

PRODUCT INFORMATION



N-Glycine 3'-sulfo Galactosylsphingosine

Item No. 27200

CAS Registry No.: 2620491-81-6
Formal Name: (2R,3R,4S,5S,6R)-2-(((2S,3R,E)-2-(2-aminoacetamido)-3-hydroxyoctadec-4-en-1-yl)oxy)-3,5-dihydroxy-6-(hydroxymethyl) tetrahydro-2H-pyran-4-yl hydrogen sulfate

Synonyms: N-Glycinated *lyso*-sulfatide, N-Glycinated 3'-sulfo Galactosylsphingosine, N-Glycine Lyso 3'-sulfo Galactosylceramide, N-Glycine *lyso*-sulfatide, N-Glycine Sphingosine-1-Galactoside-3-Sulfate, N-Glycine Sulfogalactosylsphingosine

MF: C₂₆H₅₀N₂O₁₁S

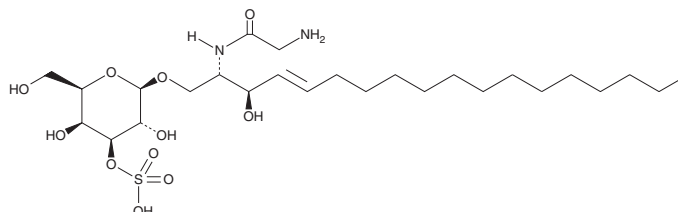
FW: 598.7

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

N-Glycine 3'-sulfo galactosylsphingosine is supplied as a solid. A stock solution may be made by dissolving the N-glycine 3'-sulfo galactosylsphingosine in the solvent of choice, which should be purged with an inert gas. N-Glycine 3'-sulfo galactosylsphingosine is soluble in ethanol and a 70:30:4 solution of chloroform:methanol:water.

Description

N-Glycine 3'-sulfo galactosylsphingosine is a glycinated form of 3'-sulfo galactosylsphingosine (Item No. 25316). 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, 02/03/2023

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