

PRODUCT INFORMATION

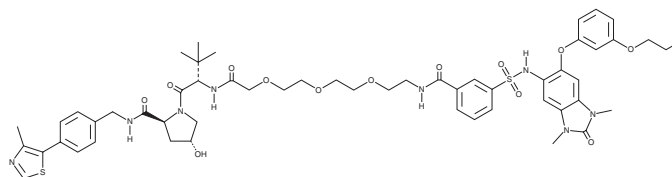


dTRIM24

Item No. 45749

CAS Registry No.: 2170695-14-2
Formal Name: (2S,4R)-1-((S)-15-(*tert*-butyl)-1-(3-(*N*-(1,3-dimethyl-2-oxo-6-(3-propoxyphenoxy)-2,3-dihydro-1H-benzo[d]imidazol-5-yl)sulfamoyl)phenyl)-1,13-dioxo-5,8,11-trioxa-2,14-diazahexadecan-16-oyl)-4-hydroxy-N-(4-(4-methylthiazol-5-yl)benzyl)pyrrolidine-2-carboxamide

MF: C₅₅H₆₈N₈O₁₃S₂
FW: 1,113.3
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

dTRIM24 is supplied as a solid. A stock solution may be made by dissolving the dTRIM24 in the solvent of choice, which should be purged with an inert gas. dTRIM24 is soluble (≥10 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

dTRIM24 is a proteolysis-targeting chimera (PROTAC) containing the ubiquitin ligase TRIM24 ligand IACS-7e conjugated to the von Hippel-Lindau (VHL) ligand VL-269 via a PEG linker.¹ It induces degradation of TRIM24 in HEK293-FT cells at 0.25-5 μM with maximum degradation at 5 μM. dTRIM24 (5 μM) reduces growth and induces apoptosis in MOLM-13 leukemia cells.

Reference

1. Gechijian, L.N., Buckley, D.L., Lawlor, M.A., et al. Functional TRIM24 degrader via conjugation of ineffectual bromodomain and VHL ligands. *Nat. Chem. Biol.* **14**(4), 405-412 (2018).

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, 05/22/2026

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