

Asymmetric Synthesis



*New chiral diene ligand for
conjugate addition*

Featuring:

Asymmetric Reduction

Asymmetric Allylation

Asymmetric Epoxidation

Introduction

Asymmetric synthesis has become one of the most important tools for the synthesis of chiral molecules. The importance of this transformation is illustrated by the nearly endless discovery of chiral molecules used in various fields such as new drugs, fragrances, pesticides, etc. Since Knowles and Horner's pioneering work with the hydrogenation of prochiral olefins,^{1,2} a plethora of new ligands and catalysts have been developed for new asymmetric reactions. These new transformations give access to molecules in fewer steps, high yields and high enantiopurity.

In this issue of *ChemFiles*, we are pleased to present new chiral ligands and catalysts for a variety of transformations; new chiral phosphines for hydrogenation, hydrosilylation and for the reduction of ketones; a new generation of ligands such as chiral dienes for conjugate addition from Carreira and Zhou, as well as the ligands developed by Walsh. Our different Ipc_2 boranes as well as Soderquist borabicyclodecanes are also highlighted.

If you are unable to find the catalyst or ligand for your research, "Please Bother Us" with your suggestions at william.sommer@sial.com.

References: (1) Knowles, W. S. et al. *Chem. Commun.* **1968**, 1445. (2) Horner, L. et al. *Angew. Chem.* **1968**, 80, 1034.



William Sommer
Product Manager



Starting a new lab?

Let Sigma-Aldrich® help you get it done on time, on the spot and on the money.

Moving to a new location?

Make a good first impression – choose high-quality products at competitive prices from Sigma-Aldrich.

Received your first research grant?

Focus on your discoveries – not your expenses. Let Sigma-Aldrich help you make the most of your research dollars.

Join the Sigma-Aldrich New Lab Start-Up program, visit sigma-aldrich.com/newlabchem

Easy, Effective, Economical Sigma-Aldrich New Lab Start-Up Program!

- Over 100,000 quality biochemical and organic chemical products
- Instant savings of up to 70%
- Additional savings as your order value increases

sigma-aldrich.com

SIGMA-ALDRICH®

About Our Cover

The cover graphic represents the three-dimensional structure of (3a*S*, 6a*S*)-3,6-Diphenyl-1,3a,4,6a-tetrahydropentalene (hydrogen is not represented for clarity). This ligand is used in the conjugate addition of aryl boronic acids to α,β -unsaturated carbonyl compounds.

ChemFiles

Vol. 8 No. 6

**Aldrich Chemical Co., Inc.
Sigma-Aldrich Corporation**
6000 N. Teutonia Ave.
Milwaukee, WI 53209, USA

To Place Orders

Telephone 800-325-3010 (USA)
FAX 800-325-5052 (USA)

Customer & Technical Services

Customer Inquiries	800-325-3010
Technical Service	800-231-8327
SAFC®	800-244-1173
Custom Synthesis	800-244-1173
Flavors & Fragrances	800-227-4563
International	414-438-3850
24-Hour Emergency	414-438-3850
Web Site	sigma-aldrich.com
Email	aldrich@sial.com

Subscriptions

To request your **FREE** subscription to *ChemFiles*, please contact us by:

Phone: 800-325-3010 (USA)

Mail: **Attn: Marketing Communications**
Aldrich Chemical Co., Inc.
Sigma-Aldrich Corporation
P.O. Box 355
Milwaukee, WI 53201-9358

Email: sams-usa@sial.com

International customers, please contact your local Sigma-Aldrich office. For worldwide contact information, please see back cover.

ChemFiles are also available in PDF format on the Internet at sigma-aldrich.com/chemfiles.

Aldrich 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 for its particular use. See reverse side of invoice or packing slip for additional terms and conditions of sale.

All prices are subject to change without notice.

ChemFiles (ISSN 1933-9658) is a publication of Aldrich Chemical Co., Inc. Aldrich is a member of the Sigma-Aldrich Group. © 2008 Sigma-Aldrich Co.

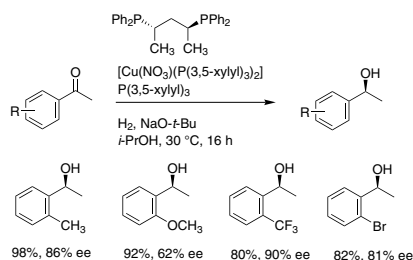
Asymmetric Reduction

Asymmetric Hydrogenation

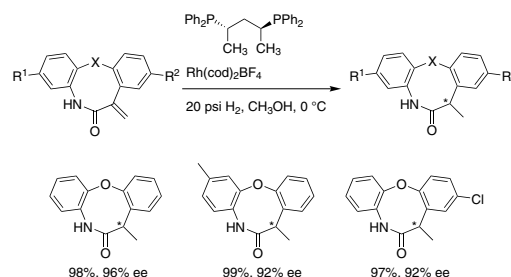
Asymmetric hydrogenation plays an important role in today's synthesis world. Knowles and Noyori pioneered this field and were rewarded with the Nobel Prize in 2001 for their work. It allows the generation of chiral centers selectively, with good yields. Furthermore, it gives access to a new pool of molecules with specific properties used in the pharmaceutical, flavors and fragrances, and agrochemical industries. However, the metal complex commonly used for these transformations is usually based on rhodium or ruthenium. The high cost and scarcity of these metal complexes make the technology prohibitive for use in large-scale synthesis. Based on the preliminary work of Stryker with copper catalysts, Shimizu et al. developed a new methodology for the hydrogenation of aryl ketones using copper complexes.¹ To achieve high enantioselectivities, (S,S)-BDPP ligand is necessary. Using a copper complex with (S,S)-BDPP, and tri(3,5-xylyl)-phosphine, a variety of aryl ketones were reduced to secondary alcohols in good to excellent yields with overall good ee's (**Scheme 1**).

The versatility of (S,S)-BDPP was demonstrated in the work done by Lu and Alper.² In their quest to synthesize chiral heterocycles, they investigated a number of catalysts and ligands for the asymmetric hydrogenation of terminal alkenes. Common ruthenium and iridium complexes were used with limited success. Furthermore, several chiral bidentate ligands such as BINAP, Monophos, DIOP, Duphos or Josiphos were tested in the transformation with little success. The best catalytic system was found to be (S,S)-BDPP with Rh(cod)₂BF₄. A variety of tricyclic lactams were synthesized with excellent enantioselectivity. The procedure typically involves the use of 1 mol% of rhodium complex and ligand in methanol at 0 °C under 20 psi of H₂. It is important to note that these heterocyclic motifs are encountered in many biologically active natural products as well as drug candidates.

References: (1) Shimizu, H. et al. *Org. Lett.* **2007**, 9, 1655. (2) Lu, S. -M. et al. *J. Am. Chem. Soc.* **2008**, 130, 6451.



Scheme 1

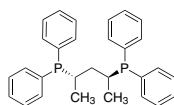


Scheme 2

(2S,4S)-2,4-Bis(diphenylphosphino)pentane, 97% NEW

[(1S,3S)-1,3-Dimethyl-1,3-propanediyl]bis(diphenylphosphine); (S,S)-BDPP
[77876-39-2]

C₂₉H₃₀P₂
FW 440.50



693618-500MG

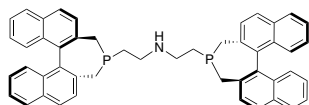
500 mg

693618-1G

1 g

2,2'-Bis[(R-1,1'-binaphthyl-2,2'-dimethyl)phosphino]diethylamine

Bis(R-2-(4,5-dihydro-3H-binaphtho[1,2-c;2',1'-e]-phosphino)ethyl)amine
C₄₈H₄₁NP₂
FW 693.79

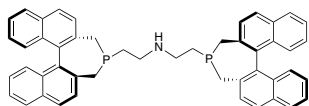


665002-250MG

250 mg

2,2'-Bis[(S-1,1'-binaphthyl-2,2'-dimethyl)phosphino]diethylamine, 97%

Bis(S-2-(4,5-dihydro-3H-binaphtho[1,2-c;2',1'-e]-phosphino)ethyl)amine
C₄₈H₄₁NP₂
FW 693.79

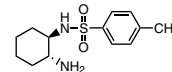


665010-250MG

250 mg

(1R,2R)-(-)-N-p-Tosyl-1,2-cyclohexanediamine, 97%

N-Toluenesulfonyl-(1R,2R)-(-)-1,2-diamino-cyclohexane; (-)-TsCYDN
C₁₃H₂₀N₂O₂S
FW 268.38

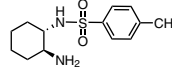


677302-1G

1 g

(1S,2S)-(+)-N-p-Tosyl-1,2-cyclohexanediamine, 97%

N-Toluenesulfonyl-(1S,2S)-(+)-1,2-diamino-cyclohexane; (+)-TsCYDN
C₁₃H₂₀N₂O₂S
FW 268.38

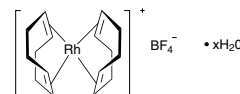


677310-1G

1 g

Bis(1,5-cyclooctadiene)rhodium(I) tetrafluoroborate hydrate NEW

[207124-65-0]
C₁₆H₂₄BF₄Rh · xH₂O
FW 406.07 (Anh)



683140-500MG

500 mg

683140-2G

2 g

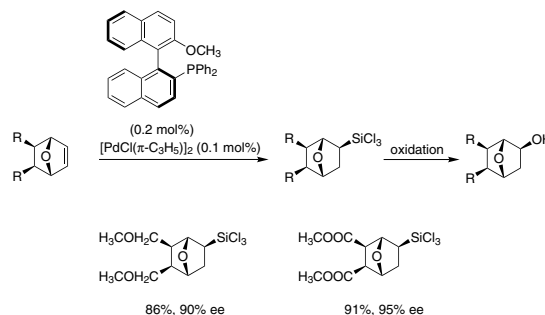
Asymmetric Hydrosilylation

The asymmetric hydrosilylation of olefins has proven to be a useful method to access a variety of chiral alcohols.¹ The introduction of the silica moiety allows further transformation of the molecule before converting it to the alcohol analogue. However, this particular reaction proved to be particularly challenging. To overcome this challenge Hayashi and co-workers synthesized a new chiral monodentate phosphine ligand using 1,1'-binaphthyl as a backbone.¹ This ligand, when used with palladium showed excellent stereoselectivities for the asymmetric hydrosilylation of 2,5-dihydrofurans.² The versatility of this ligand was showcased in the hydrosilylation of a variety of meso bicyclic olefins containing 2,5-dihydrofuran skeleton, with enantioselectivities as high as 95% (**Scheme 1**). In a typical reaction, 0.1 mol% of palladium allyl chloride dimer, with 0.2 mol% of MOP ligand is used. A reaction time of 24 hours is required to achieve good yields.

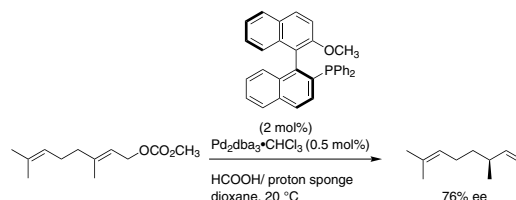
Palladium-Catalyzed Reduction of Allylic Esters

The palladium catalyzed reduction of allylic esters provides a convenient method to access chiral olefins.³ Hayashi et al. studied the activity of the MOP ligand with palladium toward the reduction of allylic esters.⁴ Mechanistic studies of the reaction have showed that a monodentate phosphine ligand is essential to achieve high selectivity. Reacting geranyl methyl carbonate with $\text{Pd}_2(\text{dba})_3$ and MOP (1 mol% respectively), quantitatively yielded (S)-3,7-dimethyl-1,6-octadiene with 76% ee (**Scheme 2**). It is important to note that a low catalyst and ligand loading with mild conditions is necessary for good selectivity.

References: (1) Hayashi, T. *Acc. Chem. Res.* **2000**, *33*, 354. (2) Uozumi, Y. et al. *Tetrahedron Lett.* **1993**, *34*, 2335. (3) Tsuji, J. *Palladium Reagents and Catalysts*; Wiley: New York, 1995. (4) Hayashi, T. et al. *J. Am. Chem. Soc.* **1994**, *116*, 775. Hayashi, T. et al. *Synthesis* **1994**, 526.



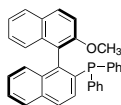
Scheme 1



Scheme 2

R-MOP

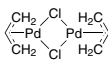
(R)-(+)-2-(Diphenylphosphino)-2'-methoxy-1,1'-binaphthyl
[145964-33-6]
 $\text{C}_{33}\text{H}_{25}\text{OP}$
FW 468.52



693537-50MG	50 mg
693537-100MG	100 mg

Allylpalladium(II) chloride dimer, 98%

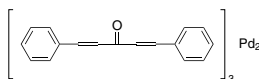
[12012-95-2]
 $\text{C}_6\text{H}_{10}\text{Cl}_2\text{Pd}_2$
FW 365.89



222380-100MG	100 mg
222380-500MG	500 mg
222380-5G	5 g

Tris(dibenzylideneacetone)dipalladium(0)

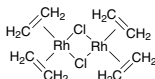
[51364-51-3]
 $\text{C}_{51}\text{H}_{42}\text{O}_3\text{Pd}_2$
FW 915.72



328774-500MG	500 mg
328774-1G	1 g
328774-5G	5 g
328774-50G	50 g
328774-100G	100 g

μ-Dichlorotetraethylene dirhodium(I)

[12081-16-2]
 $\text{C}_8\text{H}_{16}\text{Cl}_2\text{Rh}_2$
FW 388.93



656763-100MG	100 mg
656763-1G	1 g

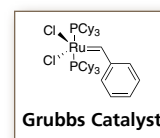
Discover Sigma-Aldrich's Newest Cheminar®

The Efficient Application of Olefin Metathesis in Pharmaceuticals and Fine Chemicals

- Featuring the latest innovative chemical synthesis technologies and products for metathesis
- Access directly via your desktop
- Convenient navigation
- Highly interactive

Contents:

- Ring Closing Metathesis
- Cross Metathesis
- Recent Catalyst Developments
- Optimizing Metathesis Reactions
- Ruthenium Catalyst Removal Techniques



To view Sigma-Aldrich's Web-based chemistry seminar series, please visit sigma-aldrich.com/cheminars

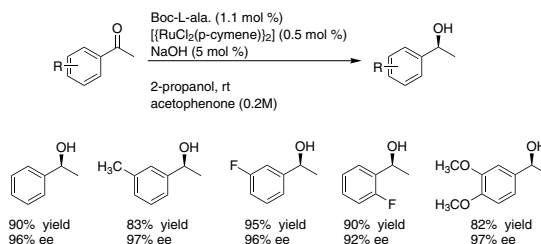
sigma-aldrich.com

SIGMA-ALDRICH®

Asymmetric Reduction of Ketones

The enantioselective reduction of ketones to alcohols gives access to a pool of chiral building blocks that can be used for the synthesis of natural products. Adolfsson and co-workers reported a novel class of ligands, based on a pseudo-dipeptide,¹ for the mild and efficient reduction of ketones using 2-propanol as a hydrogen source. The ligand is used with $[(\text{RuCl}_2(\text{p-cymene}))_2]$ in presence of NaOH for the asymmetric reduction of a variety of acetophenone derivatives. Excellent yields and ee's were reported (**Scheme 1**).

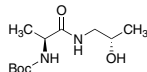
Reference: (1) Bøgevig, A. et al. *Chem. Eur. J.* **2004**, *10*, 294.; Västilä, P. et al. *Chem. Commun.* **2005**, 4039.



Scheme 1

Boc-L-alanine (2S)-2-hydroxypropylamide, 97%

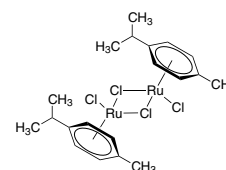
$\text{C}_{11}\text{H}_{22}\text{N}_2\text{O}_4$
FW 246.30



684414-100MG	100 mg
684414-500MG	500 mg

Dichloro(p-cymene)ruthenium(II) dimer

[52462-29-0]
 $\text{C}_{20}\text{H}_{28}\text{Cl}_4\text{Ru}_2$
FW 612.39

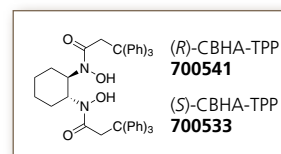
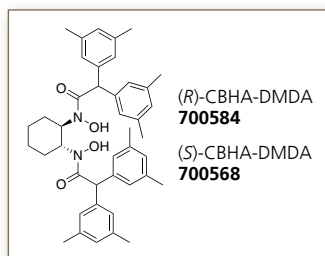
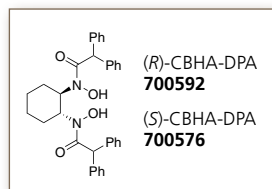
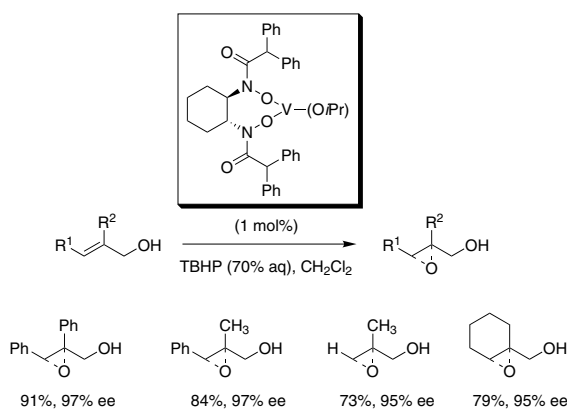


343706-1G	1 g
343706-5G	5 g

Asymmetric Epoxidation of Allylic Alcohols

Catalytic asymmetric epoxidation of olefins has become a reaction of choice to generate various chiral building blocks used in the synthesis of natural products and biologically active molecules. Yamamoto and co-workers developed bishydroxamic acid based ligands for the efficient asymmetric epoxidation of various allylic alcohols. Using only 1 mol% of catalyst (derived from vanadium and the bishydroxamic acid ligand), a variety of allylic alcohols were reacted to form enantiopure epoxides with high selectivity.

Reference: Zhang, W. et al. *Angew. Chem., Int. Ed.* **2005**, *44*, 4389.

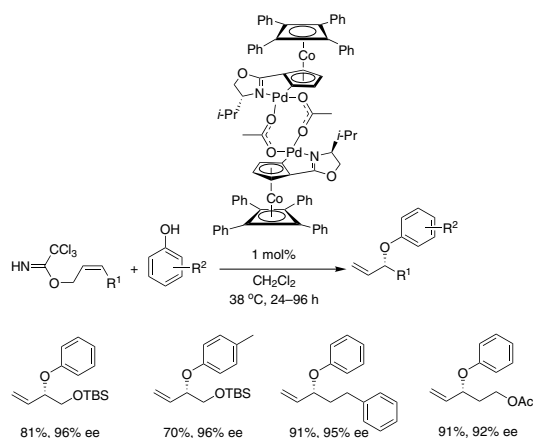


Asymmetric Allylation

Asymmetric Allylation Using COP Catalysts

Chiral allylic aryl esters are valuable building blocks for the synthesis of natural products and drug candidates. In recent years, several research groups developed new catalysts to synthesize chiral allylic esters from phenols and prochiral allylic precursors.¹ Overman and co-workers developed a new method to synthesize chiral 3-aryloxy-1-alkenes using COP-OAc catalysts.² This reaction only requires 1 mol% of catalyst producing the desired product in high yields and high ee's (**Scheme 1**). A library of allylic trichloroacetimidates and phenols were screened. This new method is compatible with the presence of base-labile functionality.

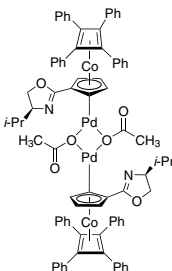
References: (1)(a) Trost, B. et al. *J. Am. Chem. Soc.* **2005**, 127, 7014. (b) Mbaye, M. D. Et al. *Chem. Commun.* **2004**, 1870. (c) Shekhar, S. et al. *J. Am. Chem. Soc.* **2006**, 128, 11770. (2) Kirsch, S. F. Et al. *Org. Lett.* **2007**, 9, 911.



Scheme 1

(S)-(+)-COP-OAc Catalyst, 95%

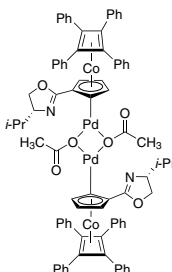
Di-μ-acetato-bis[η⁵-(S)-(*p*,*R*)-2-(2'-(4'-methylethyl)oxazolyl)cyclopentadienyl, 1-C,3'-N(η⁴-tetraphenylcyclobutadiene)cobalt]dipalladium; [Bis[μ-(acetato-κO:κO')]dipalladium] bis[1,1',1'',1'''-(η⁴-1,3-cyclobutadiene-1,2,3,4-tetrayl)tetrakis[benzene]]bis[μ-[(1-η:1,2,3,4,5-η)-2-[(4*S*)-4,5-dihydro-4-(1-methylethyl)-2-oxazolyl-κN³]-2,4-cyclopentadien-1-ylidene]]dicobalt [222400-03-5]
C₈₂H₇₂O₆N₂Co₂Pd₂
FW 1512.17



661716-250MG	250 mg
661716-1G	1 g

(R)-(-)-COP-OAc Catalyst, 95%

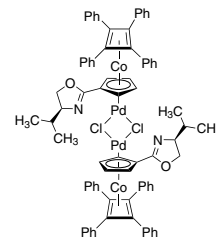
Di-μ-acetato-bis[η⁵-(*R*)-(*p*,*R*)-2-(2'-(4'-methylethyl)oxazolyl)cyclopentadienyl, 1-C,3'-N(η⁴-tetraphenylcyclobutadiene)cobalt]dipalladium [849592-74-1]
C₈₂H₇₂O₆N₂Co₂Pd₂
FW 1512.17



661708-250MG	250 mg
661708-1G	1 g

(S)-(+)-COP-Cl Catalyst

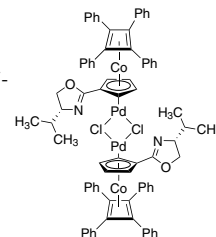
[612065-01-7]
C₇₈H₆₆Cl₂Co₂N₂O₂Pd₂
FW 1464.98



646636-250MG	250 mg
646636-1G	1 g

(R)-(-)-COP-Cl Catalyst

Di-μ-chloro-bis[η⁵-(*R*)-(*p*,*R*)-2-(2'-(4'-methylethyl)oxazolyl)cyclopentadienyl, 1-C,3'-N(η⁴-tetraphenylcyclobutadiene)cobalt]dipalladium; [Bis[μ-(chloro-κO:κO')]dipalladium]bis[1,1',1'',1'''-(η⁴-1,3-cyclobutadiene-1,2,3,4-tetrayl)tetrakis[benzene]]bis[μ-[(1-η:1,2,3,4,5-η)-2-[(4*R*)-4,5-dihydro-4-(1-methylethyl)-2-oxazolyl-κN³]-2,4-cyclopentadien-1-ylidene]]dicobalt [612065-00-6]
C₇₈H₆₆Cl₂Co₂N₂O₂Pd₂
FW 1464.98



661791-100MG	100 mg
661791-500MG	500 mg

Asymmetric Allylboration Using $\text{Ipc}_2\text{B(allyl)}$

Asymmetric allylboration of aldehydes is an extremely important method for preparation of homoallylic alcohols, as evidenced in numerous complex natural product syntheses.¹ Although a variety of reagents have been developed to execute this reaction, the most broadly adopted have been Brown's α -pinene-derived (-)- and (+)- β -allyldiisopinocampheylboranes (i.e., $\text{Ipc}_2\text{B(allyl)}$).² Typically the allylboration reagents are generated in situ and used immediately at -78 to -100 °C, since they are not stable for extended periods at ambient temperature (**Scheme 1**).

New, salt-free, $\text{Ipc}_2\text{B(allyl)}$ reagents have been developed as stable solutions in pentane or dioxane. Under refrigerator storage, no appreciable decreases in selectivity are observed after several months. They exhibit excellent reactivity and reactions can be performed at 5 °C, even in water.

References: (1)(a) Furstner, A. et al., *Angew. Chem. Int. ed.* **2006**, 45, 5506. (b) Smith, A. B. et al. *J. Am. Chem. Soc.* **2003**, 125, 350. (2) For a review of allylborane reagents, see: Ramachandran, P. V. *Aldrichimica Acta* 2002, 35, 23.

(-)- $\text{Ipc}_2\text{B(allyl)}$ borane solution

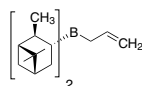
(-)- β -Allyldiisopinocampheylborane solution
[85116-38-7]

$\text{C}_{23}\text{H}_{39}\text{B}$

FW 326.37

► 1 M in dioxane

678546-5ML	5 mL
678546-25ML	25 mL



(-)- $\text{Ipc}_2\text{B(allyl)}$ borane solution

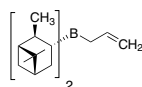
(-)- β -Allyldiisopinocampheylborane solution
[85116-38-7]

$\text{C}_{23}\text{H}_{39}\text{B}$

FW 326.37

► 1 M in pentane

678538-5ML	5 mL
678538-25ML	25 mL



(+)- $\text{Ipc}_2\text{B(allyl)}$ borane solution

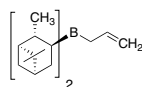
(+)- β -Allyldiisopinocampheylborane solution
[106356-53-0]

$\text{C}_{23}\text{H}_{39}\text{B}$

FW 326.37

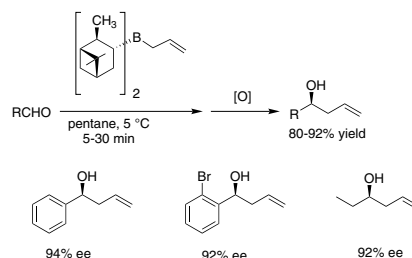
► 1 M in dioxane

678511-5ML	5 mL
678511-25ML	25 mL



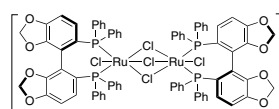
► 1 M in pentane

678503-5ML	5 mL
678503-25ML	25 mL



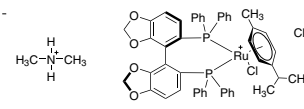
Scheme 1

Takasago Catalysts for Asymmetric Synthesis



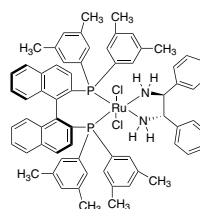
692204

JP Registration No. 3148136, 3549390



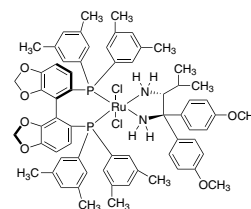
692522

JP Registration No. 3148136



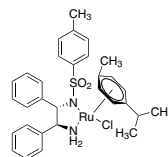
693251

JP Registration No. 2041996, 2731377, 2935453,
JP Application No. 19970359654

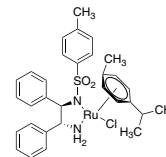


692328

JP Application No. 19970359654,
JP Registration No. 3148136



703907



703915

Sold in collaboration with Takasago for research purposes only.

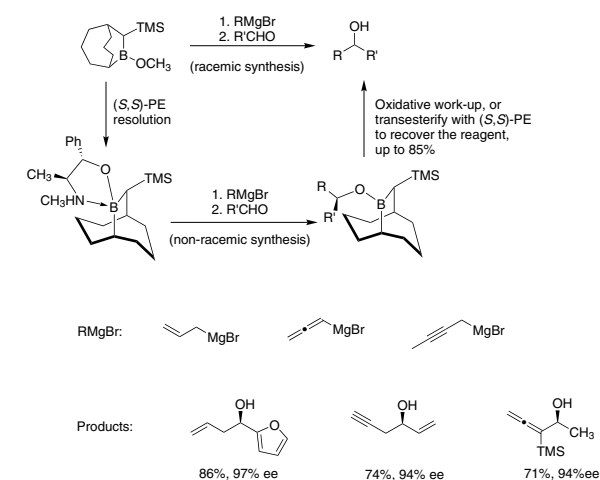
sigma-aldrich.com

SIGMA-ALDRICH
Chemistry

Soderquist Borabicyclodecanes

In the past 20 years, the asymmetric allylboration of aldehydes has become an essential reaction for the synthesis of homoallylic or homopropargylic alcohols. Soderquist et al. developed a robust, versatile and recyclable pseudoephedrine reagent improving the reaction performance and selectivity. Treatment of these air stable crystalline pseudoephedrine borinic acid complexes with the appropriate Grignard reagent gives access to homoallylic or homopropargylic alcohols and -allenyl carbinols via symmetric allyl-, propargyl-, and allenylboration reactions. Allylboration reactions occur rapidly, can be performed in the presence of Mg salts, and the degree of asymmetric induction is only minimally affected by temperature. Moreover, the reagents are easily recycled in good yield. The reagents also perform well in asymmetric crotylboration,¹ allylboration of imines,² and the alkynylboration of *N*-acylimines to give chiral *N*-propargylamides (**Scheme 1**).³

References: (1) Burgos, C. H. et al. *J. Am. Chem. Soc.* **2005**, 127, 8044. (2) Hernandez, E. et al. *Pure Appl. Chem.* **2006**, 78, 1389. (3) Gonzales, A. Z. et al. *Org. Lett.* **2006**, 8, 3331.



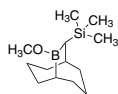
Scheme 1

B-Methoxy-10-trimethylsilyl-9-borabicyclo[3.3.2]decane, 95%

9-Methoxy-10-trimethylsilyl-9-borabicyclo[3.3.2]decane

C₁₃H₂₇BOSi

FW 238.25



676667-5G	5 g
676667-25G	25 g

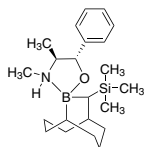
(+)-9-(1*S*, 2*S*-Pseudoephedriny)-(10*R*)-(trimethylsilyl)-9-borabicyclo[3.3.2]decane

N, α -Dimethyl- β -[[(10*R*)-10-(trimethylsilyl)-9-borabicyclo[3.3.2]dec-9-yl]oxy]-(α *S*, β *S*)-benzeneethanamine

[848618-13-3]

C₂₂H₃₈BNOSi

FW 371.44

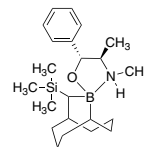


676675-1G	1 g
676675-5G	5 g

(-)-9-(1*R*,2*R*-Pseudoephedriny)-(10*S*)-(trimethylsilyl)-9-borabicyclo[3.3.2]decane

C₂₂H₃₈BNOSi

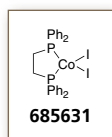
FW 371.44



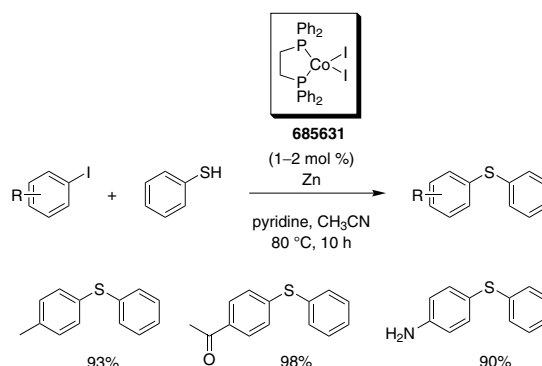
676683-1G	1 g
676683-5G	5 g

Aryl-Sulfur Bond Formation

Wong et al. developed a new cobalt catalyst for the cross coupling of aryl halides with thiophenols. A growing interest in this coupling reaction has been witnessed in the past few years due to the occurrence of carbon-sulfur bond motifs in natural products and drug candidates. Utilizing 1 to 2 mol % of CoI₂(dppe), various aryl halides were coupled to thiophenols with excellent yields.



Reference: Wong, Y.-C. et al. *Org. Lett.* **2006**, 8, 5613.



sigma-aldrich.com

SIGMA-ALDRICH®

Asymmetric Aldol Reaction Using Bis-ProPhenol Ligands

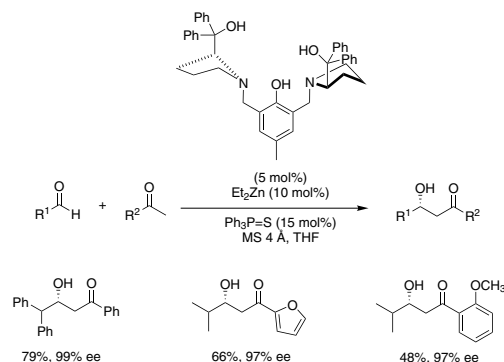
The aldol reaction is an important tool for the synthesis of complex molecules. However, the chemo- and regioselectivity of this reaction is usually poor. In view of these challenges, Trost and co-workers developed a new ligand capable of performing aldol reactions enantioselectively.¹ The approach taken by the researchers was to look at the different requirements for a good catalyst for asymmetric transformations. The final structure of the catalyst is a semi-crown complex which allows good chiral recognition as well as a path for easy exchange of product. The ligand structure is based on a *p*-cresol backbone. To probe the activity of the ligand, various aldehydes were reacted with various aryl ketones with overall good yields and excellent selectivity (**Scheme 1**).

Asymmetric Mannich-Type Reaction Using Bis-ProPhenol Ligands

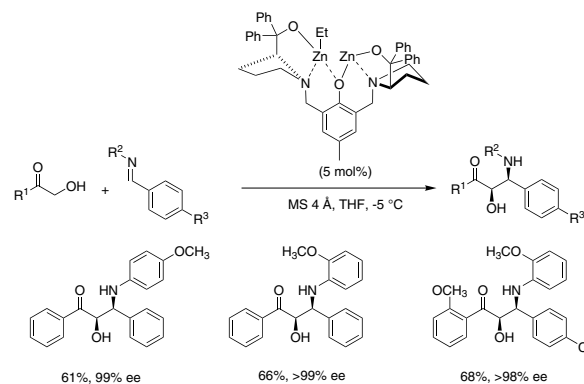
The Mannich reaction has become an essential tool for the synthesis of nitrogen-containing compounds. Trost and co-workers utilized the semi-crown ligand that they developed for asymmetric aldol reaction in an asymmetric imine-aldol reaction generating syn 1,2-amino alcohols.² A series of hydroxyacetophenone derivatives were reacted with different glyoxalate imines in the presence of a zinc complex in THF (**Scheme 2**). It was noted that more electron-rich aromatic hydroxy ketones required higher catalyst loading. Good yields and selectivity were reported for the synthesis of a variety of amino alcohols. It is important to note that these imine alcohols are important synthons that can be reacted further to yield different α -hydroxyl- β -amino acids.

A recent example of the use of Bis-ProPhenol ligands is the Friedel-Crafts alkylation of pyrroles with nitroalkenes. The Friedel-Crafts reaction has been extensively studied and used for the past 100 years. Recently, several research groups focused their attention on asymmetric Friedel-Crafts alkylation with very good success.³ Trost et al. reported a dinuclear zinc-bis-ProPhenol complex to efficiently catalyze the asymmetric Friedel-Crafts alkylation of unprotected pyrroles with nitroalkenes.⁴ The resulting products are 2-alkyl-substituted and 2,5-dialkyl-substituted pyrroles containing a chiral center at C-1 of the alkyl substituent. The significance of this transformation lies in the fact that pyrroles are relatively unstable in the acidic environment of the classical Friedel-Crafts reaction and that unprotected pyrroles can be used directly in the reaction. Furthermore, pyrroles are important motifs present in numerous natural products and medicinal reagents. The current alkylation is carried out at room temperature in THF and gives rise to moderate-to-high yields and enantioselectivities (**Scheme 3**).

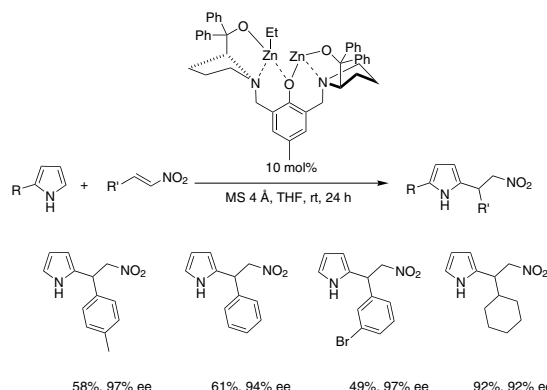
References: (1) Trost, B. M. et al. *J. Am. Chem. Soc.* **2000**, 122, 12003. (2) Trost, B. M. et al. *J. Am. Chem. Soc.* **2003**, 125, 338. (3)(a) Bandini, M. et al. *Eur. J. Org. Chem.* **2006**, 3527. (b) Jia, Y. X. et al. *J. Org. Chem.* **2006**, 71, 75. (4) Trost, B. M. et al. *J. Am. Chem. Soc.* **2008**, 130, 2438.



Scheme 1



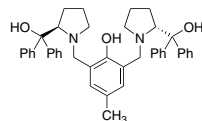
Scheme 2



Scheme 3

(*R,R*)-(-)-2,6-Bis[2-(hydroxydiphenylmethyl)-1-pyrrolidinyl-methyl]-4-methylphenol, 95%

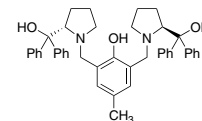
(*R,R*)-2,6-Bis[2-(2-(hydroxydiphenylmethyl)pyrrolidin-1-yl)methyl]-4-methylphenol;
Trost (*R,R*)-Bis-ProPhenol Ligand
 $C_{43}H_{46}N_2O_3$
FW 638.84



667625-1G	1 g
667625-5G	5 g

(*S,S*)-(+)-2,6-Bis[2-(hydroxydiphenylmethyl)-1-pyrrolidinyl-methyl]-4-methylphenol, 95%

(*S,S*)-2,6-Bis[2-(2-(hydroxydiphenylmethyl)pyrrolidin-1-yl)methyl]-4-methylphenol;
Trost (*S,S*)-Bis-ProPhenol Ligand
 $C_{43}H_{46}N_2O_3$
FW 638.84



668370-1G	1 g
668370-5G	5 g

Asymmetric Conjugate Addition

In the past 5 years, chiral diene ligands have surfaced for a variety of asymmetric transformations. Hayashi¹ and Carreira² pioneered this field by synthesizing chiral diene ligands that formed stable complexes with metals and exerted high catalytic activity as well as high enantioselectivity. Carreira and co-workers reported the asymmetric synthesis of 3,3-diarylpropanals using a chiral bicyclo[2.2.2]octadiene as a ligand with a rhodium complex.³ This new synthesis provides access to important building blocks that are otherwise not readily available. The reaction offers a general synthesis of a wide variety of diaryl propanals. This rhodium catalyzed conjugate addition of aryl boronic acids is both chemo- and regio-selective wherein conjugate addition is favored over 1,2-addition. In a typical reaction, 3.3 mol% of the ligand and 3 mol% of the rhodium complex in a mixture of methanol and water is required. Good yields and selectivity were reported with both electron rich and electron poor boronic acids and various cinnamaldehydes (**Scheme 1**).

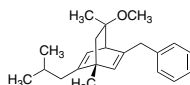
Following on the work of Carreira and Hayashi, Feng et al. developed a new type of diene ligand based on a nonbridged bicyclo[3.3.0]octadiene framework.⁴ The two cis-fused cyclopentadiene rings generate a wedge structure providing a chiral environment giving excellent enantiocontrol in the reaction when coordinated to the metal. The activity of the ligand was evaluated in the asymmetric arylation of *N*-tosylarylimines with aryl boronic acids. Using 3 mol% of [RhCl(C₂H₄)₂]₂ and 3.3 mol% of the ligand, a variety of aryl boronic acids and *N*-tosylarylimines were screened. It is important to note that the reactants screened had diverse steric and electronic properties and showed little influence on the yield and enantioselectivity of the reaction (**Scheme 2**).

To expand the scope of this new ligand, Feng et al. investigated the asymmetric 1,4-addition of aryl boronic acids to α,β -unsaturated carbonyl compounds.⁵ Using 2.5 mol% of [RhCl(C₂H₄)₂]₂ and 5.5 mol% of the ligand, a variety of sterically hindered, electron donating and electron withdrawing substituted aryl boronic acids were screened (**Scheme 3**). The nature of the substituent on the aryl boronic acid had no influence on the yield and the enantioselectivity.

References: (1) Hayashi, T. et al. *J. Am. Chem. Soc.* **2003**, 125, 11508. (2) Fischer, C. et al. *J. Am. Chem. Soc.* **2004**, 126, 1628. (3) Paquin, J. -F. et al. *J. Am. Chem. Soc.* **2005**, 127, 10850. (4) Wang, Z., -Q. Et al. *J. Am. Chem. Soc.* **2007**, 129, 5336. (5) Feng, C. -G. Et al. *Chem. Asian. J.* **2008**, ASAP

(1*R*,4*R*,8*R*)-5-Benzyl-8-methoxy-1,8-dimethyl-2-(2'-methylpropyl)-bicyclo[2.2.2]octa-2,5-diene, $\geq 98\%$

(1*R*,4*R*,8*R*)-5-Benzyl-2-isobutyl-8-methoxy-1,8-dimethyl-2-bicyclo[2.2.2]octa-2,5-diene; (1*R*,4*R*,8*R*) Carreira DOLEFIN Ligand [948594-95-4]



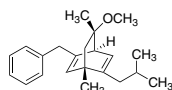
C₂₂H₃₀O
FW 310.47

669490-100MG

100 mg

(1*S*,4*S*,8*S*)-5-Benzyl-8-methoxy-1,8-dimethyl-2-(2'-methylpropyl)-bicyclo[2.2.2]octa-2,5-diene, $\geq 98\%$

(1*S*,4*S*,8*S*)-5-Benzyl-2-isobutyl-8-methoxy-1,8-dimethyl-2-bicyclo[2.2.2]octa-2,5-diene; (1*S*,4*S*,8*S*) Carreira DOLEFIN Ligand [862499-50-1]



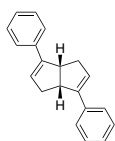
C₂₂H₃₀O
FW 310.47

672254

(3*aR*,6*aR*)-3,6-Diphenyl-1,3*a*,4,6*a*-tetrahydropentalene, 97%

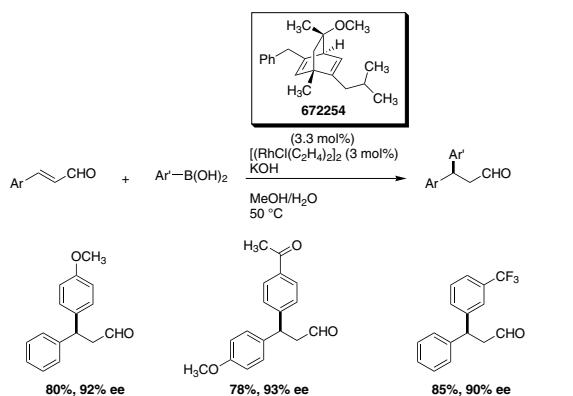
(3*aR*,6*aR*)-1,3*a*,4,6*a*-tetrahydro-3,6-diphenylpentalene [947503-81-3]

C₂₀H₁₈
FW 258.36

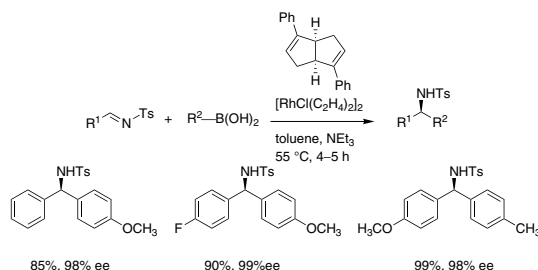


698725-50MG

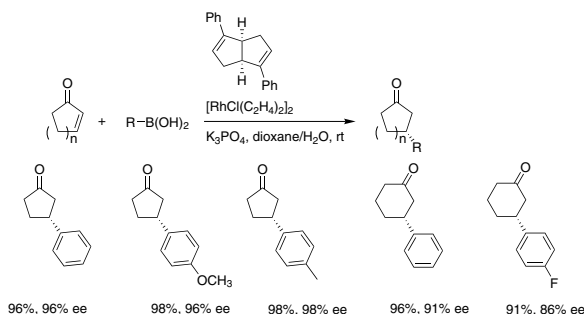
50 mg



Scheme 1



Scheme 2

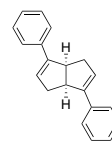


Scheme 3

(3*aS*,6*aS*)-3,6-Diphenyl-1,3*a*,4,6*a*-tetrahydropentalene, 97%

(3*aS*,6*aS*)-1,3*a*,4,6*a*-tetrahydro-3,6-diphenylpentalene [940280-80-8]

C₂₀H₁₈
FW 258.36



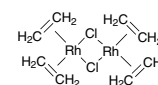
698733-50MG

50 mg

μ -Dichlorotetraethylene dirhodium(I)

[12081-16-2]

C₈H₁₆Cl₂Rh₂
FW 388.93



656763-100MG

100 mg

656763-1G

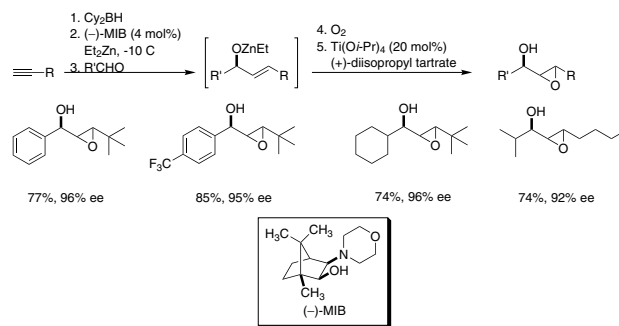
1 g

Asymmetric Epoxidation

Asymmetric Epoxidation Reaction Using MIB

Chiral epoxy alcohols are essential building blocks extensively used in the synthesis of natural products. However, the synthesis of these molecules proved to be challenging, especially allylic alcohols containing (*E*)-olefins. To access a plethora of chiral epoxy alcohols Professor Walsh and coworkers have invented a novel methodology,^{1,2} which complements the Sharpless asymmetric epoxidation. This process is mediated by a zinc catalyst built around Nugent's (–)-MIB ligand. Starting with a substituted terminal alkyne, the one pot reaction yields the desired epoxy alcohols with excellent levels of enantiocontrol, as well as yielding moderate to excellent diastereomeric ratios for a wide array of substrates.

References: (1) Jeon, S. -J. et al. *Org. Lett.* **2005**, 7, 1729. (2) Lurain, A. E. et al. *J. Org. Chem.* **2005**, 70, 1262.



Scheme 1

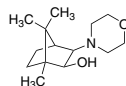
(2S)-3-exo-(Morpholino)isoborneol, 96 %

[(-)-MIB]

[287105-48-0]

C₁₄H₂₅NO₂

FW 239.35



676942-1G

1 g

676942-5G

5 g

CRC International Symposium in Stockholm

JSPS–Hokkaido University—Stockholm University

Cross-Coupling & Organometallics

November 3, 2008

**The Royal Swedish
Academy of Sciences
Stockholm**

For more information, visit
www.cat.hokudai.ac.jp

Opening Address:

Prof. Tamotsu Takahashi
Hokkaido University

Invited Speakers:

Prof. Jan-Erling Bäckvall
Stockholm University

Prof. Irina P. Beletskaya
Moscow State University

Prof. Tamejiro Hiyama
Kyoto University

Prof. Mats Larhed
Uppsala University

Prof. Shengming Ma
Shanghai Institute of Organic Chemistry

Prof. Ei-ichi Negishi
Purdue University

Prof. Michael G. Organ
York University

Prof. Akira Suzuki
Hokkaido University

Prof. Kohei Tamao
Riken

sigma-aldrich.com

SIGMA-ALDRICH®

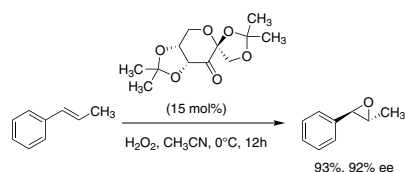
Asymmetric Epoxidation Using Shi Catalyst

Catalytic asymmetric epoxidation of alkenes has been the focus of many research efforts over the past two decades, the best known methods probably being those developed by Sharpless¹ and Jacobsen-Katsuki.² Shi has also developed a very efficient method for asymmetric epoxidation, using a ketone-derived organocatalyst based on D-fructose.³ This organocatalyst is able to epoxidize *trans* alkenes and certain *cis* alkenes with good to excellent yields and selectivities. More recently, Shi has achieved excellent results using hydrogen peroxide as an oxidant instead of Oxone, which allows a significant reduction in the amount of additional salts introduced and solvent used in the reaction (**Scheme 1**).⁴

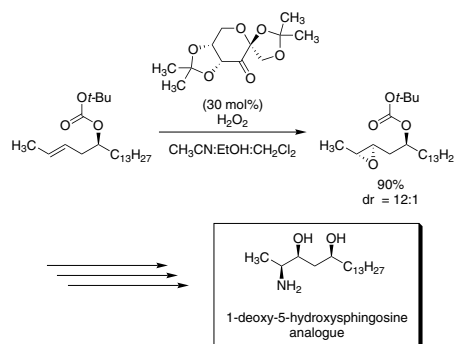
Other groups have used of Shi's methodology in pursuit of a number of unique structural moieties. For example, McDonald and co-workers recently reported a robust and selective synthesis of 2-amino-3,5-diols that employs the Shi epoxidation in a key step (**Scheme 2**).⁵ These 2-amino-3,5-diols are 1-deoxy-5-hydroxysphingosine analogues, which show promise as anticancer agents.

A recent example of the use of the Shi catalyst is the synthesis of thiosulfonate. Chiral sulfinyl derivatives have been of interest for the past three decades.⁶ This interest stemmed from the use of these molecules as chiral controllers in asymmetric synthesis, as chiral ligands and as Lewis bases. Khair et al. reported a new method to synthesize enantiomerically pure, functionalized, sterically demanding thiosulfonates.⁷ Investigating several catalysts for the oxidation of disulfides diesters and diethers, Khair et al. determined that a particular chiral pyranone derived from D-glucose, the Shi catalyst, yielded the best yields and stereoselectivities. Using the Shi catalyst in combination with oxone in acetonitrile and dimethoxymethane, the authors reported up to 89% yield and up to 96% ee's for the oxidation of disulfides (**Scheme 3**).

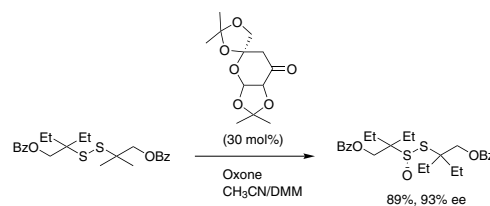
References: (1) Johnson, R. A.; Sharpless, K. B. In *Catalytic Asymmetric Synthesis* (2nd Ed.); Ojima, I., Ed.; Wiley-VCH: New York, 2000; Chapter 6A. (2) Jacobsen, E. N. In *Catalytic Asymmetric Synthesis* (1st Ed.); Ojima, I., Ed.; VCH: New York, 1993; Chapter 4.2. (3) Shi, Y. *Acc. Chem. Res.* **2004**, *37*, 488. (b) Frohn, M.; Shi, Y. *Synthesis* **2000**, 1979. (c) Wang, Z.-X. et al. *J. Am. Chem. Soc.* **1997**, *119*, 11224. (d) Tu, Y. et al. *J. Am. Chem. Soc.* **1996**, *118*, 9806. (4) Shu, L.; Shi, Y. *Tetrahedron* **2001**, *57*, 5213. (5) Wiseman, J. M. et al. *Org. Lett.* **2005**, *7*, 3155. (6) Fernandez, I.; Khair, N. *Chem. Rev.* **2003**, *95*, 1717. (7) Khair, N. et al. *Org. Lett.* **2007**, *9*, 1255.



Scheme 1



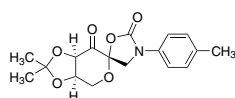
Scheme 2



Scheme 3

Shi Epoxidation Oxazolidinone Methyl Catalyst, 95%

(3a*R*,5'*S*,7a*R*)-Dihydro-2,2-dimethyl-3'-(4-methylphenyl)-spiro[6*H*-1,3-dioxolo[4,5-*c*]pyran-6,5'-oxazolidine]-2',7(4*H*)-dione [403501-30-4]

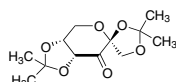


C₁₇H₁₉NO₆
FW 333.34

693456-100MG 100 mg
693456-500MG 500 mg

Shi Epoxidation Diketal Catalyst, 98%

1,2:4,5-Di-*O*-isopropylidene-β-D-*erythro*-2,3-hexodiulo-2,6-pyranose [18422-53-2]



C₁₂H₁₈O₆
FW 258.27

520160-5G 5 g

Got ChemNews?



A Monthly Chemistry E-Newsletter
from Sigma-Aldrich®

Visit sigma-aldrich.com/chemnews

sigma-aldrich.com

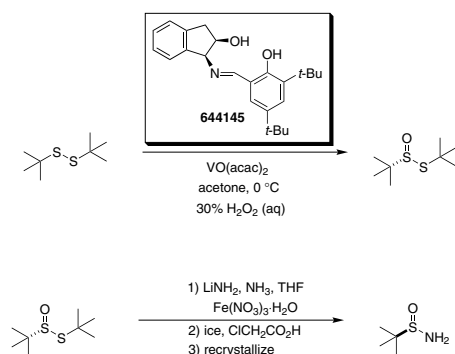
SIGMA-ALDRICH®

Asymmetric Oxidation

Asymmetric Oxidation of Di-*tert*-butyl Disulfide

The seminal work of Ellman and coworkers provided the first example of the catalytic asymmetric oxidation of *tert*-butyl disulfide. The desymmetrization involves reaction of *tert*-butyl disulfide with catalytic VO(acac)₂ and a chiral indanol ligand in the presence of hydrogen peroxide as the stoichiometric oxidant (**Scheme 1**). Upon monooxidation to provide *tert*-Butyl *tert*-butanethiosulfinate, the sulfur-sulfur bond is cleaved via displacement with LiNH₂ to afford the sulfonamide. The sulfonamides are often reacted with aldehydes to provide upon addition of organometallic reagents and deprotection, chiral amines (one of many uses of chiral sulfinimines). This sequence has been demonstrated to be effective on a large scale, utilizes relatively non-toxic solvents such as acetone (first step), and is economically feasible.

References: (1)(a) Weix, D. J.; Ellman, J. A. *Org. Lett.* **2003**, 5, 1317. (b) Weix, D. J.; Ellman, J. A. *Org. Syn.* **2005**, 82, 157. (c) Cogan, D. A.; Liu, G.; Kim, K.; Backes, B. J.; Ellman, J. A. *J. Am. Chem. Soc.* **1998**, 120, 8011. (d) Morton, D.; Stockman, R. A. *Tetrahedron* **2006**, 62, 8869.

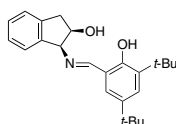


Scheme 1

(1*S*,2*R*)-1-[(3,5-Di-*tert*-butyl-2-hydroxybenzylidene)amino]-2-indanol, 97%

(1*S*,2*R*)-1-[(2-Hydroxy-3,5-di-*tert*-butylbenzylidene)amino]indan-2-ol
[212378-89-7]

C₂₄H₃₁NO₂
FW 365.51

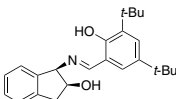


644145-1G	1 g
644145-5G	5 g

(1*R*,2*S*)-1-[(3,5-Di-*tert*-butyl-2-hydroxybenzylidene)amino]-2-indanol, 97%

(1*R*,2*S*)-1-[(2-Hydroxy-3,5-di-*tert*-butylbenzylidene)amino]indan-2-ol
[275374-67-9]

C₂₄H₃₁NO₂
FW 365.51

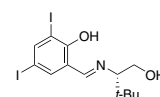


642959-1G	1 g
642959-5G	5 g

(*S*)-(-)-2-[(1-Hydroxy-3,3-dimethylbutan-2-ylimino)methyl]-4,6-diiodophenol, 97%

2-[[[(1*S*)-1-(Hydroxymethyl)-2,2-dimethylpropyl]imino]methyl]-4,6-diiodophenol
[477339-39-2]

C₁₃H₁₇I₂NO₂
FW 473.09



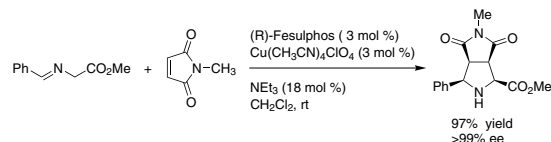
677558-100MG	100 mg
677558-500MG	500 mg

Asymmetric Cycloaddition

Asymmetric 1,3-dipolar Cycloaddition Using Fesulphos

The asymmetric 1,3-dipolar cycloaddition reaction is of the utmost importance for the enantioselective synthesis of five-membered heterocycles. Cabrera et al. introduced a new family of ligands consisting of a planar chiral *P,S*-ligand, named Fesulphos, for the 1,3-dipolar cycloaddition of azomethine halides.¹ The catalytic reaction is carried out with the Fesulphos ligand, a copper salt, and triethylamine in methylene chloride. This new catalytic system demonstrated complete enantiocontrol (ee >99%) with conversion up to 97% (**Scheme 1**).

Reference: (1) Cabrera, S. et al. *J. Am. Chem. Soc.* **2005**, 127, 16394.

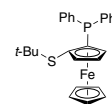


Scheme 1

(*R*_p)-2-(*tert*-Butylthio)-1-(diphenylphosphino)ferrocene, 98%

(1*S*)-1-(*tert*-Butylthio)-2-(diphenylphosphino)ferrocene;
(*R*)-FeSulPhos
[503859-61-8]

C₂₆H₂₇FePS
FW 458.38



687561-100MG	100 mg
687561-500MG	500 mg

Introducing Aldrich's **NEW** Chemical Synthesis Website

Check out our new website for synthetic chemistry. Designed and produced by chemists, the site is packed with the latest product/service information and productivity tools to accelerate your research success.

Featuring:

- **Chem Product Central**—40,000+ products categorized based on compound class and application area
 - Complete libraries of **Aldrichimica Acta** and **ChemFiles**
 - **Cheminars**®—Open access to seminal presentations on innovative synthetic technologies
 - **ChemBlogs**—Weblogs on latest hot papers
- ... and so much more!



If you love the Aldrich Handbook, you will love **sigma-aldrich.com/gochem**

Join our community and take our **Did You Know? Challenge**

Website registrants will receive a free T-shirt and be entered into our October drawings for a chance to win either a \$250 Apple® gift card or a \$250 Sigma-Aldrich® voucher. One winner will be randomly selected on each business day in October 2008.



To take the Challenge and register for the October drawings, visit **sigma-aldrich.com/gochem**

Terms and Conditions:

No purchase necessary. You must be over 18 to participate in this contest. The Did You Know? Challenge is limited to one entry per registration to our Community. Registrant must score 100% to be entered into the daily drawing for the Apple Gift Card or Aldrich voucher. Winners' names will be posted on the website each business day. Employees, officers and directors of Sigma-Aldrich and their respective family and/or household members are not eligible. Additional restrictions may apply. Void where prohibited. T-shirts, heterocyclic ring systems, numbering posters and ionic liquid charts are available while supplies last. Please allow 6 to 8 weeks for delivery. Apple Gift Cards can be applied only to qualified purchases directly from Apple at an Apple Store, the Apple Store online, or Apple Telesales (1-800-MY-APPLE) in the United States. Apple is a registered trademark of Apple Computer, Inc. All rights reserved. Apple is not a participant or sponsor of this promotion.

Argentina

SIGMA-ALDRICH DE ARGENTINA S.A.
Free Tel: 0810 888 7446
Tel: (+54) 11 4556 1472
Fax: (+54) 11 4552 1698

Australia

SIGMA-ALDRICH PTY LTD.
Free Tel: 1800 800 097
Free Fax: 1800 800 096
Tel: (+61) 2 9841 0555
Fax: (+61) 2 9841 0500

Austria

SIGMA-ALDRICH HANDELS GmbH
Tel: (+43) 1 605 81 10
Fax: (+43) 1 605 81 20

Belgium

SIGMA-ALDRICH NV/S.A.
Free Tel: 0800 14747
Free Fax: 0800 14745
Tel: (+32) 3 899 13 01
Fax: (+32) 3 899 13 11

Brazil

SIGMA-ALDRICH BRASIL LTDA.
Free Tel: 0800 701 7425
Tel: (+55) 11 3732 3100
Fax: (+55) 11 5522 9895

Canada

SIGMA-ALDRICH CANADA LTD.
Free Tel: 1800 565 1400
Free Fax: 1800 265 3858
Tel: (+1) 905 829 9500
Fax: (+1) 905 829 9292

China

SIGMA-ALDRICH (SHANGHAI)
TRADING CO. LTD.
Free Tel: 800 819 3336
Tel: (+86) 21 6141 5566
Fax: (+86) 21 6141 5567

Czech Republic

SIGMA-ALDRICH spol. s r. o.
Tel: (+420) 246 003 200
Fax: (+420) 246 003 291

Denmark

SIGMA-ALDRICH DENMARK A/S
Tel: (+45) 43 56 59 10
Fax: (+45) 43 56 59 05

Finland

SIGMA-ALDRICH FINLAND OY
Tel: (+358) 9 350 9250
Fax: (+358) 9 350 92555

France

SIGMA-ALDRICH CHIMIE S.à.r.l.
Free Tel: 0800 211 408
Free Fax: 0800 031 052
Tel: (+33) 474 82 28 00
Fax: (+33) 474 95 68 08

Germany

SIGMA-ALDRICH CHEMIE GmbH
Free Tel: 0800 51 55 000
Free Fax: 0800 64 90 000
Tel: (+49) 89 6513 0
Fax: (+49) 89 6513 1160

Greece

SIGMA-ALDRICH (O.M.) LTD.
Tel: (+30) 210 994 8010
Fax: (+30) 210 994 3831

Hungary

SIGMA-ALDRICH Kft
Ingyenes telefonszám: 06 80 355 355
Ingyenes fax szám: 06 80 344 344
Tel: (+36) 1 235 9055
Fax: (+36) 1 235 9050

India

SIGMA-ALDRICH CHEMICALS
PRIVATE LIMITED
Telephone
Bangalore: (+91) 80 6621 9600
New Delhi: (+91) 11 4358 8000
Mumbai: (+91) 22 2570 2364
Hyderabad: (+91) 40 4015 5488
Fax
Bangalore: (+91) 80 6621 9650
New Delhi: (+91) 11 4358 8001
Mumbai: (+91) 22 2579 7589
Hyderabad: (+91) 40 4015 5466

Ireland

SIGMA-ALDRICH IRELAND LTD.
Free Tel: 1800 200 888
Free Fax: 1800 600 222
Tel: +353 (0) 402 20370
Fax: + 353 (0) 402 20375

Israel

SIGMA-ALDRICH ISRAEL LTD.
Free Tel: 1 800 70 2222
Tel: (+972) 8 948 4100
Fax: (+972) 8 948 4200

Italy

SIGMA-ALDRICH S.r.l.
Numero Verde: 800 827018
Tel: (+39) 02 3341 7310
Fax: (+39) 02 3801 0737

Japan

SIGMA-ALDRICH JAPAN K.K.
Tel: (+81) 3 5796 7300
Fax: (+81) 3 5796 7315

Korea

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

Malaysia

SIGMA-ALDRICH (M) SDN. BHD
Tel: (+60) 3 5635 3321
Fax: (+60) 3 5635 4116

Mexico

SIGMA-ALDRICH QUÍMICA, S.A. de C.V.
Free Tel: 01 800 007 5300
Free Fax: 01 800 712 9920
Tel: 52 722 276 1600
Fax: 52 722 276 1601

The Netherlands

SIGMA-ALDRICH CHEMIE BV
Free Tel: 0800 022 9088
Free Fax: 0800 022 9089
Tel: (+31) 78 620 5411
Fax: (+31) 78 620 5421

New Zealand

SIGMA-ALDRICH NEW ZEALAND LTD.
Free Tel: 0800 936 666
Free Fax: 0800 937 777
Tel: (+61) 2 9841 0555
Fax: (+61) 2 9841 0500

Norway

SIGMA-ALDRICH NORWAY AS
Tel: (+47) 23 17 60 60
Fax: (+47) 23 17 60 50

Poland

SIGMA-ALDRICH Sp. z o.o.
Tel: (+48) 61 829 01 00
Fax: (+48) 61 829 01 20

Portugal

SIGMA-ALDRICH QUÍMICA, S.A.
Free Tel: 800 202 180
Free Fax: 800 202 178
Tel: (+351) 21 924 2555
Fax: (+351) 21 924 2610

Russia

SIGMA-ALDRICH RU, LLC
Tel: +7 (495) 621 6037
+7 (495) 621 5828
Fax: +7 (495) 621 5923

Singapore

SIGMA-ALDRICH PTE. LTD.
Tel: (+65) 6779 1200
Fax: (+65) 6779 1822

Slovakia

SIGMA-ALDRICH spol. s r. o.
Tel: (+421) 255 571 562
Fax: (+421) 255 571 564

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

Sweden

SIGMA-ALDRICH SWEDEN AB
Tel: (+46) 8 742 4200
Fax: (+46) 8 742 4243

Switzerland

SIGMA-ALDRICH CHEMIE GmbH
Free Tel: 0800 80 00 80
Free Fax: 0800 80 00 81
Tel: (+41) 81 755 2828
Fax: (+41) 81 755 2815

United Kingdom

SIGMA-ALDRICH COMPANY LTD.
Free Tel: 0800 717 181
Free Fax: 0800 378 785
Tel: (+44) 1747 833 000
Fax: (+44) 1747 833 313
SAFC (UK) Tel: 01202 712305

United States

SIGMA-ALDRICH
P.O. Box 14508
St. Louis, Missouri 63178
Toll-Free: 800 325 3010
Toll-Free Fax: 800 325 5052
Call Collect: (+1) 314 771 5750
Tel: (+1) 314 771 5765
Fax: (+1) 314 771 5757

Internet

sigma-aldrich.com



Mixed Sources
Product group from well-managed
forests, controlled sources and
recycled wood or fiber
www.fsc.org Cert no. SGS-COC-003338
© 1996 Forest Stewardship Council

**World Headquarters**

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

Order/Customer Service (800) 325-3010 • Fax (800) 325-5052

Technical Service (800) 325-5832 • sigma-aldrich.com/techservice

Development/Bulk Manufacturing Inquiries SAFC™ (800) 244-1173

©2008 Sigma-Aldrich Co. All rights reserved. SIGMA, SAFC, SAFC™, SIGMA-ALDRICH, ALDRICH, FLUKA, SUPELCO, are trademarks belonging to Sigma-Aldrich Co. and its affiliate Sigma-Aldrich Biotechnology, L.P. 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. Eppendorf is a registered trademark of Eppendorf-Netheler-Hinz GmbH. Sepharose is a registered trademark of GE Healthcare. Coomassie is a registered trademark of Imperial Chemical Industries Ltd. ProteoSilver is a trademark of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co. Cheminars and TraceSELECT are registered trademarks of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co.

LAH
70636-504740
0108

*Accelerating Customers'
Success through Leadership
in Life Science, High
Technology and Service*

SIGMA-ALDRICH®

Sigma-Aldrich
3050 Spruce Street
St. Louis, MO 63103 USA
sigma-aldrich.com