

PRODUCT INFORMATION

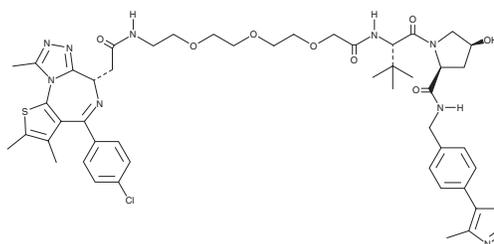


cis-MZ1

Item No. 45053

CAS Registry No.: 1797406-72-4
Formal Name: (4S)-N-[14-[(6S)-4-(4-chlorophenyl)-2,3,9-trimethyl-6H-thieno[3,2-f][1,2,4]triazolo[4,3-a][1,4]diazepin-6-yl]-1,13-dioxo-3,6,9-trioxa-12-azatetradec-1-yl]-3-methyl-L-valyl-4-hydroxy-N-[[4-(4-methyl-5-thiazolyl)phenyl]methyl]-L-prolinamide

MF: C₄₉H₆₀ClN₉O₈S₂
FW: 1,002.6
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

cis-MZ1 is supplied as a solid. A stock solution may be made by dissolving the cis-MZ1 in the solvent of choice, which should be purged with an inert gas. cis-MZ1 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

cis-MZ1 is an inactive diastereomer of the proteolysis-targeting chimera (PROTAC) MZ1 (Item No. 21622) that has been used as a negative control for MZ1-induced degradation of bromodomain-containing protein 4 (BRD4).¹

Reference

1. Zengerle, M., Chan, K.-H., and Ciulli, A. Selective small molecule induced degradation of the BET bromodomain protein BRD4. *ACS Chem. Biol.* **10**(8), 1770-1777 (2015).

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, 02/06/2026

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