

PRODUCT INFORMATION



C6 Biotin Ganglioside G_{T1b} (d18:1/6:0) (ammonium salt)

Item No. 40501

Synonyms: Biotin-C6:0-G_{T1b},
N-hexanoyl-Biotin-Ganglioside G1,
N-hexanoyl-Biotin-Trisialoganglioside G_{T1b}

MF: C₉₃H₁₆₅N₁₁O₄₉S

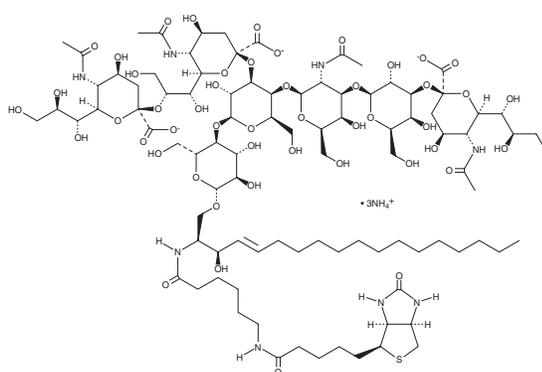
FW: 2,253.4

Purity: ≥95% (mixture of sphingoid chains)

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.

Description

C6 Biotin Ganglioside G_{T1b} (d18:1/6:0) is an affinity probe that allows ganglioside G_{T1b} (Item Nos. 31712 | 31592 | 15888) 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.

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