

PRODUCT INFORMATION



C6 Glucosylceramide (d18:2/6:0)

Item No. 24474

Formal Name: N-((2S,3R,4E,8E)-3-hydroxy-1-(((2R,3R,4S,5S,6R)-3,4,5-trihydroxy-6-(hydroxymethyl)tetrahydro-2H-pyran-2-yl)oxy)octadeca-4,8-dien-2-yl)hexanamide

Synonyms: GluCer(d18:2/6:0), Glucosylceramide (d18:2/6:0), N-hexanoyl Glucosylceramide

MF: C₃₀H₅₅NO₈

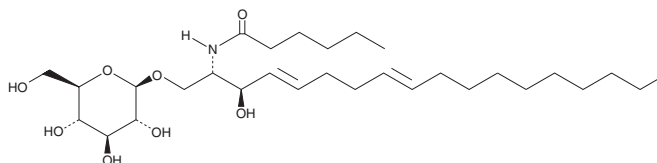
FW: 557.8

Purity: ≥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

C6 Glucosylceramide (d18:2/6:0) is supplied as a solid. A stock solution may be made by dissolving the C6 glucosylceramide (d18:2/6:0) in the solvent of choice, which should be purged with an inert gas. C6 Glucosylceramide (d18:2/6:0) is soluble in organic solvents such as chloroform, methanol, DMSO, and dimethyl formamide.

Description

C6 Glucosylceramide is a glucosylated form of C6 ceramide (Item No. 62525). As this product is derived from a natural source, there may be variations in the sphingoid backbone.

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, 11/10/2022

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