

PRODUCT INFORMATION



1-Palmitoyl-*rac*-glycerol

Item No. 26993

CAS Registry No.: 542-44-9
Formal Name: hexadecanoic acid,
2,3-dihydroxypropyl ester
Synonyms: 1-Hexadecanoyl-*rac*-glycerol,
1-Monohexadecanoin,
1-Monopalmitin,
MG(16:0/0:0/0:0)

MF: C₁₉H₃₈O₄

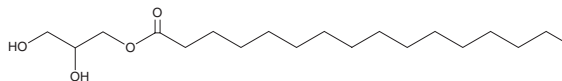
FW: 330.5

Purity: ≥95%

Supplied as: A crystalline 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-Palmitoyl-*rac*-glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1-palmitoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1-Palmitoyl-*rac*-glycerol is soluble in the organic solvents ethanol and dimethyl formamide. The solubility of 1-palmitoyl-*rac*-glycerol in these solvents is approximately 15 mg/ml.

Description

1-Palmitoyl-*rac*-glycerol is a monoacylglycerol that contains palmitic acid (Item No. 10006627) at the *sn*-1 position. It is an emulsifying agent and has been used to increase the viscoelasticity of processed cheese.¹

Reference

1. Bunka, F., Pavlínek, V., Hrabě, J., *et al.* Effect of 1-monoglycerides on viscoelastic properties of processed cheese. *Int. J. Food Prop.* **10(4)**, 819-828 (2007).

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, 10/24/2022

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