

## WACKER—Company Profile

WACKER is a globally operating chemical company headquartered in Munich, Germany. Its portfolio focuses on semiconductor technology, silicone chemistry, polymers and fine chemicals, as well as polysilicon. In line with these core areas of expertise, WACKER organized its activities into a chemical and a semiconductor segment. The former comprises WACKER SILICONES, WACKER POLYMERS, WACKER FINE CHEMICALS and WACKER POLYSILICON; the latter comprises Siltronic.



The WACKER Group posted sales of around €2.5 billion in 2004. Germany accounted for 23 percent of these sales, Europe (excluding Germany) for 33 percent, and the Americas and Asia Pacific (plus the rest of the world) for 22 percent each. WACKER employs roughly 14,700 people at 20 production sites in Europe, the Americas and Asia, and at over 100 sales offices around the globe. Founded in 1914 by Dr. Alexander Wacker, the WACKER Group is 100 percent controlled by the Wacker family.

WACKER's business strategy is geared to long-term profitable growth and, in turn, to securing a sustained increase in corporate value. The Group's portfolio strategy is designed to ensure a well-balanced mix of mature, profitable sectors and newer, high-growth activities and also to reinforce WACKER's core expertise in the chemical sector.

Fourteen percent of WACKER's sales revenue is earmarked for investment, with six percent of this sum being spent on R&D. Corporate strategy is geared to enhancing the Group's global presence and strengthening its position as a technology leader and innovation partner for its customers. Thus, WACKER ranks among the world's most research-intensive chemical companies. WACKER's corporate research facility, the "Consortium für elektrochemische Industrie," is an indispensable part of the continuous innovation process.

At WACKER, health, safety, and the environment are top priorities. Every year, the Group invests over €300 million in the safety and environmental compatibility of its products and plants, in employee health and safety, as well as in basic and advanced training. A substantial reduction in emissions and resource consumption was documented since 2000, despite rising production volumes. WACKER's 2003 accident rate of less than five per million working hours was well below the German chemical industry's average.

Members of the Group Executive Board are Dr. Peter-Alexander Wacker (President & CEO), Dr. Rudolf Staudigl, and Dr. Joachim Rauhut.

## Sigma-Aldrich Worldwide Locations

### 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: +420 246 003 200  
Fax: +420 246 003 291

### Denmark

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

### Finland

SIGMA-ALDRICH FINLAND  
Tel: (09) 350 9250  
Fax: (09) 350 92555

### 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  
New Delhi: 91-11-5165 4255  
Mumbai: 91-22-2570 2364  
Hyderabad: 91-40-5584 5488  
Fax  
Bangalore: 91-80-5112-7473  
New Delhi: 91-11-5165 4266  
Mumbai: 91-22-2579 7589  
Hyderabad: 91-40-5584 5466

### Ireland

SIGMA-ALDRICH IRELAND LTD.  
Free Tel: 1800 200 888  
Free Fax: 1800 600 222  
Tel: 353 1 4041900  
Fax: 353 1 4041910

### Israel

SIGMA-ALDRICH ISRAEL LTD.  
Free Tel: 1-800-70-2222  
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  
Tel: 61 2 9841 0500  
Fax: 61 2 9841 0500

### Norway

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

### Poland

SIGMA-ALDRICH Sp. z o.o.  
Tel: 061 829 01 00  
Fax: 061 829 01 20

### Portugal

SIGMA-ALDRICH QUÍMICA, S.A.  
Free Tel: 800 202180  
Free Fax: 800 202178  
Tel: 21 9242555  
Fax: 21 9242610

### Russia

SIGMA-ALDRICH RUS, LLC  
Tel: +7 (095) 975-1917/3321  
Fax: +7 (095) 975-4792

### Singapore

SIGMA-ALDRICH PTE. LTD.  
Tel: 65-67791200  
Fax: 65-67791822

### South Africa

SIGMA-ALDRICH SOUTH AFRICA (PTY) LTD.  
Free Tel: 0800 1100 75  
Free Fax: 0800 1100 79  
Tel: 27 11 979 1188  
Fax: 27 11 979 1119

### Spain

SIGMA-ALDRICH QUÍMICA S.A.  
Free Tel: 900 101376  
Free Fax: 900 102028  
Tel: 91 661 99 77  
Fax: 91 661 96 42

### Sweden

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

### Switzerland

SIGMA-ALDRICH CHEMIE LLC  
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  
SAFC (UK): 01202 712305

### 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

 **ChemFiles**

Chemical Synthesis  
**Sigma-Aldrich Minibulk Concept**

Vol. 5  
Supplement

**WACKER**



Development/Bulk Manufacturing Inquiries **SAFC**  
World Headquarters • 3050 Spruce St., St. Louis, MO 63103



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-de Haën  SUPELCO  ISOTEC

©2005 Sigma-Aldrich Co. All rights reserved.  
SIGMA, SAFC, SAFC, SIGMA-ALDRICH, ISOTEC, ALDRICH, FLUKA, and SUPELCO are trademarks belonging to Sigma-Aldrich Co. and its affiliate Sigma-Aldrich Biotechnology LP. Riedel-de Haën, trademark under license from Riedel-de Haën GmbH. Sigma 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. Synthesis Power is a registered trademark of WACKER FINE CHEMICALS.



To order: Contact your local Sigma-Aldrich office (see back cover),  
or visit [sigma-aldrich.com](http://sigma-aldrich.com)

IJX

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



SIGMA-ALDRICH

## The Sigma-Aldrich Minibulk Concept

Sigma-Aldrich is the exclusive worldwide distribution partner for scale-up laboratory quantities of WACKER FINE CHEMICALS. Selected products are available in the following package sizes:

- gram samples
- returnable containers
- original WACKER drums
- multidrum pallets

Larger quantities for customers are supplied directly from WACKER. This partnership allows our customers the seamless availability of products from development to manufacturing scale, eliminating the need to qualify a new source as the project moves along the development pathway.

For each WACKER product, a Sigma-Aldrich catalog number has been created with the note "WACKER quality." The current portfolio consists of more than 40 organic intermediates and 20 organosilanes.

## WACKER FINE CHEMICALS—A Partner for the Life Sciences and Consumer Care Industries

As an expert in organic synthesis, silane chemistry, and biotechnology, WACKER FINE CHEMICALS provides tailored solutions for its customers in the life sciences and consumer care industries.

WACKER FINE CHEMICALS is a preferred partner for challenging chemical and biotech custom manufacturing projects that serve as the basis for novel pharmaceutical active ingredients. Our TRIPLE SYNTHESIS POWER®, based on the unique combination of three technology platforms—organic synthesis, silane chemistry, and biotechnology—gives you access to a broad range of complex organic intermediates, innovative silanes, and chiral products and also enables us to develop cost-efficient, innovative routes for the most demanding custom synthesis projects.

The range of products further includes cyclodextrins, cysteines, and unnatural amino acids, as well as various chemical-starting materials. Major customers include the pharmaceutical, agrochemical, household goods, and personal care, food and textile industries.



WACKER FINE CHEMICALS has production sites in Germany and the United States. The business division generated sales of some €95 million in 2004 and employs a workforce of about 300.

Wacker-Chemie GmbH  
WACKER FINE CHEMICALS  
Hanns-Seidel-Platz 4  
81737 München, Germany  
Tel: +49 89 6279 1484  
Fax: +49 89 6279 2921  
info.finechemicals@wacker.com

**WACKER** FINE CHEMICALS

## Available WACKER Products

Cat. No.	Product Name	WACKER Abbrev.	Min. Assay	CAS No.	Mol. Weight	Sample	Minibulk	Original Drum	Pallet	WACKER Direct Supply
<b>Organic Intermediates</b>										
<b>Ethers</b>										
10917	Dibutyl ether	DBE	≥99.3% GC	142-96-1	130.23	1 kg	25 kg	155 kg	4*155 kg	>3 pallets
80103	Dibutyl ether CD-R grade	DBE CD-R Grade	≥99.3% GC	142-96-1	130.23	1 kg	25 kg	155 kg	4*155 kg	>3 pallets
<b>Aldehydes and Acetals</b>										
10891	3-Oxobutylaldehyde dimethylacetal	AADMA	≥95.0% GC	5436-21-5	132.16	250 g	25 kg	-	-	>1 drum
41585	Ketene dimethyl acetal	Ketendi-methylacetal	≥96.0% GC	922-69-0	88.11	100 g	5 kg	-	-	>1 drum
<b>Ketones</b>										
10916	Acetylacetone	ACAC	≥99.3% GC	123-54-6	100.12	1 kg	25 kg	185 kg	4*185 kg	>3 pallets
10890	2,5-Hexanedione	ACAN	≥97.0% GC	110-13-4	114.14	250 g	25 kg	-	-	>1 drum
<b>Esters</b>										
10914	Isopropenyl acetate	IPA	≥98.8% GC	108-22-5	100.12	1 kg	25 kg	170 kg	4*170 kg	>3 pallets
10913	Methyl acetoacetate	ACM	≥99.0% GC	105-45-3	116.12	1 kg	25 kg	220 kg	4*220 kg	>3 pallets
10912	Ethyl acetoacetate	ACE	≥99.0% GC	141-97-9	130.14	1 kg	25 kg	210 kg	4*210 kg	>3 pallets
10893	Allyl acetoacetate	ACALLYL	≥98.5% GC	1118-84-9	142.15	1 kg	25 kg	210 kg	-	>1 pallet
10869	2-Methoxyethyl acetoacetate	ACMDE	≥96.0% GC	22502-03-0	160.17	1 kg	25 kg	200 kg	-	>1 pallet
10911	tert-Butyl acetoacetate	ACTB	≥98.0% GC	1694-31-1	158.19	1 kg	25 kg	190 kg	-	>1 pallet
10902	Isobutyl acetoacetate	ACIB	≥96.0% GC	7779-75-1	158.19	1 kg	25 kg	210 kg	-	>1 pallet
10903	Isopropyl acetoacetate	ACIP	≥98.0% GC	542-08-5	144.17	1 kg	25 kg	200 kg	-	>1 pallet
10898	Benzyl acetoacetate	AC-benzyl	≥96.0% GC	5396-89-4	192.21	1 kg	25 kg	210 kg	-	>1 pallet
10899	Methyl 3-oxoheptanoate	VEM	≥96.0% GC	39815-78-6	158.19	100 g	5 kg	-	-	>1 drum
10910	Methyl 3-oxovalerate	PEM	≥98.5% GC	30414-53-0	130.14	250 g	25 kg	-	-	>1 drum
10900	Methyl 4-methyl-3-oxovalerate	IBEM	≥98.0% GC	42558-54-3	144.17	100 g	25 kg	-	-	>1 drum
10901	Methyl 3-oxohexanoate	BEM	≥98.0% GC	30414-54-1	144.17	250 g	25 kg	-	-	>1 drum
10892	Ethyl 3-(4-methoxyphenyl)-3-oxopropionate	MOBEE	≥97.0% GC	28821-83-6	222.24	250 g	25 kg	-	-	>1 drum
10896	Vinyl laurate	VL	≥98.0% GC	2146-71-6	226.36	1 kg	25 kg	167 kg	-	>1 pallet
<b>Nitrogen Compounds</b>										
10908	N,N-Diethylacet-oacetamide	DMAA	≥97.0% GC	2235-46-3	157.21	1 kg	50 kg	205 kg	-	>1 pallet
10909	N,N-Dimethyl-aceto-acetamide solution	DEAA	80% in water	2044-64-6	129.16	1 kg	50 kg	220 kg	-	>1 pallet
<b>Halogenated Compounds</b>										
10907	Chloroacetaldehyde solution	CLACH	45% in water	107-20-0	78.50	1 kg	25 kg	230 kg	-	>1 pallet
10906	Chloroacetaldehyde dimethyl acetal	CADMA	≥98.0% GC	97-97-2	124.57	1 kg	25 kg	210 kg	-	>1 pallet
11927	2-Chloropropionaldehyde dimethyl acetal	CL-PADMA	≥96.0% GC	87894-32-4	138.59	250 g	25 kg	-	-	>1 drum
10905	Chloroacetone	MCA	≥96.0% GC	78-95-5	92.52	1 kg	25 kg	225 kg	-	>1 pallet
10877	1,1-Dichloro-2-propanone	1,1-DCA	≥96.0% GC	513-88-2	126.97	1 kg	25 kg	-	-	>1 drum
10878	1,3-Dichloroacetone solution	1,3-DCA	50% in acetone	534-07-6	126.97	250 g	25 kg	-	-	>1 drum
10879	1,1,3-Trichloroacetone	TCA	≥86.5% GC	921-03-9	161.41	1 kg	25 kg	-	-	>1 drum
10894	Hexachloro-2-propanone	HCA	≥99.0% GC	116-16-5	264.75	1 kg	25 kg	330 kg	-	>1 pallet
10904	3-Chloro-2-butanone	Chlor-MEK	≥96.0% GC	4091-39-8	106.55	1 kg	25 kg	-	-	>1 drum
10872	3-Chloro-2,4-pentanedione	3-CL-ACAC	≥97.0% GC	1694-29-7	134.56	250 g	25 kg	-	-	>1 drum
10895	Ethyl 2-chloroacetoacetate	2-CL-ACE	≥96.0% GC	609-15-4	164.59	1 kg	25 kg	210 kg	-	>1 pallet
10876	Methyl 2-chloro-3-oxovalerate	2-CL-PEM	≥93.0% GC	114192-09-5	164.59	250 g	25 kg	-	-	>1 drum
10882	N,N-Dimethyl-2-chloroacetoacetamide	DMCAA	≥95.0% GC	5810-11-7	163.60	1 kg	25 kg	-	-	>1 drum

Cat. No.	Product Name	WACKER Abbrev.	Min. Assay	CAS No.	Mol. Weight	Sample	Minibulk	Original Drum	Pallet	WACKER Direct Supply
<b>Sulfur Compounds</b>										
10883	1,4-Dithiane-2,5-diol	Dithian	≥96.0% GC	40018-26-6	152.24	100 g	5 kg	40 kg	-	>1 pallet
<b>5-Ring Heterocycles</b>										
05091	3,5-Dimethylpyrazole	3,5-DMP	≥99.0% GC	67-51-6	96.13	100 g	1 kg; 25 kg	100 kg	-	>1 pallet
79705	3,5-Dimethylisoxazole	Dimexazol	≥98.0% GC	300-87-8	97.12	100 g	25 kg	-	-	>1 drum
<b>Silanes</b>										
<b>Silylating Agents</b>										
95541	Chlorotrimethylsilane	Silan M3	≥99.0% GC	75-77-4	108.64	500 g	5 kg; 25 kg	170 kg	-	>1 pallet
80430	Dichlorodimethylsilane	Silan M2	≥99.6% GC	75-78-5	129.06	500 g	5 kg; 25 kg	-	-	>1 drum
75986	Chlorotriethylsilane	Silan E3	≥99.0% GC	994-30-9	150.72	500 g	5 kg	-	-	>1 can
90174	Chlorotrisopropylsilane	Silan IP3	≥97.0% GC	13154-24-0	192.80	500 g	5 kg	-	-	>1 can
86944	Hexamethyldisilazane	Silazan HMN	≥97.0% GC	999-97-3	161.39	500 g	5 kg; 20 kg	-	-	>1 can
07048	Triisopropylsilyltrifluoromethane sulfonate	Silan IP3-Triflat	≥97.0%	80522-42-5	306.42	250 g	1 kg	-	-	>1 can
29383	Trimethylsilyl trifluoromethane sulfonate	Silan TMS-Triflat	≥97.0%	27607-77-8	222.26	250 g	1 kg	-	-	>1 can
76016	Dichloro-methyl-octadecylsilane	Silan ODM	≥92.0% GC	5157-75-5	367.51	500 g	5 kg	-	-	>1 can
71034	Chloro-dimethyl-octadecylsilane solution	Silan ODM-tol. Lsg.	70% in toluene	18643-08-8	347.09	500 g	5 kg	-	-	>1 drum
94394	Dimethoxy-methyl-octadecylsilane	Silan ODM-	≥92.0%	70851-50-2	358.67	500 g	5 kg	-	-	>1 can
80631	Sodium bis(trimethylsilyl) amide solution	HMN-Na-Salzlg. 40%/THF	40% in THF	1070-89-9	183.37	500 g	5 kg	-	-	>1 can
<b>Reducing Agents</b>										
89706	Triethylsilane	Silan E3H	≥98.0% GC	617-86-7	116.28	500 g	5 kg	-	-	>1 can
92231	1,1,3,3-Tetramethyl-disiloxane	Silan HSi2	≥98.0% GC	3277-26-7	134.32	500 g	5 kg	-	-	>1 can
<b>Building Blocks and Synthons</b>										
87406	Chloro(chloromethyl) dimethylsilane	Silan CMM2	≥98.0% GC	1719-57-9	143.09	250 g	2.5 kg	-	-	>1 can
70750	Dichloro(chloromethyl) methylsilane	Silan CMM1	≥98.0% GC	1558-33-4	163.51	500 g	5 kg	-	-	>1 can
18146	(Chloromethyl) trimethylsilane	Silan CMM3	≥98.0%	2344-80-1	122.67	250 g	2.5 kg	-	-	>1 can
77829	Trichloro(chloromethyl)silane	Silan CM	≥95.0%	1558-25-4	183.92	500 g	5 kg	-	-	>1 can
73314	Ethynyltrimethylsilane	TMSA	≥98.0% GC	1066-54-2	98.22	500 g	5 kg	-	-	>1 can
89787	(Trimethylsilyloxy)ethylene	TMSO-Ethen	≥96.0% GC	6213-94-1	116.23	500 g	5 kg	-	-	>1 can
<b>Amino Acids</b>										
79256	S-(4-Tolyl)-L-cysteine	-	purum ≥95.0% HPLC	-	211.28	1 g; 5 g	-	-	-	-
78904	S-(2-Thiazolyl)-L-cysteine	-	purum ≥95.0% HPLC	405150-20-1	204.27	1 g	-	-	-	-
95631	S-(2-Thienyl)-L-cysteine	-	purum ≥95.0% HPLC	405150-23-4	203.28	1 g	-	-	-	-
06993	3-(1-Pyrazolyl)-L-alanine	-	purum ≥95.0% HPLC	2734-48-7	155.15	1 g	-	-	-	-
12227	3-(1,2,4-Triazole-1-yl)-L-alanine	-	purum ≥95.0% HPLC	4819-36-7	156.14	1 g	-	-	-	-
67387	N-Boc-pyrazolyl-L-alanine	-	purum ≥97.0% HPLC	21012-18-0	255.27	0.5 g	-	-	-	-
50996	N-Boc-triazolyl-L-alanine	-	purum	-	256.26	0.5 g	-	-	-	-
51916	N-Fmoc-pyrazolyl-L-alanine	-	purum ≥97.0% HPLC	-	377.39	0.5 g	-	-	-	-
53229	N-Fmoc-triazolyl-L-alanine	-	purum ≥97.0% HPLC	-	378.38	0.5 g	-	-	-	-