

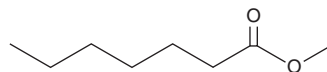
# PRODUCT INFORMATION



## Heptanoic Acid methyl ester

Item No. 33673

CAS Registry No.: 106-73-0  
Synonyms: C7:0 methyl ester, Methyl Heptanoate, SFE 8:0  
MF:  $C_8H_{16}O_2$   
FW: 144.2  
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

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

### Description

Heptanoic acid methyl ester is an esterified form of heptanoic acid (Item No. 33672). It is a target odor compound in Cascade and Willamette hops.<sup>1</sup>

### Reference

1. Sharp, D.C., Townsend, M.S., Qian, Y., *et al.* Effect of harvest maturity on the chemical composition of Cascade and Willamette hops. *J. Am. Soc. Brew. Chem.* **72**(4), 231-238 (2014).

#### 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