

PRODUCT INFORMATION



5-Hexynoic Acid

Item No. 35507

CAS Registry No.: 53293-00-8

Synonyms: ALK-4, 5-Hexynoate

MF: $C_6H_8O_2$

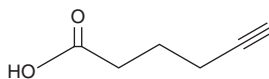
FW: 112.1

Purity: $\geq 95\%$

Supplied as: A neat liquid

Storage: -20°C

Stability: ≥ 4 years



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

Laboratory Procedures

5-Hexynoic acid is supplied as a neat liquid. A stock solution may be made by dissolving the 5-hexynoic acid in the solvent of choice, which should be purged with an inert gas. 5-Hexynoic acid is soluble (≥ 10 mg/ml) in acetonitrile and methanol.

Description

5-Hexynoic acid is a protein labeling compound.¹ It has been used to label proteins at acylation sites in a fatty acid synthase-dependent manner. 5-Hexynoic acid can be further conjugated to fluorescent probes or biotin *via* click chemistry for protein identification.

Reference

1. Karthigeyan, K.P., Zhang, L., Loiselle, D.R., *et al.* A bioorthogonal chemical reporter for fatty acid synthase-dependent protein acylation. *J. Biol. Chem.* **297**(5), 101272 (2021).

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, 12/03/2024

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