

PRODUCT INFORMATION



N-Glycine Glucosylsphingosine (d18:1)

Item No. 24473

Formal Name: 2-amino-N-((2S,3R,E)-3-hydroxy-1-(((2R,3R,4S,5S,6R)-3,4,5-trihydroxy-6-(hydroxymethyl)tetrahydro-2H-pyran-2-yl)oxy)octadec-4-en-2-yl)acetamide

Synonyms: N-Glycinated Glucosylsphingosine, N-Glycinated 1-β-D-lyso-Glucosylceramide

MF: C₂₆H₅₀N₂O₈

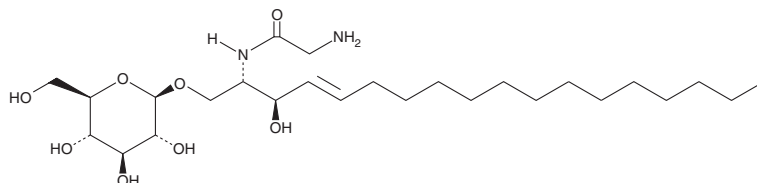
FW: 518.7

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

N-Glycine glucosylsphingosine (d18:1) is supplied as a solid. A stock solution may be made by dissolving the N-glycine glucosylsphingosine (d18:1) in the solvent of choice. N-Glycine glucosylsphingosine (d18:1) is soluble in chloroform, methanol, and an 8:2 solution of DMSO:DMF.

Description

N-glycine Glucosylsphingosine is a glycinated form of 1-β-D-glucosylsphingosine (Item No. 23211).

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, 04/23/2018

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