

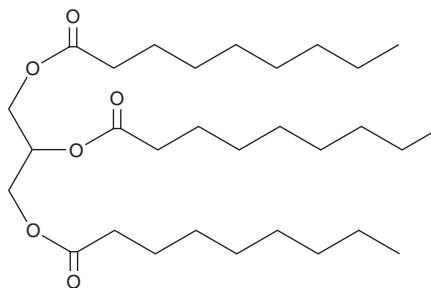
PRODUCT INFORMATION



1,2,3-Trinonanoyl Glycerol

Item No. 28177

CAS Registry No.: 126-53-4
Formal Name: nonanoic acid, 1,1',1''-(1,2,3-propanetriyl) ester
Synonyms: Glyceryl Trinonanoate, NSC 5647, Trinonanoin, Tripelargonin
MF: C₃₀H₅₆O₆
FW: 512.8
Purity: ≥95%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

1,2,3-Trinonanoyl glycerol is supplied as a liquid. A stock solution may be made by dissolving the 1,2,3-trinonanoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Trinonanoyl glycerol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 1,2,3-trinonanoyl glycerol in ethanol is approximately 25 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Description

1,2,3-Trinonanoyl glycerol is a triacylglycerol that contains nonanoic acid at the *sn*-1, *sn*-2, and *sn*-3 positions. It has been found in *Schizochytrium* biomass.¹ 1,2,3-Trinonanoyl-*rac*-glycerol increases blood levels of the ketone body D-(-)-3-hydroxybutyrate in newborn rhesus monkeys.²

References

1. Mioso, R., Toledo Marante, F.J., González, J.E., *et al.* Metabolite profiling of *Schizochytrium* sp. by GC-MS, an oleaginous microbial source of biodiesel. *Braz. J. Microbiol.* **45(2)**, 403-409 (2014).
2. Tetrick, M.A., Greer, F.R., and Benevenga, N.J. Blood D-(-)-3-hydroxybutyrate concentrations after oral administration of trioctanoin, trinonanoin, or tridecanoin to newborn rhesus monkeys (*Macaca mulatta*). *Comp. Med.* **60(6)**, 486-490 (2010).

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, 06/05/2020

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