

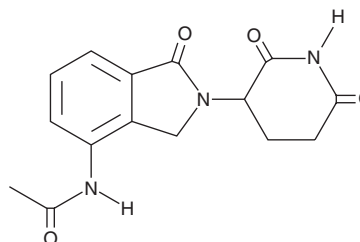
PRODUCT INFORMATION



N-acetyl Lenalidomide

Item No. 34421

CAS Registry No.: 1421593-80-7
Formal Name: N-[2-(2,6-dioxo-3-piperidiny)-2,3-dihydro-1-oxo-1H-isoindol-4-yl]-acetamide
MF: $C_{15}H_{15}N_3O_4$
FW: 301.3
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 219 nm
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

N-acetyl Lenalidomide is supplied as a solid. A stock solution may be made by dissolving the N-acetyl lenalidomide in the solvent of choice, which should be purged with an inert gas. N-acetyl Lenalidomide is soluble in the organic solvent ethanol.

Description

N-acetyl Lenalidomide is a metabolite of lenalidomide (Item No. 14643).¹ It is formed by hydrolysis of lenalidomide in plasma.

Reference

1. Chen, N., Wen, L., Lau, H., *et al.* Pharmacokinetics, metabolism and excretion of [^{14}C]-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, 11/22/2022

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