

PRODUCT INFORMATION



C12 Lissamine-rhodamine Galactosylceramide (d18:1/12:0)

Item No. 24470

Formal Name: 2-(6-(diethylamino)-3-(diethyliminio)-3H-xanthen-9-yl)-5-(N-(12-(((2S,E)-3R-hydroxy-1-((3R,4S,5R-trihydroxy-6R-(hydroxymethyl)tetrahydro-2H-pyran-2R-yl)oxy)octadec-4-en-2-yl)amino)-12-oxododecyl)sulfamoyl)benzenesulfonate

Synonyms: Lissamine-rhodamine B-dodecanoyl Galactosylceramide, Lissamine-rhodamine GalCer(d18:1/12:0)

MF: C₆₃H₉₈N₄O₁₄S₂

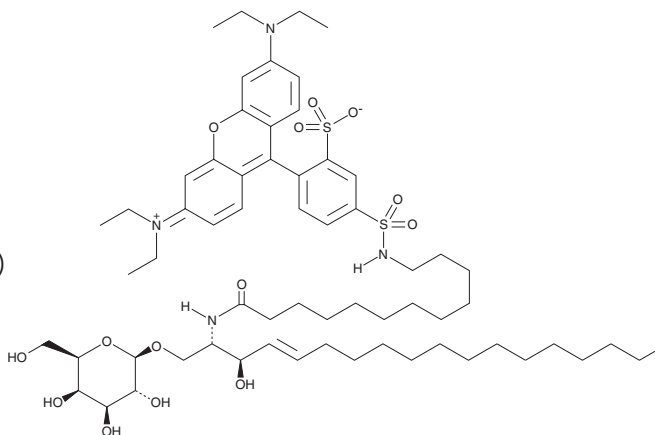
FW: 1,199.6

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

C12 Lissamine-rhodamine galactosylceramide (d18:1/12:0) is supplied as a solid. A stock solution may be made by dissolving the C12 lissamine-rhodamine galactosylceramide (d18:1/12:0) in the solvent of choice, which should be purged with an inert gas. C12 Lissamine-rhodamine galactosylceramide (d18:1/12:0) is soluble in DMSO, dimethyl formamide, and an 8:2 solution of chloroform:methanol.

Description

C12 Lissamine-rhodamine galactosylceramide is a fluorescently labeled version of C12 galactosylceramide (Item No. 24347). Lissamine-rhodamine B displays excitation and emission maxima of 560 and 575 nm, respectively.¹ As this product is derived from a natural source, there may be variations in the sphingoid backbone.

Reference

1. Glunde, K., Foss, C.A., Takagi, T., *et al.* Synthesis of 6'-O-lissamine-rhodamine B-glucosamine as a novel probe for fluorescence imaging of lysosomes in breast tumors. *Bioconjug. Chem.* **16**(4), 843-851 (2005).

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, 12/17/2022

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