

PRODUCT INFORMATION



TBS-Corey Lactone Aldehyde

Item No. 70031

CAS Registry No.: 64091-14-1
Formal Name: [3aR-(3aα,4α,5β,6aα)]-5-tert-butyl-dimethylsilyl)hexahydro-2-oxo-2H-cyclopenta[b]furan-4-carboxaldehyde

MF: C₁₄H₂₄O₄Si

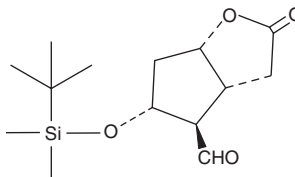
FW: 284.4

Purity: ≥95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥1 year



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

Laboratory Procedures

TBS-Corey lactone aldehyde is supplied as a crystalline solid. A stock solution may be made by dissolving the TBS-Corey lactone aldehyde in an organic solvent purged with an inert gas. TBS-Corey lactone aldehyde is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TBS-Corey lactone aldehyde in these solvents is approximately 20 mg/ml.

TBS-Corey lactone aldehyde is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TBS-Corey lactone aldehyde should first be dissolved in DMF and then diluted with the aqueous buffer of choice. TBS-Corey lactone aldehyde has a solubility of approximately 50 µg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

TBS-Corey lactone aldehyde is a versatile, hydroxyl-protected intermediate for the synthesis of prostaglandins and prostaglandin analogs. Prostaglandin derivatives with modifications in the alpha chain, the omega chain, or both, are readily accessible from TBS-Corey lactone aldehyde.¹

Reference

1. Bindra, J.S. and Bindra, R. Prostaglandin synthesis. *Academic Press, Inc., New York* (1977).

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, 12/13/2018

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