

# PRODUCT INFORMATION



## Monodocosahexaenoin

Item No. 35091

**Formal Name:** 4Z,7Z,10Z,13Z,16Z,19Z-docosahexaenoic acid, 2,3-dihydroxypropyl ester with 1,2,3-propanetriol

**MF:**  $C_{25}H_{38}O_4$

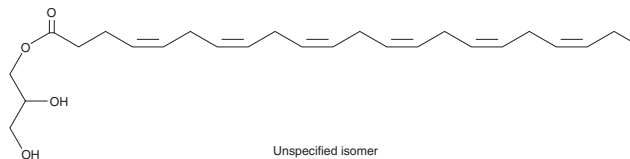
**FW:** 402.6

**Purity:**  $\geq 95\%$

**Supplied as:** A neat oil

**Storage:**  $-20^{\circ}\text{C}$

**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Monodocosahexaenoin is supplied as a neat oil. A stock solution may be made by dissolving the monodocosahexaenoin in the solvent of choice, which should be purged with an inert gas. Monodocosahexaenoin is soluble in chloroform.

### Description

Monodocosahexaenoin is a monoacylglycerol that contains docosahexaenoic acid (DHA; Item No. 90310).

#### 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](http://WWW.CAYMANCHEM.COM)