

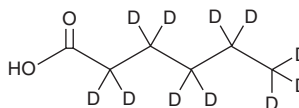
PRODUCT INFORMATION



Hexanoic Acid-d₁₁

Item No. 28082

CAS Registry No.: 95348-44-0
Synonyms: C6:0-d₁₁, Caproic Acid-d₁₁, FA 6:0-d₁₁
MF: C₆HD₁₁O₂
FW: 127.2
Chemical Purity: ≥98% (Hexanoic Acid)
Deuterium
Incorporation: ≥99% deuterated forms (d₁-d₁₁); ≤1% d₀
Supplied as: A semi-solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Low melting



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

Laboratory Procedures

Hexanoic acid-d₁₁ is intended for use as an internal standard for the quantification of hexanoic acid by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Hexanoic acid-d₁₁ is supplied as a semi-solid. A stock solution may be made by dissolving the hexanoic acid-d₁₁ in the solvent of choice, which should be purged with an inert gas. Hexanoic acid-d₁₁ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of hexanoic acid-d₁₁ in these solvents is approximately 30 mg/ml.

Description

Hexanoic acid is a short-chain saturated fatty acid that has been found in various animal fats used in biodiesel production.¹

Reference

1. Sales, E.A., Ghirardi, M.L., and Jorquera, O. Subcritical ethylic biodiesel production from wet animal fat and vegetable oils: A net energy ratio analysis. *Energy Convers. Manag.* **141**, 216-223 (2017).

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, 03/25/2024

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