

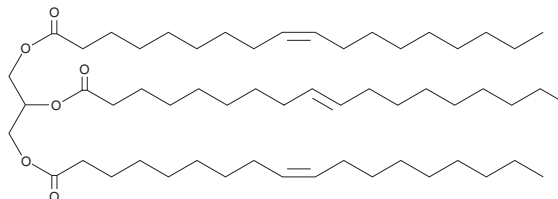
# PRODUCT INFORMATION



## 1,3-Dioleoyl-2-Elaidoyl Glycerol

Item No. 29235

**CAS Registry No.:** 110170-81-5  
**Formal Name:** (9Z,9'E,9''Z)-9-octadecenoic acid,  
1,2,3-propanetriyl ester  
**Synonyms:** 1,3-Olein-2-Elaidin,  
TG(18:1(9Z)/18:1(9E)/18:1(9Z))  
**MF:** C<sub>57</sub>H<sub>104</sub>O<sub>6</sub>  
**FW:** 885.4  
**Purity:** ≥95%  
**Supplied as:** An oil  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

1,3-Dioleoyl-2-elaidoyl glycerol is supplied as an oil. A stock solution may be made by dissolving the 1,3-dioleoyl-2-elaidoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Dioleoyl-2-elaidoyl glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 1,3-dioleoyl-2-elaidoyl glycerol in these solvents is approximately 10 mg/ml.

1,3-Dioleoyl-2-elaidoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,3-dioleoyl-2-elaidoyl glycerol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 1,3-Dioleoyl-2-elaidoyl glycerol has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

1,3-Dioleoyl-2-elaidoyl glycerol is a triacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 and *sn*-3 positions and elaidic acid (Item No. 90250) at the *sn*-2 position. It has been used to study the effects of unsaturation and *sn*-2 substitution on the polymorphic behavior of triacylglycerols.<sup>1</sup>

### Reference

1. Kodali, D.R., Atkinson, D., Redgrave, T.G., *et al.* Structure and polymorphism of 18-carbon fatty acyl triacylglycerols: Effect of unsaturation and substitution in the 2-position. *J. Lipid Res.* **28(4)**, 403-413 (1987).

#### 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, 11/15/2022

#### 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