

# PRODUCT INFORMATION



## 1,2,3-Tri-13(Z)-Docosenoyl Glycerol

Item No. 26894

**CAS Registry No.:** 2752-99-0  
**Formal Name:** (13Z,13'Z,13''Z)-13-docosenoic acid,

**Synonyms:** 1,1,1''-(1,2,3-propanetriyl) ester  
Glycerol Tri-13(Z)-Docosenoate,  
Glycerol Trierucate,  
TG(22:1/22:1/22:1),  
Tri-13(Z)-Docosenoin, Trierucin

**MF:** C<sub>69</sub>H<sub>128</sub>O<sub>6</sub>

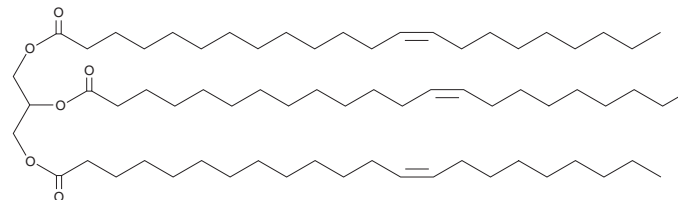
**FW:** 1,053.8

**Purity:** ≥95%

**Supplied as:** A solution in chloroform

**Storage:** -20°C

**Stability:** ≥2 years



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

### Description

1,2,3-Tri-13(Z)-docosenoyl glycerol is a triacylglycerol that contains 13(Z)-docosenoic acid (Item No. 90175) at the *sn*-1, *sn*-2, and *sn*-3 positions and has been found in nasturtium seed oil.<sup>1</sup> 1,2,3-Tri-13(Z)-docosenoyl glycerol (30% of fat calories in a 60% fat diet), in combination with sunflowerseed oil, induces severe cardiac lipidosis in rats after 1 week, with severity decreasing by 24 weeks.<sup>2</sup> 1,2,3-Tri-13(Z)-docosenoyl glycerol is a minor component of Lorenzo's oil.<sup>3</sup>

### References

1. Deineka, V.I., Deineka, L.A., Fofanov, G.M., *et al.* Reversed-phase HPLC of seed oils to establish triglyceride fatty acid composition. *Rastitel'nye Resursy* **40(1)**, 104-112 (2004).
2. Abdellatif, A.M. and Vles, R.O. Short-term and long-term pathological effects of glyceryl trierucate and of increasing levels of dietary rapeseed oil in rats. *Nutr. Metab.* **15(4)**, 219-231 (1973).
3. Moser, H.W., Moser, A.B., Hollandsworth, K., *et al.* "Lorenzo's oil" therapy for X-linked adrenoleukodystrophy: Rationale and current assessment of efficacy. *J. Mol. Neurosci.* **33(1)**, 105-113 (2007).

#### 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, 06/09/2020

#### 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