

PRODUCT INFORMATION



1,2-Dioleoyl-3-linoleoyl-*rac*-glycerol

Item No. 62610

CAS Registry No.: 2190-20-7

Formal Name: 9Z,12Z-octadecadienoic acid
2,3-bis[[[(9Z)-1-oxo-9-octadecen-1-yl]
oxy]propyl ester

Synonyms: OOL, TG(18:1/18:1/18:2),
TG(18:1(9Z)/18:1(9Z)/18:2(9Z,12Z)),
TG(54:4)

MF: $C_{57}H_{102}O_6$

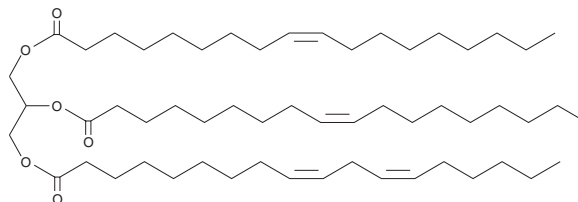
FW: 883.4

Purity: $\geq 95\%$

Supplied as: A solution in methyl acetate

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-Dioleoyl-3-linoleoyl-*rac*-glycerol (OOL) is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the OOL under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol and dimethyl formamide purged with an inert gas can be used. The solubility of OOL in these solvents is approximately 10 mg/ml.

OOL is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of OOL should be diluted with the aqueous buffer of choice. OOL has a solubility of 500 $\mu\text{g/ml}$ 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

OOL is a triacylglycerol containing oleic acid at the *sn*-1 and *sn*-2 positions and linoleic acid at the *sn*-3 position. It is found in a variety of seed and vegetable oils, including pumpkin seed, olive, and sesame oils.¹⁻³

References

1. Ghaleb, M.L., Farines, M., and Soulier, J. Chemical composition of seed oils of pumpkin, gourd and melon. *Rev. Fr. Corps Gras* **38**(1-2), 17-22 (1991).
2. Ruiz- Samblás, C., Rodríguez-Cuadros, L., González-Casado, A., *et al.* A straightforward quantification of triacylglycerols (and fatty acids) in monovarietal extra virgin olive oils by high-temperature GC. *Anal. Methods* **4**(3), 753-758 (2012).
3. U.S.Pharmacopeia and National Formulary Official Monographs NF 19; Sesame Oil. *in* USP 24-NF 19, 3rd Supplement. 19th edition, 3138-3139 (2001).

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, 10/10/2022

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