

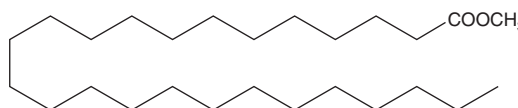
# PRODUCT INFORMATION



## Pentacosanoic Acid methyl ester

Item No. 10006451

**CAS Registry No.:** 55373-89-2  
**Formal Name:** pentacosanoic acid, methyl ester  
**Synonyms:** Methyl Pentacosanoate, SFE 26:0  
**MF:** C<sub>26</sub>H<sub>52</sub>O<sub>2</sub>  
**FW:** 396.7  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Pentacosanoic acid methyl ester is supplied as a crystalline solid. A stock solution may be made by dissolving the pentacosanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. Pentacosanoic acid methyl ester is soluble in the organic solvent ethanol at a concentration of approximately 0.5 mg/ml.

Pentacosanoic acid methyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pentacosanoic acid methyl ester should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Pentacosanoic acid methyl ester has a solubility of approximately 0.5 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

Long chain saturated fatty acids and their esters are common components of plant lipids. Pentacosanoic acid methyl ester is the methyl ester of the 25-carbon, fully saturated fatty acid pentacosanoic acid. This product is designed to be used as a standard in total fatty acid analysis.

#### 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, 03/29/2024

#### 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