

# PRODUCT INFORMATION



## Hexanoic Acid methyl ester

Item No. 33675

CAS Registry No.: 106-70-7

Synonyms: C6:0 methyl ester, Caproic Acid methyl ester, Methyl Caproate, Methyl Hexanoate, NSC 5023, SFE 7:0

MF:  $C_7H_{14}O_2$

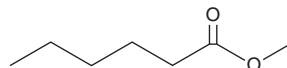
FW: 130.2

Purity:  $\geq 98\%$

Supplied as: A liquid

Storage:  $-20^{\circ}C$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

Hexanoic acid methyl ester is supplied as a liquid. A stock solution may be made by dissolving the hexanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. Hexanoic acid methyl ester is soluble in ethanol, chloroform, and ethyl ether.

### Description

Hexanoic acid methyl ester is an esterified form of hexanoic acid and a major volatile compound in strawberries.<sup>1</sup> Formulations containing hexanoic acid methyl ester have been used as cosmetic flavoring and fragrance agents.

### Reference

1. Miszczak, A., Forney, C.F., and Prange, R.K. Development of aroma volatiles and color during postharvest ripening of 'Kent' strawberries. *J. Amer. Soc. Hort. Sci.* **120**(4), 650-655 (1995).

#### 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)