

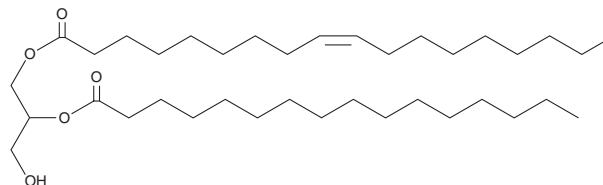
PRODUCT INFORMATION



1-Oleoyl-2-Palmitoyl-*rac*-glycerol

Item No. 27166

CAS Registry No.: 3331-34-8
Formal Name: 9Z-octadecenoic acid, 3-hydroxy-2-[(1-oxohexadecyl)oxy]propyl ester
Synonyms: 18:1/16:0-DG, DG(18:1/16:0/0:0), 1,2-OP, 1-Olein-2-Palmitin
MF: C₃₇H₇₀O₅
FW: 595.0
Purity: ≥95%
Supplied as: A neat oil
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-Oleoyl-2-palmitoyl-*rac*-glycerol (1,2-OP) is supplied as a neat oil. A stock solution may be made by dissolving the 1,2-OP in the solvent of choice, which should be purged with an inert gas. 1,2-OP is slightly soluble in chloroform and ethyl acetate.

Description

1,2-OP is a diacylglycerol containing oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 position and palmitic acid (Item No. 10006627) at the *sn*-2 position. It enhances the inhibition of superoxide anion production by cytochalasin B (Item No. 11328) when used at a concentration of 10 μM.¹ The concentration of 1,2-OP decreases from 25.7 to 11.15% during the dry-curing process of Iberian ham.²

References

1. Wagh, V.D., Korinek, M., Lo, I.-W., *et al.* Inflammation modulatory phorbol esters from the seeds of *Aquilaria malaccensis*. *J. Nat. Prod.* **80**(5), 1421-1427 (2017).
2. Narváez-Rivas, M., Vicario, I.M., Constante, E.G., *et al.* Changes in the concentrations of free fatty acid, monoacylglycerol, and diacylglycerol in the subcutaneous fat of Iberian ham during the dry-curing process. *J. Agric. Food Chem.* **55**(26), 10953-10961 (2007).

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, 11/17/2022

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