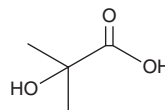


PRODUCT INFORMATION

2-Hydroxyisobutyric Acid

Item No. 43901

CAS Registry No.: 594-61-6
Formal Name: 2-hydroxy-2-methyl-propanoic acid
Synonyms: Acetonic Acid, 2-HIBA, 2-Methylactic Acid, NSC 4505, NSC 402158
MF: C₄H₈O₃
FW: 104.1
Purity: ≥95%
Supplied as: A 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

2-Hydroxyisobutyric acid is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxyisobutyric acid in the solvent of choice, which should be purged with an inert gas. 2-Hydroxyisobutyric acid is soluble (≥10 mg/ml) in ethanol and DMSO.

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 2-hydroxyisobutyric acid can be prepared by directly dissolving the solid in aqueous buffers. 2-Hydroxyisobutyric acid is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Hydroxyisobutyric acid is a metabolism modifying agent.^{1,2} It inhibits glutathione-induced decreases in NAD/NADH ratios, ATP levels, and maximal and basal oxygen consumption rates in PC12 adrenal medulla cells.¹ 2-Hydroxyisobutyric acid (10 mM) increases lipid droplet size, decreases reactive oxygen species (ROS), and increases lifespan in *C. elegans*.² It reduces blood glucose levels, decreases the latency to escape in the Barnes maze test, and increases brain Bdnf, nerve growth factor (Ngf), and glutathione (GSH) levels in a *db/db* mouse model of diabetic neuropathy when administered at doses of 15 or 30 mg/kg per day.¹

References

1. Xie, M., Gu, S., Liu, Y., et al. 2-Hydroxyisobutyric acid targeted binding to MT-ND3 boosts mitochondrial respiratory chain homeostasis in hippocampus to rescue diabetic cognitive impairment. *Redox Biol.* **79**, 103446 (2025).
2. Schifano, E., Conta, G., Preziosi, A., et al. 2-hydroxyisobutyric acid (2-HIBA) modulates ageing and fat deposition in *Caenorhabditis elegans*. *Front. Mol. Biosci.* **9**, 986022 (2022).

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, 08/18/2025

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