

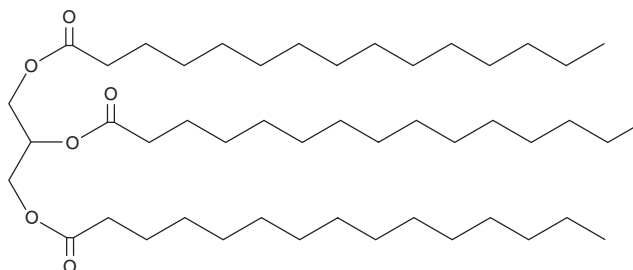
# PRODUCT INFORMATION



## 1,2,3-Tripentadecanoyl Glycerol

Item No. 26962

**CAS Registry No.:** 7370-46-9  
**Formal Name:** pentadecanoic acid, 1,1',1''-(1,2,3-propanetriyl) ester  
**Synonyms:** TG(15:0/15:0/15:0), Tripentadecanoin  
**MF:** C<sub>48</sub>H<sub>92</sub>O<sub>6</sub>  
**FW:** 765.2  
**Purity:** ≥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-Tripentadecanoyl glycerol is supplied as a solid. A stock solution may be made by dissolving the 1,2,3-tripentadecanoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Tripentadecanoyl glycerol is soluble in the organic solvent dimethyl formamide at a concentration of approximately 10 mg/ml.

### Description

1,2,3-Tripentadecanoyl glycerol is a triacylglycerol that contains pentadecanoic acid (Item No. 17399) at the *sn*-1, *sn*-2, and *sn*-3 positions. It has been found in various grapes.<sup>1</sup> 1,2,3-Tripentadecanoyl glycerol has been used as a standard for the detection of triacylglycerols in human plasma by LC-MS.<sup>2</sup>

### References

1. Della Corte, A., Chitarrini, G., Di Gangi, I.M., *et al.* A rapid LC-MS/MS method for quantitative profiling of fatty acids, sterols, glycerolipids, glycerophospholipids and sphingolipids in grapes. *Talanta* **140**, 52-61 (2015).
2. Castro-Perez, J.M., Kamphorst, J., DeGroot, J., *et al.* Comprehensive LC-MSE lipidomic analysis using a shotgun approach and its application to biomarker detection and identification in osteoarthritis patients. *J. Proteome. Res.* **9**(5), 2377-2389 (2010).

#### 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, 12/20/2022

#### 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