

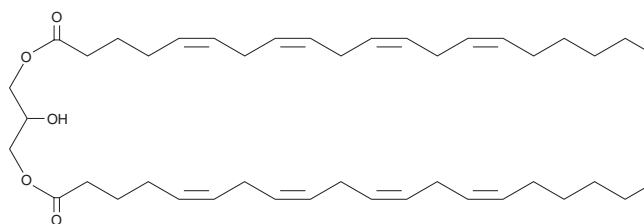
PRODUCT INFORMATION



Diarachidonin

Item No. 26928

CAS Registry No.: 82231-61-6
Formal Name: (5Z,5'Z,8Z,8'Z,11Z,11'Z,14Z,14'Z)-5,8,11,14-eicosatetraenoic acid, diester with 1,2,3-propanetriol
MF: C₄₃H₆₈O₅
FW: 665.0
Purity: ≥98%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



Unspecified isomer

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

Laboratory Procedures

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

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

Diarachidonin is a diacylglycerol that contains the ω-6 polyunsaturated 20-carbon fatty acid arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) 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, 10/01/2019

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