

# PRODUCT INFORMATION



## 1,2,3-Triheneicosanoyl-*rac*-glycerol

Item No. 10010540

**CAS Registry No.:** 26536-14-1  
**Formal Name:** heneicosanoic acid,  
1,2,3-propanetriyl ester  
**Synonyms:** Glycerol Triheneicosanoate,  
TG(21:0/21:0/21:0),  
Triheneicosanoin,  
Triheneicosanoyl Glycerol

**MF:** C<sub>66</sub>H<sub>128</sub>O<sub>6</sub>

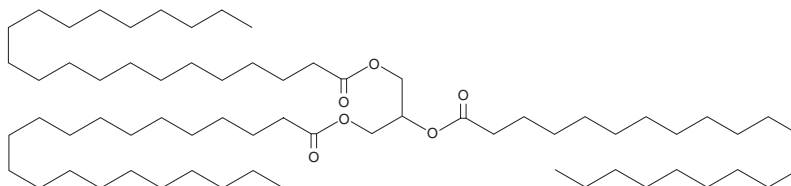
**FW:** 1,017.7

**Purity:** ≥95%

**Supplied as:** A crystalline solid

**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-Triheneicosanoyl-*rac*-glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2,3-triheneicosanoyl-*rac*-glycerol in the solvent of choice. 1,2,3-Triheneicosanoyl-*rac*-glycerol is soluble in the organic solvent dimethyl formamide, which should be purged with an inert gas, at a concentration of approximately 10 mg/ml.

### Description

1,2,3-Triheneicosanoyl-*rac*-glycerol is a triacylglycerol containing heneicosanoic acid (Item No. 22593) groups at the *sn*-1, *sn*-2, and *sn*-3 positions. It has been used as an internal standard for the quantification of fatty acids in human breast milk samples and triacylglycerols in rat heart.<sup>1,2</sup>

### References

1. Hoppu, U., Isolauri, E., Laakso, P., *et al.* Probiotics and dietary counselling targeting maternal dietary fat intake modifies breast milk fatty acids and cytokines. *Eur. J. Nutr.* **51(2)**, 211-219 (2012).
2. Han, X., Abendschein, D.R., Kelley, J.G., *et al.* Diabetes-induced changes in specific lipid molecular species in rat myocardium. *Biochem J.* **352(Pt. 1)**, 79-89 (2000).

#### 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, 05/06/2019

#### 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