

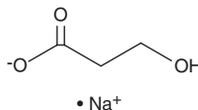
PRODUCT INFORMATION



3-Hydroxypropionic Acid (sodium salt)

Item No. 27603

CAS Registry No.: 6487-38-3
Synonym: Sodium β -hydroxypropionate
Formal Name: 3-hydroxy-propanoic acid, monosodium salt
MF: $C_3H_5O_3 \cdot Na$
FW: 112.1
Purity: $\geq 95\%$
Supplied as: A crystalline 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

3-Hydroxypropionic acid (sodium salt) is supplied as a crystalline solid. Aqueous solutions of 3-hydroxypropionic acid (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 3-hydroxypropionic acid (sodium salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

3-Hydroxypropionic acid is a building block in the synthesis of a variety of compounds with industrial applications, including acrylic acid, acrylamide, and 1,3-propanediol.^{1,2} Accumulation of 3-hydroxypropionic acid is also used as a biomarker of defects in propionic acid and branched-chain amino acid catabolism in patients with methylmalonic acidemia or propionic acidemia.³

References

1. David, Y., Oh, Y.H., Baylon, M.G., *et al.* Microbial production of 3-hydroxypropionic acid. *Industrial biotechnology: Products and processes*. Wittman, C. and Liao, J.C., editors, 1st edition, Wiley-VCH (2017).
2. Matsakas, L., Hrůzová, K., Rova, U., *et al.* Biological production of 3-hydroxypropionic acid: An update on the current status. *Fermentation* **4(13)**, 4010013 (2018).
3. Baumgartner, M.R., Hörster, F., Dionisi-Vici, C., *et al.* Proposed guidelines for the diagnosis and management of methylmalonic and propionic acidemia. *Orphanet J. Rare Dis.* **9:130**, (2014).

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, 11/11/2022

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