

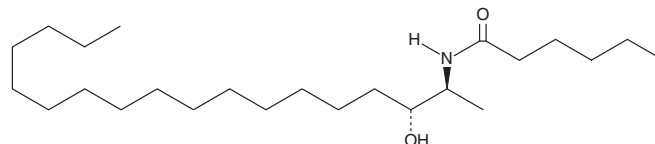
PRODUCT INFORMATION



C6 dihydro 1-Deoxyceramide (m18:0/6:0)

Item No. 25494

Formal Name: N-((2S,3R)-3-hydroxyoctadecan-2-yl)hexanamide
Synonyms: C6 dihydro Ceramide (m18:0/6:0), Cer(m18:0/6:0), N-hexanoyl-1-deoxy-Sphinganine
MF: C₂₄H₄₉NO₂
FW: 383.7
Purity: ≥98%
Supplied as: A crystalline 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

C6 dihydro 1-Deoxyceramide (m18:0/6:0) is supplied as a crystalline solid. A stock solution may be made by dissolving the C6 dihydro 1-deoxyceramide (m18:0/6:0) in the solvent of choice. C6 dihydro 1-Deoxyceramide (m18:0/6:0) is soluble in the organic solvent dimethyl formamide, which should be purged with an inert gas, at a concentration of approximately 0.5 mg/ml.

Description

C6 dihydro 1-Deoxyceramide (m18:0/6:0) is a short-chain atypical ceramide containing a 1-deoxysphinganine (m18:0) (Item No. 13511) backbone. 1-Deoxysphingolipids are formed when serine palmitoyltransferase condenses palmitoyl-CoA with alanine instead of serine during sphingolipid synthesis.^{1,2}

References

- Steiner, R., Saied, E.M., Othman, A., *et al.* Elucidating the chemical structure of native 1-deoxysphingosine. *J. Lipid Res.* **57(7)**, 1194-1203 (2016).
- Alecu, I., Othman, A., Penno, A., *et al.* Cytotoxic 1-deoxysphingolipids are metabolized by a cytochrome P450-dependent pathway. *J. Lipid Res.* **58(1)**, 60-71 (2017).

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, 02/16/2024

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