

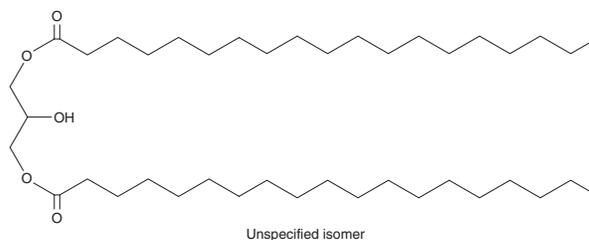
# PRODUCT INFORMATION



## Dinonadecanoin

Item No. 26981

**CAS Registry No.:** 372516-94-4  
**Formal Name:** nonadecanoic acid, diester with 1,2,3-propanetriol  
**MF:** C<sub>41</sub>H<sub>80</sub>O<sub>5</sub>  
**FW:** 653.1  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

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

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

Dinonadecanoin is a diacylglycerol that contains the saturated fatty acid nonadecanoic acid (Item No. 19723) 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, 12/08/2022

#### 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