

PRODUCT INFORMATION



1,3-Dipalmitoyl-2-Eicosapentaenoyl Glycerol

Item No. 26840

CAS Registry No.: 148691-18-3
Formal Name: 5Z,8Z,11Z,14Z,17Z-eicosapentaenoic acid, 2-[(1-oxohexadecyl)oxy]-1-[[[(1-oxohexadecyl)oxy]methyl]ethyl] ester

Synonyms: 1,3-Palmitin-2-Eicosapentaenoin, TG(16:0/20:5/16:0)

MF: C₅₅H₉₆O₆

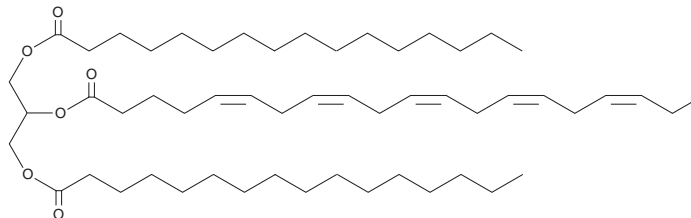
FW: 853.4

Purity: ≥98%

Supplied as: A 10 mg/ml solution in chloroform

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

1,3-Dipalmitoyl-2-eicosapentaenoyl glycerol is supplied as a solution in chloroform. To change the solvent, simply evaporate the chloroform under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol and dimethyl formamide purged with an inert gas can be used at a concentration of approximately 10 mg/ml.

1,3-Dipalmitoyl-2-eicosapentaenoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 1,3-dipalmitoyl-2-eicosapentaenoyl glycerol should be diluted with the aqueous buffer of choice. 1,3-Dipalmitoyl-2-eicosapentaenoyl glycerol has a solubility of 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

1,3-Dipalmitoyl-2-eicosapentaenoyl glycerol is a triacylglycerol that contains palmitic acid (Item No. 10006627) at the *sn*-1 and *sn*-3 positions and eicosapentaenoic acid (Item Nos. 90110 | 90110.1 | 21908) at the *sn*-2 position. Dietary administration of 1,3-dipalmitoyl-2-eicosapentaenoyl glycerol decreases hepatic triacylglycerol levels in mice.¹

Reference

1. Yoshinaga, K., Sasaki, K., Watanabe, H., *et al.* Differential effects of triacylglycerol positional isomers containing n-3 series highly unsaturated fatty acids on lipid metabolism in C57BL/6J mice. *J. Nutr. Biochem.* **26**(1), 57-63 (2015).

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, 03/25/2026

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