

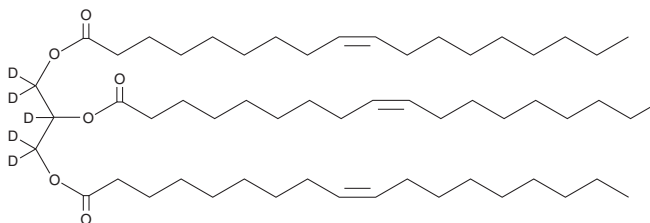
# PRODUCT INFORMATION



## 1,2,3-Trioleoyl Glycerol-d<sub>5</sub>

Item No. 41987

**CAS Registry No.:** 60892-83-3  
**Formal Name:** 9Z-octadecenoic acid, 1,2,3-propanetriyl-1,1,2,3,3-d<sub>5</sub> ester  
**Synonyms:** Glyceryl Trioleate-d<sub>5</sub>, TG(18:1/18:1/18:1)-d<sub>5</sub>, Triolein-d<sub>5</sub>  
**MF:** C<sub>57</sub>H<sub>99</sub>D<sub>5</sub>O<sub>6</sub>  
**FW:** 890.5  
**Chemical Purity:** ≥95% (1,2,3-Trioleoyl Glycerol)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>5</sub>); ≤1% d<sub>0</sub>  
**Supplied as:** A solution in ethanol  
**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-Trioleoyl glycerol-d<sub>5</sub> is intended for use as an internal standard for the quantification of 1,2,3-trioleoyl glycerol (Item No. 26871) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

1,2,3-Trioleoyl glycerol-d<sub>5</sub> is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. 1,2,3-Trioleoyl glycerol-d<sub>5</sub> is slightly soluble (0.1-1 mg/ml) in chloroform and methanol.

### Description

1,2,3-Trioleoyl glycerol is a triacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) at the *sn*-1, *sn*-2, and *sn*-3 positions. It inhibits oxidized LDL-induced decreases in cell viability and superoxide dismutase (SOD) and glutathione peroxidase (GPX) activities and increases in apoptosis in endothelial cells when used at a concentration of 10 μM.<sup>1</sup> 1,2,3-Trioleoyl glycerol has been found in sunflower, corn, and extra virgin olive oils.<sup>2,3</sup> It has been used as a standard for the quantification of triacylglycerols in extra virgin olive oil by high-temperature gas chromatography-(ionic trap)mass spectrometry (HTGC-(IT)MS).<sup>3</sup> 1,2,3-Trioleoyl glycerol is a major component of Lorenzo's oil.<sup>4</sup>

### References

1. Luo, T., Deng, Z.Y., Li, X.P., *et al.* Triolein and trilinolein ameliorate oxidized low-density lipoprotein-induced oxidative stress in endothelial cells. *Lipids* **49**(5), 495-504 (2014).
2. Gao, B., Luo, Y., Lu, W., *et al.* Triacylglycerol compositions of sunflower, corn and soybean oils examined with supercritical CO<sub>2</sub> ultra-performance convergence chromatography combined with quadrupole time-of-flight mass spectrometry. *Food Chem.* **218**, 569-574 (2017).
3. Ruiz-Samblás, C., Rodríguez-Cuadros, L., González-Casado, A., *et al.* A straightforward quantification of triacylglycerols (and fatty acids) in monovarietal extra virgin olive oils by high-temperature GC. *Anal. Methods* **4**(3), 753-758 (2012).
4. 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, 10/30/2024

#### 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