

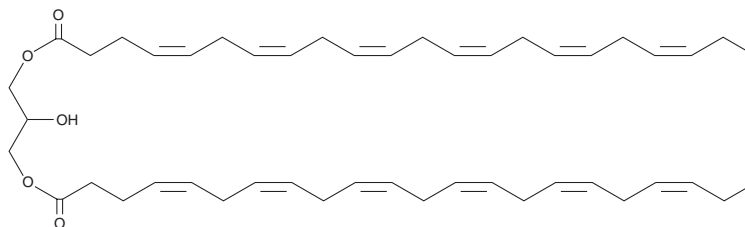
PRODUCT INFORMATION



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

1,3-Didocosahexaenoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,3-didocosahexaenoyl glycerol should first be dissolved in ethanol and then 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, 06/05/2020

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