

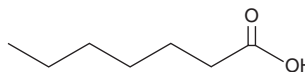
PRODUCT INFORMATION



Heptanoic Acid

Item No. 33672

CAS Registry No.: 111-14-8
Synonyms: C7:0, FA 7:0, Enanthic Acid, NSC 2192
MF: $C_7H_{14}O_2$
FW: 130.2
Purity: $\geq 98\%$
Supplied as: A liquid
Storage: $-20^\circ C$
Stability: ≥ 2 years



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

Laboratory Procedures

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

Description

Heptanoic acid is a medium-chain saturated fatty acid. It is a volatile component of the rancid odor of defective olive oil.¹ It has also been found as a volatile component in pig farm wastewater.² Heptanoic acid induces neurite outgrowth, a marker of neuronal differentiation, in PC12 cells in a concentration-dependent manner and increases the length of neurites when used at a concentration of 7.5 mM.³

References

1. Oliver-Pozo, C., Aparicio-Ruiz, R., Romero, I., *et al.* Analysis of volatile markers for virgin olive oil aroma defects by SPME-GC/FID: Possible sources of incorrect data. *J. Agric. Food Chem.* **63(48)**, 10477-10483 (2015).
2. Yo, S.-P. Analysis of volatile fatty acids in wastewater collected from a pig farm by a solid phase microextraction method. *Chemosphere* **38(4)**, 823-834 (1999).
3. Kamata, Y., Shiraga, H., Tai, A., *et al.* Induction of neurite outgrowth in PC12 cells by the medium-chain fatty acid octanoic acid. *Neuroscience* **146(3)**, 1073-1081 (2007).

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, 02/20/2024

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