

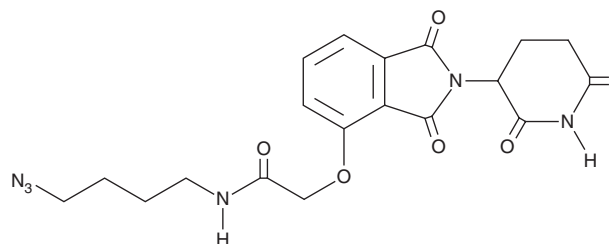
PRODUCT INFORMATION



Azido-Thalidomide

Item No. 26146

CAS Registry No.: 2098488-36-7
Formal Name: N-(4-azidobutyl)-2-[[2-(2,6-dioxo-3-piperidiny)-2,3-dihydro-1,3-dioxo-1H-isoindol-4-yl]oxy]-acetamide
Synonym: Click Tag™ Azido-Thalidomide
MF: C₁₉H₂₀N₆O₆
FW: 428.4
Purity: ≥95%
UV/Vis.: λ_{max}: 219 nm
Supplied as: A crystalline 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

Azido-thalidomide is supplied as a crystalline solid. A stock solution may be made by dissolving the azido-thalidomide in the solvent of choice, which should be purged with an inert gas. Azido-thalidomide is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of azido-thalidomide in these solvents is approximately 12 mg/ml.

Azido-thalidomide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, azido-thalidomide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Azido-thalidomide has a solubility of approximately 0.11 mg/ml in a 1:8 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Azido-thalidomide is a building block in the synthesis of proteolysis-targeting chimera (PROTAC) technologies.¹ It contains thalidomide, which is a ligand for the E3 ubiquitin ligase cereblon, and can be conjugated to alkynylated ligands for target proteins *via* click chemistry. Azido-thalidomide has been used in the synthesis of a SirReal-based PROTAC for the degradation of sirtuin 2 (Sirt2) in cells.

Reference

- Schiedel, M., Herp, D., Hammelmann, S., *et al.* Chemically induced degradation of Sirtuin 2 (Sirt2) by a proteolysis targeting chimera (PROTAC) based on sirtuin rearranging ligands (SirReals). *J. Med. Chem.* **61**(2), 482-491 (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, 08/23/2023

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