

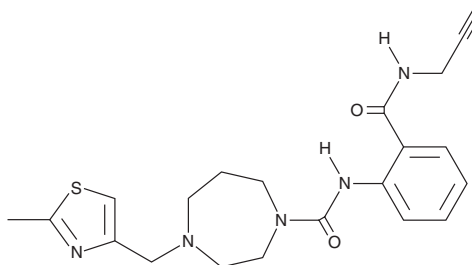
PRODUCT INFORMATION



MTDB-Alkyne

Item No. 38926

CAS Registry No.: 2923694-73-7
Formal Name: hexahydro-4-[(2-methyl-4-thiazolyl)methyl]-N-[2-[(2-propyn-1-ylamino)carbonyl]phenyl]-1H-1,4-diazepine-1-carboxamide
Synonyms: Click Tag™ MTDB-Alkyne, MTDB-Alk
MF: C₂₁H₂₅N₅O₂S
FW: 411.5
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 252 nm
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥3 years



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

Description

MTDB-alkyne is a clickable RNA pseudoknot binder.¹ It is composed of the β-coronavirus pseudoknot ligand MTDB that has been functionalized with a clickable alkyne group. MTDB-alkyne has been used in the synthesis of proximity-induced nucleic acid degrader (PINAD) molecules for the degradation of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) RNA.

Reference

1. Mikutis, S., Rebelo, M., Yankova, E., *et al.* Proximity-induced nucleic acid degrader (PINAD) approach to targeted RNA degradation using small molecules. *ACS Cent. Sci.* **9**(5), 892-904 (2023).

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, 01/26/2026

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