

PRODUCT INFORMATION

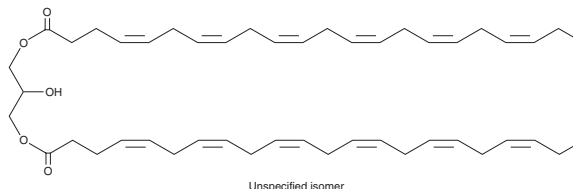


Didocosahexaenoin

Item No. 27003

CAS Registry No.: 124538-05-2
Formal Name: (4'Z,7'Z,10'Z,13'Z,16'Z,19'Z)-4Z,7Z,10Z,13Z,16Z,19Z-docosahexaenoic acid, diester with 1,2,3-propanetriol

MF: C₄₇H₆₈O₅
FW: 713.1
Purity: ≥98%
Supplied as: A liquid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

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

Didocosahexaenoin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, didocosahexaenoin should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Didocosahexaenoin 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

Didocosahexaenoin is a diacylglycerol that contains docosahexaenoic acid (DHA; Item No. 90310) at two 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