

PRODUCT INFORMATION



1-Palmitoyl-2-Oleoyl-3-Linolenoyl-*rac*-glycerol

Item No. 29061

CAS Registry No.: 78144-20-4

Formal Name: (9Z,12Z,15Z)-9,12,15-octadecatrienoic acid, 3-[(1-oxohexadecyl)oxy]-2-[[[(9Z)-1-oxo-9-octadecen-1-yl]oxy]propyl ester

Synonyms: 1-Palmitin-2-Olein-3-Linolenin, TG(16:0/18:1/18:3)

MF: $C_{55}H_{98}O_6$

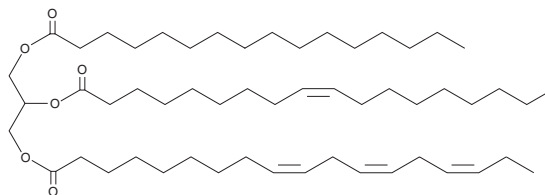
FW: 855.4

Purity: $\geq 95\%$

Supplied as: A solution in ethyl acetate

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-Palmitoyl-2-oleoyl-3-linolenoyl-*rac*-glycerol is supplied as a solution in ethyl acetate. To change the solvent, simply evaporate the 1-palmitoyl-2-oleoyl-3-linolenoyl-*rac*-glycerol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 1-palmitoyl-2-oleoyl-3-linolenoyl-*rac*-glycerol in these solvents is approximately 10 mg/ml.

1-Palmitoyl-2-oleoyl-3-linolenoyl-*rac*-glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethyl acetate solution of 1-palmitoyl-2-oleoyl-3-linolenoyl-*rac*-glycerol should be diluted with the aqueous buffer of choice. 1-Palmitoyl-2-oleoyl-3-linolenoyl-*rac*-glycerol has a solubility of 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-Palmitoyl-2-oleoyl-3-linolenoyl-*rac*-glycerol is a triacylglycerol that contains palmitic acid (Item No. 10006627), oleic acid (Item Nos. 90260 | 24659), and α -linolenic acid (Item Nos. 90210 | 21910) at the *sn*-1, *sn*-2, and *sn*-3 positions, respectively.

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, 05/17/2023

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