

PRODUCT INFORMATION



C6 Biotin Ganglioside G_{M3} (d18:1/6:0) (ammonium salt)

Item No. 31590

Formal Name: (2S,4S,5R,6R)-5-acetamido-2-(((2S,3R,4S,5S,6R)-2-(((2R,3S,4R,5R,6R)-4,5-dihydroxy-6-(((2S,3R,E)-3-hydroxy-2-(6-(5-((3aS,4S,6aR)-2-oxohexahydro-1H-thieno[3,4-d]imidazol-4-yl)pentanamido)hexanamido)octadec-4-en-1-yl)oxy)-2-(hydroxymethyl)tetrahydro-2H-pyran-3-yl)oxy)-3,5-dihydroxy-6-(hydroxymethyl)tetrahydro-2H-pyran-4-yl)oxy)-4-hydroxy-6-((1R,2R)-1,2,3-trihydroxypropyl)tetrahydro-2H-pyran-2-carboxylate, monoammonium salt

Synonyms: Biotin-C6:0-G_{M3},
N-hexanoyl-Biotin-Monosialoganglioside G_{M3}

MF: C₅₇H₉₈N₅O₂₃S • NH₄

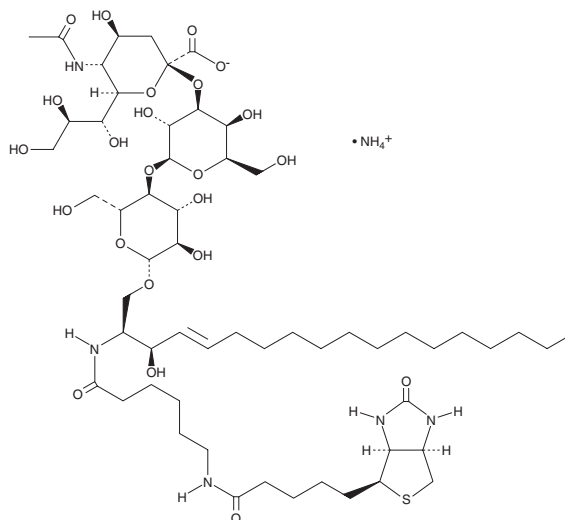
FW: 1,271.5

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 ganglioside G_{M3} (d18:1/6:0) (ammonium salt) is supplied as a solid. A stock solution may be made by dissolving the C6 biotin ganglioside G_{M3} (d18:1/6:0) (ammonium salt) in the solvent of choice, which should be purged with an inert gas. C6 Biotin ganglioside G_{M3} (d18:1/6:0) (ammonium salt) is soluble in 2:1:0.1 solution of chloroform:methanol:water. C6 Biotin ganglioside G_{M3} (d18:1/6:0) (ammonium salt) is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

C6 Biotin ganglioside G_{M3} is an affinity probe that allows ganglioside G_{M3} (Item No. 15587) to be detected or immobilized through interaction with the biotin ligand. As this product is derived from a natural source, there may also 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, 12/22/2022

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