

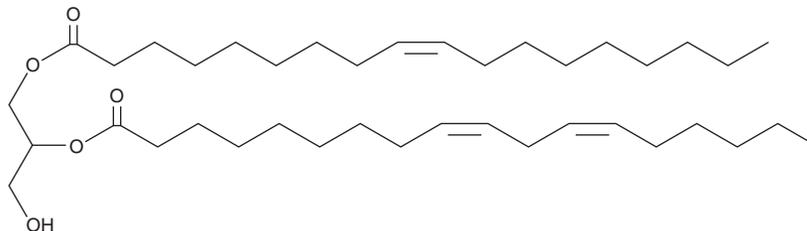
PRODUCT INFORMATION



1-Oleoyl-2-Linoleoyl-*rac*-glycerol

Item No. 37241

CAS Registry No.: 106292-55-1
Formal Name: 9Z,12Z-octadecadienoic acid, 1-(hydroxymethyl)-2-[[[(9Z)-1-oxo-9-octadecen-1-yl]oxy]ethyl ester
Synonyms: DG(18:1/18:2/0:0), 1-Olein-2-Linolein-*rac*-glycerol
MF: C₃₉H₇₀O₅
FW: 619.0
Purity: ≥95%
Supplied as: A 10 mg/ml solution in ethanol
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-Oleoyl-2-linoleoyl-*rac*-glycerol is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as chloroform, dichloromethane, DMSO, ethyl acetate, and hexane purged with an inert gas can be used.

Description

1-Oleoyl-2-linoleoyl-*rac*-glycerol is a diacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) and linoleic acid (Item Nos. 90150 | 90150.1 | 21909) at the *sn*-1 and *sn*-2 positions, respectively. It has been found in raw and dry-cured Iberian ham.¹

Reference

1. 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, 12/10/2025

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