

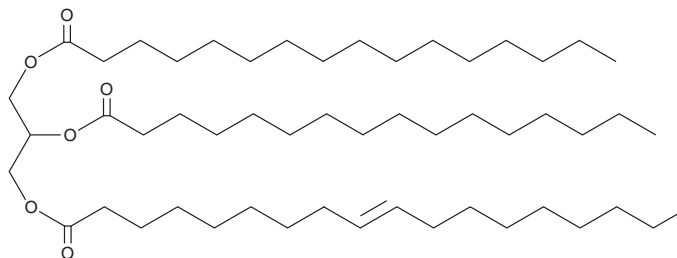
# PRODUCT INFORMATION



## 1,2-Dipalmitoyl-3-Elaidoyl-*rac*-glycerol

Item No. 28865

**CAS Registry No.:** 93452-43-8  
**Formal Name:** (9E)-9-octadecenoic acid, 2,3-bis[(1-oxohexadecyl)oxy]propyl ester  
**Synonyms:** 1,2-Dipalmitin-3-Elaidin,  
1,2-Dipalmitin-3-Elaidoyl-*rac*-glycerol,  
TG(16:0/16:0/18:1(9E))  
**MF:** C<sub>53</sub>H<sub>100</sub>O<sub>6</sub>  
**FW:** 833.4  
**Purity:** ≥95%  
**Supplied as:** A crystalline 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-Dipalmitoyl-3-elaidoyl-*rac*-glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-dipalmitoyl-3-elaidoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1,2-Dipalmitoyl-3-elaidoyl-*rac*-glycerol is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml.

### Description

1,2-Dipalmitoyl-3-elaidoyl-*rac*-glycerol is a triacylglycerol that contains palmitic acid (Item No. 10006627) at the *sn*-1 and *sn*-2 positions and elaidic acid (Item No. 90250) at the *sn*-3 position. It has been used as a standard for the identification of triacylglycerols in rapeseed and palm oils.<sup>1</sup>

### Reference

1. Macher, M.-B. and Holmqvist, A. Triacylglycerol analysis of partially hydrogenated vegetable oils by silver ion HPLC. *J. Sep. Sci.* **24**(3), 179-185 (2001).

**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, 11/07/2022

### 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