

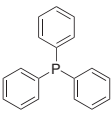
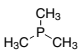
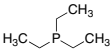
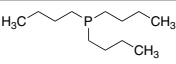
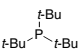
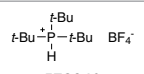
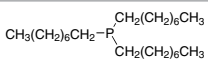
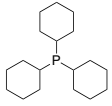
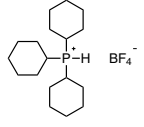
Phosphine Ligand Application Guide

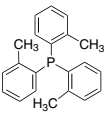
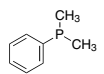
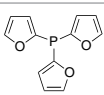
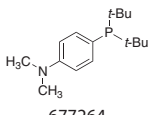
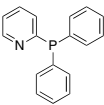
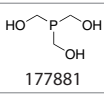
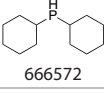
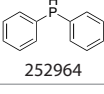
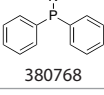
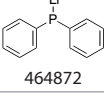
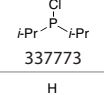
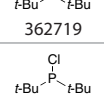
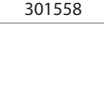
Add  Aldrich[™]

Phosphine Ligands

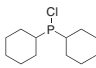
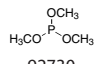
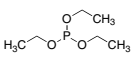
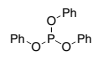
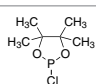
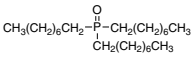
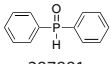
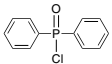
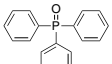
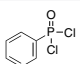
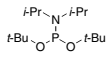
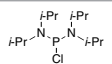
Transition-metal-catalyzed cross-coupling is an efficient tool for constructing C–C, C–N and C–O bonds in organic synthesis. Intensive interest has been focused on the development of the most effective ligands that improve catalyst performance. Ligands play a key role in stabilizing and activating the central metal atom and fine-tuning the selectivity of the transformation. Phosphines continue to be the most significant class of ligands for cross coupling. This brochure highlights the common applications and advantages of monodentate, bidentate, Buchwald, cataCXium®, and Dalphos ligands, all of which are commercially available from Aldrich Chemistry.

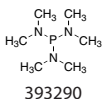
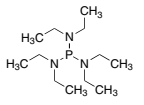
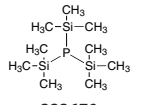
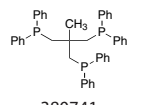
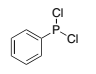
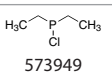
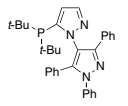
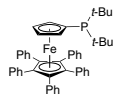
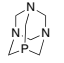
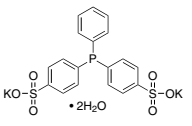
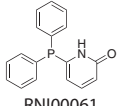
Monodentate Phosphine Ligands and Precursors

Structure and Cat. No.	Common Name	CAS No.	Mol. Wt.	Negishi	Suzuki	Stille	Heck	Sonogashira	Buchwald-Hartwig (C–N or C–O)	Additional Applications
 93090 93092 T84409	PPh ₃	603-35-0	262.29	X	X	X	X	X		<ul style="list-style-type: none"> • Morita-Baylis-Hillman reaction
 323322 324108 (1.0 M in THF) 324116 (1.0 M in toluene)	PMe ₃	594-09-2	76.08							<ul style="list-style-type: none"> • Mitsunobu reaction • Transformation of azides into carbamates • Aziridines from azidoalcohols • Reagent for iminophosphoranes • Aza-Wittig reaction
 346179 (1.0 M in THF) 245275	PEt ₃	554-70-1	118.16							<ul style="list-style-type: none"> • α-Arylation of ketones
 T49484 247049	P(<i>n</i> -Bu) ₃	998-40-3	202.32			X				<ul style="list-style-type: none"> • 1,4-addition • Acylation • Thioetherification of alcohols • Preparation of active esters
 570958 655325 (1.0 M in toluene)	P(<i>t</i> -Bu) ₃	13716-12-6	202.32	X	X	X	X	X	X	<ul style="list-style-type: none"> • Addition of organoboronic acids to aldehydes • Addition of alkynes to benzoylchlorides
 578940	TTBP · HBF ₄	131274-22-1	290.13	X	X	X	X	X	X	<ul style="list-style-type: none"> • Precursor for P(<i>t</i>-Bu)₃ • Carbonylation • Addition of boronate esters to alkynes and alkenes
 117854 718165	P(<i>n</i> -Oct) ₃	4731-53-7	370.64							<ul style="list-style-type: none"> • Surfactant for nanomaterial and nanocrystal synthesis • A general phosphorus source for the low-temperature conversion of metals into metal phosphides
 261971	PCy ₃	2622-14-2	280.43		X		X	X		<ul style="list-style-type: none"> • Hiyama coupling • Krische formal [4+2] of dienes and vicinal diols
 631493	PCy ₃ · HBF ₄	58656-04-5	368.24		X	X		X		<ul style="list-style-type: none"> • Control regiochemistry of hydrostannations of 1-alkynes • Intramolecular arylation • Synthesis of indolines from <i>N</i>-alkyl-2-bromoanilines

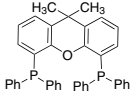
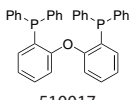
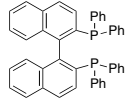
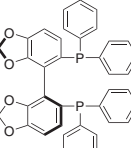
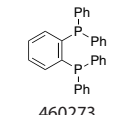
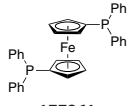
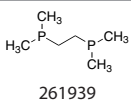
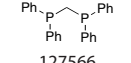
Structure and Cat. No.	Common Name	CAS No.	Mol. Wt.	Negishi	Suzuki	Stille	Heck	Sonogashira	Buchwald-Hartwig (C–NorC–O)	Additional Applications
 287822	P(o-tol) ₃	6163-58-2	304.37	X	X	X	X		X	• Silylation
 265020	Me ₂ PPh	672-66-2	138.15							• Ligand for transition metal complexes
 383767	TFP	5518-52-5	232.17	X	X	X	X	X	X	• Cycloisomerization/oxidation of alcohols
 677264	APhos	932710-63-9	265.37		X					• A-Phos ligand for cross-coupling
 392960	Diphenyl-2-pyridylphosphine	37943-90-1	263.27							• Carbonylations • Dehydrogenative coupling • Distannylation • Silylation • Carbostannylation • Hydration • Mitsunobu reaction
 177881	Tris(hydroxymethyl)phosphine	2767-80-8	124.08							• Precursor for metal complexes
 666572	HPCy ₂	829-84-5	198.28		X					• Precursor for ligands and metal complexes
 252964	HPPH ₂	829-85-6	186.19							• Precursor for the phosphorus-based alkenation reagents • Source of phosphorus-centered radical
 380768	KPPH ₂	15475-27-1	224.28							• Precursor for novel sulfur and phosphorus containing oxazolines used in asymmetric catalysis
 464872	Ph ₂ PLi	65567-06-8	192.12							• Dealkylation of alkyl aryl ethers • Dehydroxylation of α-hydroxy ketones
 337773	P(i-Pr) ₂ Cl	40244-90-4	152.60							• O-Arylation of phenols • Precursor for ligands and metal complexes
 362719	HP(t-Bu) ₂	819-19-2	146.21		X		X	X		• Precursor for ligands and metal complexes
 301558	P(t-Bu) ₂ Cl	13716-10-4	180.66		X					

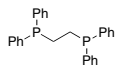
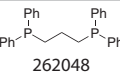
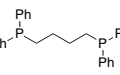
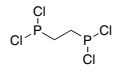
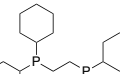
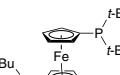
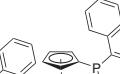
Monodentate Phosphine Ligands and Precursors – continued

Structure and Cat. No.	Common Name	CAS No.	Mol. Wt.	Negishi	Suzuki	Stille	Heck	Sonogashira	Buchwald-Hartwig (C-N or C-O)	Additional Applications
 481408	Cy ₂ P(OMe)Cl	16523-54-9	232.73							<ul style="list-style-type: none"> • Precursor for phosphine ligands
 92730 240907	P(OMe) ₃	121-45-9	124.08		X		X			<ul style="list-style-type: none"> • Mild desulfurization reagent
 T61204	P(OEt) ₃	122-52-1	166.16							<ul style="list-style-type: none"> • Reducing agent • Precursor for phosphonates or phosphates • Used with copper(I) iodide to form a stable metal complex
 T84654	P(OPh) ₃	101-02-0	310.28			X	X			<ul style="list-style-type: none"> • Conversion of alcohols to halides • Wittig-type reaction • Deoxygenation • Cycloaddition reactions • Allylic alkylations • Cyclohydrocarbonylation/CO insertion • Wittig-type olefinations
 447536	2-Chloro-4,4,5,5-tetramethyl-1,3,2-dioxaphospholane	14812-59-0	182.59							<ul style="list-style-type: none"> • Reagent for the phosphitylation of alcohols and heteroatomic nucleophiles
 223301 346187	TOPO®	78-50-2	386.63							<ul style="list-style-type: none"> • Agent for extraction of metals • Capping ligand for the production of quantum dots • Solvent for the synthesis and solubilizes the growing nanoparticles
 287881	HPOPh ₂	4559-70-0	202.19		X			X		<ul style="list-style-type: none"> • Hydrophosphinylation of terminal alkynes • Selective hydrophosphinylation of olefins • Precursor for diphenylphosphino-containing chiral ligands
 230235	Ph ₂ POCl	1499-21-4	236.63							<ul style="list-style-type: none"> • Acid activation • Alkylphosphine oxide formation • Amine protection
 T84603	TPPO	791-28-6	278.28							<ul style="list-style-type: none"> • Catalyst for allylation of aldehydes or <i>N</i>-aryl hydrazones • Promoter for coupling reaction of 1-alkenyl with sulfonyl chlorides • Cocatalyst for enantioselective cyanosylations • Tuning ligand for asymmetric epoxidation • Scavenger for enantioselective Michael reactions of enones
 389560	PhPOCl ₂	824-72-6	194.98							<ul style="list-style-type: none"> • Precursor for phosphonate ligands and metal complexes
 419362	Di- <i>tert</i> -butyl <i>N,N</i> -diisopropylphosphoramidite	137348-86-8	277.38							<ul style="list-style-type: none"> • Precursor for ligands • Reagent for the synthesis of nucleosides and nucleotides
 341347	Bis(diisopropylamino)chlorophosphine	56183-63-2	266.79							<ul style="list-style-type: none"> • Reagent for the synthesis of oligonucleotides

Structure and Cat. No.	Common Name	CAS No.	Mol. Wt.	Negishi	Suzuki	Stille	Heck	Sonogashira	Buchwald-Hartwig (C-N or C-O)	Additional Applications
 393290	HMPT	1608-26-0	163.20		X					<ul style="list-style-type: none"> Synthesis of epoxides, arene oxides, and carbonates Conversion of alcohols to alkyl chlorides Conversion of disulfides to sulfides Deoxygenation of sulfoxides and azoxyarenes Reduction of ozonides Replaces Ph_3P in the Wittig reaction
 253189	$\text{P}(\text{NEt}_3)_3$	2283-11-6	247.36							<ul style="list-style-type: none"> Precursor for phosphine ligands and metal complexes
 333670	$\text{P}(\text{SiMe}_3)_3$	15573-38-3	250.54							<ul style="list-style-type: none"> Phosphorus source Alternative to phosphine gas Precursor of $(\text{Me}_3\text{Si})_2\text{PLi}$
 380741	Triphos	22031-12-5	624.67							<ul style="list-style-type: none"> Reagent for metal complexes
 D71984	PhPCl_2	644-97-3	178.98							<ul style="list-style-type: none"> Precursor for phosphine ligands
 573949	Et_2PCl	686-69-1	124.55							<ul style="list-style-type: none"> Precursor for phosphine ligands
 676632	BippyPhos	894086-00-1	506.62					X		<ul style="list-style-type: none"> Amidation Coupling of hydroxylamines with aryl halides
 675784	QPhos	312959-24-3	710.71		X			X	X	<ul style="list-style-type: none"> α-Arylation of esters, amides and aldehydes
 695467	PTA	53597-69-6	157.15		X			X		<ul style="list-style-type: none"> Morita-Baylis-Hillman reactions Hydrogenations Olefin hydroformylations Cyclization
 698539	Bis(<i>p</i> -sulfonatophenyl) phenylphosphine dihydrate dipotassium salt	151888-20-9	534.62		X					
 RNI00061	6-DPPon	64741-28-2	279.1							<ul style="list-style-type: none"> New Breit's Regioselective Hydroformylation Ligand

Bidentate Phosphine Ligands and Precursors

Structure and Cat. No.	Common Name	CAS No.	Mol. Wt.	Negishi	Suzuki	Stille	Heck	Sonogashira	Buchwald-Hartwig (C-N or C-O)	Additional Applications
 526460	XantPhos	161265-03-8	578.62		X	X			X	<ul style="list-style-type: none"> C-P bond formation Regioselective hydroformylation of allylamines Borylation of nitriles
 510017	DPEPhos	166330-10-5	538.55	X	X		X		X	<ul style="list-style-type: none"> Hydroamination
 481084	(±)-BINAP	98327-87-8	622.67			X			X	<ul style="list-style-type: none"> Intramolecular acylation of aryl bromides Minisci reaction with alcohols Decarboxylative benzylation of diphenylglycinate Imines Regioselective hydroarylation Decarboxylative cycloaddition S-arylation
 693006	(S)-SEGPPOS®	210169-54-3	610.57							<ul style="list-style-type: none"> Stereoselective preparation of homoallylic alcohols via Ir-catalyzed stereoselective transfer hydrogenative crotylation of an allylic acetate with alcohols or aldehydes
 460273	dppbenz	13991-08-7	446.46							<ul style="list-style-type: none"> Cycloadditions Addition of arylboronic acids Baeyer-Villiger oxidation Reduction of propargyl oxiranes and hydroboration of styrene
 177261	dppe	12150-46-8	554.38	X	X	X	X	X	X	<ul style="list-style-type: none"> Tsuji-Trost allylation Carbonylation of vinyl triflates Enantioselective ene reaction Addition of organoboronic acids to aldehydes
 261939	dmpe	23936-60-9	150.14							<ul style="list-style-type: none"> Ligand for metal complexes
 127566	dppm	2071-20-7	384.39		X		X			<ul style="list-style-type: none"> Ligand for metal complexes

Structure and Cat. No.	Common Name	CAS No.	Mol. Wt.	Negishi	Suzuki	Stille	Heck	Sonogashira	Buchwald-Hartwig (C-N or C-O)	Additional Applications
 106496	dppe	1663-45-2	398.42	X	X		X			<ul style="list-style-type: none"> • Allylation of ketones • Decarboxylation of allyl esters • [3 + 2] cycloaddition • Carbonylation • Homo-Diels-Alder cycloaddition of norbornadiene with alkynes
 262048	dppp	6737-42-4	412.44	X	X			X		<ul style="list-style-type: none"> • Deoxygenation of phenols via the corresponding triflate • Substitution of aryl bromides with vinyl ethers • Carbonylation of aryl triflates • Addition of β-dicarbonyl enolates to dienes • Coupling of aryl and alkyl halides with Grignard reagents
 261947	dppb	7688-25-7	426.47	X	X	X	X	X		<ul style="list-style-type: none"> • Lactonization of unsaturated alcohol • Salkylation of allyl acetates • Hydroacylation • Alkyne isomerizations • Intramolecular allylation of alkenes • Alkyne-CO₂ coupling
 261920	1,2-Bis(dichlorophosphino)ethane	28240-69-9	231.81							<ul style="list-style-type: none"> • Precursor for ligands and transition metal complexes
 479500	dcpe	23743-26-2	422.61	X	X					<ul style="list-style-type: none"> • Decarboxylative coupling • Ligand for metal complexes
 695149	D'BPF	84680-95-5	474.42		X				X	<ul style="list-style-type: none"> • Arylation • Indolization
 L511501	dnfp	N/A	754.61				X			<ul style="list-style-type: none"> • Zhou regioselective arylation of α-olefins

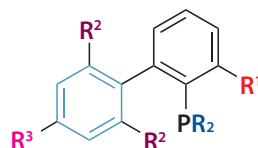
Buchwald Ligands

Buchwald ligands are bulky electron-rich dialkylbiaryl phosphines and are known to improve reactivity in palladium catalysis. These highly active reagents have been extensively applied in the synthesis of pharmaceuticals, natural products, polymers, and new materials. The structure of the dialkylbiaryl ligand is directly correlated to the efficiency of catalysts containing these ligands.

Structural features of Dialkylbiaryl Phosphine Ligands

R²: Increases catalyst stability by preventing cyclometalation, Encourages formation of L₁Pd(0)

R³: When not equal to H, usually only for ease of synthesis

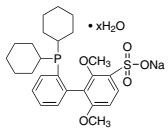
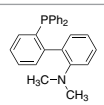
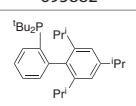
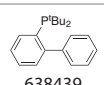
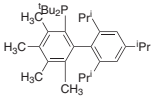
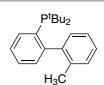
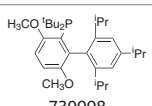
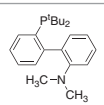
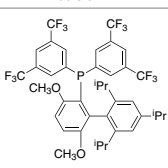


: Prevents oxidation at P by O₂, Accelerates reductive elimination

R¹: Substitution promotes reductive elimination

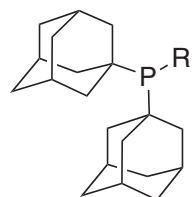
R: Electron-rich groups accelerate rate of oxidative addition

Structure and Cat. No.	Common Name	CAS No.	Mol. Wt.	Negishi	Suzuki	Stille	Heck	Sonogashira	Buchwald-Hartwig (C-N or C-O)	Additional Applications
 638099	Cyclohexyl JohnPhos	247940-06-3	350.48	X	X				X	<ul style="list-style-type: none"> • Hiyama coupling • Methylation of aryl halides • C-B bond formation • Oxidation of alcohols to ketones • α-Arylation of ketones
 638021	DavePhos	213697-53-1	393.54	X	X				X	<ul style="list-style-type: none"> • Kumada-Corriu coupling and α-Arylation of ketones • Hydroamination
 638064	XPhos	564483-18-7	476.72	X	X	X	X	X	X	<ul style="list-style-type: none"> • Hiyama coupling • Carbonyl enolate coupling
 638072	SPhos	657408-07-6	410.53	X	X		X		X	<ul style="list-style-type: none"> • Kumada-Corriu coupling
 695262	MePhos	251320-86-2	364.5		X				X	<ul style="list-style-type: none"> • α-Arylation of ketones
 663131	RuPhos	787618-22-8	466.64	X	X			X	X	<ul style="list-style-type: none"> • α-Arylation of oxindoles
 718742	BrettPhos	1070663-78-3	536.77						X	<ul style="list-style-type: none"> • Trifluoromethylation of aryl chlorides

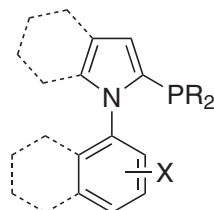
Structure and Cat. No.	Common Name	CAS No.	Mol. Wt.	Negishi	Suzuki	Stille	Heck	Sonogashira	Buchwald-Hartwig (C-N or C-O)	Additional Applications
 677280	³ SPhos	1049726-96-6	512.57		X					<ul style="list-style-type: none"> 1, 2-addition of arylboronic acids to aldehydes and ketones
 695882	PhDavePhos	240417-00-9	381.45					X		<ul style="list-style-type: none"> Heteroarene benzylation
 638080	tBuXPhos	564483-19-8	424.64					X		<ul style="list-style-type: none"> α-Arylation of tetramic acids Decarboxylative coupling of aromatic acids and aryl iodides Carboxylation of aryl bromides with CO₂
 638439	JohnPhos	224311-51-7	298.40		X		X	X		<ul style="list-style-type: none"> C-Si bond formation
 675938	Tetramethyl di-tBuXPhos	857356-94-6	480.75					X		<ul style="list-style-type: none"> Conversion of aryl halides to phenols
 695211	tBuMePhos	255837-19-5	312.43				X	X		<ul style="list-style-type: none"> α-Arylation of ketones Formate reduction of N-heterocyclic allylic acetates
 730998	tBuBrettPhos	1160861-53-9	484.69					X		<ul style="list-style-type: none"> O-Arylation of ethyl acetohydroximate C-F bond formation Conversion of aryl and vinyl triflates to bromides and chlorides Conversion of aryl chlorides, triflates, and nonaflates to nitroaromatics
 695874	tBuDavePhos	224311-49-3	341.47					X		<ul style="list-style-type: none"> α-Arylation of esters C-Si bond formation
 731013	JackiePhos	1160861-60-8	796.66					X		<ul style="list-style-type: none"> Coupling of aryl nonaflates and triflates with secondary amides, ureas, carbamates, and sulfonamides Coupling of aryl chlorides with secondary amides, carbamates, and sulfonamides

cataCXium® Ligands

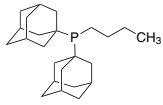
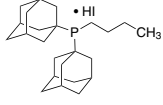
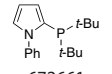
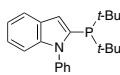
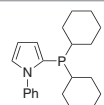
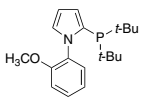
cataCXium Ligands are highly effective ligands for palladium catalyzed cross-coupling reactions. Di-adamantylalkylphosphine, known as cataCXium A, is a bulky and electron-rich phosphine ligand used for Heck and Suzuki couplings, Buchwald-Hartwig amination of arylchlorides, and α -arylation reactions of ketones. In reactions, these ligands allow for low catalyst loadings used under mild conditions. Recently, another class of basic, sterically hindered phosphines featuring phosphino-substituted *N*-aryl pyrroles (cataCXium P) has shown high catalyst turnover numbers for the Suzuki coupling of both electron-rich and electron-deficient aryl chlorides.



cataCXium® A



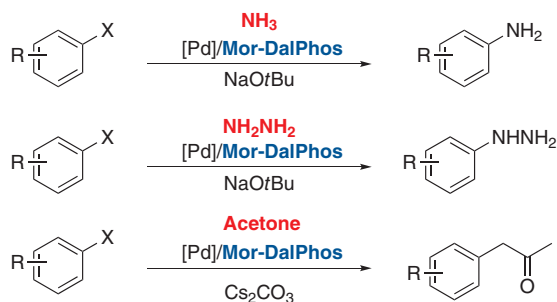
cataCXium® P

Structure and Cat. No.	Common Name	CAS No.	Mol. Wt.	Negishi	Suzuki	Stille	Heck	Sonogashira	Buchwald-Hartwig (C-N or C-O)	Additional Applications
 671479	cataCXium® A	321921-71-5	358.54		X		X	X	X	<ul style="list-style-type: none"> Formylation of aryl bromides Arylation
 671584	cataCXium® AHI	714951-87-8	486.45		X				X	<ul style="list-style-type: none"> Arylation
 672661	cataCXium® PtB	672937-61-0	287.38		X				X	<ul style="list-style-type: none"> Dehydrogenation of iso-propanol
 672343	cataCXium® PIntB	740815-37-6	337.44		X			X	X	
 672017	cataCXium® PCy	672937-60-9	339.45		X		X			<ul style="list-style-type: none"> Amination of alcohols with ammonia
 672564	cataCXium® POMetB	1053658-91-5	317.41						X	

DalPhos Ligands

The bulky di(1-adamantyl)phosphino [P(1-Ad)₂] fragment is common to the DalPhos ligand scaffold. These chelating N,P ligands are useful for Pd-catalyzed C-N and C-C bond formation. The more reactive Mor-DalPhos improves the scope and utility of ammonia coupling at room temperature, and is also effective in coupling of hydrazine and acetone.

Pd Catalyzed Monoarylation of Ammonia, Hydrazine and Acetone



Structure and Cat. No.	Common Name	CAS No.	Mol. Wt.	Negishi	Suzuki	Stille	Heck	Sonogashira	Buchwald-Hartwig (C-N or C-O)	Additional Applications
 751413	Me-DalPhos	1219080-77-9	421.6						X	
 751618	Mor-DalPhos	1237588-12-3	463.63						X	• Monoarylation of hydrazine and acetone
 759198	Di(1-adamantyl)-1-piperidinylphenylphosphine	1237588-13-4	461.66						X	• New

Sigma-Aldrich® Worldwide Offices

Argentina

Free Tel: 0810 888 7446
Tel: (+54) 11 4556 1472
Fax: (+54) 11 4552 1698

Australia

Free Tel: 1800 800 097
Free Fax: 1800 800 096
Tel: (+61) 2 9841 0555
Fax: (+61) 2 9841 0500

Austria

Tel: (+43) 1 605 81 10
Fax: (+43) 1 605 81 20

Belgium

Tel: (+32) 3 899 13 01
Fax: (+32) 3 899 13 11

Brazil

Free Tel: 0800 701 7425
Tel: (+55) 11 3732 3100
Fax: (+55) 11 5522 9895

Canada

Free Tel: 1800 565 1400
Free Fax: 1800 265 3858
Tel: (+1) 905 829 9500
Fax: (+1) 905 829 9292

Chile

Tel: (+56) 2 495 7395
Fax: (+56) 2 495 7396

People's Republic of China

Free Tel: 800 819 3336
Tel: (+86) 21 6141 5566
Fax: (+86) 21 6141 5567

Czech Republic

Tel: (+420) 246 003 200
Fax: (+420) 246 003 291

Denmark

Tel: (+45) 43 56 59 00
Fax: (+45) 43 56 59 05

Finland

Tel: (+358) 9 350 9250
Fax: (+358) 9 350 92555

France

Free Tel: 0800 211 408
Free Fax: 0800 031 052
Tel: (+33) 474 82 28 88
Fax: (+33) 474 95 68 08

Germany

Free Tel: 0800 51 55 000
Free Fax: 0800 64 90 000
Tel: (+49) 89 6513 0
Fax: (+49) 89 6513 1169

Hungary

Tel: (+36) 1 235 9055
Fax: (+36) 1 235 9068

India

Telephone

Bangalore: (+91) 80 6621 9400
New Delhi: (+91) 11 4358 8000
Mumbai: (+91) 22 4087 2364
Pune: (+91) 20 4146 4700
Hyderabad: (+91) 40 3067 7450
Kolkata: (+91) 33 4013 8000

Fax

Bangalore: (+91) 80 6621 9550
New Delhi: (+91) 11 4358 8001
Mumbai: (+91) 22 2579 7589
Pune: (+91) 20 4146 4777
Hyderabad: (+91) 40 3067 7451
Kolkata: (+91) 33 4013 8016

Ireland

Free Tel: 1800 200 888
Free Fax: 1800 600 222
Tel: +353 (0) 402 20370
Fax: + 353 (0) 402 20375

Israel

Free Tel: 1 800 70 2222
Tel: (+972) 8 948 4222
Fax: (+972) 8 948 4200

Italy

Free Tel: 800 827 018
Tel: (+39) 02 3341 7310
Fax: (+39) 02 3801 0737

Japan

Tel: (+81) 3 5796 7300
Fax: (+81) 3 5796 7315

Korea

Free Tel: (+82) 80 023 7111
Free Fax: (+82) 80 023 8111
Tel: (+82) 31 329 9000
Fax: (+82) 31 329 9090

Luxembourg

Tel: (+32) 3 899 1301
Fax: (+32) 3 899 1311

Malaysia

Tel: (+60) 3 5635 3321
Fax: (+60) 3 5635 4116

Mexico

Free Tel: 01 800 007 5300
Free Fax: 01 800 712 9920
Tel: (+52) 722 276 1600
Fax: (+52) 722 276 1601

The Netherlands

Tel: (+31) 78 620 5411
Fax: (+31) 78 620 5421

New Zealand

Free Tel: 0800 936 666
Free Fax: 0800 937 777
Tel: (+61) 2 9841 0555
Fax: (+61) 2 9841 0500

Norway

Tel: (+47) 23 17 60 00
Fax: (+47) 23 17 60 10

Poland

Tel: (+48) 61 829 01 00
Fax: (+48) 61 829 01 20

Portugal

Free Tel: 800 202 180
Free Fax: 800 202 178
Tel: (+351) 21 924 2555
Fax: (+351) 21 924 2610

Russia

Tel: (+7) 495 621 5828
Fax: (+7) 495 621 6037

Singapore

Tel: (+65) 6779 1200
Fax: (+65) 6779 1822

Slovakia

Tel: (+421) 255 571 562
Fax: (+421) 255 571 564

South Africa

Free Tel: 0800 1100 75
Free Fax: 0800 1100 79
Tel: (+27) 11 979 1188
Fax: (+27) 11 979 1119

Spain

Free Tel: 900 101 376
Free Fax: 900 102 028
Tel: (+34) 91 661 99 77
Fax: (+34) 91 661 96 42

Sweden

Tel: (+46) 8 742 4200
Fax: (+46) 8 742 4243

Switzerland

Free Tel: 0800 80 00 80
Free Fax: 0800 80 00 81
Tel: (+41) 81 755 2511
Fax: (+41) 81 756 5449

Thailand

Tel: (+66) 2 126 8141
Fax: (+66) 2 126 8080

United Kingdom

Free Tel: 0800 717 181
Free Fax: 0800 378 785
Tel: (+44) 01747 833 000
Fax: (+44) 01747 833 574

United States

Toll-Free: 800 325 3010
Toll-Free Fax: 800 325 5052
Tel: (+1) 314 771 5765
Fax: (+1) 314 771 5757

Vietnam

Tel: (+84) 8 3516 2810
Fax: (+84) 8 6258 4238

Internet

sigma-aldrich.com

*Enabling Science to
Improve the Quality of Life*

Order/Customer Service (800) 325-3010 • Fax (800) 325-5052
Technical Service (800) 325-5832 • sigma-aldrich.com/techservice
Development/Custom Manufacturing Inquiries **SAFC® (800) 244-1173**
Safety-related Information sigma-aldrich.com/safetycenter

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