

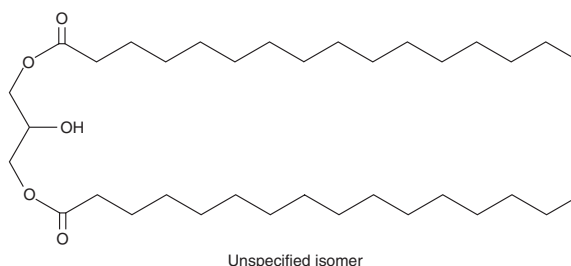
PRODUCT INFORMATION



Dipalmitin

Item No. 28172

CAS Registry No.: 26657-95-4
Formal Name: hexadecanoic acid, diester with 1,2,3-propanetriol
Synonym: Dipalmitoyl-*rac*-glycerol
MF: C₃₅H₆₈O₅
FW: 568.9
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

Dipalmitin is supplied as a solid. A stock solution may be made by dissolving the dipalmitin in the solvent of choice, which should be purged with an inert gas. Dipalmitin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of dipalmitin 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 dipalmitin can be prepared by directly dissolving the solid in aqueous buffers. The solubility of dipalmitin 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

Dipalmitin is a diacylglycerol that contains the saturated 16-carbon fatty acid palmitic acid (Item No. 10006627) at two 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, 11/18/2022

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