

# PRODUCT INFORMATION



## C6 Biotin Glucosylceramide (d18:1/6:0)

Item No. 23208

**CAS Registry No.:** 2692623-22-4  
**Formal Name:** (3aS,4S,6aR)-N-[6-[[[(1S,2R,3E,7E)-1-[(β-D-glucopyranosyloxy)methyl]-2-hydroxy-3,7-heptadecadien-1-yl]amino]-6-oxohexyl]hexahydro-2-oxo-1H-thieno[3,4-d]imidazole-4-pentanamide

**Synonyms:** Biotin GluCer(18:1/6:0), Biotin Glucosylceramide (d18:1/6:0), C6 Glucosylceramide-biotin, N-C6:0-Glucosylceramide-biotin, N-C6:0-β-D-Glucosylsphingosine-biotin, N-Hexanoyl-biotin-Glucosylceramide

**MF:** C<sub>40</sub>H<sub>70</sub>N<sub>4</sub>O<sub>10</sub>S

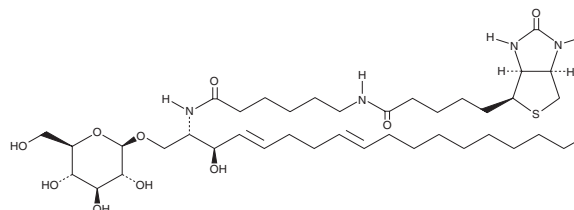
**FW:** 799.1

**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 Biotin glucosylceramide (d18:1/6:0) is supplied as a solid. A stock solution may be made by dissolving the C6 biotin glucosylceramide (d18:1/6:0) in the solvent of choice. C6 Biotin glucosylceramide (d18:1/6:0) is soluble in methanol, dimethyl formamide, and in a 2:1 solution of chloroform:methanol.

### Description

C6 Biotin glucosylceramide is an affinity probe that allows C6 glucosylceramide to be detected or immobilized through interaction with the biotin ligand. 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, 12/21/2022

#### 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