

PRODUCT INFORMATION



1,2,3-Tri-13(E)-Docosenoyl Glycerol

Item No. 26991

CAS Registry No.: 37635-44-2

Formal Name: (13E,13'E,13''E)-13-docosenoic acid,
1,2,3-propanetriyl ester

Synonyms: Brassidic Acid Triglyceride, Glycerol
Tridocosenoate, TG(22:1/22:1/22:1),
Tribrassidin, Tridocosenoyl Glycerol,
Tri-13(E)-Docosenoin

MF: $C_{69}H_{128}O_6$

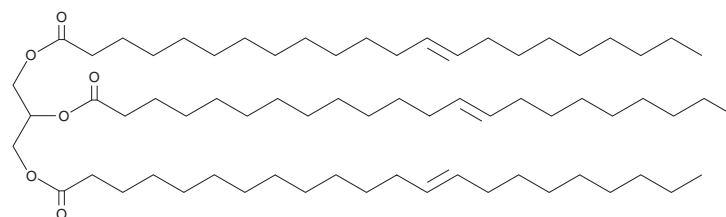
FW: 1,053.8

Purity: $\geq 98\%$

Supplied as: A 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-Tri-13(E)-docosenoyl glycerol is supplied as a solid. A stock solution may be made by dissolving the 1,2,3-tri-13(E)-docosenoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Tri-13(E)-docosenoyl glycerol is slightly soluble in chloroform and methanol.

Description

1,2,3-Tri-13(E)-docosenoyl glycerol is a triacylglycerol that contains 13(E)-docosenoic acid (Item No. 22580) at the *sn*-1, *sn*-2, and *sn*-3 positions. Dietary administration of 1,2,3-tri-13(E)-docosenoyl glycerol (15% of the diet) in weanling rats fed a calcium-deficient diet transiently increases heart triglyceride levels, which return to baseline after 28 days.¹

Reference

1. Astorg, P.O. Heart lipidosis induced by short-term feeding of *cis*- or *trans*-docosenoic acids in weanling or 7-week-old rats. *Ann. Nutr. Metab.* **25(4)**, 201-207 (1981).

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, 01/20/2026

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