

PRODUCT INFORMATION



TL12-186

Item No. 40970

CAS Registry No.: 2250025-88-6

Formal Name: N-[2-[2-[2-[4-[4-[[5-chloro-4-[[2-[(1-methylethyl)sulfonyl]phenyl]amino]-2-pyrimidinyl]amino]phenyl]-1-piperazinyl]ethoxy]ethoxy]ethyl]-2-[[2-(2,6-dioxo-3-piperidinyl)-2,3-dihydro-1,3-dioxo-1H-isoindol-4-yl]amino]-acetamide

MF: $C_{44}H_{51}ClN_{10}O_9S$

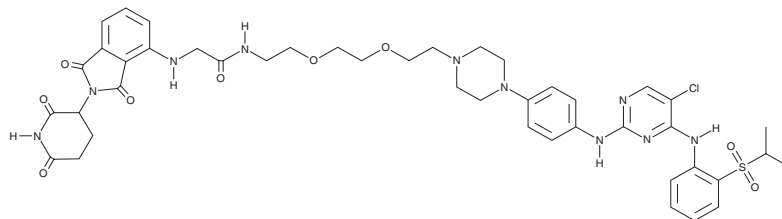
FW: 931.5

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

TL12-186 is supplied as a solid. A stock solution may be made by dissolving the TL12-186 in the solvent of choice, which should be purged with an inert gas. TL12-186 is soluble (≥ 10 mg/ml) in DMSO.

Description

TL12-186 is a proteolysis-targeting chimera (PROTAC) containing a multi-kinase inhibitor linked to a cereblon (CRBN) binding moiety.¹ It binds to CRBN ($IC_{50} = 12$ nM) and inhibits 193 kinases by greater than 90% in cell-free assays. TL12-186 (1-10,000 nM) inhibits the proliferation of wild-type CRBN-expressing but not CRBN^{-/-} MOLM-14 and MOLT-4 cells.

Reference

1. Huang, H.-T., Dobrovolsky, D., Paulk, J., et al. A chemoproteomic approach to query the degradable kinome using a multi-kinase degrader. *Cell Chem. Biol.* **25(1)**, 88-99 (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, 06/25/2024

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