

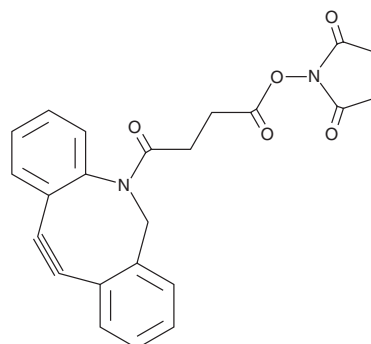
PRODUCT INFORMATION



DBCO-NHS ester

Item No. 40461

CAS Registry No.: 1353016-71-3
Formal Name: 11,12-didehydro-γ-oxo-dibenz[b,f]azocine-5(6H)-butanoic acid, 2,5-dioxo-1-pyrrolidinyl ester
Synonyms: DBCO N-hydroxysuccinimide ester, DIBAC N-hydroxysuccinimide ester, DIBAC-NHS ester, Dibenzocyclooctyne N-hydroxysuccinimide ester
MF: C₂₃H₁₈N₂O₅
FW: 402.4
Purity: ≥98%
Supplied as: A 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

DBCO-NHS ester is supplied as a solid. A stock solution may be made by dissolving the DBCO-NHS ester in the solvent of choice, which should be purged with an inert gas. DBCO-NHS ester is soluble in acetonitrile and DMSO.

Description

DBCO-NHS ester is an azido-reactive click-chemistry reagent.¹ It has been conjugated to various anticancer agents for targeting azido-labeled cancer cells in mice.

Reference

1. Wang, H., Wang, R., Cai, K., *et al.* Selective *in vivo* metabolic cell-labeling-mediated cancer targeting. *Nat. Chem. Biol.* **13**(4), 415-424 (2017).

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