

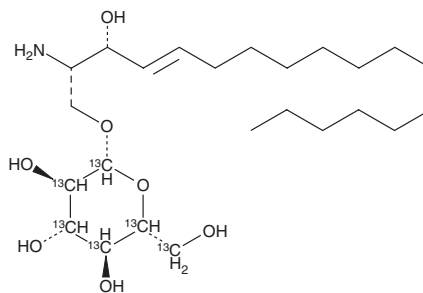
PRODUCT INFORMATION



$^{13}\text{C}_6$ Glucosylsphingosine (d18:1)

Item No. 23212

CAS Registry No.: 299172-48-8
Formal Name: 2S-amino-3R-hydroxy-4E-octadecen-1-yl,
 β -D-glucopyranoside-1,2,3,4,5,6- $^{13}\text{C}_6$
Synonyms: 1- β -D-Glucosyl-1,2,3,4,5,6- $^{13}\text{C}_6$ -sphingosine,
 $^{13}\text{C}_6$ -Glucosylsphingosine
MF: $\text{C}_{18}[^{13}\text{C}]_6\text{H}_{47}\text{NO}_7$
FW: 467.6
Purity: $\geq 98\%$
Supplied as: A 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

$^{13}\text{C}_6$ Glucosylsphingosine (d18:1) is supplied as a solid. A stock solution may be made by dissolving the $^{13}\text{C}_6$ glucosylsphingosine (d18:1) in the solvent of choice. $^{13}\text{C}_6$ Glucosylsphingosine (d18:1) is soluble in ethanol, methanol, and a 2:1 solution of chloroform:methanol.

Description

$^{13}\text{C}_6$ Glucosylsphingosine is an isotopically enriched form of 1- β -D-glucosylsphingosine (d18:1) (Item No. 23211) that is intended for use as an internal standard for the quantification of 1- β -D-glucosylsphingosine by GC- or LC-MS. 1- β -D-Glucosylsphingosine is a lysolipid derivative of glucosylcerebroside that decreases activity of glucocerebrosidase in LA-N-2 cells in a dose-dependent manner.^{1,2}

References

1. Murugesan, V., Chuang, W.-L., Liu, J., *et al.* Glucosylsphingosine is a key biomarker of Gaucher disease. *Am. J. Hematol.* **91(11)**, 1082-1089 (2016).
2. Schueler, U.H., Kolter, T., Kaneski, C.R., *et al.* Toxicity of glucosylsphingosine (glucopsychosine) to cultured neuronal cells: A model system for assessing neuronal damage in Gaucher disease type 2 and 3. *Neurobiol. Dis.* **14(3)**, 595-601 (2003).

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, 09/22/2022

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