

PRODUCT INFORMATION



PT-179

Item No. 43028

CAS Registry No.: 2924858-25-1

Formal Name: 2-(2,6-dioxo-3-piperidinyl)-5-(4-morpholinyl)-1H-isoindole-1,3(2H)-dione

MF: C₁₇H₁₇N₃O₅

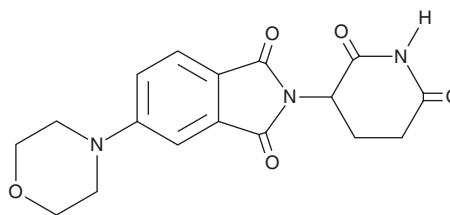
FW: 343.3

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

PT-179 is supplied as a solid. A stock solution may be made by dissolving the PT-179 in the solvent of choice, which should be purged with an inert gas. PT-179 is soluble (≥10 mg/ml) in DMSO.

Description

PT-179 is a molecular glue and ligand for the E3 ubiquitin ligase cereblon.¹ It has been used in conjunction with the zinc finger degron SD40 to induce degradation of SD40-tagged proteins.

Reference

1. Mercer, J.A.M., DeCarlo, S.J., Roy Burman, S.S., *et al.* Continuous evolution of compact protein degradation tags regulated by selective molecular glues. *Science* **383**(6688), eadk4422 (2024).

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, 04/07/2025

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