

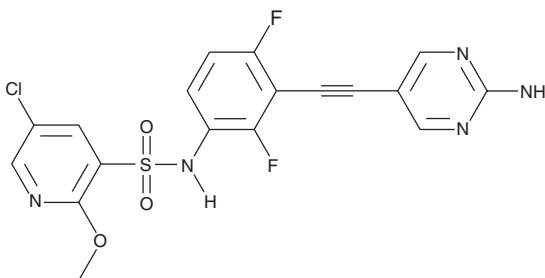
PRODUCT INFORMATION



GCN2iB

Item No. 35897

CAS Registry No.: 2183470-12-2
Formal Name: N-[3-[2-(2-amino-5-pyrimidinyl)ethynyl]-2,4-difluorophenyl]-5-chloro-2-methoxy-3-pyridinesulfonamide
MF: C₁₈H₁₂ClF₂N₅O₃S
FW: 451.8
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

GCN2iB is supplied as a solid. A stock solution may be made by dissolving the GCN2iB in the solvent of choice, which should be purged with an inert gas. GCN2iB is soluble in organic solvents such as acetonitrile and DMSO.

Description

GCN2iB is an inhibitor of general control nonderepressible 2 kinase (GCN2; IC₅₀ = 2.4 nM).¹ It is selective for GCN2 over a panel of 465 kinases but does inhibit MAP2K5, eukaryotic translation initiation factor 2α kinase 2 (EIF2AK2), and MAP3K20 at 1 μM. GCN2iB (1 μM), in combination with L-asparaginase, reduces the viability of mouse embryonic fibroblasts (MEFs). It decreases the levels of activating transcription factor 4 (ATF4) and phosphorylated GCN2 in combination with L-asparaginase in CCRF CEM leukemia cells when used at a concentration of 0.4 μM. *In vivo*, GCN2iB (10 mg/kg twice per day), in combination with L-asparaginase, reduces tumor volume in CCRF CEM and MV4-11 leukemia mouse xenograft models.

Reference

1. Nakamura, A., Nambu, T., Ebara, S., *et al.* Inhibition of GCN2 sensitizes ASNS-low cancer cells to asparaginase by disrupting the amino acid response. *Proc. Natl. Acad. Sci. USA* **115(33)**, E7776-E7785 (2018).

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, 08/20/2024

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