

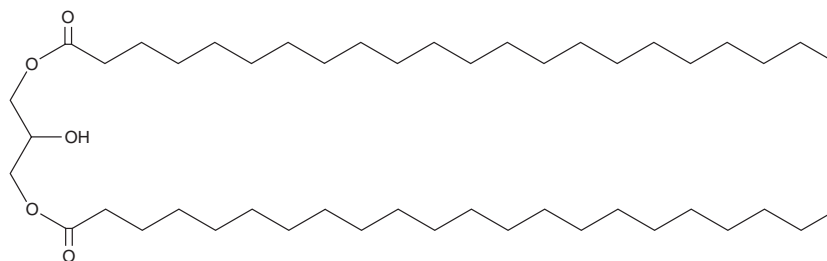
# PRODUCT INFORMATION



## Didocosanoin

Item No. 26963

**CAS Registry No.:** 99880-64-5  
**Formal Name:** docosanoic acid, diester with 1,2,3-propanetriol  
**Synonym:** Dibehenin  
**MF:** C<sub>47</sub>H<sub>92</sub>O<sub>5</sub>  
**FW:** 737.2  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Unspecified isomer

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

### Laboratory Procedures

Didocosanoin is supplied as a solid. A stock solution may be made by dissolving the didocosanoin in the solvent of choice, which should be purged with an inert gas. Didocosanoin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of didocosanoin in these solvents is approximately 0.25, 30, and 20 mg/ml, respectively.

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. Organic solvent-free aqueous solutions of didocosanoin can be prepared by directly dissolving the solid in aqueous buffers. The solubility of didocosanoin in PBS, pH 7.2, is approximately 0.7 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Didocosanoin is a diacylglycerol that contains the saturated 22-carbon fatty acid docosanoic acid (Item No. 9000338) 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, 09/19/2023

### 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