

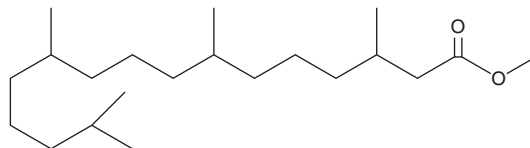
# PRODUCT INFORMATION



## Phytanic Acid methyl ester

Item No. 10007422

**CAS Registry No.:** 1118-77-0  
**Formal Name:** 3,7,11,15-tetramethyl-hexadecanoic acid, methyl ester  
**Synonyms:** Methyl Phytanate, SFE 21:0  
**MF:** C<sub>21</sub>H<sub>42</sub>O<sub>2</sub>  
**FW:** 326.6  
**Purity:** ≥95%  
**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

Phytanic acid methyl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the phytanic acid methyl ester under a gentle stream of nitrogen and immediately add the solvent of choice. Phytanic acid methyl ester is soluble in the organic solvent ethanol. The solubility of phytanic acid methyl ester in ethanol is approximately 5 mg/ml.

### Description

Phytanic acid methyl ester is a long-chain fatty acid methyl ester and esterified form of phytanic acid (Item No. 90360) that has been found in Greek tobacco.<sup>1</sup>

### Reference

1. Kimland, B., Aasen, A.J., and Enzell, C.R. Tobacco chemistry. 10. Volatile neutral constituents of Greek tobacco. *Acta Chem. Scand.* **26(6)**, 2177-2184 (1972).

#### 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, 03/27/2024

#### 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