

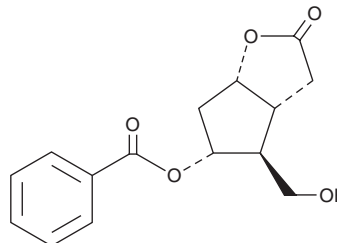
PRODUCT INFORMATION



Corey Lactone Benzoate

Item No. 10007983

CAS Registry No.: 39746-00-4
Formal Name: (3aR,4S,5R,6aS)-5-(benzoyloxy)hexahydro-4-(hydroxymethyl)-2H-cyclopenta[b]furan-2-one
Synonyms: (-)-Corey Lactone Benzoate,
(-)-Corey Lactone Benzoate Alcohol
MF: C₁₅H₁₆O₅
FW: 276.3
Purity: ≥98%
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

Corey lactone benzoate is supplied as a crystalline solid. A stock solution may be made by dissolving the corey lactone benzoate in the solvent of choice, which should be purged with an inert gas. Corey lactone benzoate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of corey lactone benzoate in these solvents is approximately 39.7, 25.3, and 40.3 mg/ml, respectively.

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. Organic solvent-free aqueous solutions of corey lactone benzoate can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of corey lactone benzoate in PBS (pH 7.2) is approximately 60 µg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Corey lactone benzoate is a precursor that has been used in the synthesis of prostaglandins (PGs), as well as the agonist of the PGE₂ receptor subtype EP₄ rivenprost (Item No. 13618).^{1,2}

References

1. Fleming, I. and Winter, S.B.D. Stereocontrol in organic synthesis using silicon-containing compounds. A formal synthesis of prostaglandins controlling the stereochemistry at C-15 using a silyl-to-hydroxy conversion following a stereochemically convergent synthesis of an allylsilane. *J. Chem. Soc. Perkin Trans. 1* **17**, 2687-2700 (1998).
2. Ohta, C., Kuwabe, S.i., Shiraishi, T., et al. An improved synthesis of the selective EP₄ receptor agonist ONO-4819. *J. Org. Chem.* **74(21)**, 8298-8308 (2009).

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, 05/15/2023

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