

PRODUCT INFORMATION



Bio-Ben

Item No. 39299

Formal Name: N-(1-hydroxy-1,3-dihydrobenzo[c][1,2]oxaborol-6-yl)-5-((3aS,4S,6aR)-2-oxohexahydro-1H-thieno[3,4-d]imidazol-4-yl)pentanamide

Synonym: Biotin-Benzoboroxole

MF: C₁₇H₂₂BN₃O₄S

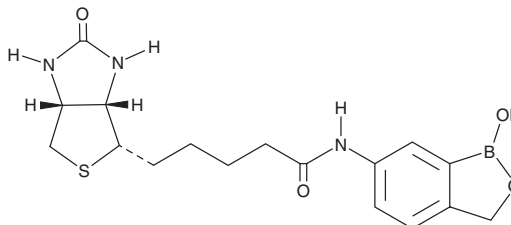
FW: 375.2

Purity: ≥98%

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

Bio-ben is supplied as a crystalline solid. A stock solution may be made by dissolving the bio-ben in the solvent of choice, which should be purged with an inert gas. Bio-ben is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of bio-ben in these solvents is approximately 5 mg/ml.

Description

Bio-ben is a probe for sulfenic acid. It has been used to label free sulfenic acid, as well as cysteine sulfenic acid in proteins and peptides.¹

Reference

1. Niu Y, Chen Z, Jiang Z, et al. Detection of cysteine sulfenic acid on *E. coli* proteins with a biotin-benzoboroxole probe. *ACS Chem. Biol.* **18**(6), 1351-1359 (2023).

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, 10/10/2025

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