA ID 204	960	2	MCE filter PVC filter	----- -----	23381 23387	225-5 225-8-01	XRF
Quartz	OSHA ID 142	816	1.7	PVC filter 10mm nylon cyclone	----- -----	23387 NA	225-8-01	XRD
Quartz, Respirable, in Coal Dust	NIOSH 7603	300-1000	1.7 or 2.2	PVC filter	-----	23387	225-8-01	IR
Quicklime (Calcium)	NIOSH 7020	20-400	1-3	MCE filter	-----	23381	225-5	AAS
Quinone	NIOSH 4/S181	24	0.01-0.2	XAD-2	ORBO-605	20049	NA	HPLC-UV
Reactive Acidic and Basic Gases and Particulate Matter	EPA IP-9			annular denuder system				IC
Resorcinol	NIOSH 5701	5-160	0.2-1	XAD-7/glass fiber filter (OVS)	ORBO-47P	custom	226-57	GC-FID
Rhodium, Metal, Fumes & Dust	NIOSH 3/S188	720	1.5	MCE filter	-----	23381	225-5	AAS
Rhodium, Soluble Salts	NIOSH 3/S189	370	1.5	MCE filter	-----	23381	225-5	AAGF
Ribavirin	NIOSH 5027	5-1000	1-4	glass fiber filter	-----	23378	225-7	HPLC-UV
Ronnel	NIOSH 5600	12-60	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Rotenone	NIOSH 5007	8-400	1-4	PTFE filter	-----	23383	225-1710	HPLC-UV
Rubidium	NIOSH 1/173	180	1.5	MCE filter	-----	23381	225-5	AAS
Selenium	NIOSH 7300	5-2000	1-4	MCE filter	-----	23381	225-5	ICP-AES
Selenium	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Semivolatiles	EPA 0010			glass fiber filter XAD-2	----- Supelpak 2	custom 20279	NA	GC-MS
Semivolatiles (SASS)	EPA 0020			glass fiber filter XAD-2	----- Supelpak 2	custom 20279	NA	GC-MS
Sevin	NIOSH 5006	20-400	1-3	glass fiber filter	-----	23378	225-7	VIS
Sevin	OSHA 63	60	1	XAD-2/glass fiber filter (OVS)	ORBO-49P	20350	226-30-16	HPLC-UV
Silica, Amorphous	NIOSH 7501	50-400	1.7 or 2.2	PVC filter	-----	23387	225-8-01	XPD
Silica, Crystalline	NIOSH 7601	400-800	1.7 or 2.2	MCE filter or PVC filter + 10mm nylon cyclone	----- ----- -----	23381 23387 NA	225-5 225-8-01	VIS

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equip	Analytical Technique
Silica, Crystalline	NIOSH 7602	400-800	1.7 or 2.2	MCE filter or PVC filter + 10mm nylon cyclone	----- ----- -----	23381 23387 NA	225-5 225-8-01	IR
Silica, Crystalline, Respirable	NIOSH 7500	400-1000	1.7 or 2.2	PVC filter	-----	23387	225-8-01	XPD
Silica (Quartz, Respirable) also see Quartz	OSHA ID 142	816	1.7	PVC filter	-----	23387	225-8-01	XRD
Silica, Quartz, in Coal Dust	NIOSH 7603	300-1000	1.7 or 2.2	PVC filter	-----	23387	225-8-01	IR
Silicon	NIOSH 1/173	180	1.5	MCE filter	-----	23381	225-5	AAS
Silver	NIOSH 7300	250-2000	1-4	MCE filter	-----	23381	225-5	ICP-AES
Silver (as Ag)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Sodium Azide	OSHA ID 211	>5	1.0	silica gel c/w NaOH PVC filter	ORBO-560 -----	custom 23387	226-55 225-8-01	IC-UV
Sodium Dichloroisocyanurate Dihydrate	NIOSH 5/314	400	1.7	PVC filter	-----	23387	225-8-01	titration
Sodium Fluoride	NIOSH 7902	12-800	1-2	MCE filter + support pad c/w Na <sub>2</sub> CO <sub>3</sub>	845	custom	225-9005 225-9001	ISE
Sodium Fluoroacetate	NIOSH 5/S301	480	1.5-2	Nucleopore cellulosic membrane	-----	special order		IC-ELCD
Sodium Hexafluoroaluminate	NIOSH 7902	12-800	1-2	MCE filter + support pad c/w Na <sub>2</sub> CO <sub>3</sub>	845	custom	225-9005 225-9001	ISE
Sodium Hexafluoroaluminate	NIOSH 7906	120-800	1-2	MCE filter + support pad c/w Na <sub>2</sub> CO <sub>3</sub>	845	custom	225-9005 225-9001	IC
Sodium Hydroxide	NIOSH 7401	70-1000	1-4	PTFE filter	-----	23383	225-1710	titration
Sodium Hydroxide	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Sodium Polyacrylate	NIOSH 5035	50-1500	1-3	PVC filter	-----	23387	225-8-01	ICP-AES or AAS
Solder Metals	OSHA ID 206	480	2	MCE filter	-----	23381	225-5	ICP-AES
Stibine	NIOSH 6008	4-50	0.01-0.2	silica gel c/w mercuric chloride	ORBO-551	custom	226-10-02	VIS
Stoddard Solvent	NIOSH 1550	1.3-20	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Stoddard Solvent	OSHA 48	3	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Strychnine	NIOSH 5016	70-1000	1-3	glass fiber filter	-----	23378	225-7	HPLC-UV
Styrene (Phenylethylene)	OSHA 9	10-15	0.2-1.0	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Styrene	NIOSH 1501	1-14	0.01-1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Styrene	OSHA 89	12	0.05	charcoal c/w TBC	ORBO-356	20047	NA	GC-FID
Styrene Oxide	NIOSH 5/303	13	0.01-0.2	Tenax	ORBO-406	custom	226-35-06	GC-FID
Sulfur Dioxide	NIOSH 6004	4-200	0.5-1.5	MCE filter + support pad c/w Na <sub>2</sub> CO <sub>3</sub>	845	custom	225-9005 225-9001	IC
Sulfur Dioxide	OSHA ID 104	15-60	1	MCE filter fritted bubbler	----- -----	23381 64835-U	225-5 225-36-2	IC
Sulfur Dioxide	OSHA ID 200	1.5-12	0.1	carbon c/w KOH	ORBO-1010	custom	226-80	IC
Sulfur Hexafluoride	NIOSH 5/S244	3-4	0.01-0.05	5L sampling bag	-----	24655	232-05	GC-TCD
Sulfur Hexafluoride	NIOSH 6602	80% vol	0.01-0.05	5L sampling bag	-----	24655	232-05	GC-ECD
Sulfuric Acid	NIOSH 7903	3-100	0.2-0.5	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Sulfuric Acid	OSHA ID 113	480	2	MCE filter	-----	23381	225-5	IC
Sulfuric Acid	OSHA ID 165SG	96	0.20	silica gel/glass fiber filter	ORBO-53	20265	226-10-03	IC
Sulfuric Acid Mist	ASTM D4856	40	1	MCE filter	-----	23381	225-5	IC
Sulfuryl Fluoride	NIOSH 6012	1.3-10	0.05-0.1	coconut charcoal	ORBO-32J	custom	226-16	IC
Sulprofos	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Super Absorbent Polymers	NIOSH 5035	50-1500	1-3	PVC filter	-----	23387	225-8-01	ICP-AES or AAS
2,4,5-T	NIOSH 5001	15-200	1-3	glass fiber filter	-----	23378	225-7	HPLC-UV
Talc, Respirable	NIOSH 355 2nd ed. update	200	1.7	PVC filter 10mm nylon cyclone	----- -----	23387 NA	225-8-01	XRD
Tellurium	NIOSH 7300	25-2000	1-4	MCE filter	-----	23381	225-5	ICP-AES
Tellurium	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Tellurium Hexafluoride	NIOSH 3/S187	400	1	coconut charcoal	ORBO-32S	20267-U	226-01	AAS
Temik	OSHA 74	480	1	XAD-2/glass fiber filter (OVS)	ORBO-49P	20350	226-30-16	GC-NPD

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Terbufos	NIOSH 5600	12-240	0.2-1	XAD-2/quartz filter (OVS)	ORBO-49Q	custom	226-58	GC-FPD
Terpenes	NIOSH 1552	2-30	0.01-2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
o-Terphenyl	NIOSH 5021	2-30	1-3	PTFE filter	-----	23390-U	225-17-07	GC-FID
1,1,2,2-Tetrabromoethane	NIOSH 2003	50-100	0.2-1	silica gel	ORBO-52S	20229	226-10	GC-FID
Tetrabutyltin	NIOSH 5504	50-500	1-1.5	XAD-2 glass fiber filter	ORBO-614 -----	custom 23378	226-30 225-7	HPLC-AAGF
1,2,4,5-Tetrachlorobenzene	NIOSH 5517	3-12	0.01-0.2	XAD-2 PTFE filter	ORBO-44 -----	20260-U special order	226-30-04	GC-ECD
1,1,1,2-Tetrachloro-2,2-difluoroethane	NIOSH 1016	0.5-2	0.01-0.035	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,1,1,2-Tetrachloro-2,2-difluoroethane	OSHA 7	2	0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,1,2,2-Tetrachloro-1,2-difluoroethane	NIOSH 1016	0.5-2	0.01-0.035	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,1,2,2-Tetrachloro-1,2-difluoroethane	OSHA 7	2	0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,1,2,2-Tetrachloroethane	NIOSH 1019	3-30	0.01-0.2	petroleum charcoal	ORBO-303	20040-U	226-38	GC-FID
1,1,2,2-Tetrachloroethane	OSHA 7	10	0.2	petroleum charcoal	ORBO-303	20040-U	226-38	GC-FID
Tetrachloroethylene (Perchloroethylene)	NIOSH 1003	0.2-40	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Tetrachloroethylene	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Tetrachloronaphthalene	NIOSH 2/S130	100	1.3	glass fiber filter fritted bubbler	----- -----	23378 64835-U	225-7 225-36-2	GC-FID
2,3,4,6-Tetrachlorophenol	OSHA 45	48	0.2	XAD-7 XAD-7/glass fiber filter (set of 3)	ORBO-602	custom	226-97	HPLC-UV
Tetraethyl Lead	NIOSH 2533	30-200	0.01-1	XAD-2	ORBO-605	20049	NA	GC-PID
Tetraethyl Pyrophosphate	NIOSH 2504	20-48	0.01-0.2	(2) Chromosorb 102	ORBO-707	custom	226-105	GC-FPD
Tetraethyltin	OSHA 110	3-48	0.2	XAD-7 (Supelpak 70)	ORBO-47	20349	226-95	GC-FID
Tetraethylene Glycol	NIOSH 5523	5-60	0.5-2	XAD-7/glass fiber filter (OVS)	ORBO-47P	custom	226-57	GC-FID
Tetrahydrofuran	NIOSH 1609	1-9	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Tetrahydrofuran	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Tetramethyl Lead (as Pb)	NIOSH 2534	15-100	0.01-0.2	XAD-2	ORBO-609	20051	226-30-06	GC-PID
Tetramethylperoxidicarbonyldiamide (Thiram)	NIOSH 5005	10-400	1-4	PTFE filter	-----	23383	225-1710	HPLC-UV
Tetramethyl Succinonitrile	NIOSH 3/S155	50	1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Tetramethyl Succinonitrile	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Tetramethyl Thiourea	NIOSH 3505	50-250	0.2-1	impinger	-----	20270-U	225-36-1	VIS
Tetranitromethane	NIOSH 3/S224	250	1	impinger	-----	20270-U	225-36-1	GC-FID
Tetranitromethane	NIOSH 3513	20-250	0.5-1	impinger	-----	20270-U	225-36-1	GC-FID
Tetryl (2,4,6-trinitrophenylmethylatramine)	NIOSH 3/S225	100	1.5	MCE filter	-----	23381	225-5	VIS
Thallium	NIOSH 7300	25-2000	1-4	MCE filter	-----	23381	225-5	ICP-AES
Thallium (as Tl)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Thiophene	NIOSH 1/255	70	1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FPD
Thiram	NIOSH 5005	10-400	1-4	PTFE filter	-----	23383	225-1710	HPLC-UV
Tin (as Sn), Tin Oxide	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Tin (Organic Compounds)	NIOSH 5504	50-500	1-1.5	XAD-2 glass fiber filter	ORBO-614 -----	custom 23378	226-30 225-7	HPLC-AAGF
Titanium	NIOSH 7300	5-100	1-4	MCE filter	-----	23381	225-5	ICP-AES

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equip	Analytical Technique
Titanium Diboride	NIOSH 6/323	500	1.7	PVC filter	-----	23387	225-8-01	XRD
Titanium Dioxide	NIOSH 3/S385	100	1.5	MCE filter	-----	23381	225-5	AAS
o-Tolidine	NIOSH 5013	150-500	1-3	PTFE filter	-----	23389	225-17	HPLC-UV
o-Tolidine	OSHA 71	100	1	glass fiber filter c/w H2SO4	822	20065	225-9004	GC-ECD
Toluene	NIOSH 1500	2-8	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Toluene	NIOSH 1501	1-8	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Toluene	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Toluene	NIOSH 4000	15m-8h		3M #3500 diffusive sampler	-----	20249-U		GC-FID
Toluene	OSHA 111	12	0.05	coconut charcoal or Carbosieve SIII (Anasorb 747) or 3M #3520 diffusive sampler	ORBO-32S ORBO-91 -----	20267-U 20360 207121	226-01 226-81A 575-002	GC-FID
Toluene	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
2,4-Toluenediamine	NIOSH 5516	30-500	1	impinger	-----	20270-U	225-36-1	HPLC-UV
2,4-Toluenediamine	OSHA 65	100	1	glass fiber filter c/w H2SO4	822	20065	225-9004	GC-ECD
2,6-Toluenediamine	NIOSH 5516	30-500	1	impinger	-----	20270-U	225-36-1	HPLC-UV
2,6-Toluenediamine	OSHA 65	100	1	glass fiber filter c/w H2SO4	822	20065	225-9004	GC-ECD
2,4-Toluene Diisocyanate (TDI)	ASTM D5836	15	1	glass fiber filter c/w 1-2PP	ORBO-80	20811	225-9002	HPLC-UV
2,4-Toluene Diisocyanate (TDI)	OSHA 42	15	1	glass fiber filter c/w 1-2PP	ORBO-80	20811	225-9002	HPLC-UV
Toluene-2,4-diisocyanate	NIOSH 5522	15-380	1-2	impinger	-----	20270-U	225-36-1	HPLC-FL
Toluene-2,4-diisocyanate (TDI)	NIOSH 2535	2-170	0.2-1	glass wool c/w NPMPA	ORBO-1007	custom	special order	HPLC-UV
Toluene-2,4-diisocyanate (TDI)	NIOSH 5521	5-500	1	impinger	-----	20270-U	225-36-1	HPLC-UV ELCD
2,6-Toluene Diisocyanate (TDI)	ASTM D5836	15	1	glass fiber filter c/w 1-2PP	ORBO-80	20811	225-9002	HPLC-UV
2,6-Toluene Diisocyanate (TDI)	OSHA 42	15	1	glass fiber filter c/w 1-2PP	ORBO-80	20811	225-9002	HPLC-UV
Toluene-2,6-diisocyanate	NIOSH 5521	5-500	1	impinger	-----	20270-U	225-36-1	HPLC-UV ELCD
Toluene-2,6-diisocyanate	NIOSH 5522	15-380	1-2	impinger	-----	20270-U	225-36-1	HPLC-FL
p-Toluenesulfonic Acid	NIOSH 5043	10-1000	1-3	glass fiber filter	-----	23376	225-16	HPLC-UV
m-Toluidine	OSHA 73	100	1	glass fiber filter c/w H2SO4	822	20065	225-9004	GC-ECD
o-Toluidine	NIOSH 2002	10-150	0.02-1	silica gel	ORBO-52S	20229	226-10	GC-FID
o-Toluidine	NIOSH 2017	5-50	0.2	silica gel + glass fiber filter c/w H2SO4	ORBO-507 822	20870-U 20065	226-15 225-9004	GC-FID
o-Toluidine	OSHA 73	100	1	glass fiber filter c/w H2SO4	822	20065	225-9004	GC-ECD
p-Toluidine	OSHA 73	100	1	glass fiber filter c/w H2SO4	822	20065	225-9004	GC-ECD
Tremolite Fibers	NIOSH 7400	400-	0.5-16	MCE filter	-----	23380-U	225-19	PCM
Tribromomethane	NIOSH 1003	4-70	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Tributyl Phosphate	NIOSH 5034	2-100	1-3	MCE filter	-----	23381	225-5	GC-FPD
Tributyltin Chloride	NIOSH 5504	50-500	1-1.5	XAD-2 glass fiber filter	ORBO-614 -----	custom 23378	226-30 225-7	HPLC-AAGF
1,2,4-Trichlorobenzene	NIOSH 5517	3-12	0.01-0.2	XAD-2 PTFE filter	ORBO-44 -----	20260-U special order	226-30-04	GC-ECD
1,1,1-Trichloroethane	NIOSH 1003	0.1-8	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,1,1-Trichloroethane	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
1,1,1-Trichloroethane	OSHA 14	3	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,1,2-Trichloroethane	NIOSH 1003	2-60	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,1,2-Trichloroethane	OSHA 11	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Trichloroethylene	NIOSH 1022	1-30	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Trichloroethylene	NIOSH 3701	<80% vol	0.02-0.05	5L sampling bag	-----	24655	232-05	GC-PID
Trichloroethylene	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Trichlorofluoromethane	NIOSH 1006	0.3-7	0.01-0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Trichloroisocyanuric Acid	NIOSH 5/314	400	1.7	PVC filter	-----	23387	225-8-01	titration
Trichloronaphthalene	NIOSH 2/S128	100	1.3	glass fiber filter fritted bubbler	----- -----	23378 64835-U	225-7 225-36-2	GC-FID
1,2,3-Trichloropropane	NIOSH 1003	0.6-60	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
1,2,3-Trichloropropane	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,1,2-Trichloro-1,2,2-trifluoroethane (Freon113)	NIOSH 1020	0.1-3	0.01-0.05	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
1,1,2-Trichloro-1,2,2-trifluoroethane (Freon113)	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
1,1,2-Trichloro-1,2,2-trifluoroethane (Freon 113)	OSHA 113	1	0.05	(2) Carbosieve S-III (Anasorb CMS)	ORBO-91	20360	226-121	GC-FID
Tricyclohexyltin Hydroxide	NIOSH 5504	50-500	1-1.5	XAD-2 glass fiber filter	ORBO-614 -----	custom 23378	226-30 225-7	HPLC-AAGF
Triethanolamine	NIOSH 3509	5-300	0.5-1	impinger	-----	20270-U	225-36-1	IC
Triethylamine	NIOSH 3/S152	100	1	fritted bubbler	-----	64835-U	225-36-2	GC-FID
Triethylene Glycol	NIOSH 5523	5-60	0.5-2	XAD-7/glass fiber filter (OVS)	ORBO-47P	custom	226-57	GC-FID
Triethylenetetramine (TETA)	NIOSH 2540	1-20	0.01-0.1	XAD-2 c/w 10% NITC	ORBO-654	custom	226-30-18	HPLC-UV
Triethylenetetramine (TETA)	OSHA 60	10	0.1	XAD-2 c/w 10% NITC	ORBO-654	custom	226-30-18	HPLC-UV
Trifluorobromomethane	NIOSH 1017	0.1-1	0.01-0.05	(2) coconut charcoal	ORBO-32S + ORBO-32L	20267-U 20228	226-01 226-09	GC-FID
Trimellitic Anhydride	NIOSH 5036	400-1000	1.5-2	37mm PVC-copolymer filter 0.8um	-----	special order		GC-FID
Trimellitic Anhydride	OSHA 98	480	2	glass fiber filter c/w DMBA & DOP	894	custom	225-9010	HPLC-UV
2,4,7-Trinitrofluorene-9-one	NIOSH 5018	100-500	1-3	PTFE filter	-----	special order	225-17-04	HPLC-UV
2,4,6-Trinitrotoluene (TNT)	OSHA 44	60	1	Tenax/glass fiber filter	ORBO-79	custom	NA	GC-TEA
Triorthocresyl Phosphate	NIOSH 5037	2-100	1-3	MCE filter	-----	23381	225-5	GC-FPD
Triphenyl Phosphate	NIOSH 5038	10-400	1-3	MCE filter	-----	23381	225-5	GC-FPD
Tungsten	OSHA ID 213	30-480	2	MCE filter	-----	23381	225-5	ICP-AES
Tungsten, Insoluble	NIOSH 7074	200-1000	1-4	MCE filter	-----	23381	225-5	AAS
Tungsten, Soluble	NIOSH 7074	200-1000	1-4	MCE filter	-----	23381	225-5	AAS
Turpentine	NIOSH 1551	1-10	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Turpentine	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
V M & P Naphtha	NIOSH 1550	1.30-20	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Valeraldehyde	NIOSH 2536	0.5-10	0.01-0.04	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-FID
Valeraldehyde	NIOSH 2539	5	0.01-0.05	XAD-2 c/w 2-HMP	ORBO-23	20257-U	226-118	GC-FID
Valeraldehyde	OSHA 85	3	0.05	glass fiber filter c/w DNPH	827	20069	225-9003	HPLC-UV
Vanadic Anhydride	NIOSH 7504	200-1000	1.7 or 2.2	PVC filter	-----	23387	225-8-01	XPD
Vanadic Oxide	NIOSH 7504	200-1000	1.7 or 2.2	PVC filter	-----	23387	225-8-01	XPD
Vanadium	NIOSH 7300	5-2000	1-4	MCE filter	-----	23381	225-5	ICP-AES
Vanadium (V2O5) Dust (as V)	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Vanadium (V2O5) Fume (as V)	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Vanadium Oxides	NIOSH 7504	200-1000	1.7 or 2.2	PVC filter	-----	23387	225-8-01	XPD
Vanadium Pentoxide	OSHA ID 185	816	1.7	PVC filter	-----	23387	225-8-01	XRF/XRD
Vinyl Acetate	NIOSH 1453	0.085-24	0.1-0.2	Carboxen 564	ORBO-92	20362	NA	GC-FID
Vinyl Acetate	OSHA 51	24	0.1	Carboxen 564	ORBO-92	20362	NA	GC-FID
Vinyl Benzene (Styrene)	NIOSH 1501	1-14	0.01-1	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Vinyl Bromide	NIOSH 1009	2-10	0.01-0.2	coconut charcoal	ORBO-32L	20228	226-09	GC-FID
Vinyl Bromide	OSHA 8	5	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Vinyl Chloride	ASTM D4766	25	0.05-0.1	coconut charcoal	ORBO-32J	custom	226-16	GC-FID
Vinyl Chloride	NIOSH 1007	0.7-5	0.05	(2) coconut charcoal	ORBO-301	20039	NA	GC-FID
Vinyl Chloride	OSHA 75	3	0.05	Carbosieve S-III	ORBO-91	20360	226-121	GC-FID

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Vinyl Toluene	NIOSH 1501	24	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Vinyl Toluene	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Vinylidene Chloride (1,1-Dichloroethene)	NIOSH 1015	2.5-7	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Vinylidene Chloride	OSHA 19	3	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
Volatile Organic Compounds (VOST)	EPA 0030	20	1	Tenax + Tenax/charcoal	VOST 100 + VOST 200	20074-U 20075-U	226-134 226-135	GC-MS
Volatile Organic Compounds (VOST)	EPA 0031	20	1	(2) Tenax + (1) Carbosieve S-III (Anasorb 747)	VOST 100 + VOST 150	20074-U custom	226-134 226-133	GC-MS
Volatile Organic Compounds (VOCs)	EPA IP-1A			canister	-----	NA	NA	GC-MS
Volatile Organic Compounds (VOCs)	EPA IP-1B	17-23	0.02-0.06	3-bed thermal desorption tube	Carbotrap 300	20379		TD-GC/MS
Volatile Organic Compounds (VOCs)	EPA TO-1	varies		3-bed thermal desorption tube	Carbotrap 300	20379		GC-MS
Volatile Organic Compounds (VOCs)	EPA TO-2	20-100	0.020-0.5	3-bed thermal desorption tube	Carbotrap 300	20379		GC-MS
Volatile Organic Compounds (VOCs)	EPA TO-3	0.1-1	0.01-0.1	3-bed thermal desorption tube	Carbotrap 300	20379		GC/ECD/FID
Volatile Organic Compounds (VOCs)	EPA TO-14			canister	-----	NA	NA	GC-MS or NPD/FID/ECD/ PID
Volatile Organic Compounds (VOCs)	EPA TO-15		0.0035-0.007	canister focusing trap	----- Purge Trap H	NA 20321	NA NA	GC-MS
Volatile Organic Compounds (VOCs)	EPA TO-16			FTIR and other remote sensing instruments				FTIR
Volatile Organic Compounds (VOCs)	EPA TO-17	1 & 4	0.01-0.2	multi-bed thermal desorption tube	Carbotrap 217 Carbotrap 317 Air Toxics Air Toxics Carbotrap 300 Carbotrap 300	20895-U 20877 25086 25051 25085 25050		TD-GC/MS
Volatile Organic Compounds (VOCs)	ASTM D6196	<10	<0.2	multi-bed thermal desorption tube	Carbotrap 217 Carbotrap 317 Air Toxics Air Toxics Carbotrap 300 Carbotrap 300	20895-U 20877 25086 25051 25085 25050		TD-GC/MS
Volatile Organic Compounds (VOCs)	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243		TD-GC/MS
Warfarin	NIOSH 5002	200-1000	1-4	PTFE filter	-----	23383	225-1710	HPLC-UV
Welding Fumes	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Xylene	NIOSH 2549	1-6	0.01-0.05	3-bed thermal desorption tube	Carbotrap 349	20243	NA	TD-GC/MS
Xylene	OSHA 7	10	0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
m-Xylene	NIOSH 1501	2-23	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
o-Xylene	NIOSH 1501	2-23	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
p-Xylene	NIOSH 1501	2-23	0.01-0.2	coconut charcoal	ORBO-32S	20267-U	226-01	GC-FID
m-Xylene- $\alpha,\alpha'$ -diamine	OSHA 105	15	1	glass fiber filter c/w H2SO4	822	20065	225-9004	HPLC-UV
p-Xylene- $\alpha,\alpha'$ -diamine	OSHA 105	15	1	glass fiber filter c/w H2SO4	822	20065	225-9004	HPLC-UV
2,4-Xylidine	NIOSH 2002	3-20	0.02-0.2	silica gel	ORBO-52S	20229	226-10	GC-FID
Yttrium	NIOSH 7300	5-1000	1-4	MCE filter	-----	23381	225-5	ICP-AES
Zinc, Zinc Compounds (as Zn)	NIOSH 7030	2-400	1-3	MCE filter	-----	23381	225-5	AAS
Zinc, Zinc Compounds (as Zn)	NIOSH 7300	5-200	1-4	MCE filter	-----	23381	225-5	ICP-AES

Compound	Method	Volume Liters	Rate Liters/Min	Medium and/or Sample Collection Device	Supelco Model No.	Supelco Cat. No.	SKC Equiv	Analytical Technique
Zinc, Zinc Compounds (as Zn)	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Zinc, Zinc Compounds (as Zn)	OSHA ID 125G	480	2	MCE filter	-----	23381	225-5	ICP-AES
Zinc Chloride Fume	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Zinc Oxide	NIOSH 7502	10-400	1-3	PVC filter	-----	special order		XPD
Zinc Oxide	OSHA ID 143	30-960	2	PVC filter	-----	23387	225-8-01	XRD
Zineb	OSHA 107	500	2	MCE filter	-----	23380-U	225-19	HPLC-UV
Zirconium	NIOSH 7300	5-200	1-4	MCE filter	-----	23381	225-5	ICP-AES
Zirconium	OSHA ID 121	480-960	2	MCE filter	-----	23381	225-5	AAS
Zirconium Oxides	NIOSH 4/250	90	1.7	PVC filter	-----	23387	225-8-01	XRD

If you have questions concerning this guide, please contact our Technical Service specialists. For additional method information, visit the agency websites listed below.



Supelco Park  
Bellefonte, PA 16823-0048 USA

### Technical Service

Telephone 800-359-3041  
814-359-3041  
Fax 800-359-3044  
814-359-5468  
email [techservice@sial.com](mailto:techservice@sial.com)

### Ordering

To place an order or to inquire about an existing order.

Telephone 800-247-6628  
800-325-3010  
Fax 800-325-5052  
email [supelco@sial.com](mailto:supelco@sial.com)

### Custom Products/Quotations

Gas chromatography, liquid chromatography, and sample preparation

Telephone 800-247-6628  
814-359-3441  
Fax 800-447-3044

Chemical standards and reagents

Telephone 814-359-5419  
Fax 814-359-5750

### Free Literature

Web Site [www.sigma-aldrich.com](http://www.sigma-aldrich.com)  
Telephone 800-447-3044  
814-359-3441

### Agency Websites

NIOSH [www.cdc.gov/niosh/homepage.html](http://www.cdc.gov/niosh/homepage.html)  
OSHA [www.osha.gov](http://www.osha.gov)  
EPA [www.epa.gov](http://www.epa.gov)  
ASTM [www.astm.org](http://www.astm.org)

Order/Customer Service 800-247-6628, 800-325-3010  
Fax 800-325-5052 • E-mail [supelco@sial.com](mailto:supelco@sial.com)

Technical Service 800-359-3041, 814-359-3041  
Fax 800-359-3044, 814-359-5468 • E-mail [techservice@sial.com](mailto:techservice@sial.com)

SUPELCO • 595 North Harrison Road, Bellefonte, PA 16823-0048 • 814-359-3441

We are committed to the success of our Customers, Employees and Shareholders through leadership in Life Science, High Technology and Service.

The SIGMA-ALDRICH Family  SIGMA  ALDRICH  Fluka  Riedel-deHaen  SUPELCO  ISOTEC

©1999 Sigma-Aldrich Co. Printed in USA. Supelco brand products are sold through Sigma-Aldrich, Inc. Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip. Supelco is a registered trademark of Sigma-Aldrich Co.

BUO T411091D



[sigma-aldrich.com/supelco](http://sigma-aldrich.com/supelco)