

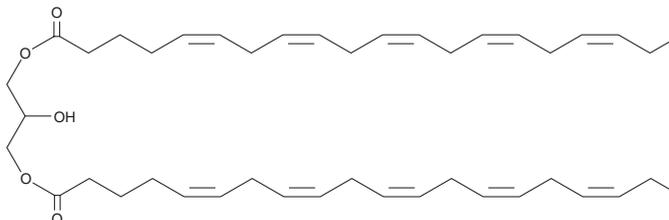
PRODUCT INFORMATION



1,3-Dieicosapentaenoyl Glycerol

Item No. 27010

CAS Registry No.: 140670-41-3
Formal Name: (5Z,5'Z,8Z,8'Z,11Z,11'Z,14Z,14'Z,17Z,17'Z)-5,8,11,14,17-eicosapentaenoic acid, 2-hydroxy-1,3-propanediyl ester
Synonyms: DG(20:5/0:0/20:5), 1,3-Dieicosapentaenoin
MF: C₄₃H₆₄O₅
FW: 661.0
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

1,3-Dieicosapentaenoyl glycerol is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. 1,3-Dieicosapentaenoyl glycerol is soluble in the organic solvent dimethyl formamide at a concentration of approximately 10 mg/ml. 1,3-Dieicosapentaenoyl glycerol is slightly soluble in chloroform.

1,3-Dieicosapentaenoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 1,3-dieicosapentaenoyl glycerol should be diluted with the aqueous buffer of choice. 1,3-Dieicosapentaenoyl 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-Dieicosapentaenoyl glycerol is a diacylglycerol that contains eicosapentaenoic acid (Item Nos. 90110 | 90110.1 | 21908) at the *sn*-1 and *sn*-3 positions.

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, 01/20/2026

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