

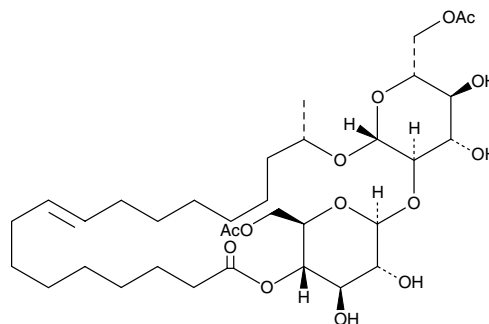
Product Information



1',4''-Sophorolactone 6',6''-diacetate

Item No. 14718

CAS Registry No.: 148409-20-5
Formal Name: 17S-[[[6-O-acetyl-2-O-(6-O-acetyl-β-D-glucopyranosyl)-β-D-glucopyranosyl]oxy]-9Z-octadecenoic acid, 1,4''-ester
MF: C₃₄H₅₆O₁₄
FW: 688.8
Purity: ≥80%
Stability: ≥1 year at room temperature
Supplied as: A white to slightly brownish powder



Laboratory Procedures

For long term storage, we suggest that 1',4''-sophorolactone 6',6''-diacetate be stored as supplied at room temperature. It should be stable for at least one year.

1',4''-Sophorolactone 6',6''-diacetate is supplied as a white to slightly brownish powder. A stock solution may be made by dissolving the 1',4''-sophorolactone 6',6''-diacetate in the solvent of choice. 1',4''-Sophorolactone 6',6''-diacetate is soluble in organic solvents such as ethanol, methanol, and ethyl acetate, which should be purged with an inert gas.

1',4''-Sophorolactone 6',6''-diacetate is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

1',4''-Sophorolactone 6',6''-diacetate is a major glycolipid produced from fermentation of sugar by yeasts including *C. bombicola*. This lactone form of sophorolipid has found use as a biosurfactant, possessing lower surface tension compared to the acid form of sophorolipids.¹

Reference

1. Nguyen, T.T. and Sabatini, D.A. Characterization and emulsification properties of rhamnolipid and sophorolipid biosurfactants and their applications. *Int. J. Mol. Sci.* **12(2)**, 1232-1244 (2011).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/14718

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 08/21/2013

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com