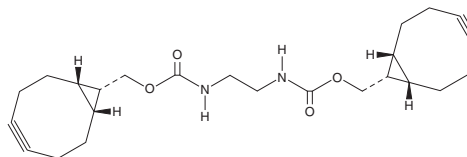


# PRODUCT INFORMATION



## BCN-E-BCN Item No. 21287

**Formal Name:** bis(((1R,8S,9s)-bicyclo[6.1.0]non-4-yn-9-yl)methyl) ethane-1,2-diyl dicarbamate  
**Synonym:** Click Tag™ BCN-E-BCN  
**MF:** C<sub>24</sub>H<sub>32</sub>N<sub>2</sub>O<sub>4</sub>  
**FW:** 412.5  
**Purity:** ≥95%  
**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

BCN-E-BCN is supplied as a crystalline solid. A stock solution may be made by dissolving the BCN-E-BCN in the solvent of choice. BCN-E-BCN is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of BCN-E-BCN in these solvents is approximately 3 and 1 mg/ml, respectively.

BCN-E-BCN is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BCN-E-BCN should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BCN-E-BCN has a solubility of approximately 25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

BCN-E-BCN is a strained cycloalkyne probe for detecting proteins that have been sulfenylated, the first intermediate step in protein oxidation.<sup>1</sup> BCN-E-BCN is selective for the sulfenic acid form of thiol oxidation and does not label free thiol, sulfinic, or sulfonic forms of proteins. The BCN-E-BCN protein conjugate can be detected using copper-free click chemistry to attach azide-bearing biotin or fluorescent tags. It is highly reactive and faster-acting than other sulfenic acid probes.

### Reference

1. McGarry, D.J., Shchepinova, M.M., Lilla, S., *et al.* A cell-permeable bicyclooctyne as a novel probe for the identification of protein sulfenic acids. *ACS Chem. Biol.* **11(12)**, 3300-3304 (2016).

#### 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, 02/15/2024

#### 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