

PRODUCT INFORMATION



Citrulline-specific Probe-biotin

Item No. 17450

CAS Registry No.: 2468149-70-2
Formal Name: (3aR,4R,6aS)-rel-hexahydro-N-[2-[[4-[1-[1-methyl-2-oxo-2-[[3-(2-oxoacetyl)phenyl]amino]ethyl]-1H-1,2,3-triazol-4-yl]-1-oxobutyl]amino]ethyl]-2-oxo-1H-thieno[3,4-d]imidazole-4-pentanamide

MF: C₂₉H₃₈N₈O₆S

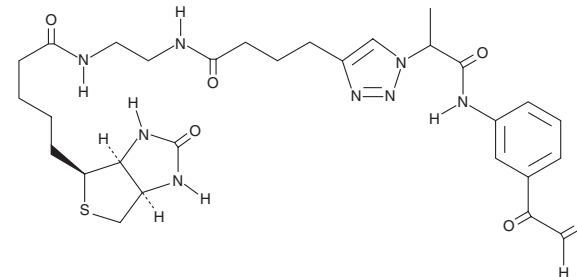
FW: 626.7

Purity: ≥75%

Supplied as: A crystalline solid

Storage: -80°C

Stability: ≥2 years



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

Description

Citrulline-specific probe-biotin is an affinity probe that allows for detection or immobilization of citrullinated proteins through interaction with the biotin ligand. It contains a phenylglyoxal group that selectively labels citrulline over arginine at acidic pH.¹ Citrulline-specific probe-biotin is a biotinylated form of citrulline-specific probe-rhodamine (Item No. 16172) that contains biotin in place of the rhodamine fluorophore.

Reference

1. Bicker, K.L., Subramanian, V., Chumanevich, A.A., *et al.* Seeing citrulline: Development of a phenylglyoxal-based probe to visualize protein citrullination. *J. Am. Chem. Soc.* **134**(41), 17015-17018 (2012).

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, 11/18/2020

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