

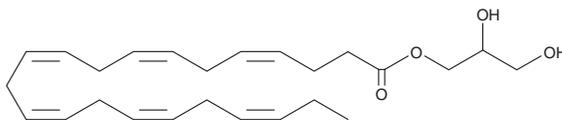
# PRODUCT INFORMATION



## 1-Docosahexaenoyl-*rac*-glycerol

Item No. 35871

**CAS Registry No.:** 140670-40-2  
**Formal Name:** 4Z,7Z,10Z,13Z,16Z,19Z-docosahexaenoic acid, 2,3-dihydroxypropyl ester  
**Synonyms:** 1-Docosahexaenoin, MAG-DHA, MG(22:6/0:0/0:0)  
**MF:** C<sub>25</sub>H<sub>38</sub>O<sub>4</sub>  
**FW:** 402.6  
**Purity:** ≥95% (mixture of isomers)  
**Supplied as:** A 10 mg/ml solution in acetonitrile  
**Storage:** -80°C  
**Stability:** ≥1 year



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

### Laboratory Procedures

1-Docosahexaenoyl-*rac*-glycerol is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile under a gentle stream of nitrogen and immediately add the solvent of choice. 1-Docosahexaenoyl-*rac*-glycerol is soluble in chloroform.

### Description

1-Docosahexaenoyl-*rac*-glycerol is a monoacylglycerol that contains docosahexaenoic acid (DHA; Item Nos. 90310 | 17950) at the *sn*-1 position. It reduces smooth muscle tension induced by endothelin-1 (ET-1) in isolated human pulmonary arteries (HPAs) when used at concentrations ranging from 1 to 30 μM.<sup>1</sup> 1-Docosahexaenoyl-*rac*-glycerol also reduces ET-1-induced increases in HPA cell proliferation and VEGF levels.

### Reference

1. Morin, C., Fortin, S., and Rousseau, E. Docosahexaenoic acid monoacylglyceride decreases endothelin-1 induced Ca<sup>2+</sup> sensitivity and proliferation in human pulmonary arteries. *Am. J. Hypertens.* **25(7)**, 756-763 (2012).

#### 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, 12/01/2025

#### 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