

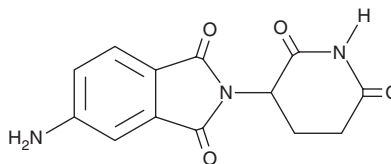
# PRODUCT INFORMATION



## 5-Aminothalidomide

Item No. 39514

**CAS Registry No.:** 191732-76-0  
**Formal Name:** 5-amino-2-(2,6-dioxo-3-piperidinyl)-1H-isoindole-1,3(2H)-dione  
**MF:** C<sub>13</sub>H<sub>11</sub>N<sub>3</sub>O<sub>4</sub>  
**FW:** 273.2  
**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

5-Aminothalidomide is supplied as a solid. A stock solution may be made by dissolving the 5-aminothalidomide in the solvent of choice, which should be purged with an inert gas. 5-Aminothalidomide is soluble in acetonitrile and DMSO.

### Description

5-Aminothalidomide is a derivative of the immunomodulatory compound (±)-thalidomide (Item No. 14610). It inhibits the proliferation of EA.hy926 endothelial cells (IC<sub>50</sub> = 200 μM).<sup>1</sup> 5-Aminothalidomide is an intermediate in the synthesis of a PROTAC c-MET degrader with antiproliferative activity in c-MET-sensitive EBC-1 and Hs 746T cancer cells.<sup>2</sup> It is also an intermediate in the synthesis of a thalidomide derivative that reduces clinical scores in a mouse model of experimental autoimmune encephalomyelitis (EAE).<sup>1</sup>

### References

1. Contino-Pépin, C., Parat, A., Périno, S., *et al.* Preliminary biological evaluations of new thalidomide analogues for multiple sclerosis application. *Bioorg. Med. Chem. Lett.* **19(3)**, 878-881 (2009).
2. Li, P., Jia, C., Fan, Z., *et al.* Discovery of novel exceptionally potent and orally active c-MET PROTACs for the treatment of tumors with MET alterations. *Acta Pharm. Sin. B* **13(6)**, 2715-2735 (2023).

#### 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, 02/19/2024

#### 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