

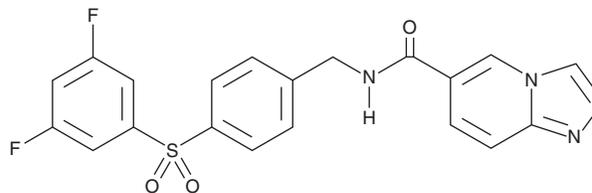
PRODUCT INFORMATION



GNE-617

Item No. 22463

CAS Registry No.: 1362154-70-8
Formal Name: N-[[4-[(3,5-difluorophenyl)sulfonyl]phenyl]methyl]-imidazo[1,2-a]pyridine-6-carboxamide
MF: C₂₁H₁₅F₂N₃O₃S
FW: 427.4
Purity: ≥98%
UV/Vis.: λ_{max}: 238 nm
Supplied as: A crystalline 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

GNE-617 is supplied as a crystalline solid. A stock solution may be made by dissolving the GNE-617 in the solvent of choice, which should be purged with an inert gas. GNE-617 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GNE-617 in these solvents is approximately 20 mg/ml.

GNE-617 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GNE-617 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GNE-617 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GNE-617 is a potent inhibitor of nicotinamide phosphoribosyltransferase (Namt; IC₅₀ = 5 nM).¹ It has antiproliferative activity in U251 glioblastoma, HT-1080 fibrosarcoma, PC3 prostate, MiaPaCa2 pancreatic, and HCT116 colon cancer cell lines (IC₅₀s = 1.8, 2.1, 2.7, 7.4, and 2.0 nM, respectively). In a U251 mouse xenograft model, GNE-617 (25 mg/kg once per day) led to rapid tumor regression.

Reference

1. Zheng, X., Bauer, P., Baumeister, T., *et al.* Structure-based discovery of novel amide-containing nicotinamide phosphoribosyltransferase (nampt) inhibitors. *J. Med. Chem.* **56(16)**, 6413-6433 (2013).

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, 01/03/2023

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