

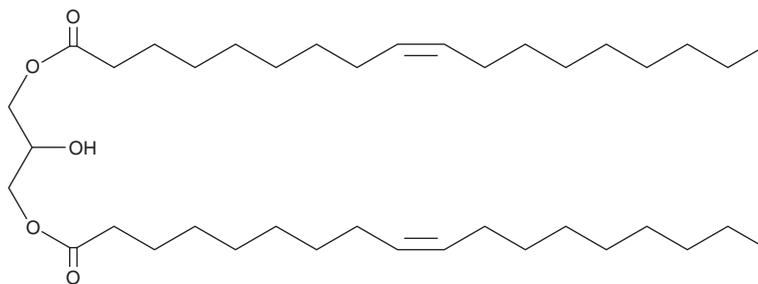
# PRODUCT INFORMATION



## Diolein

Item No. 26976

**CAS Registry No.:** 25637-84-7  
**Formal Name:** 9Z-octadecenoic acid, diester with 1,2,3-propanetriol  
**Synonym:** Dioleoylglycerol  
**MF:** C<sub>39</sub>H<sub>72</sub>O<sub>5</sub>  
**FW:** 621.0  
**Purity:** ≥98%  
**Supplied as:** A liquid  
**Storage:** -20°C  
**Stability:** ≥2 years



Unspecified isomer

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Diolein is supplied as a liquid. A stock solution may be made by dissolving the diolein in the solvent of choice, which should be purged with an inert gas. Diolein is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of diolein in these solvents is approximately 10 mg/ml.

Diolein is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, diolein should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Diolein 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

Diolein is a diacylglycerol that contains the monounsaturated 18-carbon fatty acid oleic acid (Item Nos. 90260 | 24659) at two positions.

#### 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, 08/08/2019

#### 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