

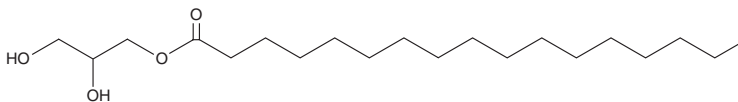
# PRODUCT INFORMATION



## 1-Heptadecanoyl-*rac*-glycerol

Item No. 28180

**CAS Registry No.:** 5638-14-2  
**Formal Name:** heptadecanoic acid,  
2,3-dihydroxypropyl ester  
**Synonyms:** MG(17:0/0:0/0:0),  
1-Heptadecanoin  
**MF:** C<sub>20</sub>H<sub>40</sub>O<sub>4</sub>  
**FW:** 344.5  
**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-Heptadecanoyl-*rac*-glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1-heptadecanoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1-Heptadecanoyl-*rac*-glycerol is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 1-heptadecanoyl-*rac*-glycerol in these solvents is approximately 1 and 3 mg/ml, respectively.

### Description

1-Heptadecanoyl-*rac*-glycerol is a monoacylglycerol that contains heptadecanoic acid (Item No. 19722) at the *sn*-1 position. It is active against the bacteria *E. aerogens*, *E. cloacae*, *P. mirabilis*, and *S. faecalis* (MIC = 78 µg/ml for all).<sup>1</sup> 1-Heptadecanoyl-*rac*-glycerol has been found in *T. africana*, *I. sonorae*, and wheat bran.<sup>1-3</sup>

### References

1. Kuete, V., Metuno, R., Ngameni, B., *et al.* Antimicrobial activity of the methanolic extracts and compounds from *Treculia africana* and *Treculia acuminata* (Moraceae). *S. Afr. J. Bot.* **74**(1), 111-115 (2008).
2. Fernández-Galicia, E., Calada, F., Roman-Romos, R., *et al.* Monoglycerides and fatty acids from *Ibervillea sonorae* root: Isolation and hypoglycemic activity. *Planta Med.* **73**(3), 236-240 (2007).
3. Prinsen, P., Gutiérrez, A., Faulds, C.B., *et al.* Comprehensive study of valuable lipophilic phytochemicals in wheat bran. *J. Agric. Food Chem.* **62**(7), 1664-1673 (2014).

#### 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, 12/16/2022

#### 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