

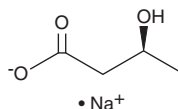
PRODUCT INFORMATION



(S)-3-Hydroxybutyrate (sodium salt)

Item No. 21822

CAS Registry No.: 127604-16-4
Formal Name: 3S-hydroxy-butanoic acid, monosodium salt
Synonyms: (S)-3-Hydroxybutanoate, (S)-3-Hydroxybutyric Acid
MF: C₄H₇O₃ • Na
FW: 126.1
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

(S)-3-Hydroxybutyrate (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the (S)-3-hydroxybutyrate (sodium salt) in the solvent of choice, which should be purged with an inert gas. (S)-3-Hydroxybutyrate (sodium salt) is soluble in the organic solvent ethanol at a concentration of approximately 5 mg/ml.

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 (S)-3-hydroxybutyrate (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of (S)-3-hydroxybutyrate (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

(S)-3-Hydroxybutyrate is a building block used in the synthesis of optically active fine chemicals, such as vitamins, antibiotics, pheromones, and flavor compounds.¹ (S)-3-Hydroxybutyrate is an enantiomer of the naturally occurring ketone body β-hydroxybutyrate, which is the (R) isomer.

Reference

1. Tseng, H.-C., Martin, C.H., Nielsen, D.R., *et al.* Metabolic engineering of *Escherichia coli* for enhanced production of (R)- and (S)-3-hydroxybutyrate. *Appl. Environ. Microbiol.* **75(10)**, 3137-3145 (2009).

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, 10/24/2022

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