

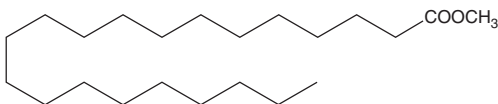
# PRODUCT INFORMATION



## Heneicosanoic Acid methyl ester

Item No. 26865

**CAS Registry No.:** 6064-90-0  
**Formal Name:** heneicosanoic acid, methyl ester  
**Synonyms:** C21:0 methyl ester,  
Heneicosylic Acid methyl ester,  
Methyl Heneicosanoate,  
Methyl Heneicosylate, SFE 22:0  
**MF:**  $C_{22}H_{44}O_2$   
**FW:** 340.6  
**Purity:**  $\geq 98\%$   
**Supplied as:** A solid  
**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

Heneicosanoic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the heneicosanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. Heneicosanoic acid methyl ester is soluble in organic solvents such as ethanol, chloroform, DMSO, and dimethyl formamide (DMF). The solubility of heneicosanoic acid methyl ester in ethanol and chloroform is approximately 10 and 50 mg/ml, respectively. Heneicosanoic acid methyl ester is slightly soluble in DMSO and DMF.

### Description

Heneicosanoic acid methyl ester is an ester form of heneicosanoic acid (Item No. 22593). It has been used as an internal standard in the quantification of fatty acids in dairy products and fatty acid methyl esters in fish by GC-MS.<sup>1,2</sup>

### References

1. Lin, H., Boylston, T.D., Chang, M.J., *et al.* Survey of the conjugated linoleic acid contents of dairy products. *J. Dairy Sci.* **78**(11), 2358-2365 (1995).
2. Spiric, A., Trbovic, D., Vranic, D., *et al.* Statistical evaluation of fatty acid profile and cholesterol content in fish (common carp) lipids obtained by different sample preparation procedures. *Anal. Chim. Acta* **672**(1-2), 66-71 (2010).

#### 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, 02/14/2024

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