

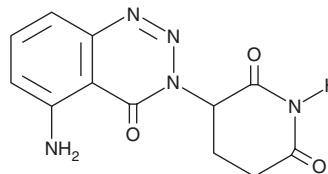
PRODUCT INFORMATION



TD-106

Item No. 44836

CAS Registry No.: 2250288-69-6
Formal Name: 3-(5-amino-4-oxo-1,2,3-benzotriazin-3(4H)-yl)-2,6-piperidinedione
MF: C₁₂H₁₁N₅O₂
FW: 273.3
Purity: ≥95%
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

TD-106 is supplied as a solid. A stock solution may be made by dissolving the TD-106 in the solvent of choice, which should be purged with an inert gas. TD-106 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

TD-106 is a modulator of cereblon.¹ It binds to cereblon in a cellular thermal shift assay (CETSA) using HEK293 cells. TD-106 (1-1,000 nM) induces degradation of the cereblon targets Ikaros family zinc finger protein 1 (IKZF1) and IKZF3 in NCI H929 multiple myeloma cells and is cytotoxic to NCI H929 cells (CC₅₀ = 0.039 μM). It also reduces tumor growth in a TMD8 diffuse large B cell lymphoma (DLBCL) mouse xenograft model when administered at a dose of 50 mg/kg per day. TD-106 has been used in the synthesis of proteolysis-targeting chimeras (PROTACs) with anticancer activity.

Reference

1. Kim, S.A., Go, A., Jo, S.H., *et al.* A novel cereblon modulator for targeted protein degradation. *Eur. J. Med. Chem.* **166**, 65-74 (2019).

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, 01/22/2026

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