

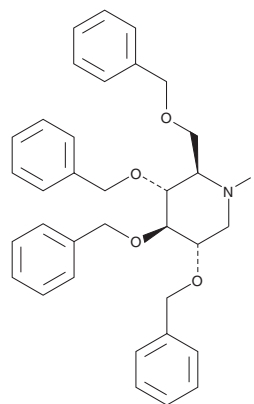
PRODUCT INFORMATION



Deoxynojirimycin Tetrabenzyl Ether

Item No. 10011914

CAS Registry No.: 69567-11-9
Formal Name: 3R,4R,5S-tris(phenylmethoxy-2R-[(phenylmethoxy)methyl]-piperidine
Synonyms: DNJ Tetrabenzyl Ether, dNM Tetrabenzyl Ether
MF: C₃₄H₃₇NO₄
FW: 523.7
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

Deoxynojirimycin (dNM) tetrabenzyl ether is supplied as a crystalline solid. A stock solution may be made by dissolving the dNM tetrabenzyl ether in the solvent of choice, which should be purged with an inert gas. dNM tetrabenzyl ether is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of dNM tetrabenzyl ether in these solvents is approximately 0.25, 20, and 15 mg/ml, respectively.

dNM tetrabenzyl ether is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, dNM tetrabenzyl ether should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. dNM tetrabenzyl ether has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

dNM tetrabenzyl ether is an intermediate for the synthesis of glucosylceramide synthase inhibitors such as 1-dNM, a glucose analog that potently inhibits α -glucosidase I and II.¹

Reference

1. Matos, C.R.R., Lopes, R.S.C., and Lopes, C.C. Synthesis of 1-deoxynojirimycin and N-butyl-1-deoxynojirimycin. *Synthesis* 4, 571-573 (1999).

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, 04/05/2024

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