

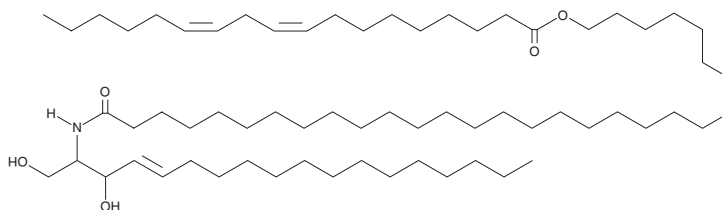
# PRODUCT INFORMATION



## EOS (d18:1/30:0/18:2)

Item No. 24394

**CAS Registry No.:** 97040-38-5  
**Formal Name:** 30-[[2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]amino]-30-oxotriacontyl ester, 9,12-octadecadienoic acid  
**Synonyms:** N-(30-linoleoyloxy-triacontanoyl)-Sphingosine, C48:2 EOS (d18:1/30:0w18:2 (9Z,12Z))  
**MF:** C<sub>66</sub>H<sub>125</sub>NO<sub>5</sub>  
**FW:** 1,012.7  
**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

EOS (d18:1/30:0/18:2) is supplied as a solid. A stock solution may be made by dissolving the EOS (d18:1/30:0/18:2) in the solvent of choice, which should be purged with an inert gas. EOS (d18:1/30:0/18:2) is soluble in chloroform, methanol, and dimethyl formamide.

### Description

EOS is a ceramide found in the outer layer of the epidermis in mammals.<sup>1</sup> It is comprised of an ω-hydroxy very long-chain ceramide (C28-36) esterified to the essential fatty acid linoleic acid (Item No. 90150). The consecutive regio- and stereospecific oxygenation of the linoleate portion of EOS by 12(R)-lipoxygenase (12(R)-LO) and eLOX3 is essential for the maintenance of the epidermal barrier to prevent water loss. Following oxygenation, the oxidized linoleate is hydrolyzed, leaving the ω-hydroxy end of the very long-chain fatty acid to covalently bind the protein layer, forming the corneocyte lipid envelope and sealing the gap between the extracellular lipid lamellae and the cornified cell envelope of the corneocyte. EOS (d18:1/30:0/18:2) is a sphingolipid that has been found in intact and desquamated human stratum corneum as well as porcine epidermis.<sup>2,3</sup>

### References

1. Zheng, Y., Yun, H., Boeglin, W.E., *et al.* Lipoxygenases mediate the effect of essential fatty acid in skin barrier formation: A proposed role in releasing omega-hydroxyceramide for construction of the corneocyte lipid envelope. *J. Biol. Chem.* **286**(27), 24046-24056 (2011).
2. Long, S.A., Wertz, P.W., Strauss, J.S., *et al.* Human stratum corneum polar lipids and desquamation. *Arch. Dermatol. Res.* **277**(4), 284-287 (1985).
3. Wertz, P.W., Abraham, W., Cho, E.S., *et al.* Linoleate-rich O-acylsphingolipids of mammalian epidermis: Structures and effects of essential fatty acid deficiency. *Prog. Lipid Res.* **25**, 383-389 (1986).

#### 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, 12/06/2022

#### 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