

PRODUCT INFORMATION



1,3-Diarachidonoyl-2-Oleoyl Glycerol

Item No. 29232

CAS Registry No.: 869990-37-4
Formal Name: (5Z,5'Z,8Z,8'Z,11Z,11'Z,14Z,14'Z)-5,8,11,14-eicosatetraenoic acid, 1,1'-[2-[[[(9Z)-1-oxo-9-octadecen-1-yl]oxy]-1,3-propanediyl] ester

Synonyms: 1,3-Diarachidonin-2-Olein, TG(20:4/18:1/20:4)

MF: C₆₁H₁₀₀O₆

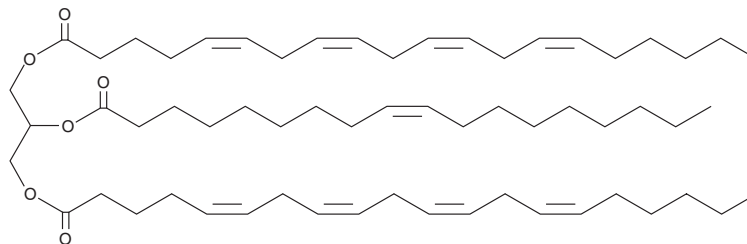
FW: 929.4

Purity: ≥95%

Supplied as: An oil

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-Diarachidonoyl-2-oleoyl glycerol is supplied as an oil. A stock solution may be made by dissolving the 1,3-diarachidonoyl-2-oleoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Diarachidonoyl-2-oleoyl glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 1,3-diarachidonoyl-2-oleoyl glycerol in these solvents is approximately 10 mg/ml.

1,3-Diarachidonoyl-2-oleoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,3-diarachidonoyl-2-oleoyl glycerol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 1,3-Diarachidonoyl-2-oleoyl 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-Diarachidonoyl-2-oleoyl glycerol is a triacylglycerol that contains arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) at the *sn*-1 and *sn*-3 positions and oleic acid (Item Nos. 90260 | 24659) at the *sn*-2 position.

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, 08/18/2025

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