

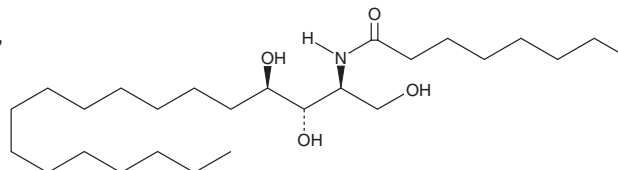
PRODUCT INFORMATION



C8 Phytoceramide (t18:0/8:0)

Item No. 24357

CAS Registry No.: 249728-93-6
Formal Name: N-[(1S,2S,3R)-2,3-dihydroxy-1-(hydroxymethyl)heptadecyl]-octanamide
Synonyms: Cer(t18:0/8:0), Ceramide (t18:0/8:0), N-Octanoyl Phytosphingosine, C8 Phytosphingosine (t18:0/8:0)
MF: C₂₆H₅₃NO₄
FW: 443.7
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

C8 Phytoceramide (t18:0/8:0) is supplied as a solid. A stock solution may be made by dissolving the C8 phytoceramide (t18:0/8:0) in the solvent of choice. C8 Phytoceramide (t18:0/8:0) is soluble in ethanol and methanol and a 1:1 (warm) solution of chloroform:methanol.

Description

C8 Phytoceramide is a cell-permeable analog of phytosphingosine (Item No. 20217). It is composed of a phytosphingosine backbone amine-linked to a C8 fatty acid chain. C8 Phytoceramide has been used as a standard for the determination of ceramides in commercial cosmetics using LC-MS.¹

Reference

1. Lee, M.H., Lee, G.H., and Yoo, J.S. Analysis of ceramides in cosmetics by reversed-phase liquid chromatography/electrospray ionization mass spectrometry with collision-induced dissociation. *Rapid Commun. Mass Spectrom.* **17**(1), 64-75 (2003).

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, 12/19/2022

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