

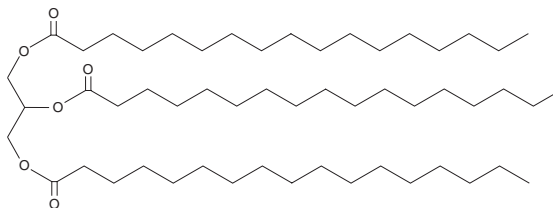
PRODUCT INFORMATION



1,2,3-Triheptadecanoyl Glycerol

Item No. 16489

CAS Registry No.: 2438-40-6
Formal Name: heptadecanoic acid,
1,1',1''-(1,2,3-propanetriyl) ester
Synonyms: Glycerol Triheptadecanoate,
TG(17:0/17:0/17:0)
MF: C₅₄H₁₀₄O₆
FW: 849.4
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

1,2,3-Triheptadecanoyl glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2,3-triheptadecanoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Triheptadecanoyl glycerol is soluble in the organic solvent dimethyl formamide at a concentration of approximately 10 mg/ml.

Description

1,2,3-Triheptadecanoyl glycerol is a triacylglycerol that contains heptadecanoic acid (Item No. 19722) at the *sn*-1, *sn*-2, 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, 09/29/2022

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