

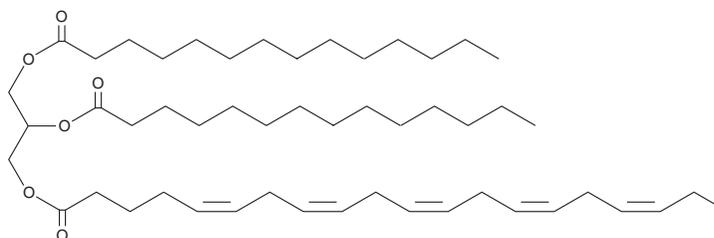
# PRODUCT INFORMATION



## 1,2-Dimyristoyl-3-Eicosapentaenoyl-*rac*-glycerol

Item No. 26826

**CAS Registry No.:** 116198-40-4  
**Formal Name:** (all-*Z*)-5,8,11,14,17-eicosapentaenoic acid, 2,3-*bis*[(1-oxotetradecyl)oxy]propyl ester  
**Synonyms:** MME, 1,2-Myristin-3-Eicosapentaenoin, 14:0/14:0/20:5-TG, TG(14:0/14:0/20:5)  
**MF:** C<sub>51</sub>H<sub>88</sub>O<sub>6</sub>  
**FW:** 797.2  
**Purity:** ≥98%  
**Supplied as:** A liquid  
**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-Dimyristoyl-3-eicosapentaenoyl-*rac*-glycerol is supplied as a liquid. A stock solution may be made by dissolving the 1,2-dimyristoyl-3-eicosapentaenoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1,2-Dimyristoyl-3-eicosapentaenoyl-*rac*-glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 1,2-dimyristoyl-3-eicosapentaenoyl-*rac*-glycerol in these solvents is approximately 10 mg/ml. 1,2-Dimyristoyl-3-eicosapentaenoyl-*rac*-glycerol is slightly soluble in chloroform.

1,2-Dimyristoyl-3-eicosapentaenoyl-*rac*-glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,2-dimyristoyl-3-eicosapentaenoyl-*rac*-glycerol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 1,2-Dimyristoyl-3-eicosapentaenoyl-*rac*-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

1,2-Dimyristoyl-3-eicosapentaenoyl-*rac*-glycerol is a triacylglycerol that contains myristic acid (Item No. 13351) at the *sn*-1 and *sn*-2 positions and eicosapentaenoic acid (Item Nos. 90110 | 90110.1 | 21908) at the *sn*-3 position. Relative quantities of 1,2-dimyristoyl-3-eicosapentaenoyl-*rac*-glycerol in the alga *T. minutus* positively correlate with increases in temperature.<sup>1</sup>

### Reference

- Řezanka, T., Lukavský, J., Sigler, K., *et al.* Temperature dependence of production of structured triacylglycerols in the alga *Trachydiscus minutus*. *Phytochemistry* **110**, 37-45 (2015).

**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, 01/16/2026

### 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