

PRODUCT INFORMATION



1,2,3-Tri-12(Z)-Heneicosanoyl Glycerol

Item No. 26844

CAS Registry No.: 169054-26-6
Formal Name: (12Z,12'Z,12''Z)-12-heneicosenoic acid, 1,1',1''-(1,2,3-propanetriyl) ester
Synonyms: Glycerol 12(Z)-Triheneicosanoate, 21:1(12Z)/21:1(12Z)/21:1(12Z)-TG, TG(21:1(12Z)/21:1(12Z)/21:1(12Z)), 12(Z)-Triheneicosanoylglycerol, cis-12-Triheneicosenoic Acid, 12(Z)-Triheneicosenoin, 1,2,3-Triheneicosenoylglycerol

MF: C₆₆H₁₂₂O₆

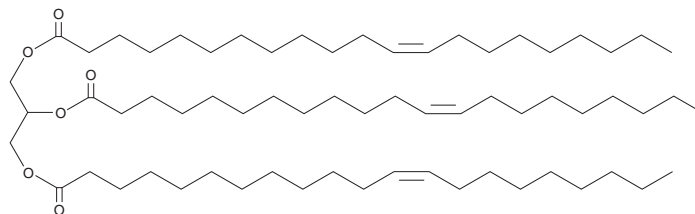
FW: 1,011.7

Purity: ≥98%

Supplied as: An oil

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-Tri-12(Z)-heneicosanoyl glycerol is supplied as an oil. A stock solution may be made by dissolving the 1,2,3-tri-12(Z)-heneicosanoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Tri-12(Z)-heneicosanoyl glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 1,2,3-tri-12(Z)-heneicosanoyl glycerol in these solvents is approximately 10 mg/ml.

Description

1,2,3-Tri-12(Z)-heneicosanoyl glycerol is an unsaturated triacylglycerol containing heneicosenoic acid 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, 12/21/2022

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