

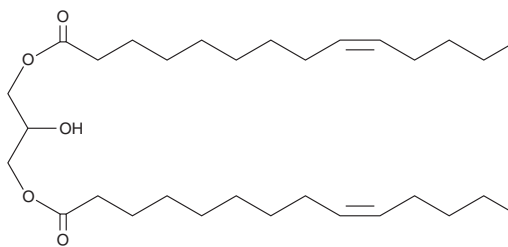
# PRODUCT INFORMATION



## Dimyristolein

Item No. 26977

**CAS Registry No.:** 372516-89-7  
**Formal Name:** (9Z,9'Z)-9-tetradecenoic acid, diester with 1,2,3-propanetriol  
**MF:** C<sub>31</sub>H<sub>56</sub>O<sub>5</sub>  
**FW:** 508.8  
**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

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

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

Dimyristolein is a diacylglycerol that contains the monounsaturated 14-carbon fatty acid myristoleic acid (Item No. 9002461) 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, 11/26/2019

### 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