

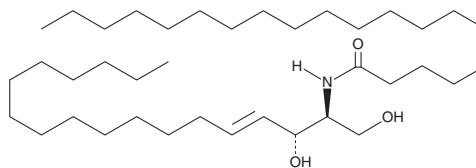
PRODUCT INFORMATION



C20 Ceramide (d18:1/20:0)

Item No. 10724

CAS Registry No.: 7344-02-7
Formal Name: N-[2R-hydroxy-1S-(hydroxymethyl)-3E-heptadecen-1-yl]-eicosanamide
Synonyms: Cer(d18:1/20:0), Ceramide (d18:1/20:0), N-arachidoyl-D-erythro-Sphingosine
MF: C₃₈H₇₅NO₃
FW: 594.0
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

C20 Ceramide (d18:1/20:0) is supplied as a crystalline solid. A stock solution may be made by dissolving the C20 ceramide (d18:1/20:0) in the solvent of choice, which should be purged with an inert gas. C20 Ceramide (d18:1/20:0) is soluble in the organic solvent dimethyl formamide at a concentration of approximately 0.15 mg/ml.

Description

Ceramides are generated from sphingomyelin through activation of sphingomyelinases or through the *de novo* synthesis pathway. Certain forms of ceramide have been shown to mediate cellular responses such as apoptosis, growth arrest, and differentiation in certain cell types.^{1,2} C20 Ceramide is a natural 20:0 ceramide that is abundant in the brain. It is synthesized *de novo* by ceramide synthases 1 and 2.^{3,4}

References

- Ogretmen, B., Pettus, B.J., Rossi, M.J., *et al.* Biochemical mechanisms of the generation of endogenous long chain ceramide in response to exogenous short chain ceramide in the A547 human lung adenocarcinoma cell line. *J. Biol. Chem.* **277**(15), 12960-9 (2002).
- Pandey, S., Murphy, R.F., and Agrawal, D.K. Recent advances in the immunobiology of ceramide. *Exp. Mol. Pathol.* **82**(3), 298-309 (2007).
- Cowart, L.A. Sphingolipids: Players in the pathology of metabolic disease. *Trends Endocrinol. Metab.* **20**(1), 34-42 (2008).
- Pewzner-Jung, Y., Ben-Dor, S., and Futerman, A.H. When do lasses (longevity assurance genes) become CerS (ceramide synthases)? *J. Biol. Chem.* **281**(35), 25001-5 (2006).

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