



ChemFiles

Volatile Precursors for NanoFabrication

Vol. 4 No. 3

ALD Precursors for
High κ and Barrier
Materials

Precursors for Low κ
Materials

Metal Amidinates

High Purity Inorganics
and Gases

III-V Material Sources

Manufacturing
Capabilities

Custom Packaging

Equipment, Books and
Software

Precursors for Nanofabrication: For Your Lab or for Your Fab

Sigma-Aldrich is pleased to present our complete line of high-purity organometallics and inorganics for micro and nanoelectronic applications. In this brochure you will find products for all aspects of atomic layer (ALD) and chemical vapor deposition (CVD). We feature precursors for high and low (κ) dielectric materials, transition metals for barrier and metallization applications, III-V materials, and epitaxy gases. This brochure features a discussion of the ALD process and technical

synopses on key application areas, equipment, books and software of interest to researchers and engineers.

Ready to begin? Look through our product index below by end material type and then go to the page to find the products you need. Do you have a precursor in mind but can't find it in this brochure? Need a precursor for a rare-earth oxide or a single source nitride precursor? Contact our Organometallics Product Manager at matsci@sial.com.

Material Deposited	Product Name	Page
Al ₂ O ₃ , Al, AlN, AlP AlAs, LaAlO ₃ , Aluminates	Aluminum <i>sec</i> -butoxide	2
	Aluminum tribromide	6
	Aluminum trichloride	6
	Diethylaluminum ethoxide	2
	Tris(ethylmethylamido)aluminum	2
	Triethylaluminum	2
	Triisobutylaluminum	2
	Trimethylaluminum	2
	Tris(diethylamido)aluminum	2
	Tris(ethylmethylamido)aluminum	2
AlAs, GaAs, InAs	Trimethylarsine	8
MgB ₂ , BN, B, B ₄ C, B ₂ O ₃ B doping	Diborane (10% in Hydrogen)	10
	Diborane-d ₆ (10% in D ₂ or He)	10
	Trimethylboron	10
	Trimethylboron-d ₉	10
Co, CoO, CoSi ₂	Bis(N,N'-Diisopropylacetamidinato)cobalt(II)	4
	Dicarbonyl(cyclopentadienyl)cobalt(I)	9
Cu, YBaCuO _x , CuO	(N,N'-Diisopropylacetamidinato)copper(I)	4
Fe, FeO	Bis(N,N'-di- <i>tert</i> -butylacetamidinato)iron(II)	4
Ga ₂ O ₃ , Ga, GaN, GaP, GaAs	Gallium tribromide	6
	Gallium trichloride (bead or slug)	6
	Triethylgallium	7
	Triisopropylgallium	7
	Trimethylgallium	7
	Tris(dimethylamido)gallium	7
	Tri- <i>tert</i> -butylgallium	7
Ge, GeO ₂ , GeSi	Digermane (10% in H ₂)	10
	Germane	10
	Tetramethylgermanium	10
HfO ₂ , Hf ₃ N ₄	Hafnium(IV) chloride	6
	Hafnium(IV) <i>tert</i> -butoxide	2
	Tetrakis(diethylamido)hafnium(IV)	2
	Tetrakis(dimethylamido)hafnium(IV)	2
	Tetrakis(ethylmethylamido)hafnium(IV)	2
In ₂ O ₃ , InN, InP, InAs In containing Solar Cells Indium Tin Oxide	Indium trichloride	6
	Indium(I) iodide (Anhydrous beads)	6
	Indium acetylacetonate	7
	Triethylindium	7
La ₂ O ₃ , LaAlO ₃	Tris(N,N'-Di- <i>tert</i> -butylacetamidinato)lanthanum(III)	4
Mg dopant in III-V	Bis(pentanethylcyclopentadienyl)magnesium	9
MoS ₂ , MoO ₂ , Mo	Molybdenum hexacarbonyl	6
	Molybdenum(V) chloride	6
	Molybdenum(VI) fluoride	6
GaN, InGaN, AlGaN, Si ₃ N ₄	N,N-dimethylhydrazine	8
	Ammonia	8
	Azidotrimethylsilane	8
Nb ₂ O ₅	Niobium(V) chloride	6
	Niobium(V) ethoxide	9

Material Deposited	Product Name	Page
Ni, NiO	Bis(methylcyclopentadienyl)nickel(II)	9
P doping, InP, GaP	Phosphine	10
	Phosphine-d ₃	10
	<i>tert</i> -Butylphosphine	8
	Tris(trimethylsilyl)phosphine	8
Pt, PtO ₂	Cyclopentadienyl(trimethyl)platinum(IV)	9
Ru, RuO ₂	Bis(ethylcyclopentadienyl)ruthenium(II)	9
Sb Source	Trimethylantimony	8
	Tris(dimethylamido)antimony	8
SiO ₂ /Si ₃ N ₄ /SiC	2,4,6,8-Tetramethylcyclotetrasiloxane	4
	Dimethoxydimethylsilane	4
	Disilane	10
	Methylsilane	10
	Octamethylcyclotetrasiloxane	4
	Silane	10
	Silane-d ₄	10
	Tris(isopropoxy)silanol	4
	Tris(<i>tert</i> -butoxy)silanol	4
	Tris(<i>tert</i> -pentoxy)silanol	4
Ta ₂ O ₅ /TaN	Pentakis(dimethylamido)tantalum(V)	9
	Tantalum(V) chloride	6
	Tantalum(V) ethoxide	9
	Tris(diethylamino)(<i>tert</i> -butylimido)-tantalum(V)	9
TiN/TiO ₂	Bis(diethylamido)bis(dimethylamido)-titanium(IV)	3
	Tetrakis(diethylamido)titanium(IV)	3
	Tetrakis(dimethylamido)titanium(IV)	3
	Tetrakis(ethylmethylamido)titanium(IV)	3
	Titanium(IV) bromide	6
	Titanium(IV) chloride	6
	Titanium(IV) <i>tert</i> -butoxide	3
V ₂ O ₅	Vanadium(V) oxytriisopropoxide	9
W, WO ₂ , WO ₃ , WC	Bis(<i>tert</i> -butylimido)bis(dimethylamido)-tungsten(VI)	9
	Tungsten hexacarbonyl	6
	Tungsten(VI) chloride	6
	Tungsten(VI) fluoride	6
Y ₂ O ₃ , YBaCuOx	Tris(N,N-bis(trimethylsilyl)amide)yttrium(III)	9
	Yttrium(III) butoxide solution 0.5M in toluene	9
ZnO	Diethylzinc	9
Zr ₃ N ₄ , ZrO ₂	Tetrakis(diethylamido)zirconium(IV)	3
	Tetrakis(dimethylamido)zirconium(IV)	3
	Tetrakis(ethylmethylamido)zirconium(IV)	3
	Zirconium(IV) bromide	6
	Zirconium(IV) chloride	6
	Zirconium(IV) <i>tert</i> -butoxide	3

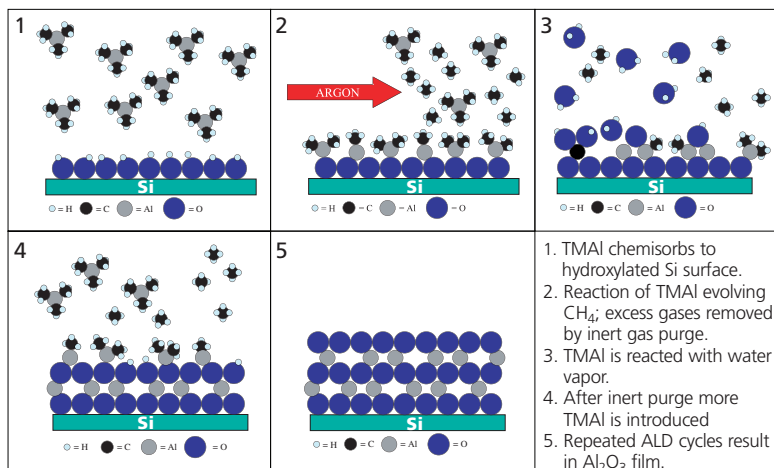
TO ORDER: Contact your local Sigma-Aldrich office (see back cover), call 1-800-558-9160 (USA), or visit sigma-aldrich.com.

Atomic Layer Deposition

Atomic Layer Deposition (ALD) is recognized as the key technology for the semiconductor industry to break below the 90 nm device node barrier, keeping device miniaturization on track with Moore's Law.¹ ALD processes can be used for the deposition of barrier materials,² nucleation layers, metallization,³ and high κ dielectrics.⁴ The ALD process is able to meet the demands of microelectronics miniaturization because it is a layer-by-layer build up of materials starting with the chemisorption of molecular precursors (Figure). These precursors are reacted at the substrate surface and the resulting layer is purged of by-products with inert gas.⁵ The process is repeated as needed to attain desired film characteristics. Because ALD is a self-limiting deposition process, it offers a high degree of control over film composition, thickness and 100% step-coverage, even over high aspect ratio steps.⁶ Contrast this with traditional CVD processes where gas-phase chemical reactions produce the final material to be deposited on the surface. While effective for traditional thick and thin-film applications, CVD does not offer the layer-by-layer control and step coverage of ALD. Furthermore, recent advances in processing technology are allowing for ALD processes to be used in high throughput.

The ALD mechanism is composed of two separate self-limiting steps: the substrate saturation process and the reactivity between the surface groups and precursor molecules. These requirements place several demands on the ALD precursor molecule. Not every molecule is ideal but we can help you find the material ideal for your application.

Sigma-Aldrich, a leader in *High Technology* products is proud to present our line of Volatile Precursors for Nanofabrication. These products encompass a wide variety of products to get you the final material phases you need.



Ideal ALD Precursor Characteristics

- High vapor pressure
- Thermal stability prior to deposition
- Ease of handling and transfer
- Chemisorbs to substrate
- Aggressive reaction with complementary precursors
- Noncorrosive to substrate
- High purity
- Low hazard by-products

High κ Materials

The miniaturization of integrated devices brings serious challenges to the continued use of current materials. The capacitors used in traditional microelectronic circuits, composed of $\text{Si}/\text{SiO}_2/\text{metal}$, will not be able to function due to the physical limits of the SiO_2 dielectric. At the nanometer dimension, the SiO_2 dielectric constant (κ) is not large enough to prevent leakage currents, leading to the unwanted discharge of the capacitor. New higher κ materials are under consideration and are listed in the table. 1.5-10 nm thick layers of Zr, Hf and Al oxides on silicon, grown by ALD processes exhibit much lower gate leakage than SiO_2 of equivalent thickness. Sigma-Aldrich offers a wide array of volatile Al, Hf and Zr precursors for your high κ deposition processes. We can tailor ligand designed to give you the physical properties and purity you demand for your systems.

Candidate High κ Materials

Material	Dielectric constant (κ)
$\text{Si}_3\text{N}_4/\text{SiO}_2$	5-6
Sc_2O_3	>10
Al_2O_3	8-9
Y_2O_3	15
HfO_2	21
ZrO_2	22
LaAlO_3	25
Ta_2O_5	25
La_2O_3	25-30
SrTiO_3	200
$\text{Ba}_x\text{Sr}_{1-x}\text{O}_3$	300



Ready to scale up? For competitive quotes on larger quantities or custom synthesis, contact Sigma-Aldrich Fine Chemicals at 1-800-336-9719 (USA), or visit www.sigma-aldrich.com/safc.

Precursors for High κ Materials

Aluminum

Trimethylaluminum

J1,0001-3

Purity: 99.9999%

CAS No. 75-24-1

MF: (C₃H₉)Al

FW: 72.09

mp: 15 °C

bp: 125–126 °C

Density: 0.752 g/mL



Typical ICP-MS results for J1,0001-3 in ppb

Ag	<1	Ge	3	Pt	<1
As	<1	Hf	8	Sb	<1
Au	<1	Hg	<1	Sn	<1
B	<1	In	<1	Sr	<1
Ba	4	K	<1	Ta	<1
Be	<1	La	<1	Th	<1
Bi	<1	Li	<1	Ti	4
Ca	6	Mg	1	Tl	<1
Cd	<1	Mn	<1	U	<1
Co	<1	Mo	1	V	<1
Cr	3	Na	44	W	<1
Cu	1	Nb	<1	Zn	2
Fe	4	Ni	2	Zr	<1
Ga	<1	Pb	<1		

Aluminum sec-butoxide

51,160-9

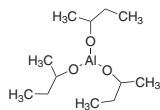
Purity: 99.99%

CAS No. 2269-22-9

FW: 246.33

bp: 200–206 °C @ 30 mmHg

Density: 0.967 g/mL



Hafnium

Tetrakis(dimethylamido)hafnium(IV)

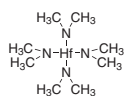
45,519-9

Purity: 99.99%

CAS No. 19962-11-9

FW: 354.80

mp: 26–29 °C



Tetrakis(ethylmethylamido)hafnium(IV)

55,312-3

Purity: 99.99%

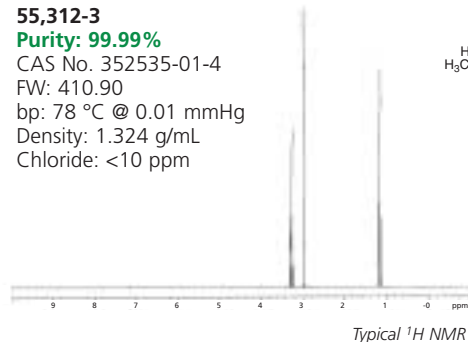
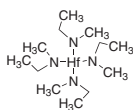
CAS No. 352535-01-4

FW: 410.90

bp: 78 °C @ 0.01 mmHg

Density: 1.324 g/mL

Chloride: <10 ppm



Triethylaluminum

J1,0001-8

Purity: 99.99%

CAS No. 97-93-8

MF: (C₂H₅)₃Al

FW: 114.17

mp: -50 °C

bp: 128–130 °C @ 50 mmHg

Density: 0.835 g/mL



Tris(diethylamido)aluminum

J1,0003-0

CAS No. 352546-72-6

FW: 486.74

mp: 28–31 °C

bp: 50 °C @ 20 mmHg

Density: 0.915 g/mL

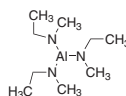


Also Available

J1,0002-9 Tris(ethylmethylamido)aluminum

25,674-9 Diethylaluminum ethoxide 97%

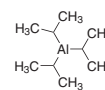
25,720-6 Triisobutylaluminum



J1,0002-9



25,674-9



25,720-6

Tetrakis(diethylamido)hafnium(IV)

45,520-2

Purity: 99.99%

CAS No. 19962-12-0

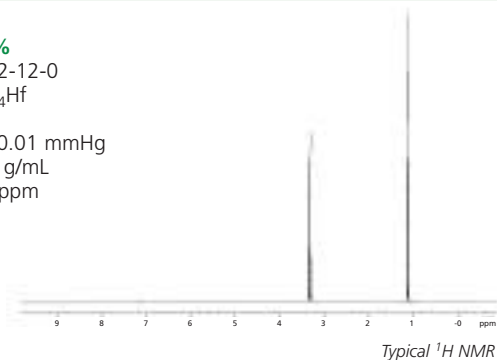
MF: [(C₂H₅)₂N]₄Hf

FW: 467.01

bp: 130 °C @ 0.01 mmHg

Density: 1.249 g/mL

Chloride: <10 ppm



Hafnium(IV) tert-butoxide

44,554-1

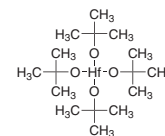
Purity: 99.99%

CAS No. 2172-02-03

FW: 473.95

bp: 90 °C @ 5 mmHg

Density: 1.166 g/mL



Precursors for High κ Materials—cont'd

Zirconium

Zirconium(IV) *tert*-butoxide

56,003-0

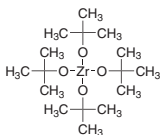
Purity: 99.999%

CAS No. 2081-12-1

FW: 383.68

bp: 81 °C @ 3 mmHg

Density: 0.985 g/mL



Tetrakis(dimethylamido)zirconium(IV)

57,921-1

Purity: 99.99%

CAS No. 19756-04-8

MF: $[(CH_3)_2N]_4Zr$

FW: 267.53

mp: 57–60 °C

Tetrakis(ethylmethylamido)zirconium(IV)

55,313-1

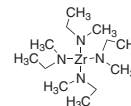
Purity: 99.99%

CAS No. 175923-04-3

FW: 323.63

bp: 81 °C @ 0.1 mmHg

Density: 1.049 g/mL



Tetrakis(diethylamido)zirconium(IV)

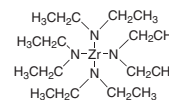
J1,0002-3

Purity: 99.99%

CAS No. 13801-49-5

FW: 379.74

bp: 128 °C @ 0.05 mmHg



Titanium Precursors for Barrier Materials

Tetrakis(diethylamido)titanium(IV)

46,986-6

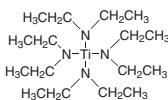
Purity: 99.99%

CAS No. 4419-47-0

FW: 336.42

bp: 112 °C @ 0.1 mmHg

Density: 0.931 g/mL



Tetrakis(dimethylamido)titanium(IV)

46,985-8

Purity: 99.999%

CAS No. 3275-24-9

MF: $[(CH_3)_2N]_4Ti$

FW: 224.21

bp: 50 °C @ 05 mmHg

Density: 0.947 g/mL

Tetrakis(ethylmethylamido)titanium(IV)

47,353-7

Purity: 99.99%

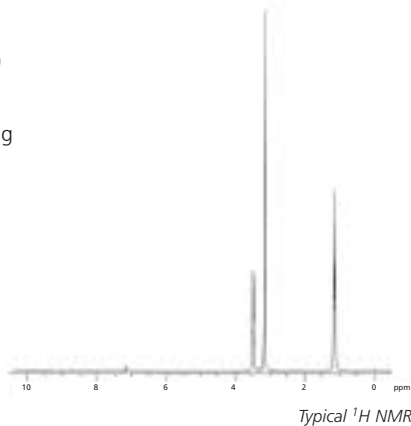
CAS No. 308103-54-0

MF: $[(CH_3)(C_2H_5)N]_4Ti$

FW: 280.32

bp: 80 °C @ 0.1 mmHg

Density: 0.923 g/mL



Bis(diethylamido)bis(dimethylamino)titanium(IV)

J1,0002-6

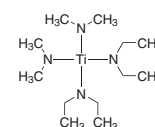
Purity: 99.99%

CAS No. 123798-13-0

FW: 280.31

bp: 80 °C @ 0.1 mmHg

Density: 0.931 g/mL



Titanium(IV) *tert*-butoxide

46,255-1

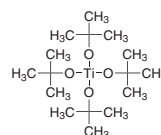
CAS No. 3087-39-6

MF: $[(CH_3)_3CO]_4Ti$

FW: 340.36

bp: 70 °C @ 0.2 mmHg

Density: 0.881 g/mL



Ready to scale up? For competitive quotes on larger quantities or custom synthesis, contact Sigma-Aldrich Fine Chemicals at 1-800-336-9719 (USA), or visit www.sigma-aldrich.com/safc.

Low κ Precursors

Low κ materials address technical challenges of increased chip performance by allowing for the reduction of spacing between metals in multi-level interconnects, interlayer dielectrics and passivation layers.⁸ Shrinking of device sizes leads to issues with resistance capacitance (RC) coupling, resulting in power dissipation, propagation delay and noise between circuits called crosstalk. Current low κ strategies include organosilicate glass (OSG) films for the 90-100 nm node ($\kappa < 3$), carbon doped

silicon oxide films ($\kappa < 3$), and porous silicate films for the 65 nm node ($\kappa < 2.6$). Porous films have some copper compatibility issues and reducing the κ of such films is accompanied by a loss of mechanical properties. Future low κ materials will be an extension of current CVD/ALD deposited non-porous films. Sigma-Aldrich has a variety of high quality low κ precursors. If you can't find what you want check the website or send a question to matsci@sial.com.

2,4,6,8-Tetramethylcyclotetrasiloxane

51,299-0

Purity: 99.999%

CAS No. 2370-88-9

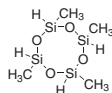
FW: 240.51

mp: -69 °C

bp: 134 mmHg

Density: 0.986 g/mL

Chloride: <10 ppm



Dimethoxydimethylsilane

55,668-8

Purity: 99.999%

CAS No. 1112-39-6

FW: 120.23

bp: 81.4 °C

Density: 0.880 g/mL



Octamethylcyclotetrasiloxane

23,569-5

Purity: 98+%

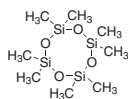
CAS No. 556-67-2

FW: 296.62

mp: 17-18 °C

bp: 175-175 °C

Density: 0.956 g/mL



Amidinate Precursors

Developed with Professor Roy Gordon at Harvard, Sigma-Aldrich presents these exciting new precursors for Atomic Layer Deposition. The amidinates are volatile, homoleptic transition metal precursors, which meet the rigorous standards for ALD applications. When employed in ALD processes, these precursors allow for the deposition of pure metals when reacted

with H₂ gas or oxides when co-reacted with water vapor.⁹ The precursors can be vaporized appreciably at temperatures between 50-100 °C for the transition metal compounds and at 130 °C for the lanthanum compound. Deposition temperatures range between 150-350 °C.

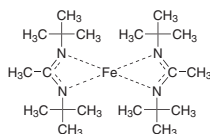
Bis(*N,N'*-di-*tert*-butylacetamidinato)iron(II)

J1,0000-6

Purity: 99%

CAS No. 635680-56-7

FW: 394.42



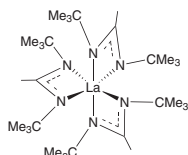
Tris(*N,N'*-Di-*tert*-butylacetamidinato)lanthanum(III)

J1,0000-8

Purity: 99.999%

FW: 646.77

mp: 165 °C



(*N,N'*-Diisopropylacetamidinato)copper(I)

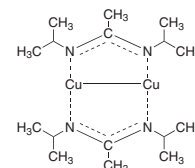
J1,0000-9

Purity: 99%

CAS No. 635680-64-7

FW: 409.56

mp: 152-154 °C



Bis(*N,N'*-Diisopropylacetamidinato)cobalt(II)

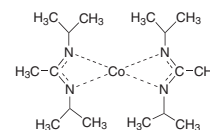
J1,0001-0

Purity: 99.98%

CAS No. 635680-58-9

FW: 341.40

mp: 80-83 °C



TO ORDER: Contact your local Sigma-Aldrich office (see back cover), call 1-800-558-9160 (USA), or visit sigma-aldrich.com.

Sigma-Aldrich Fine Chemicals Capabilities:

Manufacturing Capabilities for Volatile Precursors

Sigma Aldrich Fine Chemicals (SAFC) is known as a worldwide leader in the manufacture of high purity chemicals and biotechnology products. We have decades of experience in the production, purification and handling of air-sensitive organometallics at our ISO9001 certified Sheboygan, Wisconsin facility, manufacturing high-purity inorganics through our Aldrich-APL facility in Urbana, Illinois and stable isotope materials at our ISOTEC™ plant in Miamisburg, Ohio. We have translated this expertise into a world-class capability to manufacture a broad range of high technology products. No molecule is beyond our grasp, whether you need a few grams for proof-of-concept or you are ready to go to production, SAFC can provide you with the quality materials you need to get the job done.



SAFC Process Development Lab

As the microelectronics industry continues to change, we have the knowledge and facilities to meet your needs through our dedicated R&D and production sites. Along with these resources and our strong collaborations with key researchers in the field, Sigma-Aldrich continues to develop new products for micro and nanofabrication. We can provide these products with the highest purities (5N, 6N, 7N and higher) up to multi-kilogram scales. We back up our quality with a full range of analytical tests such as ICP-MS. Our volatile precursor products are packaged to specification using our state-of-the-art liquid filling station, to ensure you get what you order every time. Need custom packaging? We have a variety of custom packaging options as shown on pages 10 and 11.



Liquid filling station

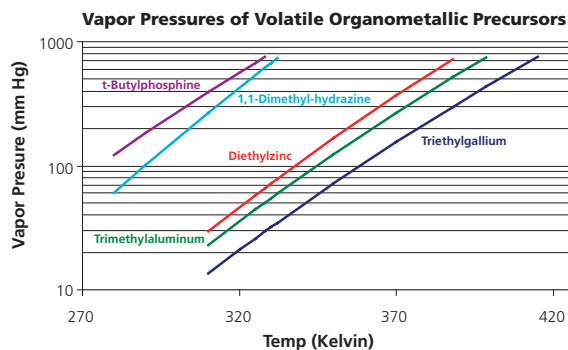


Sheboygan, WI Manufacturing

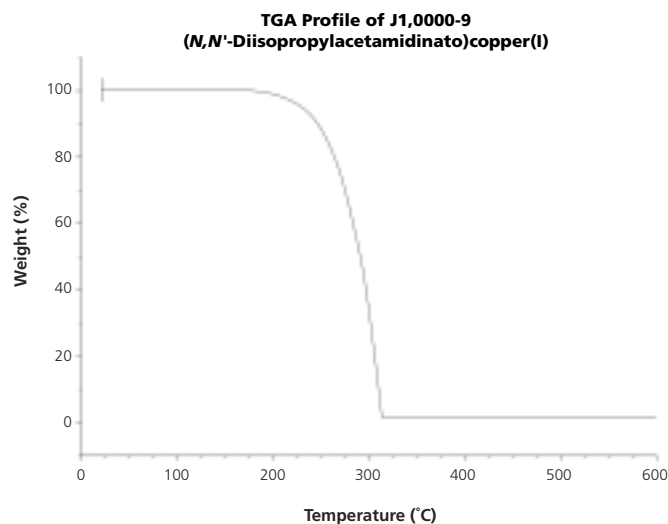
Analytical Capabilities for Volatile Precursors

ALD and MOCVD processes require key volatility, decomposition, product quality and safety and handling data. Sigma-Aldrich can provide the data you need to make the best choice of precursor including:

- Vapor Pressure Equation or Graph
- DSC and TGA for decomposition/thermal properties
- ICP-MS for ultra trace elements
- MSDS on all products for safety and handling



For product data send an inquiry to matsci@sial.com; for MSDS and safety go to the website www.sigma-aldrich.com and search by product number.



Courtesy of Roy Gordon at Harvard University.



Ready to scale up? For competitive quotes on larger quantities or custom synthesis, contact Sigma-Aldrich Fine Chemicals at 1-800-336-9719 (USA), or visit www.sigma-aldrich.com/safc.

High Purity Inorganics

Aldrich-APL (AAPL) offers a complete line of the highest purity inorganic salts and materials for the micro and nanoelectronics market. Through proprietary processes, AAPL manufactures metal halide salts as anhydrous beads or powders. These ultra-high purity materials have total metals impurities 4N-5N or higher and



water contents less than 100 ppm. AAPL can bring its expertise in reactive gas processing, high-temperature synthesis and metal distillations to meet your materials demands. AAPL products are carefully packaged under moisture and air-free conditions, ensuring the quality you expect. Whatever your scale, your demands can be met. If you are looking for a special material and cannot find it here, contact us at matsci@sial.com.

Gallium trichloride (beads)

42,712-8

Purity: 99.999+%

CAS No. 13450-90-3

MF: GaCl₃

FW: 176.08

Density: 2.47 g/mL

Also available as powder or slugs.
Call for your product configuration

Typical ICP-MS Results for 42,712-8 in ppm

Ag	>0.04	Eu	>0.04	Na	<0.04	Sm	<0.04
Al	<0.04	Fe	0.58	Nb	<0.04	Sn	<0.04
As	<0.04	Ga	Major	Nd	<0.04	Sr	<0.04
Au	<0.04	Gd	<0.04	Ni	0.15	Ta	<0.04
B	0.17	Ge	<0.04	Os	<0.04	Tb	<0.04
Ba	0.18	Hf	<0.04	Pb	<0.04	Te	<0.04
Be	<0.04	Hg	<0.04	Pd	<0.04	Th	<0.04
Bi	<0.04	Ho	<0.04	Pr	0.16	Ti	<0.04
Ca	1.42	In	<0.04	Pt	<0.04	Tl	<0.04
Cd	<0.04	Ir	<0.04	Rb	<0.04	Tm	<0.04
Ce	0.25	K	1.12	Re	<0.04	U	<0.04
Co	<0.04	La	<0.04	Rh	<0.04	V	<0.04
Cr	<0.04	Li	<0.04	Ru	<0.04	W	<0.04
Cs	<0.04	Lu	<0.04	Sb	<0.04	Y	<0.04
Cu	<0.04	Mg	<0.04	Sc	<0.04	Yb	<0.04
Dy	<0.04	Mn	<0.04	Se	<0.04	Zn	<0.04
Er	<0.04	Mo	<0.04	Si	3.72	Zr	<0.04

Hafnium(IV) chloride

59,059-2

Purity: 99.9%

CAS No. 13499-05-3

MF: HfCl₄

FW: 320.30

Aluminum trichloride (powder)

56,391-9

Purity: 99.999%

CAS No. 7446-70-0

MF: AlCl₃

FW: 133.34

Sublimes at 181 °C

Typical ICP-MS Results for 56,391-9 in ppm

Ag	<0.04	Eu	0.07	Na	1.55	Sm	0.7
Al	<0.04	Fe	0.38	Nb	<0.04	Sn	<0.04
As	<0.04	Ga	1.01	Nd	<0.04	Sr	<0.04
Au	<0.04	Gd	<0.04	Ni	0.13	Ta	<0.04
B	0.23	Ge	<0.04	Os	<0.04	Tb	<0.04
Ba	<0.04	Hf	0.08	Pb	0.5	Te	0.05
Be	<0.04	Hg	<0.04	Pd	<0.04	Th	<0.04
Bi	<0.04	Ho	<0.04	Pr	<0.04	Ti	0.2
Ca	1.03	In	<0.04	Pt	0.05	Tl	<0.04
Cd	<0.04	Ir	<0.04	Rb	<0.04	Tm	<0.04
Ce	<0.04	K	<0.04	Re	<0.04	U	<0.04
Co	<0.04	La	<0.04	Rh	<0.04	V	<0.04
Cr	.023	Li	<0.04	Ru	<0.04	W	<0.04
Cs	0.05	Lu	<0.04	Sb	<0.04	Y	<0.04
Cu	0.12	Mg	1.29	Sc	<0.04	Yb	<0.04
Dy	<0.04	Mn	<0.04	Se	0.11	Zn	1.05
Er	<0.04	Mo	<0.04	Si	1.07	Zr	<0.04

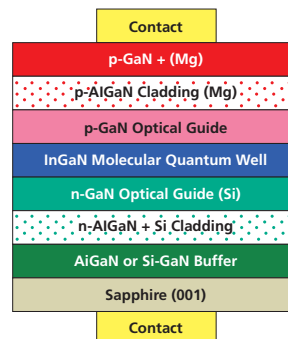
Catalog #	Name	Formula	Metals Purity %	mp °C	bp °C/mmHg*	Sublime °C
44,960-1	Aluminum tribromide	AlBr ₃	99.999	97.5	263.3/747	
45,086-3	Gallium tribromide	GaBr ₃	99.999	121.5	279	
42,941-4	Indium trichloride	InCl ₃	99.999			300
41,365-8	Indium(I) iodide	InI	99.999 beads	351		
57,776-6	Molybdenum hexacarbonyl	Mo(CO) ₆	99.9		156	
64,245-2	Molybdenum(V) chloride	MoCl ₅	99.99 powder	194	268	
39,909-4	Molybdenum(VI) fluoride	MoF ₅	99.9	17.5	37	
51,069-6	Niobium(V) chloride	NbCl ₅	99.995	204.7	254	
51,068-8	Tantalum(V) chloride	TaCl ₅	99.999	216	242	
45,160-6	Titanium(IV) bromide	TiBr ₄	99.99	39	230	
25,431-2	Titanium(IV) chloride	TiCl ₄	99.995	-24	136.4	
39,910-8	Tungsten(VI) fluoride	WF ₆	99.9	2.3	17.5	
47,295-6	Tungsten hexacarbonyl	W(CO) ₆	99.9	150		
64,516-8	Tungsten(VI) chloride	WCl ₆	99.99 powder	275	346.7	
40,056-4	Zirconium(IV) bromide	ZrBr ₄	98.0			357
64,764-0	Zirconium(IV) chloride	ZrCl ₄	99.99			331

*@760 mmHg unless otherwise stated.

TO ORDER: Contact your local Sigma-Aldrich office (see back cover), call 1-800-558-9160 (USA), or visit sigma-aldrich.com.

III-V Material Sources

At the forefront of optoelectronic, lighting and data storage applications, III-V semiconductor materials continue to generate exciting new applications. III-V quantum dot heterostructures can address the technical challenges of data transfer and storage issues in telecom and datacom systems. Blue laser GaN LEDs are expected to make a significant impact on HD-DVD storage in the coming years.¹⁰ The figure shows the necessary material layers to construct the blue-violet lasers.¹¹ GaAs and InP are primary components of low noise and power amplifiers. High brightness LEDs for mobile phones and automotive display lighting are made from a variety of III-V semiconductors. Sigma-Aldrich can provide the highest quality III-V precursors for research and production applications. For aluminum precursors see page 2.



Gallium Precursors

Trimethylgallium

J1,0001-5

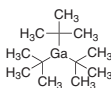
CAS No. 1445-79-0
FW: 114.83
mp: -15.8 °C
bp: 55.7 °C
Density: 1.10 g/mL



Tri-tert-butylgallium

J1,0003-4

CAS No. 55681-99-7
FW: 241.07
bp: 72 °C @ 60 mmHg

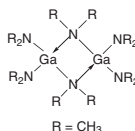


Tris(dimethylamido)gallium

54,653-4

Purity: 98%

CAS No. 57731-40-5
FW: 403.90
mp: 104–105.5 °C



Triethylgallium

J1,0001-7

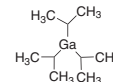
CAS No. 1115-99-7
FW: 156.91
mp: -82 °C
bp: 109 °C @ 300 mmHg



Triisopropylgallium

J1,0001-6

CAS No. 54514-59-9
FW: 198.99
bp: 95-98 °C @ 60 mmHg



Indium Precursors

Triethylindium

57,091-5

CAS No. 923-34-2
FW: 202.00
mp: -32 °C
bp: 184 °C
Density: 1.260 g/mL

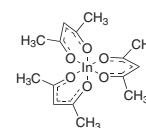


Indium Acetylacetonate

I,330-0

Purity: 99.99%

CAS No. 14405-45-9
FW: 412.15
mp: 187-189 °C



References

- Ritala, M.; Leskelä M.; Handbook of Thin Film Materials, Vol. 1; Nalwa, H.S. (ed); Academic Press: New York, **2001**, p 103-159.
- Hoyas, A.M.; Schumacher, J.; Shamiryan, D.; Waeterloos, J.; Besling, W.; Celis, J.P.; Maex, K.; *J. App. Phys.*, **2004**, 95, 381.
- Becker, J.S.; Suh, S.; Wang, S.; Gordon, R.G.; *Chem Mater.* **2003**, 15, 2969.
- Kim, S.H.; Oh, S.; Kim, K.B.; *Appl. Phys. Lett.*, **2003**, 82, 25.
- Park, J.S.; Park, H.S.; Kang, S.W.; *J. Electrochem. Soc.*, **2002**, 149, C28.
- Klaus, J.W.; Ferro, S.J.; George, S.M.; *J. Electrochem. Soc.*, **2001**, 147, 1175.
- Aaltonen, T.; Ritala, M.; Sajavaara, T.; Keinonen, J.; Leskelä, M.; *Chem. Mater.*, **2003**, 15, 1924.
- Lim, B.S.; Rahtu, A.; Gordon, R.G.; *Nature Materials*, **2003**, 2, 749.
- Elam, J.W.; Nelson, C.E.; Grobbs, R.K.; George, S.M.; *Surf. Sci.*, **2001**, 479, 121.
- Putkonen, M.; Niinistö, J.; Kukli, K.; Sajavaara, T.; Karppinen, M.; Yamauchi, H.; Niinistö, L. *Chem. Vap. Dep.*, **2003**, 9, 207.
- Lee, S.Y.; Kim, H.; McIntyre, P.; Saraswat K.C.; Byun J.S.; *Appl. Phys. Lett.* **2003**, 82, 2874.
- Kukli, K.; Ritala, M.; Sajavaara, T.; Keinonen, J.; Leskelä, M.; *Chem. Vap. Dep.*, **2002**, 8, 199.
- Sneh, O.; Clark-Phelps, R.B.; Londergan, A.R.; Winkler, J.L.; Seidel, T.E.; *Thin Solid Films*, **2002**, 402, 248.
- Crowell, J.E.; *J. Vac. Sci. Technol.* **2003**, A21, S88.
- Ritala, M.; Kukli, K.; Rahtu, A.; Raisanen, P.I.; Sajavaara, T.; Keinonen, J.; Leskelä, M.; *Science*, **2000**, 288, 319.
- George, S.M.; Klaus, J.W.; Ferguson, J.D.; *Mater. Res., Soc. Symp. Proc.*, **2000**, 616, 93.
- Gordon, R.G.; Hausmann, D.; Shepard, J.; *Chem. Vap. Dep.*, **2003**, 9, 73.
- Becker, J.S.; Gordon, R.G.; *Appl. Phys. Lett.* **2003**, 82, 2239.
- Hand, A. *Semiconductor International*, **2003**, On-line edition. <http://www.reedelectronics.com/semiconductor/article/CA293949>.
- Shamiryan, D.; Abell, T.; Iacopi, F.; Maex, K.; *Materials Today*, **2004**, 7, 34.
- Lim, B.S.; Rahtu, A.; Park, J.S.; Gordon, R.G.; *Inorg. Chem.*, **2003**, 42, 7951.
- Compound Semiconductor*, **2004**, April, p13-15.
- Monemar, B.; Pozina G., *Prog. Quantum Electronics*, **2000**, 24, 239-290.



Ready to scale up? For competitive quotes on larger quantities or custom synthesis, contact Sigma-Aldrich Fine Chemicals at 1-800-336-9719 (USA), or visit www.sigma-aldrich.com/safc.

III-V Material Sources—cont'd

N Sources

Ammonia

29,499-3

Purity: 99.99%

CAS No. 7664-41-7

MF: NH₃

FW: 17.03

mp: -78 °C

bp: -33 °C

Density: 0.6 g/mL

Azidotrimethylsilane

15,507-1

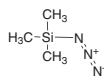
Purity: 95%

CAS No. 4648-54-8

FW: 115.21

bp: 52-53 °C @ 175 mmHg

Density: 0.868 g/mL



N,N-dimethylhydrazine

40,680-F

CAS No. 57-14-7

FW: 60.10

bp: 60-62 °C

Density: 0.791 g/mL



P,As and Sb Sources

tert-Butylphosphine

49,547-6

Purity: 99.999%

CAS No. 2501-94-2

FW: 90.11

bp: 54-55 °C

Density: 0.738g/mL



Tris(trimethylsilyl)phosphine

33,367-0

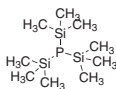
Purity: 95%

CAS No. 15573-38-3

FW: 250.55

bp: 243-244 °C

Density: 0.863 g/mL



Trimethyl Antimony

J1,0003-3

CAS No. 594-10-5

FW: 166.86



Tris(dimethylamido)antimony

55,397-2

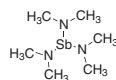
Purity: 99.99%

CAS No. 7289-92-1

FW: 253.98

bp: 32-34 °C @ 0.45 mmHg

Density: 1.325 g/mL



Trimethylarsine

55,131-7

Purity: 99.999%

CAS No. 593-88-4

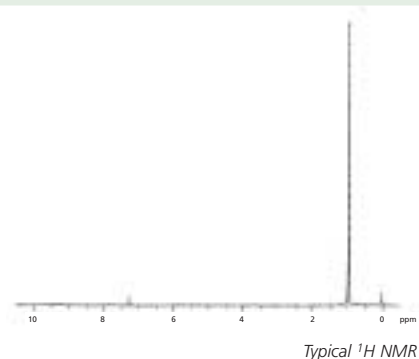
MF: (CH₃)₃As

FW: 120.03

bp: 51 °C

Density: 1.133 g/mL

Chloride: <100ppm



Typical ICP-MS Results for 55,131-7 in ppb

Ag	<1	Fe	<1	Nb	<1	Sm	<1
Al	1	Ga	<1	Nd	<1	Sn	<1
Au	3	Gd	<1	Ni	<1	Sr	<1
B	2	Ge	<1	Os	<1	Ta	<1
Ba	<1	Hf	<1	P	<1	Tb	<1
Be	<1	Hg	<1	Pb	<1	Te	<1
Bi	<1	Ho	<1	Pd	<1	Th	<1
Ca	<1	In	<1	Pr	<1	Ti	<1
Cd	<1	Ir	<1	Pt	<1	Tl	<1
Ce	<1	K	<1	Rb	<1	Tm	<1
Co	<1	La	<1	Re	<1	U	<1
Cr	<1	Li	<1	Rh	<1	V	<1
Cs	<1	Lu	<1	Ru	<1	W	<1
Cu	4	Mg	<1	Sb	3	Y	<1
Dy	<1	Mn	<1	Sc	<1	Yb	<1
Er	<1	Mo	<1	Se	<1	Zn	1
Eu	<1	Na	2	Si	<1	Zr	<1

Precursors for Metallization and Barrier Materials*

Dicarbonyl(cyclopentadienyl)cobalt(I)

24,525-9

CAS No. 12078-25-0

FW: 180.05

bp: 139–140 °C @ 710 mmHg



Bis(pentamethylcyclopentadienyl)magnesium

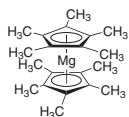
51,254-0

Purity: 99.999%

CAS No. 74507-64-5

FW: 294.77

mp: 230 °C (D)



Niobium(V) ethoxide

J1,0003-6

Purity: 99.999%

CAS No. 3236-82-6

MF: (C₂H₅O)₅Nb

FW: 318.21

Bis(methylcyclopentadienyl)nickel(II)

51,047-5

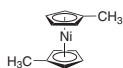
Purity: 97%

CAS No. 1293-95-4

FW: 216.96

mp: 34–36 °C

bp: 85–90 °C @ 1 mmHg



Bis(ethylcyclopentadienyl)ruthenium(II)

64,866-3

CAS No. 32992-96-4

FW: 235.29

mp: 6 °C

bp: 100 °C @ 0.01 mmHg



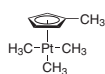
Cyclopentadienyl(trimethyl)platinum(IV)

64,560-5

CAS No. 94442-22-5

FW: 319.31

mp: 30–31 °C



Tantalum(V) ethoxide

51,855-7

Purity: 99.999%

CAS No. 6074-84-6

MF: (C₂H₅O)₅Ta

FW: 406.26

mp: 21 °C

bp: 155 °C @ 0.01 mmHg

Density: 1.566 g/mL

Pentakis(dimethylamido)tantalum(V)

49,686-3

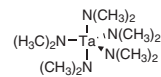
Purity: 99.99%

CAS No. 19824-59-0

FW: 401.33

mp: 100 °C (D)

Density: 1.252 g/mL



Tris(diethylamido)(tert-butylimido)tantalum(V)

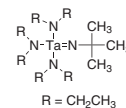
52,128-0

Purity: 99.99%

CAS No. 169896-41-7

FW: 468.46

Density: 1.252 g/mL



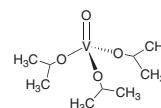
Vanadium(V) oxytriisopropoxide

40,492-6

CAS No. 5588-84-1

FW: 244.21

bp: 80–82 °C @ 2 mmHg



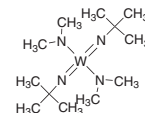
Bis(tert-butylimido)bis(dimethylamido)tungsten(VI)

57,793-6

CAS No. 406462-43-9

MF: C₁₂H₃₀N₄W

FW: 414.23



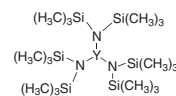
Tris(N,N-bis(trimethylsilyl)amide)yttrium(III)

52,451-4

CAS No. 41836-28-6

FW: 570.07

mp: 161–166 °C



Yttrium(III) butoxide solution 0.5M in toluene

51,066-1

CAS No. 111941-71-0

MF: (C₄H₉O)₃Y

FW: 308.25

bp: 109 °C

Diethylzinc

25,678-1

CAS No. 557-20-0

MF: (C₂H₅)₂Zn

FW: 123.49

mp: -28 °C

bp: 117 °C

Density: 1.205 g/mL

*For titanium precursors see page 3.



Ready to scale up? For competitive quotes on larger quantities or custom synthesis, contact Sigma-Aldrich Fine Chemicals at 1-800-336-9719 (USA), or visit www.sigma-aldrich.com/safc.

Order: 1.800.558.9160 Technical Service: 1.800.231.8327

ALDRICH

High Purity Deposition Gases

High Purity CVD Gases are used in the deposition of amorphous and epitaxial silicon, SiGe alloys, and dielectrics as well as the manufacture of integrated circuits and photovoltaics. The gases are also used as dopants in CVD and ion-implantation processes. Deuterated gases have been shown to improve device performance and reliability. ISOTECH™, a key member of the Sigma-Aldrich family, produces our stable

isotope product lines. Led by an impressive group of expert scientists, ISOTECH™ routinely engages in multiple step syntheses of complex labeled molecules including the stable isotope labeling patterns ¹³C, ¹⁵N, ¹⁷O, ¹⁸O, ²H. ISOTECH™ has a large inventory of stable isotope labeled starting material to accommodate any researcher's need. Contact ISOTECH™ at www.sigma-aldrich.com/isotec.

Catalog Number	Product Name	Formula	Purity (%)	bp °C	Vapor density (vs. Air)	Critical Temp °C	Water	H ₂	N ₂	O ₂	CH ₄	Other
							(ppm)					
46,305-1	Diborane (10% in H ₂)	B ₂ H ₆	99.99*	-92.5	–	–	–	–	–	–	–	
46,313-2	Diborane-d ₆ (10% in D ₂)	B ₂ D ₆	99.99*	-92.5	–	–	–	–	–	–	–	
46,314-0	Diborane-d ₆ (10% in He)	B ₂ D ₆	99.99*	-92.5	–	–	–	–	–	–	–	
46,307-8	Digermene (10% in H ₂)	Ge ₂ H ₆	99.99	31.5	–	150.9	–	–	–	–	–	
46,304-3	Disilane	Si ₂ H ₆	99.998	-14.5	–	150.9	<1	–	<1	<1	–	<50 higher silanes
46,302-7	Germane	GeH ₄	99.997	-88.4	1.53 (-142 °C)	35	<1	<50	<2	<0.5	–	<20 digermene
46,299-3	Methylsilane	CH ₃ SiH ₃	99.9	-57	0.628 (-58 °C)		–	–	<10	<2	<50	<100 higher silanes
29,564-7	Phosphine	PH ₃	99.9995	-87.5	1.15	51.3	<1	–	<1	<1	<1	<1 AsH ₃
46,312-4	Phosphine-d ₃	PD ₃	98 atom %	-87.5	–	–	–	–	–	–	–	
33,389-1	Silane	SiH ₄	99.998	-112	1.1	-3.4	<1	<100	<2	–	–	<5 higher silanes
46,311-6	Silane-d ₄	SiD ₄	98 atom%	-112	–	–	–	–	–	–	–	
39,635-4	Tetramethylgermanium	(CH ₃) ₄ Ge	98%	43.5	0.978	–	–	–	–	–	–	
46,309-4	Trimethylboron	(CH ₃) ₃ B	98.35	-20.2	2.3	–	–	<50	<10	<1	<100	<1000 N (CH ₃) ₃
46,342-6	Trimethylboron-d ₉	(CD ₃) ₃ B	99 atom %	-20.2	2.3	–	–	–	–	–	–	

*As diborane

Custom Packaging



Sigma-Aldrich has a variety of packaging options to ensure the quality and safe handling of all of our volatile precursors. Most of our products come in standard packaging depending on size, quantity and material state. Custom packing is optional for all of our products. If a product catalog number starts with a "J", then packaging configuration is up to you at the time you order.

Borosilicate or Quartz Ampoules

For handling small quantities in a dry-box. Shipped in a clam-shell.

Sure/Seal™ Bottles

The standard container for air sensitive materials.

Gas Lecture Bottles

Electronic grade gases packed in clean room conditions.

Sure/Pac™ Cylinder

Equipped with brass ball valve and treated plug.

Stainless Steel Bubblers

Standard and custom sizes for your deposition systems.

TO ORDER: Contact your local Sigma-Aldrich office (see back cover), call 1-800-558-9160 (USA), or visit sigma-aldrich.com.

Stainless Steel Bubblers

For Deposition Precursors

Sigma-Aldrich offers stainless steel bubblers designed to deliver our ultra high-purity organometallic precursors to the reaction zone during the fabrication process. All of our volatile precursors can be prepackaged in these bubblers to ensure you are getting the highest quality material available.

In addition to our standard sizes, Sigma-Aldrich can custom build bubblers to your equipment specifications. Please inquire about custom bubbler configurations by using the drawing below, calculate your dimensions then contact Sigma-Aldrich at matsci@sial.com.

Stock Stainless Steel Bubblers

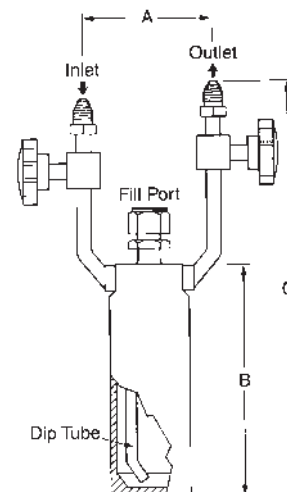
- Vertical, dual-valve bubblers equipped with a 1/2 in. fill port and 1/4 in. VCR™ inlet and outlet valves.
- Electropolished 316L stainless steel that is free of mechanical defects and foreign chemical contaminants.
- Sealed in a clean-room environment to ensure product quality and integrity.
- Pneumatically tested to ensure leak-free service.
- Available from stock in the sizes listed below.

Size	Dimensions (A x B x C) in.	Cat. No.
300 mL	3.25 x 7.36 x 12.81	Z52,706-8
1.2 L	2.77 x 8.00 x 10.28	Z55,336-0

Custom Stainless Steel Bubblers

- Full range of sizes available.
- Specify the dimensional requirements (A, B, C) as shown in the diagram.
- Select inlet and outlet valve connections.

VCR is a trademark of Swagelok Co.



Aldrich 3-Stage Oil Diffusion Pump

Aldrich 3-stage oil diffusion pumps provide fast 375 L/s pumping speed that outperforms traditional glass diffusion pumps. They feature all glass construction with a removable jet assembly for easy cleaning. 3-stage oil diffusion pumps safely reduce sublimation and distill action temperature and are chemically resistant to acidic solvents. Sigma-Aldrich also offers all the necessary equipment to run and maintain your pump.

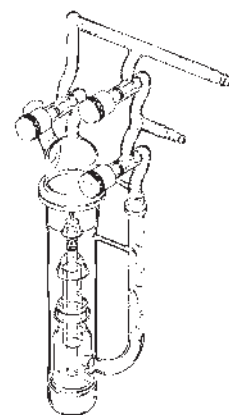
Model	Overall dim. (in.)	Pumping speed (L/s)	Max. vacuum (torr)	Catalog Number
With Gauge Port	32 x 12	375	10	Z54,478-7
Without Gauge Port	32 x 12	375	10	Z55,537-1

Equipment Required for Diffusion Pump Operation

Description

Dow Corning Diffusion Pump Oil (500 mL)	Type 704* 44,597-5	Type 705** 44,598-3
Leybold Trivac B, D8B Mechanical Vac. Pump, 197 Lpm speed	115 V Z28,459-9	220 V Z28,282-0
Glas-Col 500 mL Heating Mantle For heating mantle	115 V Z28,485-8	230 V Z28,514-5
Digitrol II Heat Controller	115 V Z28,549-8	240 V Z28,550-1
Corning 5x7" Stir Plate	115 V Z26,206-4	230 V Z26,207-2
Dewar flask, 10 L capacity	Z10,565-1	
Safety shield	Z23,140-1	
Vacuum Trap 1 L bulb capacity	F 9776	

*For Pressures 10⁻⁵ to 10⁻⁶ torr **For Pressures 10⁻⁹ to 10⁻¹⁰ torr



Ready to scale up? For competitive quotes on larger quantities or custom synthesis, contact Sigma-Aldrich Fine Chemicals at 1-800-336-9719 (USA), or visit www.sigma-aldrich.com/safc.

Deposition Substrates and Silicon Wafers

Successful deposition of materials often depends on the proper substrates, with a good lattice match and a clean, even surface. Aldrich is proud to offer a variety of new single crystal substrates and silicon wafers in multiple orientations to meet your deposition needs. All these materials undergo

rigorous analytical testing including trace metal analysis and x-ray diffraction in order to insure the highest purity standards. If there is an orientation or material you cannot find please contact us at matsci@sial.com.

Single Crystal Substrates (10 x 10 x 0.5 mm, single side polished)

Composition	Formula	mp (°C)	Unit Cell (Å)	Orientation							
				Hex (0001)	Cubic (100)	Cubic (110)	Cubic (111)	Rhomb (100) ¹	Tetrag (100)	Tetrag (001)	Tetrag (110)
Aluminum oxide	Al ₂ O ₃	2040	a=4.758 c=12.992	63,487-5	–	–	–	–	–	–	–
Lanthanum aluminum oxide	LaAlO ₃	2080	a=5.357 c=13.22	–	–	–	–	63,473-5	–	–	–
LSAT Lanthanum aluminum oxide-strontium aluminum tantalum oxide	(LaAlO ₃) _{0.3} (Sr ₂ AlTaO ₆) _{0.7}	1840	a=3.868	–	63,505-0	–	–	–	–	–	–
Magnesium aluminate	MgAl ₂ O ₄	2130	a=8.083	–	63,507-3	63,484-0	63,483-2	–	–	–	–
Magnesium oxide	MgO	2852	a=4.216	–	63,464-6	63,470-0	63,469-7	–	–	–	–
Silicon dioxide, optical grade	SiO ₂	1610	a=4.914 c=5.405	63,486-7	–	–	–	–	–	–	–
Strontium lanthanum aluminate	SrLaAlO ₄	1650	a=3.756 c=12.63	–	–	–	–	–	63,511-1	63,489-1	–
Strontium titanate	SrTiO ₃	2080	a=3.905	–	63,468-9	63,467-0	63,816-1	–	–	–	–
Titanium(IV) oxide rutile	TiO ₂	1840	a=4.5936 c=2.9582	–	–	–	–	–	63,504-9	63,505-7	63,506-5
Yttrium vanadate	YVO ₄	1825	a=7.12 c=6.29	–	–	–	–	–	–	–	63,510-3

1) Cubic @>435 °C, a=3.821 Å

Silicon Wafers (0.5 mm thick, single side polished)

Catalog Number	Orientation	Radius (in.)	Semiconductor Type	Dopant	Resistivity (Ω cm)
64,668-7	(100)	2	–	none	10 ² –10 ⁴
64,710-1	(111)	2	–	none	10 ² –10 ⁴
64,753-5	(100)	3	–	none	10 ² –10 ⁴
64,754-3	(111)	3	–	none	10 ² –10 ⁴
64,767-5	(100)	2	P-type	boron	10 ³ –40
64,770-5	(111)	2	P-type	boron	10 ³ –40
64,776-4	(100)	3	P-type	boron	10 ³ –40
64,777-2	(111)	3	P-type	boron	10 ³ –40
64,778-0	(100)	2	N-type	phosphorous	10 ³ –40
64,779-9	(111)	2	N-type	phosphorous	10 ³ –40
64,780-2	(100)	3	N-type	phosphorous	10 ³ –40
64,781-0	(111)	3	N-type	phosphorous	10 ³ –40

Silicon wafers have a density of 2.33 g/mL; unit cell: cubic, a = 5.4037 Å

Gold-Coated Substrates

We now offer gold-coated substrates with 100 to 2000 Å gold coatings on: Glass microscope slides, round and square

coverslips, silicon wafers, or mica. Please visit us at www.sigma-aldrich.com for more information.

TO ORDER: Contact your local Sigma-Aldrich office (see back cover), call 1-800-558-9160 (USA), or visit sigma-aldrich.com.

Books and Software

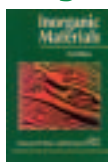
CVD of Nonmetals

Z28,630-3

W.S. Rees, Jr., Ed., John Wiley & Sons, 1992, pp. 392, Hardcover.

The basics necessary for any CVD process are discussed. Precursor requirements, with an emphasis on materials chemistry, common structures of reactants and substrates, as well as reaction control are discussed for a broad range of compositions. Technological issues, such as reactor geometries and operation parameters, are assessed and the viability of the method is compared with other techniques for the preparation of thin films.

Inorganic Materials



Z40,867-0

D.W. Bruce and D. O'Hare, Eds., John Wiley & Sons, 1997, pp. 607, Softcover.

This new edition of the acclaimed Inorganic Materials reflects the expansion in both knowledge and interest that materials chemistry has experienced in recent years. It provides a well-referenced introduction to the most exciting aspects of materials chemistry, followed by an overview with examples chosen to best emphasize the materials in question.

Characterization of Materials, 2 volume set



Z55,373-5

Elton N. Kaufmann, John Wiley & Sons, 2003, pp. 1392, Hardcover.

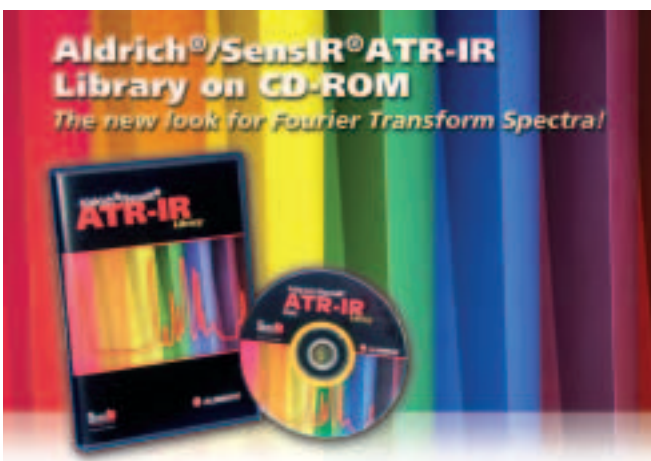
Provides comprehensive coverage of materials characterization techniques including computational and theoretical methods, crystallography, mechanical testing, thermal analysis, optical imaging and spectroscopy.

Molecular Nanoelectronics

Z55,011-6

Mark Reed and Tahkee Lee, American Scientific Publishers, 2003, pp. 400, Hardcover.

This book covers topics on transport theory through nanostructures, processing, self-assembly, device fabrication and architecture for nanoelectronic device applications. Contains about 2,000 bibliographic citations and hundreds of illustrations, figures, tables, chemical structures and equations.



Diamond ATR spectra without Nujol or KBr interference.

The Aldrich®/SensIR® ATR-IR Library (Attenuated Total Reflectance) contains high resolution FT-IR spectra for use with Thermo Electron SpecID® (Z40,748-8) spectral search software. Suitable for doing searches using spectral subtraction.

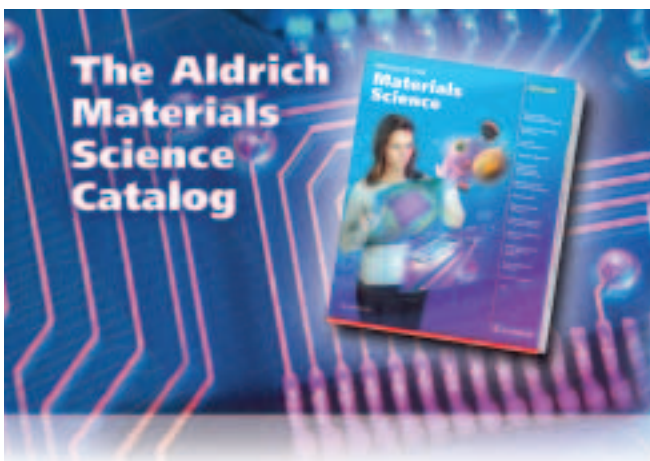
- Contains nearly 18,000 pure compounds of the most common chemicals found in the Aldrich Handbook
- Spectra archived using 2cm⁻¹ data resolution and 16-bit ordinate precision
- Searchable by ATR and standard FT-IR spectra*
- Frequency range: 650-4000cm⁻¹

Z54,751-4

Multi-unit Library discounts available.

Please inquire by phone at **800-231-8327** or email us at aldrich@sial.com.

The Library includes a wide variety of chemical products.



This "Materials by Application" Catalog Contains:

- Biocompatible/Biodegradable Materials
- Display & Photovoltaic Materials
- Fuel Cell/Battery Materials
- Magentic Materials
- Materials for Thin films/Microelectronics
- Materials for Lithography/Nanopatterning
- Nanomaterials
- Photonic/Optical Materials
- Structural materials/Electronic Ceramics
- Additional Electronic Grade Solvent/Reagents
- Reference/Calibration Standards
- Books, Equipment, and Software for Materials Scientists

Request your **FREE** copy of the Aldrich Materials Science catalog by visiting us at www.sigma-aldrich.com/matsci.



Ready to scale up? For competitive quotes on larger quantities or custom synthesis, contact Sigma-Aldrich Fine Chemicals at 1-800-336-9719 (USA), or visit www.sigma-aldrich.com/safc.

Argentina
SIGMA-ALDRICH DE ARGENTINA, S.A.
Tel: 54 11 4556 1472
Fax: 54 11 4552 1698

Australia
SIGMA-ALDRICH PTY., LIMITED
Free Tel: 1800 800 097
Free Fax: 1800 800 096
Tel: 612 9841 0555
Fax: 612 9841 0500

Austria
SIGMA-ALDRICH HANDELS GmbH
Tel: 43 1 605 81 10
Fax: 43 1 605 81 20

Belgium
SIGMA-ALDRICH NV/SA.
Free Tel: 0800-14747
Free Fax: 0800-14745
Tel: 03 899 13 01
Fax: 03 899 13 11

Brazil
SIGMA-ALDRICH BRASIL LTDA.
Tel: 55 11 3732 3100
Fax: 55 11 3733-5151

Canada
SIGMA-ALDRICH CANADA LTD.
Free Tel: 800-565-1400
Free Fax: 800-265-3858
Tel: 905-829-9500
Fax: 905-829-9292

China
SIGMA-ALDRICH CHINA INC.
Tel: 86-21-6386 2766
Fax: 86-21-6386 3966

Czech Republic
SIGMA-ALDRICH s.r.o.
Tel: 246 003 200
Fax: 246 003 291

Denmark
SIGMA-ALDRICH DENMARK A/S
Tel: 43 56 59 10
Fax: 43 56 59 05

Finland
SIGMA-ALDRICH FINLAND
Tel: 358-9-350-92 50
Fax: 358-9-350-92 555

France
SIGMA-ALDRICH CHIMIE S.à.r.l.
Tel appel gratuit: 0800 211 408
Fax appel gratuit: 0800 031 052

Germany
SIGMA-ALDRICH CHEMIE GmbH
Free Tel: 0800-51 55 000
Free Fax: 0800-649 00 00

Greece
SIGMA-ALDRICH (O.M.) LTD
Tel: 30 210 9948010
Fax: 30 210 9943831

Hungary
SIGMA-ALDRICH Kft
Tel: 06-1-235-9054
Fax: 06-1-269-6470
Ingyenes zöld telefon: 06-80-355-355
Ingyenes zöld fax: 06-80-344-344

India
SIGMA-ALDRICH CHEMICALS
PRIVATE LIMITED
Telephone
Bangalore: 91-80-5112-7272
Hyderabad:
91-40-5531 5548 / 2784 2378
Mumbai: 91-22-2579 7588 / 2570 2364
New Delhi:
91-11-2616 5477 / 2619 5360
Fax
Bangalore: 91-80-5112-7473
Hyderabad: 91-40-5531 5466
Mumbai: 91-22-2579 7589
New Delhi: 91-11-2616 5611

Ireland
SIGMA-ALDRICH IRELAND LTD.
Free Tel: 1800 200 888
Free Fax: 1800 600 222

Israel
SIGMA-ALDRICH ISRAEL LTD.
Tel: 08-948-4100
Fax: 08-948-4200

Italy
SIGMA-ALDRICH S.r.l.
Telefono: 02 33417310
Fax: 02 38010737
Numero Verde: 800-827018

Japan
SIGMA-ALDRICH JAPAN K.K.
Tokyo Tel: 03 5796 7300
Tokyo Fax: 03 5796 7315

Korea
SIGMA-ALDRICH KOREA
Tel: 031-329-9000
Fax: 031-329-9090

Malaysia
SIGMA-ALDRICH (M) SDN. BHD
Tel: 603-56353321
Fax: 603-56354116

Mexico
SIGMA-ALDRICH QUÍMICA, S.A. de C.V.
Free Tel: 01-800-007-5300
Free Fax: 01-800-712-9920

The Netherlands
SIGMA-ALDRICH CHEMIE BV
Tel Gratis: 0800-0229088
Fax Gratis: 0800-0229089
Tel: 078-6205411
Fax: 078-6205421

New Zealand
SIGMA-ALDRICH PTY., LIMITED
Free Tel: 0800 936 666
Free Fax: 0800 937 777

Norway
SIGMA-ALDRICH NORWAY AS
Tel: 23 17 60 60
Fax: 23 17 60 50

Poland
SIGMA-ALDRICH Sp. z o.o.
Tel: (+61) 829 01 00
Fax: (+61) 829 01 20

Portugal
SIGMA-ALDRICH QUÍMICA, S.A.
Free Tel: 800 20 21 80
Free Fax: 800 20 21 78

Russia
SIGMA-ALDRICH RUSSIA
OOO SAF-LAB
Tel: 095 975-1917/3321
Fax: 095 975-4792

Singapore
SIGMA-ALDRICH PTE. LTD.
Tel: 65-6271 1089
Fax: 65-6271 1571

South Africa
SIGMA-ALDRICH
SOUTH AFRICA (PTY) LTD.
Tel: 27 11 979 1188
Fax: 27 11 979 1119

Spain
SIGMA-ALDRICH QUÍMICA S.A.
Free Tel: 900101376
Free Fax: 900102028

Sweden
SIGMA-ALDRICH SWEDEN AB
Tel: 020-350510
Fax: 020-352522
Outside Sweden Tel: +46 8 7424200
Outside Sweden Fax: +46 8 7424243

Switzerland
FLUKA CHEMIE GmbH
Swiss Free Call: 0800 80 00 80
Tel: +41 81 755 2828
Fax: +41 81 755 2815

United Kingdom
SIGMA-ALDRICH COMPANY LTD.
Free Tel: 0800 717181
Free Fax: 0800 378785
Tel: 01747 833000
Fax: 01747 833313

United States
SIGMA-ALDRICH
P.O. Box 14508
St. Louis, Missouri 63178
Toll-free: 800-325-3010
Call Collect: 314-771-5750
Toll-Free Fax: 800-325-5052
Tel: 314-771-5765
Fax: 314-771-5757

Internet:
sigma-aldrich.com



P.O. Box 355
Milwaukee, WI 53201
USA

Return Service Requested

sigma-aldrich.com

Order/Customer Service 1-800-325-3010 • Fax 1-800-325-5052
Technical Service 1-800-325-5832 • sigma-aldrich.com/techservice

Development/Bulk Manufacturing Inquiries Sigma-Aldrich Fine Chemicals 1-800-336-9719

World Headquarters • 3050 Spruce St., St. Louis, MO 63103 • (314) 771-5765

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



GUN
01217-040520
0164

©2004 Sigma-Aldrich Co. Printed in USA Sigma 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. SIGMA and are registered trademarks of Sigma-Aldrich Co. and its division Sigma-Aldrich Biotechnology LP. Riedel-de Haën®: trademark under license from Riedel-de Haën GmbH.