

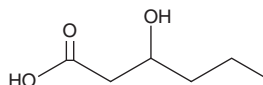
PRODUCT INFORMATION



3-hydroxy Hexanoic Acid

Item No. 24607

CAS Registry No.: 10191-24-9
Formal Name: 3-hydroxy-hexanoic acid
Synonyms: 3-hydroxy Caproic Acid, FA 6:0;O
MF: C₆H₁₂O₃
FW: 132.2
Purity: ≥95%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

3-hydroxy Hexanoic acid is supplied as a liquid. A stock solution may be made by dissolving the 3-hydroxy hexanoic acid in the solvent of choice. 3-hydroxy Hexanoic acid is soluble in organic solvents such as chloroform, ethanol, and methanol, which should be purged with an inert gas.

Description

3-hydroxy Hexanoic acid is a hydroxylated fatty acid that has been found in methyl-branched poly(3-hydroxyalkanoate) (PHA) polymers produced by *P. oleovorans*, *P. putida*, and *D. tsuruhatensis*.^{1,2} It is also found in the urine and serum of diabetic patients with ketoacidosis but not healthy patients or diabetic patients without ketoacidosis.³

References

1. Hazer, B., Lenz, R.W., and Fuller, R.C. Biosynthesis of methyl-branched poly(β-hydroxyalkanoate)s by *Pseudomonas oleovorans*. *Macromol.* **27(1)**, 45-49 (1994).
2. Razaif-Mazinah, M.R.M., Abis, S.N.S., Harun, H.I., et al. Unusual poly(3-hydroxyalkanoate) (PHA) biosynthesis behavior of *Pseudomonas putida* Bet001 and *Delftia tsuruhatensis* Bet002 isolated from palm oil mill effluent. *Biotechnol. Appl. Biochem.* **64(2)**, 259-269 (2017).
3. Niwa, T., Yamada, K., Ohki, T., et al. 3-Hydroxyhexanoic acid: An abnormal metabolite in urine and serum of diabetic ketoacidotic patients. *J. Chromatogr.* **337(1)**, 1-7 (1985).

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, 02/27/2027

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