

PRODUCT INFORMATION



DCP-Bio3

Item No. 28897

CAS Registry No.: 1001575-95-6
Formal Name: N-[4-[[5-[(3aS,4S,6aR)-hexahydro-2-oxo-1H-thieno[3,4-d]imidazol-4-yl]-1-oxopentyl]amino]butyl]-carbamic acid, 3-(2,4-dioxocyclohexyl)propyl ester

MF: C₂₄H₃₈N₄O₆S

FW: 510.6

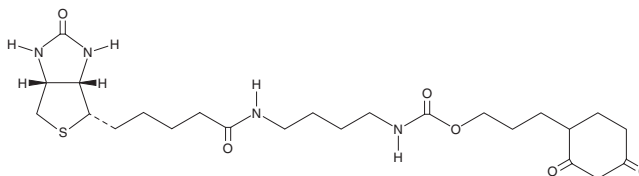
Purity: ≥90%

UV/Vis.: λ_{max}: 254 nm

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

DCP-Bio3 is supplied as a crystalline solid. A stock solution may be made by dissolving the DCP-Bio3 in the solvent of choice, which should be purged with an inert gas. DCP-Bio3 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of DCP-Bio3 in these solvents is approximately 20 mg/ml.

Description

DCP-Bio3 is a biotinylated affinity probe for the isolation of cysteine sulfenic acid-containing proteins.¹ It selectively forms a covalent adduct with a cysteine sulfenic acid form of the bacterial peroxidase alkyl hydroperoxidase (AhpC) over thiol, disulfide, and hyperoxidized forms of AhpC.

Reference

1. Poole, L.B., Klomsiri, C., Knaggs, S.A., *et al.* Fluorescent and affinity-based tools to detect cysteine sulfenic acid formation in proteins. *Bioconjug. Chem.* **18(6)**, 2004-2017 (2007).

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, 03/19/2024

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