

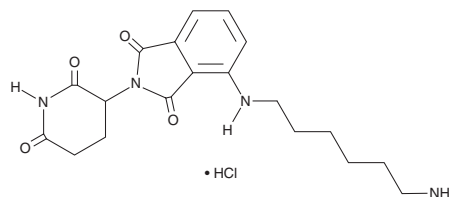
PRODUCT INFORMATION



Pomalidomide 4'-alkylC6-amine (hydrochloride)

Item No. 39985

CAS Registry No.: 2375194-37-7
Formal Name: 4-[(6-aminohexyl)amino]-2-(2,6-dioxo-3-piperidinyl)-1H-isoindole-1,3(2H)-dione, monohydrochloride
MF: $C_{19}H_{24}N_4O_4 \cdot HCl$
FW: 408.9
Purity: $\geq 98\%$
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

Pomalidomide 4'-alkylC6-amine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the pomalidomide 4'-alkylC6-amine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Pomalidomide 4'-alkylC6-amine (hydrochloride) is soluble in methanol and DMSO.

Description

Pomalidomide 4'-alkylC6-amine is a building block in the synthesis of proteolysis-targeting chimeras (PROTACs) and a derivative of the cereblon inhibitor pomalidomide (Item No. 19877).¹ It has been used in the synthesis of PROTACs containing BRD73954 (Item No. 16919) that drive histone deacetylase 8 (HDAC8) degradation.

Reference

1. Sun, Z., Deng, B., Yang, Z., *et al.* Discovery of pomalidomide-based PROTACs for selective degradation of histone deacetylase 8. *Eur. J. Med. Chem.* **239**, 114544 (2022).

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, 03/18/2024

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