

# PRODUCT INFORMATION

## 4(Z),10(Z),13(Z),16(Z)-Docosatetraenoic Acid methyl ester

Item No. 10006866

**Formal Name:** 4Z,10Z,13Z,16Z-docosatetraenoic acid, methyl ester

**Synonyms:** *cis*-4,10,13,16-Docosatetraenoic Acid methyl ester, SFE 23:4

**MF:** C<sub>23</sub>H<sub>38</sub>O<sub>2</sub>

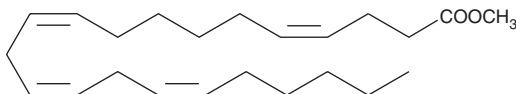
**FW:** 346.6

**Purity:** ≥98%

**Supplied as:** A solution in ethanol

**Storage:** -20°C

**Stability:** ≥1 year



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

### Laboratory Procedures

4(Z),10(Z),13(Z),16(Z)-Docosatetraenoic 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 4(Z),10(Z),13(Z),16(Z)-docosatetraenoic acid methyl ester in these solvents is approximately 30 mg/ml.

4(Z),10(Z),13(Z),16(Z)-Docosatetraenoic acid methyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 4(Z),10(Z),13(Z),16(Z)-docosatetraenoic acid methyl ester should be diluted with the aqueous buffer of choice. 4(Z),10(Z),13(Z),16(Z)-Docosatetraenoic acid methyl ester has a solubility of 0.1 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

4(Z),10(Z),13(Z),16(Z)-Docosatetraenoic acid methyl ester is a long chain polyunsaturated fatty acid. It is a minor fatty acid component of rat testis lipids.<sup>1</sup> The methyl ester is a neutral, more lipophilic form of the free acid.

### Reference

1. Yu, Q.T., Liu, B.N., Zhang, J.Y., *et al.* Location of double bonds in fatty acids of fish oil and rat testis lipids. Gas chromatography-mass spectrometry of the oxazoline derivatives. *Lipids* **24**(1), 79-83 (1989).

#### 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