

PRODUCT INFORMATION



MEL-B

Item No. 34659

Formal Name: (2R,3S,4S,5R,6R)-6-(acetoxymethyl)-5-hydroxy-2-((2S,3R)-2,3,4-trihydroxybutoxy) tetrahydro-2H-pyran-3,4-diyl dipropionate

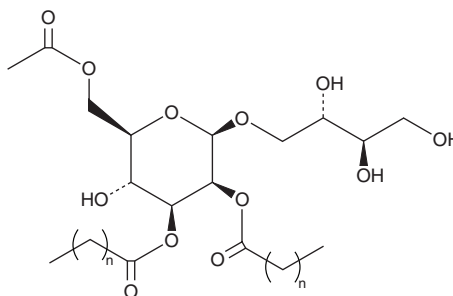
Synonym: Mannosylerythritol Lipid B

Purity: ≥95% (mixture of congeners)

Supplied as: A neat oil

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

MEL-B is supplied as a neat oil. A stock solution may be made by dissolving the MEL-B in the solvent of choice, which should be purged with an inert gas. MEL-B is soluble (≥10 mg/ml) in ethanol.

Description

MEL-B is a glycolipid biosurfactant that is produced from vegetable oils by *Pseudozyma* yeasts.¹ Mixtures of MEL-B isoforms have been used as moisturizing agents in cosmetics and are active against *S. aureus*. This product is a mixture of MEL-B isoforms with variable fatty acid chain lengths.

Reference

1. Yamauchi, S., Furukawa, M., Kawahara, A., *et al.* Roles of mannosylerythritol lipid-B components in antimicrobial activity against bovine mastitis-causing *Staphylococcus aureus*. *World J. Microbiol. Biotechnol.* **38**(3), 54 (2022).

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, 05/12/2025

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