

PRODUCT INFORMATION



Nampt-IN-1

Item No. 26066

CAS Registry No.: 1698878-14-6
Formal Name: 2-hydroxy-2-methyl-N-[1,2,3,4-tetrahydro-2-[2-(3-pyridinyloxy)acetyl]-6-isoquinoliny]-1-propanesulfonamide

Synonyms: LSN3154567,
Nicotinamide Phosphoribosyltransferase-IN-1

MF: C₂₀H₂₅N₃O₅S

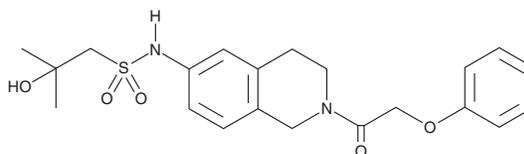
FW: 419.5

Purity: ≥98%

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

Nampt-IN-1 is supplied as a solid. A stock solution may be made by dissolving the Nampt-IN-1 in the solvent of choice, which should be purged with an inert gas. Nampt-IN-1 is soluble in DMSO.

Description

Nampt-IN-1 is a potent inhibitor of nicotinamide phosphoribosyltransferase (Nampt; IC₅₀ = 3.1 nM).¹ It is selective for Nampt over a panel of greater than 100 human kinases at 1 μM. Nampt-IN-1 inhibits NAD⁺ formation in and proliferation of HCT116 cells (IC₅₀s = 1.8 and 8.9 nM, respectively). It inhibits cell growth in a panel of cancer cell lines (IC₅₀s = <0.2-2 μM). Nampt-IN-1 inhibits NAD⁺ formation in an A2780 ovarian carcinoma mouse xenograft model (ED₅₀ = 2 mg/kg).

Reference

1. Zhao, G., Green, C.F., Hui, Y.H., *et al.* Discovery of a highly selective NAMPT inhibitor that demonstrates robust efficacy and improved retinal toxicity with nicotinic acid coadministration. *Mol. Cancer Ther.* **16**(12), 2677-2688 (2017).

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, 10/04/2022

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