

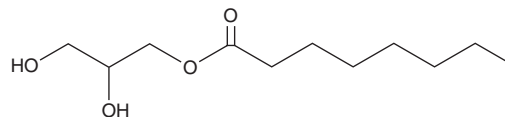
PRODUCT INFORMATION



1-Octanoyl-*rac*-glycerol

Item No. 26945

CAS Registry No.: 502-54-5
Formal Name: octanoic acid, 2,3-dihydroxypropyl ester
Synonyms: 1-Monocaprylin, 1-Monooctanoin,
1-Monooctanoyl-*rac*-glycerol,
Glycerol-1-Octanoate, MG(8:0/0:0/0:0)
MF: C₁₁H₂₂O₄
FW: 218.3
Purity: ≥95%
Supplied as: A 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-Octanoyl-*rac*-glycerol is supplied as a solid. A stock solution may be made by dissolving the 1-octanoyl-*rac*-glycerol in the solvent of choice. 1-Octanoyl-*rac*-glycerol is soluble in the organic solvent chloroform, which should be purged with an inert gas, at a concentration of approximately 100 mg/ml.

Description

1-Octanoyl-*rac*-glycerol is a monoacylglycerol that contains octanoic acid at the *sn*-1 position. It reduces the growth of a variety of bacteria and pathogenic fungi, including *C. albicans* and *C. parapsilosis*, on textiles when applied as a 15 wt% solution in ethanol.¹ 1-Octanoyl-*rac*-glycerol, in combination with ethanol or 2-mercaptoethanol, dissolves gallstones *in vitro*.²

References

1. Vltavská, P., Kašpárková, V., Janiš, R., *et al.* Antifungal and antibacterial effects of 1-monocaprylin on textile materials. *Eur. J. Lipid Sci. Technol.* **114(7)**, 849-856 (2012).
2. Smith, B.F. Dissolution of cholesterol gallstones *in vitro*: Gallstone matrix content and diameter, not cholesterol content, predict gallstone dissolution in monooctanoin. *Gastroenterology* **93(1)**, 98-105 (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, 12/08/2022

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