

# PRODUCT INFORMATION



## 3-[2-N-(Biotinyl)aminoethylthio]propanoic Acid

Item No. 16461

**CAS Registry No.:** 104582-29-8  
**Formal Name:** 3-[[2-[[5-[(3aS,4S,6aR)-hexahydro-2-oxo-1H-thieno[3,4-d]imidazol-4-yl]-1-oxopentyl]amino]ethyl]dithio]-propanoic acid

**Synonym:** SS-Biotin  
**MF:** C<sub>15</sub>H<sub>25</sub>N<sub>3</sub>O<sub>4</sub>S<sub>3</sub>

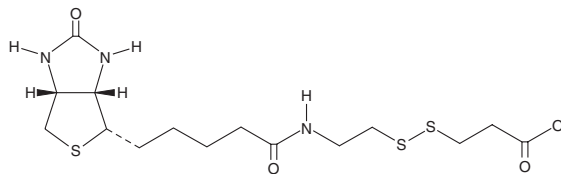
**FW:** 407.6

**Purity:** ≥98%

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥2 years



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

### Laboratory Procedures

3-[2-N-(Biotinyl)aminoethylthio]propanoic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-[2-N-(biotinyl)aminoethylthio]propanoic acid in the solvent of choice, which should be purged with an inert gas. 3-[2-N-(Biotinyl)aminoethylthio]propanoic acid is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 3-[2-N-(biotinyl)aminoethylthio]propanoic acid in these solvents is approximately 25 and 2 mg/ml, respectively.

3-[2-N-(Biotinyl)aminoethylthio]propanoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 3-[2-N-(biotinyl)aminoethylthio]propanoic acid should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 3-[2-N-(Biotinyl)aminoethylthio]propanoic acid has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

3-[2-N-(Biotinyl)aminoethylthio]propanoic acid features a biotin group and propanoic acid separated by a disulfide linkage. The carboxyl group of propanoic acid allows biotinylation of targets through either a chemical or enzymatic acylation reaction. The disulfide bond permits removal of the biotin group with reducing agents.

#### 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, 04/30/2021

#### 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