

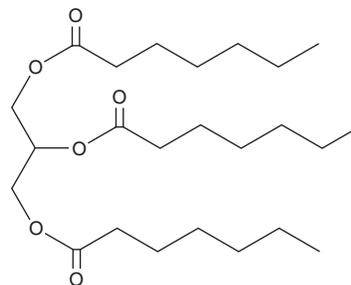
PRODUCT INFORMATION



1,2,3-Triheptanoyl Glycerol

Item No. 28174

CAS Registry No.: 620-67-7
Formal Name: heptanoic acid, 1,1',1''-(1,2,3-propanetriyl) ester
Synonyms: Glycerol Triheptanoate, TG(7:0/7:0/7:0), Triheptanoin
MF: C₂₄H₄₄O₆
FW: 428.6
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-Triheptanoyl glycerol is supplied as a liquid. A stock solution may be made by dissolving the 1,2,3-triheptanoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Triheptanoyl glycerol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 1,2,3-triheptanoyl glycerol in these solvents is approximately 30 mg/ml.

Description

1,2,3-Triheptanoyl glycerol is a triacylglycerol that contains heptanoic acid at the *sn*-1, *sn*-2, and *sn*-3 positions. It increases levels of propionyl carnitine (Item No. 9001873), pentanoyl carnitine (valeryl-L-carnitine; Item No. 26563), and heptanoyl carnitine (Item No. 26551) in mouse blood when administered as 35% of calories in the diet.¹ 1,2,3-Triheptanoyl glycerol (35% of calories in the diet) also delays the development of corneal kindling-induced seizures in mice and increases the pentylenetetrazole (PTZ; Item No. 18682) tonic extension threshold during the chronically epileptic stage in a mouse model of pilocarpine-induced status epilepticus.

Reference

1. Willis, S., Stoll, J., Sweetman, L., *et al.* Anticonvulsant effects of a triheptanoin diet in two mouse chronic seizure models. *Neurobiol. Dis.* **40(3)**, 565-572 (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, 11/29/2022

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