

PRODUCT INFORMATION



Lyso-Monosialoganglioside G_{M3} (synthetic)

Item No. 42762

CAS Registry No.: 94458-63-6

Formal Name: 2S-amino-3R-hydroxy-4E-octadecen-1-yl O-(N-acetyl- α -neuraminosyl)-(2 $\rightarrow$ 3)-O- β -D-galactopyranosyl-(1 $\rightarrow$ 4)- β -D-glucopyranoside

Synonyms: Lyso-G_{M3}, Lysoganglioside G_{M3}, lyso-Monosialoganglioside G_{M3}

MF: C₄₁H₇₄N₂O₂₀

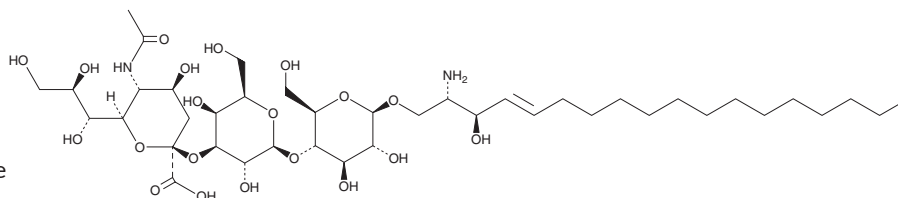
FW: 915.0

Purity: $\geq$ 98%

Supplied as: A solution in methanol

Storage: -20°C

Stability: $\geq$ 1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Lyso-monosialoganglioside G_{M3} (Lyso-G_{M3}) is a form of ganglioside G_{M3} (Item Nos. 31711 | 15587) that is lacking the fatty acyl group.¹ It inhibits EGF-induced increases in EGFR kinase activity in A431 epithelial cancer cells when used at a concentration of 500 μ M. This product is fully synthetic and has no variations in the sphingoid backbone.

Reference

1. Hanai, N., Nores, G., Torres-Méndez, C.R., *et al.* Modified ganglioside as a possible modulator of transmembrane signaling mechanism through growth factor receptors: A preliminary note. *Biochem. Biophys. Res. Commun.* **147**(1), 127-134 (1987).

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, 01/23/2025

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