

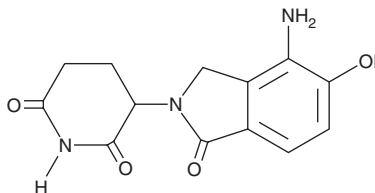
PRODUCT INFORMATION



Hydroxy Lenalidomide

Item No. 34491

CAS Registry No.: 1421593-78-3
Formal Name: 3-(4-amino-1,3-dihydro-5-hydroxy-1-oxo-2H-isoindol-2-yl)-2,6-piperidinedione
Synonym: 5-hydroxy Lenalidomide
MF: C₁₃H₁₃N₃O₄
FW: 275.3
Purity: ≥90%
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

Hydroxy lenalidomide is supplied as a solid. A stock solution may be made by dissolving the hydroxy lenalidomide in the solvent of choice, which should be purged with an inert gas. Hydroxy lenalidomide is soluble in a 1:1 solution of acetonitrile:methanol. Hydroxy lenalidomide is also soluble in DMSO.

Description

Hydroxy lenalidomide is a metabolite of the thalidomide analog and anticancer agent lenalidomide (Item No. 14643).¹

Reference

1. Chen, N., Wen, L., Lau, H., *et al.* Pharmacokinetics, metabolism and excretion of [¹⁴C]-lenalidomide following oral administration in healthy male subjects. *Cancer Chemother. Pharmacol.* **69**(3), 789-797 (2012).

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, 11/09/2022

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