

PRODUCT INFORMATION



1,3-Didocosahexaenoyl Glycerol

Item No. 27011

CAS Registry No.: 140670-42-4
Formal Name: (4Z,4'Z,7Z,7'Z,10Z,10'Z,13Z,13'Z,16Z,16'Z,19Z,19'Z)-4,7,10,13,16,19-docosahexaenoic acid, 2-hydroxy-1,3-propanediyl ester

Synonyms: DG(22:6/0:0/22:6), 1,3-Didocosahexaenoin

MF: C₄₇H₆₈O₅

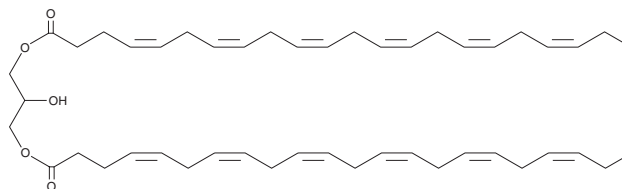
FW: 713.0

Purity: ≥98%

Supplied as: A 10 mg/ml 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-Didocosahexaenoyl glycerol is supplied as a solution in ethanol. To change the solvent, simply evaporate the 1,3-didocosahexaenoyl glycerol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as dimethyl formamide (DMF) purged with an inert gas can be used. 1,3-Didocosahexaenoyl glycerol is soluble in DMF at a concentration of approximately 10 mg/ml. 1,3-Didocosahexaenoyl glycerol is slightly soluble in chloroform.

1,3-Didocosahexaenoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 1,3-didocosahexaenoyl glycerol should be diluted with the aqueous buffer of choice. 1,3-Didocosahexaenoyl glycerol 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

1,3-Didocosahexaenoyl glycerol is a diacylglycerol that contains docosahexaenoic acid (DHA; Item No. 90310) at the *sn*-1 and *sn*-3 positions. It has been used in the synthesis of triacylglycerols that contain DHA at their *sn*-1 and *sn*-3 positions.¹

Reference

1. Gudmundsdottir, A.V., Hansen, K.-A., Magnusson, C.D., *et al.* Synthesis of reversed structured triacylglycerols possessing EPA and DHA at their terminal positions. *Tetrahedron* **71(45)**, 8544-8550 (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