

PRODUCT INFORMATION

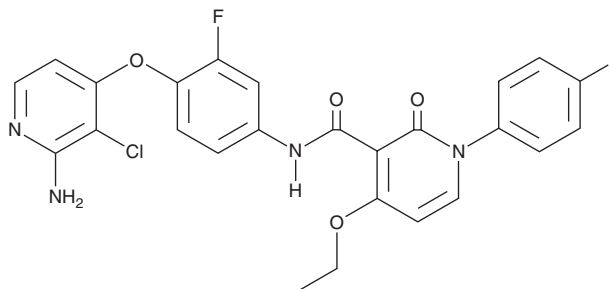


BMS-777607

Item No. 18517

CAS Registry No.: 1025720-94-8
Formal Name: N-[4-[(2-amino-3-chloro-4-pyridinyl)oxy]-3-fluorophenyl]-4-ethoxy-1-(4-fluorophenyl)-1,2-dihydro-2-oxo-3-pyridinecarboxamide

Synonym: BMS-817378
MF: C₂₅H₁₉ClF₂N₄O₄
FW: 512.9
Purity: ≥98%
UV/Vis.: λ_{max}: 204, 289 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

BMS-777607 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMS-777607 in the solvent of choice, which should be purged with an inert gas. BMS-777607 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of BMS-777607 in these solvents is approximately 5 and 20 mg/ml, respectively.

BMS-777607 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BMS-777607 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BMS-777607 has a solubility of approximately 0.1 mg/ml in a 1:7 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BMS-777607 is an inhibitor of the Met kinase family. It inhibits Ron, Met, Tyro-3, and Axl with IC₅₀ values of 1.8, 3.9, 4.3, and 1.1 nM, respectively.¹ At higher concentrations it is reported to inhibit Mer, FLT3, Aurora B, Lck, and VEGFR2 (IC₅₀s = 14, 16, 78, 120, and 180 nM, respectively).¹ In cancer cells, BMS-777607 induces polyploidy with multiple sets of chromosomes, as well as suppresses metastasis.^{1,2} In tumor xenograft models, BMS-777607 at 50 mg/kg/mouse inhibits the growth of tumor cells expressing Met/Ron.¹

References

1. Sharma, S., Zeng, J.-Y., Zhuang, C.-M., et al. Small-molecule inhibitor BMS-777607 induces breast cancer cell polyploidy with increased resistance to cytotoxic chemotherapy agents. *Mol. Cancer Ther.* **12**(5), 725-736 (2013).
2. Dai, Y. and Siemann, D.W. BMS-777607, a small-molecule met kinase inhibitor, suppresses hepatocyte growth factor-stimulated prostate cancer metastatic phenotype *in vitro*. *Mol. Cancer Ther.* **9**(6), 1554-1561 (2010).

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, 11/21/2022

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