

PRODUCT INFORMATION



1,3-Didocosahexaenoyl-2-Palmitoyl Glycerol

Item No. 26987

CAS Registry No.: 197016-86-7

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, 1,1'-[2-[(1-oxohexadecyl)oxy]-1,3-propanediyl] ester

Synonyms: 1,3-Didocosahexaenoin-2-Palmitin, DPD, TG(22:6/16:0/22:6)

MF: $C_{63}H_{98}O_6$

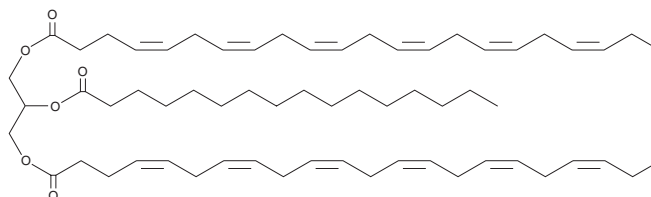
FW: 951.5

Purity: $\geq 98\%$

Supplied as: A 10 mg/ml solution in methyl acetate

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-2-palmitoyl glycerol is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. 1,3-Didocosahexaenoyl-2-palmitoyl glycerol is miscible in the organic solvent chloroform, which should be purged with an inert gas.

Description

1,3-Didocosahexaenoyl-2-palmitoyl glycerol is a triacylglycerol that contains docosahexaenoic acid (Item No. 90310) at the *sn*-1 and *sn*-3 positions and palmitic acid (Item No. 10006627) at the *sn*-2 position. It has been found in higher amounts than 1-palmitoyl-2,3-didocosahexaenoyl-*rac*-glycerol (PDD) in Steller sea lion meat and harp seal oil and at lower amounts than PDD in the meat of a variety of fishes.¹

Reference

1. Gotoh, N., Matsumoto, Y., Nagai, T., *et al.* Actual ratios of triacylglycerol positional isomers consisting of saturated and highly unsaturated fatty acids in fishes and marine mammals. *Food Chem.* **127**(2), 467-472 (2011).

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, 06/25/2026

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