

PRODUCT INFORMATION



C26 Ceramide (d18:1/26:0) (synthetic)

Item No. 42978

CAS Registry No.: 121459-09-4

Formal Name: N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-hexacosanamide

Synonyms: Cer(d18:1/26:0),
Ceramide (d18:1/26:0),
N-Hexacosanoyl-D-erythro-Sphingosine

MF: $C_{44}H_{87}NO_3$

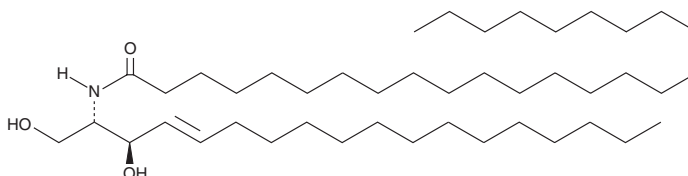
FW: 678.2

Purity: $\geq 98\%$

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

C26 Ceramide (d18:1/26:0) (synthetic) is supplied as a solid. A stock solution may be made by dissolving the C26 ceramide (d18:1/26:0) (synthetic) in the solvent of choice, which should be purged with an inert gas. C26 Ceramide (d18:1/26:0) (synthetic) is soluble in chloroform (warmed).

Description

C26 Ceramide is an endogenous ceramide.¹ C26 Ceramide levels are increased in tumor tissues from patients with hepatocellular carcinoma (HCC) compared to normal liver tissues.² C26 ceramide levels are also increased in acid ceramidase deficiency, a lysosomal disorder characterized by developmental delay, myoclonic seizures, and hematological abnormalities.³ This product is fully synthetic and has no variations in the fatty acyl chain or sphingoid backbone.

References

1. Shin, K.-O., Kim, S., Park, B.D., *et al.* N-Palmitoyl serinol stimulates ceramide production through a CB1-dependent mechanism in in vitro model of skin inflammation. *Int. J. Mol. Sci.* **22**(15), 8302 (2021).
2. Miura, K., Nagahashi, M., Prasoon, P., *et al.* Dysregulation of sphingolipid metabolic enzymes leads to high levels of sphingosine-1-phosphate and ceramide in human hepatocellular carcinoma. *Hepatol. Res.* **51**(5), 614-626 (2021).
3. Mahmoud, I.G., Elmonem, M.A., Zaki, M.S., *et al.* *ASAHI*-related disorders: Description of 15 novel pediatric patients and expansion of the clinical phenotype. *Clin. Genet.* **98**(6), 598-605 (2020).

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, 01/21/2025

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