

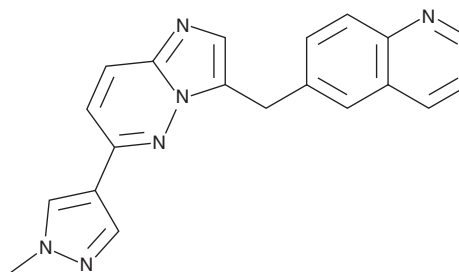
PRODUCT INFORMATION



NVP-BVU972

Item No. 34495

CAS Registry No.: 1185763-69-2
Formal Name: 6-[[6-(1-methyl-1H-pyrazol-4-yl)imidazo[1,2-b]pyridazin-3-yl]methyl]-quinoline
MF: C₂₀H₁₆N₆
FW: 340.4
Purity: ≥98%
UV/Vis.: λ_{max}: 231, 254 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

NVP-BVU972 is supplied as a solid. A stock solution may be made by dissolving the NVP-BVU972 in the solvent of choice, which should be purged with an inert gas. NVP-BVU972 is soluble in DMSO.

Description

NVP-BVU972 is an inhibitor of Met kinase (IC₅₀ = 14 nM).¹ It is selective for Met over a panel of 62 additional kinases (IC₅₀s = >1 μM for all). It inhibits proliferation of Met-amplified GTL-16 gastric, MKN45 gastric, and EBC-1 lung cancer cells (IC₅₀s = 66, 32, and 82 nM, respectively). NVP-BVU972 inhibits Met autophosphorylation in GTL-16 cells and Met phosphorylation induced by hepatocyte growth factor (HGF) in A549 cells (IC₅₀s = 7.3 and 22 nM, respectively).

Reference

1. Tiedt, R., Degenkolbe, E., Furet, P., *et al.* A drug resistance screen using a selective MET inhibitor reveals a spectrum of mutations that partially overlap with activating mutations found in cancer patients. *Cancer Res.* **71**(15), 5255-5264 (2011).

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