

PRODUCT INFORMATION

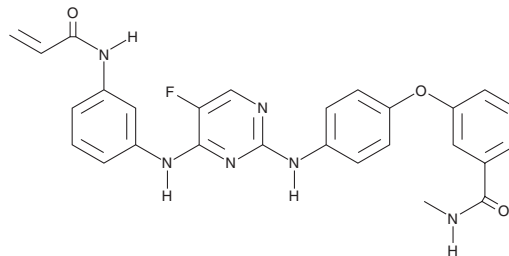


CNX-774

Item No. 21475

CAS Registry No.: 1202759-32-7
Formal Name: 4-[4-[[5-fluoro-4-[[3-[(1-oxo-2-propen-1-yl)amino]phenyl]amino]-2-pyrimidinyl]amino]phenoxy]-N-methyl-2-pyridinecarboxamide

MF: C₂₆H₂₂FN₇O₃
FW: 499.5
Purity: ≥98%
UV/Vis.: λ_{max}: 272 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

CNX-774 is supplied as a crystalline solid. A stock solution may be made by dissolving the CNX-774 in the solvent of choice. CNX-774 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of CNX-774 in these solvents is approximately 25 and 30 mg/ml, respectively.

CNX-774 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CNX-774 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CNX-774 has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CNX-774 is a potent, selective, and irreversible inhibitor of Bruton's tyrosine kinase (BTK), an important kinase in the B cell antigen receptor pathway, with IC₅₀ values of <1 nM in biochemical assays and 1-10 nM in cellular assays.¹ It inhibits proliferation in HMC-1.1 and HMC-1.2 cells (IC₅₀ = 2.82 and 0.38 μM) but has no effect on cell cycle progression.²

References

1. Akinleye, A., Chen, Y., Mukhi, N., *et al.* Ibrutinib and novel BTK inhibitors in clinical development. *J. Hematol. Oncol.* **6(59)** (2013).
2. Smiljkovic, D., Blatt, K., Stefanzi, G., *et al.* BTK inhibition is a potent approach to block IgE-mediated histamine release in human basophils. *Allergy* **1-11** (2017).

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, 12/06/2022

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