

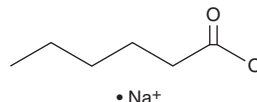
PRODUCT INFORMATION



Hexanoic Acid (sodium salt)

Item No. 45443

CAS Registry No.: 10051-44-2
Formal Name: hexanoic acid, monosodium salt
Synonyms: C6:0, Caproic Acid, FA 6:0, Sodium Hexanoate
MF: $C_6H_{11}O_2 \cdot Na$
FW: 138.1
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

Hexanoic acid (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the hexanoic acid (sodium salt) in the solvent of choice, which should be purged with an inert gas. Hexanoic acid (sodium salt) is soluble (≥ 10 mg/ml) in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of hexanoic acid (sodium salt) can be prepared by directly dissolving the solid in aqueous buffers. Hexanoic acid (sodium salt) is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Hexanoic acid is a short-chain saturated fatty acid. It has been found in various animal fats used in biodiesel production.¹ Spray application of hexanoic acid reduces *B. cinerea* lesion diameter in infected tomato plants.² It increases intratumoral numbers of CD8⁺ T cells and potentiates a reduction in tumor growth induced by an anti-programmed cell death protein 1 (PD-1) monoclonal antibody in a murine CT26 colorectal carcinoma model.³ Hexanoic acid (5% in the diet) prevents weight and fat mass gains and increases in hepatic triglyceride levels in mice fed a high-fat diet.⁴ It also increases plasma glucagon-like peptide 1 (GLP-1) levels and reduces blood glucose and plasma insulin levels in the same model.

References

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).
2. Leyva, M.O., Vicedo, B., Finiti, I., et al. Preventive and post-infection control of *Botrytis cinerea* in tomato plants by hexanoic acid. *Plant Pathol.* **57**(6), 1038-1046 (2008).
3. Seta, D., Nagahori, K., Katoh, Y., et al. Hexanoic acid intake enhances anti-tumor immune responses in colorectal cancer by reducing the immunosuppressive function of Tregs. *Mol. Nutr. Food Res.* **69**(22), e70229 (2025).
4. Ikeda, T., Takii, K., Omichi, Y., et al. Hexanoic acid improves metabolic health in mice fed high-fat diet. *Nutrients* **17**(17), 2868 (2025).

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, 03/24/2026

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