

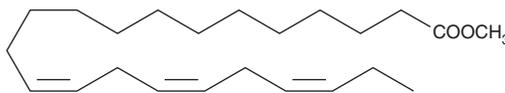
PRODUCT INFORMATION



Docosatrienoic Acid methyl ester

Item No. 20567

CAS Registry No.: 108698-01-7
Formal Name: 13Z,16Z,19Z-docosatrienoic acid, methyl ester
Synonyms: cis-13,16,19-Docosatrienoic Acid methyl ester, SFE 23:3
MF: C₂₃H₄₀O₂
FW: 348.6
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Docosatrienoic acid methyl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the docosatrienoic acid methyl ester under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of docosatrienoic acid methyl ester in these solvents is approximately 30 mg/ml.

Description

Docosatrienoic acid methyl ester is an ester version of docosatrienoic acid (Item No. 90170), a rare ω-3 fatty acid not readily detected in normal phospholipid polyunsaturated fatty acid pools that inhibits [³H]-LTB₄ binding to pig neutrophils at a concentration of 5 μM.¹ Docosatrienoic acid methyl ester is more lipid soluble than the free acid and may be more desirable than the free acid in certain formulations.

Reference

1. Yagaloff, K.A., Franco, L., Simko, B., *et al.* Essential fatty acids are antagonists of the leukotriene B₄ receptor. *Prostaglandins, Leukot. Essent. Fatty Acids* **52(5)**, 293-297 (1995).

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, 12/18/2025

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