

PRODUCT INFORMATION

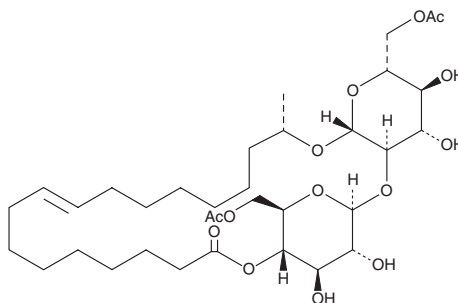


1',4''-Sophorolactone 6',6''-diacetate

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%
Supplied as: A powder
Storage: Room temperature
Stability: ≥4 years



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

Laboratory Procedures

1',4''-Sophorolactone 6',6''-diacetate is supplied as a powder. A stock solution may be made by dissolving the 1',4''-sophorolactone 6',6''-diacetate in the solvent of choice, which should be purged with an inert gas. 1',4''-Sophorolactone 6',6''-diacetate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 1',4''-sophorolactone 6',6''-diacetate in these solvents is approximately 30 mg/ml.

1',4''-Sophorolactone 6',6''-diacetate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1',4''-sophorolactone 6',6''-diacetate should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 1',4''-Sophorolactone 6',6''-diacetate has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

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).

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, 12/15/2022

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