

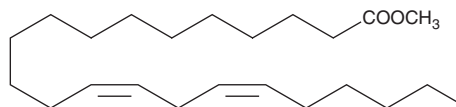
# PRODUCT INFORMATION



## Docosadienoic Acid methyl ester

Item No. 20602

**CAS Registry No.:** 61012-47-3  
**Formal Name:** 13Z,16Z-docosadienoic acid, methyl ester  
**Synonyms:** cis-13,16-Docosadienoic Acid methyl ester, SFE 23:2  
**MF:** C<sub>23</sub>H<sub>42</sub>O<sub>2</sub>  
**FW:** 350.6  
**Purity:** ≥98%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

Docosadienoic acid methyl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of docosadienoic acid methyl ester in these solvents is approximately 30 mg/ml.

### Description

Docosadienoic acid methyl ester is an esterified version of 13Z,16Z-docosadienoic acid (Item No. 20749), which is a natural ω-6 polyunsaturated fatty acid (PUFA). Docosadienoic acid is an agonist of free fatty acid receptor 4 (FFAR4, also known as GPR120) and strongly inhibits the secretion of ghrelin by isolated mouse gastric cells.<sup>1</sup>

### Reference

1. Lu, X., Zhao, X., Feng, J., *et al.* Postprandial inhibition of gastric ghrelin secretion by long-chain fatty acid through GPR120 in isolated gastric ghrelin cells and mice. *Am. J. Physiol. Gastrointest. Liver Physiol.* **303(3)**, G367-G376 (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, 02/13/2024

#### 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