

# PRODUCT INFORMATION



## ent-Corey PG-Lactone Diol

Item No. 10653

CAS Registry No.: 53110-06-8

Formal Name: (6aR)-5S-hydroxy-4S-((R,E)-3aS-hydroxyoct-1-en-1-yl)hexahydro-2H-cyclopenta[b]furan-2-one

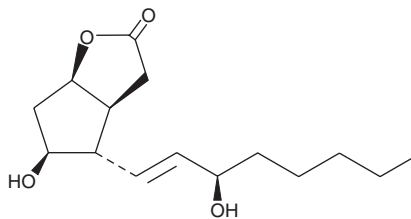
MF: C<sub>15</sub>H<sub>24</sub>O<sub>4</sub>

FW: 268.4

Purity: ≥98%

Stability: ≥1 year at -20°C

Supplied as: A solution in methyl acetate



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

### Laboratory Procedures

ent-Corey PG-lactone diol is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of ent-corey PG-lactone diol in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of ent-corey PG-lactone diol is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of ent-corey PG-lactone diol in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Corey PG-lactone diol is a versatile, chiral intermediate used in the preparation of PGs and PG analogs. ent-Corey PG-lactone diol is the opposite enantiomer of corey PG-lactone diol.

#### 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, 01/07/2021

#### 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