

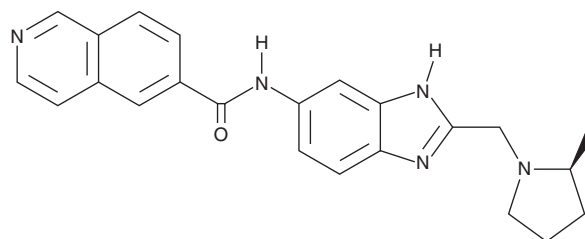
PRODUCT INFORMATION



NVS-MLLT-1

Item No. 41318

CAS Registry No.: 2561471-13-2
Formal Name: N-[2-[[[(2S)-2-methyl-1-pyrrolidinyl]methyl]-1H-benzimidazol-6-yl]-6-isoquinolinecarboxamide
MF: C₂₃H₂₃N₅O
FW: 385.5
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

NVS-MLLT-1 is supplied as a solid. A stock solution may be made by dissolving the NVS-MLLT-1 in the solvent of choice, which should be purged with an inert gas. NVS-MLLT-1 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in methanol.

Description

NVS-MLLT-1 is an inhibitor of the YEATS domain-containing proteins MLLT1 and MLLT3 (IC₅₀s = 0.15 and 0.254 μM, respectively).¹ It is selective for MLLT1 and MLLT3 over the YEATS domain-containing proteins YEATS2 and YEATS4 (IC₅₀s = >20 μM for both), as well as over the cat eye syndrome critical region protein 2 (CECR2) bromodomain (IC₅₀ = >40 μM). NVS-MLLT-1 binds to wild-type MLLT1 (IC₅₀ = 0.18 μM) and MLLT1 containing an in-frame insertion (T1 mutant) or deletion (T2 mutant) mutation (K_ds = 0.24 and 0.22 μM, respectively).² See the Structural Genomics Consortium (SGC) website for more information.

References

1. NVS-MLLT-1 a potent and selective inhibitor of YEATS proteins. Retrieved July 29, 2024, from <https://www.thesgc.org/chemical-probes/nvs-mlt-1>.
2. Ni, X., Londregan, A.T., Owen, D.R., et al. Structure and inhibitor binding characterization of oncogenic MLLT1 mutants. *ACS Chem. Biol.* **16**(4), 571-578 (2021).

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, 07/29/2024

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