

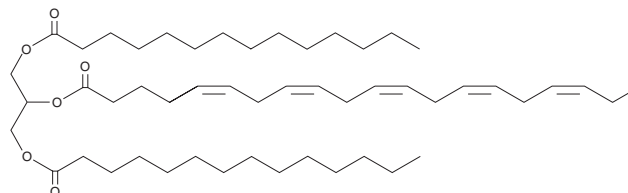
PRODUCT INFORMATION



1,3-Dimyristoyl-2-Eicosapentaenoyl Glycerol

Item No. 26986

CAS Registry No.: 191786-55-7
Formal Name: 5Z,8Z,11Z,14Z,17Z-eicosapentaenoic acid, 2-[[[1-oxotetradecyl)oxy]-1-[[[1-oxotetradecyl)oxy] methyl]ethyl ester
Synonyms: MEM, 1,3-Myristin-2-Eicosapentaenoin, 14:0/20:5/14:0-TG, TG(14:0/20:5/14:0)
MF: C₅₁H₈₈O₆
FW: 797.2
Purity: ≥98%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

1,3-Dimyristoyl-2-eicosapentaenoyl glycerol (MEM) is supplied as a liquid. A stock solution may be made by dissolving the MEM in the solvent of choice. MEM is soluble in organic solvents such as ethanol and dimethyl formamide, which should be purged with an inert gas. The solubility of MEM in these solvents is approximately 10 mg/ml. MEM is also slightly soluble in chloroform.

MEM is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MEM should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. MEM has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

MEM is a triacylglycerol that contains myristic acid (Item No. 13351) at the *sn*-1 and *sn*-3 positions and eicosapentaenoic acid (Item Nos. 90110 | 90110.1 | 21908) at the *sn*-2 position. The myristoyl groups in MEM do not affect the oxidative stability of eicosapentaenoic acid (EPA) during long-term storage at 25°C.¹

Reference

1. Endo, Y., Hoshizaki, S., and Fujimoto, K. Autoxidation of synthetic isomers of triacylglycerol containing eicosapentaenoic acid. *J. Am. Oil Chem. Soc.* **74**(5), 543-548 (1997).

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, 06/08/2020

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