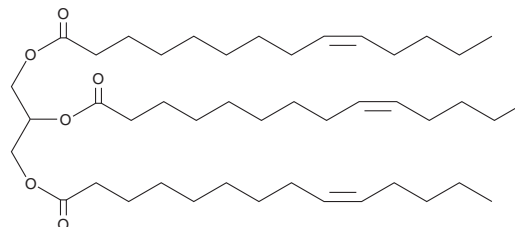


PRODUCT INFORMATION

1,2,3-Trimyristoleoyl Glycerol

Item No. 26943

CAS Registry No.: 99483-10-0
Formal Name: (9Z,9'Z,9''Z)-9-tetradecenoic acid, 1,1',1''-(1,2,3-propanetriyl) ester
Synonyms: TG(14:1/14:1/14:1), Trimyristolein
MF: C₄₅H₈₀O₆
FW: 717.1
Purity: ≥98%
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-Trimyristoleoyl glycerol is supplied as a liquid. A stock solution may be made by dissolving the 1,2,3-trimyristoleoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Trimyristoleoyl glycerol is slightly soluble in methanol and chloroform.

Description

1,2,3-Trimyristoleoyl glycerol is a triacylglycerol that contains myristoleic acid (Item No. 9002461) at the *sn*-1, *sn*-2, and *sn*-3 positions. It has been found in Idiazabal cheese.¹

Reference

1. Nájera, A.I., Perea, S., Barcina, Y., *et al.* A reverse-phase high-performance liquid chromatographic method for analyzing complex mixtures of triglycerides. Application to the fat fraction of an aged cheese. *J. Am. Oil Chem. Soc.* **76**(3), 399-407 (1999).

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, 06/08/2020

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