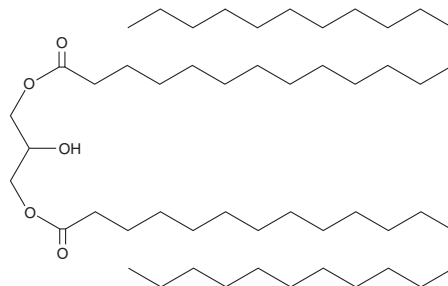


PRODUCT INFORMATION

1,3-Dilignoceroyl Glycerol

Item No. 26796

CAS Registry No.: 120639-92-1
Formal Name: tetracosanoic acid, 2-hydroxy-1,3-propanediyl ester
Synonyms: DG(24:0/0:0/24:0), 1,3-Ditetracosanoin, 1,3-Ditetracosanoyl Glycerol
MF: C₅₁H₁₀₀O₅
FW: 793.3
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-Dilignoceroyl glycerol is supplied as a solid. A stock solution may be made by dissolving the 1,3-dilignoceroyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Dilignoceroyl glycerol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 1,3-dilignoceroyl glycerol in these solvents is approximately 0.25, 30, and 20 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 1,3-dilignoceroyl glycerol can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 1,3-dilignoceroyl glycerol in PBS, pH 7.2, is approximately 0.7 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

1,3-Dilignoceroyl glycerol is a diacylglycerol that contains lignoceric acid (Item No. 13353) 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, 12/05/2022

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