

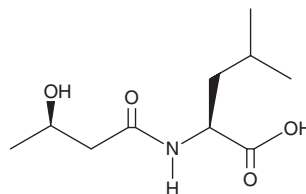
PRODUCT INFORMATION



β-Hydroxybutyrate leucine

Item No. 43377

Formal Name: ((R)-3-hydroxybutanoyl)-L-leucine
Synonym: BNB-Leu
MF: C₁₀H₁₉NO₄
FW: 217.3
Purity: ≥98%
Supplied as: A 25 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Description

β-Hydroxybutyrate leucine is a secondary ketone body metabolite composed of β-hydroxybutyrate (BHB) conjugated to the essential amino acid L-leucine (Item No. 34342).¹ It is formed endogenously by cytosolic non-specific dipeptidase 2 (CDNP2) following ketogenesis. β-Hydroxybutyrate leucine (50 mg/kg per day) reduces food intake and body weight in diet-induced obese (DIO) mice.

Reference

1. Moya-Garzon, M.D., Wang, M., Li, V.L., *et al.* A β-hydroxybutyrate shunt pathway generates anti-obesity ketone metabolites. *Cell* **188**(1), 175-186 (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, 11/21/2025

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