

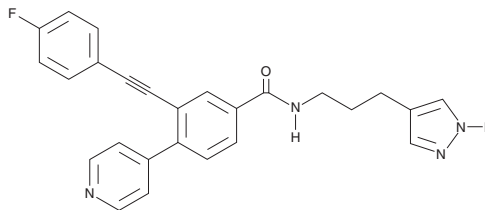
PRODUCT INFORMATION



OT-82

Item No. 38782

CAS Registry No.: 1800487-55-1
Formal Name: 3-[2-(4-fluorophenyl)ethynyl]-N-[3-(1H-pyrazol-4-yl)propyl]-4-(4-pyridinyl)-benzamide
MF: C₂₆H₂₁FN₄O
FW: 424.5
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

OT-82 is supplied as a solid. A stock solution may be made by dissolving the OT-82 in the solvent of choice, which should be purged with an inert gas. OT-82 is soluble in methanol and DMSO.

Description

OT-82 is a nicotinamide phosphoribosyltransferase (Nampt) inhibitor.¹ It inhibits recombinant human and recombinant mouse Nampt in a cell-free assay when used at concentrations ranging from 0.1 to 10 μM. OT-82 is cytotoxic against a panel of 12 hematological cancer cells (IC₅₀s = 1.01-5.64 nM). *In vivo*, OT-82 (25 and 50 mg/kg) increases survival and decreases tumor volume in MV4-11 and HEL 92.1.7 mouse xenograft models.

Reference

1. Korotchkina, L., Kazyulkin, D., Komarov, P.G., *et al.* OT-82, a novel anticancer drug candidate that targets the strong dependence of hematological malignancies on NAD biosynthesis. *Leukemia* **34(7)**, 1828-1839 (2020).

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, 03/12/2024

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