

PRODUCT INFORMATION



GNE-618

Item No. 37223

CAS Registry No.: 1362151-42-5

Formal Name: N-[[4-[[3-(trifluoromethyl)phenyl]sulfonyl]phenyl]methyl]-1H-pyrazolo[3,4-b]pyridine-5-carboxamide

MF: $C_{21}H_{15}F_3N_4O_3S$

FW: 460.4

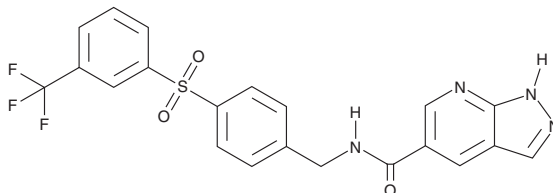
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 227 nm

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

GNE-618 is supplied as a solid. A stock solution may be made by dissolving the GNE-618 in the solvent of choice, which should be purged with an inert gas. GNE-618 is soluble in DMSO.

Description

GNE-618 is an inhibitor of nicotinamide phosphoribosyltransferase (Nampt; $IC_{50} = 6.1$ nM for the human enzyme).¹ It reduces NAD^+ levels in PC3 prostate cancer cells ($EC_{50} = 1.2$ nM) and inhibits the proliferation of A2780 ovarian cancer cells ($IC_{50} = 4.3$ nM). GNE-618 (100 mg/kg) induces tumor regression in an A2780 mouse xenograft model.

Reference

1. Zheng, X., Bair, K.W., Bauer, P., et al. Identification of amides derived from 1H-pyrazolo[3,4-b]pyridine-5-carboxylic acid as potent inhibitors of human nicotinamide phosphoribosyltransferase (NAMPT). *Bioorg. Med. Chem. Lett.* **23**(20), 5488-5497 (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, 04/26/2023

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