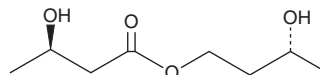


PRODUCT INFORMATION



BD-AcAc₂
Item No. 9003598

CAS Registry No.: 1208313-97-6
Formal Name: (R)-3-hydroxybutyl (R)-3-hydroxybutanoate
Synonyms: (R)-3-Hydroxybutyl (R)-3-hydroxybutyrate, ketone ester
MF: C₈H₁₆O₄
FW: 176.2
Purity: ≥95%
Supplied as: A 50 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

BD-AcAc₂ is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice.

Description

BD-AcAc₂ is a ketone ester with anti-obesity and anticonvulsant activities.^{1,2} Oral administration of BD-AcAc₂ (30% of total kcals) reduces body weight, food intake, and whole animal adiposity in mice fed an isocaloric diet.¹ BD-AcAc₂ (4 g/kg) increases the seizure threshold and blood levels of β-hydroxybutyrate in a rat model of seizures induced by pentylenetetrazole (PTZ; Item No. 18682).²

References

1. Deemer, S.E., Davis, R.A.H., Gower, B.A., *et al.* Concentration-dependent effects of a dietary ketone ester on components of energy balance in mice. *Front. Nutr.* **6**, 56 (2019).
2. Viggiano, A., Pilla, R., Arnold, P.I., *et al.* Anticonvulsant properties of an oral ketone ester in a pentylenetetrazole-model of seizure. *Brain Res.* **1618**, 50-54 (2015).

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, 01/13/2026

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