

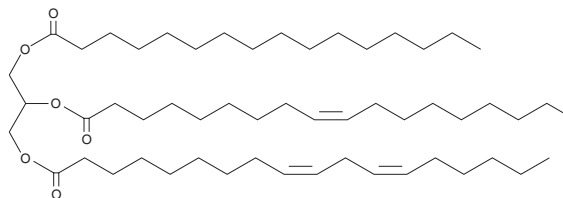
# PRODUCT INFORMATION



## 1-Palmitoyl-2-oleoyl-3-linoleoyl-*rac*-glycerol

Item No. 62660

**CAS Registry No.:** 2680-59-3  
**Formal Name:** 3-[[1-oxohexadecyl]oxy]-2-[[[(9Z)-1-oxo-9-octadecenyl]oxy]-9Z,12Z-octadecadienoic acid propyl ester  
**MF:** C<sub>55</sub>H<sub>100</sub>O<sub>6</sub>  
**FW:** 957.4  
**Purity:** ≥85%  
**Supplied as:** A solution in hexane  
**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-Palmitoyl-2-oleoyl-3-linoleoyl-*rac*-glycerol (POL) is supplied as a solution in hexane. To change the solvent, simply evaporate the hexane 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 POL in these solvents is approximately 10 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. POL is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, POL should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. POL has a solubility of approximately 0.5 mg/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

POL is a common triglyceride component of seed and vegetable oils including olive, sesame, soybean, canola, corn, hazelnut, and many others. POL is one of the standard triglyceride components used for the USP analysis of sesame oil for pharmaceutical applications.<sup>1</sup>

### Reference

1. U.S.Pharmacopeia and National Formulary Official Monographs NF 19; Sesame Oil, in USP 24-NF 19, 3rd Supplement. 19th edition, Rockville, 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, 01/04/2023

#### 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