

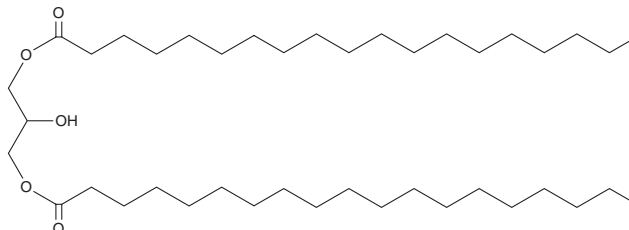
PRODUCT INFORMATION



1,3-Dinonadecanoyl Glycerol

Item No. 26770

CAS Registry No.: 372490-74-9
Formal Name: nonadecanoic acid, 1,1'-(2-hydroxy-1,3-propanediyl) ester
Synonyms: C38 DAG, DG(19:0/0:0/19:0), 9:0/0:0/19:0-DG, 1,3-Dinonadecanoin
MF: C₄₁H₈₀O₅
FW: 653.1
Purity: ≥98%
Supplied as: A solid
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-Dinonadecanoyl glycerol is supplied as a solid. A stock solution may be made by dissolving the 1,3-dinonadecanoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Dinonadecanoyl glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 1,3-dinonadecanoyl glycerol in these solvents is approximately 10 mg/ml.

1,3-Dinonadecanoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,3-dinonadecanoyl glycerol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 1,3-Dinonadecanoyl 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-Dinonadecanoyl glycerol is a diacylglycerol that contains nonadecanoic acid (Item No. 19723) at the *sn*-1 and *sn*-3 positions. It has been used as a standard for the quantification of diacylglycerols in synthetic milk fat by solid phase extraction (SPE) HPLC using an evaporative light scattering detector (ELSD).¹

Reference

1. Avalli, A. and Contarini, G. Determination of phospholipids in dairy products by SPE/HPLC/ELSD. *J. Chromatogr. A.* **1071(1-2)**, 185-190 (2005).

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, 11/17/2022

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