

PRODUCT INFORMATION



1,3-Dieicosenoyl Glycerol

Item No. 26795

CAS Registry No.: 128230-18-2
Formal Name: (11Z,11'Z)-11-eicosenoic acid, 2-hydroxy-1,3-propanediyl ester
Synonyms: 20:1/20:1-DG, DG(20:1/0:0/20:1), 11(Z)-Dieicosenoin, cis-11-Dieicosenoin, di-cis-11-Eicosenoyl-glycerol

MF: C₄₃H₈₀O₅

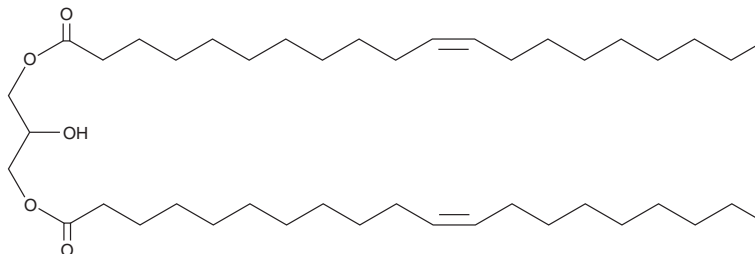
FW: 677.1

Purity: ≥98%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

1,3-Dieicosenoyl glycerol is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as dimethyl formamide, purged with an inert gas, can be used at a concentration of approximately 10 mg/ml.

1,3-Dieicosenoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 1,3-dieicosenoyl glycerol should be diluted with the aqueous buffer of choice. 1,3-Dieicosenoyl glycerol has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method.

Description

1,3-Dieicosenoyl glycerol is a diacylglycerol containing 11(Z)-eicosenoic acid (Item No. 20606) at the sn-1 and sn-3 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, 11/28/2022

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