

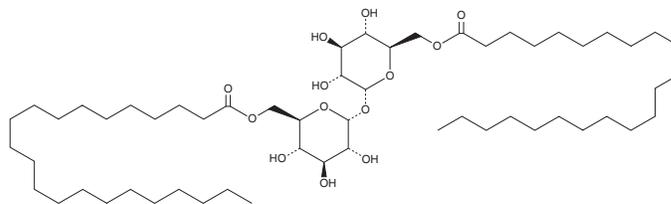
PRODUCT INFORMATION



Trehalose 6,6'-Dibehenate

Item No. 37911

CAS Registry No.: 66758-35-8
Formal Name: 6-docosanoate 6-O-(1-oxodocosyl)- α -D-glucopyranosyl α -D-glucopyranoside
Synonyms: TDB, Trehalose Dibehenate, 22:0 Trehalose
MF: C₅₆H₁₀₆O₁₃
FW: 987.4
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Trehalose 6,6'-dibehenate (TDB) is supplied as a solid. A stock solution may be made by dissolving the TDB in the solvent of choice, which should be purged with an inert gas. TDB is slightly soluble in DMSO and dimethyl formamide.

Description

TDB is a synthetic glycolipid and derivative of trehalose 6,6'-dimycolate (TDM), also known as mycobacterial cord factor.¹ It binds to a fusion protein of mouse macrophage-inducible C-type lectin (Mincle) and human IgG1 Fc in a cell-free assay when used at a concentration of 10 μ g/ml. TDB (5 μ g/ml) induces nitric oxide (NO) and G-CSF production in mouse bone marrow macrophages in a Mincle-dependent manner. Liposomes containing TDB and dimethyldioctadecylammonium (DDA; Item No. 26120) as adjuvants and the mycobacterial lipid antigens diacylated sulfoglycolipid (Ac₂SGL) and phosphatidyl-*myo*-inositol dimannoside (PIM₂) decrease spleen bacterial load and the number of lung and spleen lesions in a guinea pig model of *M. tuberculosis* infection.²

References

1. Huber, A., Kallerup, R.S., Korsholm, K.S., *et al.* Trehalose diester glycolipids are superior to the monoesters in binding to Mincle, activation of macrophages *in vitro* and adjuvant activity *in vivo*. *Innate Immun.* **22(6)**, 405-418 (2016).
2. Larrouy-Maumus, G., Layre, F., Clark, S., *et al.* Protective efficacy of a lipid antigen vaccine in a guinea pig model of tuberculosis. *Vaccine* **35(10)**, 1395-1402 (2017).

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, 05/12/2023

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