

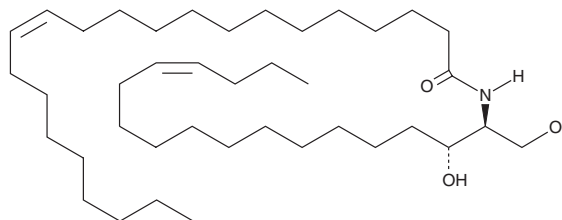
PRODUCT INFORMATION



C22:1 Ceramide (d18:1(14Z)/22:1(13Z))

Item No. 28603

Formal Name: (Z)-N-((2S,3R,Z)-1,3-dihydroxyoctadec-14-en-2-yl)docos-13-enamide
Synonyms: Cer(d18:1(14Z)/22:1(13Z)), C22:1-Cer, Ceramide (d18:1(14Z)/22:1(13Z)), N-erucoyl-4,5-dihydro-14Z-Sphingosine
MF: C₄₀H₇₇NO₃
FW: 620.1
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

C22:1 Ceramide (d18:1(14Z)/22:1(13Z)) is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. The solvent dimethyl formamide purged with an inert gas can be used. The solubility of C22:1 ceramide (d18:1(14Z)/22:1(13Z)) in this solvent is approximately 0.5 mg/ml.

Description

C22:1 Ceramide (d18:1(14Z)/22:1(13Z)) is a long-chain atypical ceramide containing a sphingosine (d18:1(14Z)) (Item No. 27019) backbone with a *cis* double bond at the 14-15 position rather than a *trans* double bond at the 4-5 position.

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, 01/30/2024

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