

PRODUCT INFORMATION

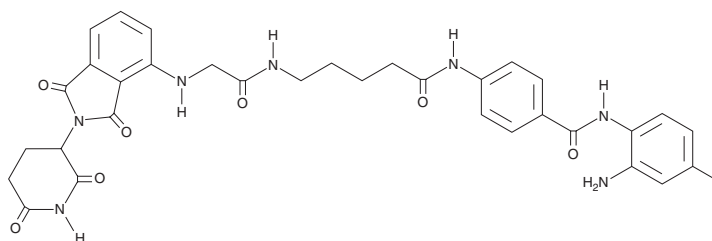


HD-TAC7

Item No. 45613

CAS Registry No.: 2978763-95-8
Formal Name: N-(2-amino-4-fluorophenyl)-4-[[5-[[2-[[2-(2,6-dioxo-3-piperidiny)]-2,3-dihydro-1,3-dioxo-1H-isoindol-4-yl]amino]acetyl]-amino]-1-oxopentyl]amino]-benzamide

MF: C₃₃H₃₂FN₇O₇
FW: 657.7
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

HD-TAC7 is supplied as a solid. A stock solution may be made by dissolving the HD-TAC7 in the solvent of choice, which should be purged with an inert gas. HD-TAC7 is slightly soluble (0.1-1 mg/ml) in methanol and DMSO.

Description

HD-TAC7 is a proteolysis-targeting chimera (PROTAC) composed of a class I histone deacetylase (HDAC) inhibitor linked to the cereblon ligand pomalidomide (Item No. 19877).¹ It binds to HDAC1, HDAC2, and HDAC3 (IC₅₀s = 3.6, 4.2, and 1.1 μM, respectively) and induces degradation of HDAC3 in RAW 264.7 cells with a half-maximal degradation concentration (DC₅₀) value of 0.32 μM, without inducing degradation of HDAC1 or HDAC2. HD-TAC7 (10 μM) increases acetylation of histone H3 lysine 27 (H3K27Ac) in RAW 264.7 macrophages but reduces the levels of NF-κB p65 and does not increase the expression of IL10 in LPS- and IFN-γ-stimulated RAW 264.7 macrophages, effects attributed to the pomalidomide moiety.

Reference

1. Cao, F., de Weerd, S., Chen, D., et al. Induced protein degradation of histone deacetylases 3 (HDAC3) by proteolysis targeting chimera (PROTAC). *Eur. J. Med. Chem.* **208**, 112800 (2020).

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, 05/06/2026

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