

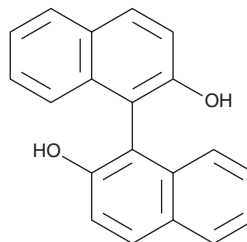
# PRODUCT INFORMATION



## (S)-(-)-1,1'-Bi-2-naphthol

Item No. 70052

**CAS Registry No.:** 18531-99-2  
**Formal Name:** (S)-(1,1')-binaphthalene)-2,2'-diol  
**Synonym:** (S)-BINOL  
**MF:** C<sub>20</sub>H<sub>14</sub>O<sub>3</sub>  
**FW:** 286.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 227 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

(S)-(-)-1,1'-bi-2-naphthol (S-BINOL) is supplied as a crystalline solid. A stock solution may be made by dissolving the S-BINOL in an organic solvent purged with an inert gas. S-BINOL is soluble in organic solvents such as ethanol (22 mg/ml), dichloromethane (19 mg/ml), dioxane (29 mg/ml), and tetrahydrofuran (30 mg/ml).

### Description

S-BINOL is a chiral auxiliary for the asymmetric reduction of 15-keto prostaglandin intermediates to 15(S)-alcohols. S-BINOL forms a complex with lithium aluminum hydride solution of ethanol. This is a versatile reagent for the asymmetric reduction of α,β-unsaturated enones at lower temperatures to the corresponding S-alcohols in ≥90% ee.

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/21/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM