

# PRODUCT INFORMATION



## 1,2,3-Tri-11(Z)-Eicosenoyl Glycerol

Item No. 26926

CAS Registry No.: 80380-39-8

Formal Name: (11Z,11'Z,11''Z)-11-eicosenoic acid,  
1,1',1''-(1,2,3-propanetriyl) ester

Synonyms: *cis*-11-Trieicosenoic Acid, *cis*-Trigondoic Acid,  
Glycerol Trieicosenoate, Glycerol Trigondosenoate,  
TG(20:1/20:1/20:1), 11(Z)-Trieicosenoin,  
Trieicosenoylglycerol, Trigondoylglycerol

MF:  $C_{63}H_{116}O_6$

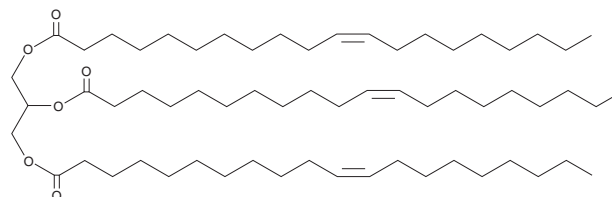
FW: 969.6

Purity:  $\geq 98\%$

Supplied as: A liquid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

1,2,3-Tri-11(Z)-eicosenoyl glycerol is supplied as a liquid. A stock solution may be made by dissolving the 1,2,3-tri-11(Z)-eicosenoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Tri-11(Z)-eicosenoyl glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 1,2,3-tri-11(Z)-eicosenoyl glycerol in these solvents is approximately 10 mg/ml.

### Description

1,2,3-Tri-11(Z)-eicosenoyl glycerol is a triacylglycerol that contains 11(Z)-eicosenoic acid (Item No. 20606) at the *sn*-1, *sn*-2, and *sn*-3 positions. It has been used as an internal standard for the quantification of triacylglycerols in seed and olive oils.<sup>1</sup>

### Reference

1. Mannina, L., Luchinat, C., Emanuele, M.C., *et al.* Acyl positional distribution of glycerol tri-esters in vegetable oils: A  $^{13}\text{C}$  NMR study. *Chem. Phys. Lipids* **103(1-2)**, 47-55 (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, 01/16/2026

#### 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