

PRODUCT INFORMATION



EOP (t18:0/30:0/18:2)

Item No. 29919

Formal Name: 30-oxo-30-((1,3,4-trihydroxyoctadecan-2-yl)amino)triacontyl (9Z,12Z)-octadeca-9,12-dienoate

Synonyms: C48:2 EOP (t18:0/30:0w18:2 (9Z,12Z)), N-(30-linoleoyloxy-triacontanoyl)-Phytosphingosine

MF: $C_{66}H_{127}NO_6$

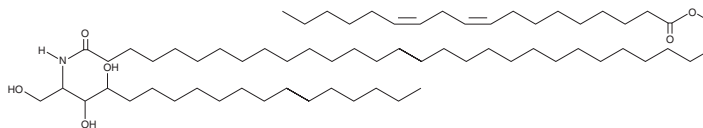
FW: 1,030.7

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

EOP (t18:0/30:0/18:2) is supplied as a solid. A stock solution may be made by dissolving the EOP (t18:0/30:0/18:2) in the solvent of choice, which should be purged with an inert gas. EOP (t18:0/30:0/18:2) is soluble in a 4:1 solution of chloroform:methanol (warmed).

Description

EOP is a phytoceramide-based analog of EOS (Item No. 24394). It is composed of an ω -hydroxy very long-chain ceramide (C28-36) esterified to the essential fatty acid linoleic acid (Item Nos. 90150 | 90150.1 | 21909). EOP has been used as a component in a lipid mixture model of the stratum corneum.¹

Reference

1. de Jager, M., Gooris, G., Ponc, M., *et al.* Acylceramide head group architecture affects lipid organization in synthetic ceramide mixtures. *J. Invest. Dermatol.* **123**(5), 911-916 (2004).

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, 12/12/2022

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