

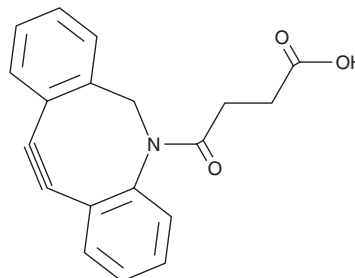
PRODUCT INFORMATION



DBCO acid

Item No. 41894

CAS Registry No.: 1353016-70-2
Formal Name: 11,12-didehydro- γ -oxo-dibenz[b,f]azocine-5(6H)-butanoic acid
Synonym: Dibenzocyclooctyne Acid
MF: C₁₉H₁₅NO₃
FW: 305.3
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

DBCO acid is supplied as a solid. A stock solution may be made by dissolving the DBCO acid in the solvent of choice, which should be purged with an inert gas. DBCO acid is sparingly soluble (1-10 mg/ml) in DMSO.

Description

DBCO acid is a copper-free click chemistry reagent.^{1,2} It has been used in the study of glycoprotein charge effects on the migration and invasion of 9-azido sialic acid-labeled cancer cells *in vitro*.¹ DBCO has also been used in the synthesis of DBCO-modified RNase A to deliver RNase A into cancer cells labeled with tetraacetyl-N-azidoacetylmannosamine (AAM) *in vitro* and *in vivo* to induce cytotoxicity and reduce tumor growth, respectively.²

References

1. Wang, M., Kejian, S., Ye, L., *et al.* Transforms of cell surface glycoproteins charge influences tumor cell metastasis via atypically inhibiting epithelial-mesenchymal transition including matrix metalloproteinases and cell junctions. *Bioconjug. Chem.* **34(8)**, 1498-1507 (2023).
2. Zhao, Z., Zhang, Z., Duan, S., *et al.* Cytosolic protein delivery via metabolic glycoengineering and bioorthogonal click reactions. *Biomater. Sci.* **9(13)**, 4639-4647 (2021).

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, 10/31/2024

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