

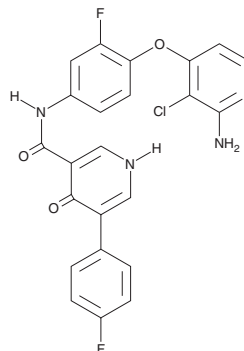
PRODUCT INFORMATION



BMS 794833

Item No. 26180

CAS Registry No.: 1174046-72-0
Formal Name: N-[4-[(2-amino-3-chloro-4-pyridinyl)oxy]-3-fluorophenyl]-5-(4-fluorophenyl)-1,4-dihydro-4-oxo-3-pyridinecarboxamide
MF: C₂₃H₁₅ClF₂N₄O₃
FW: 468.8
Purity: ≥98%
UV/Vis.: λ_{max}: 298 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 794833 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMS 794833 in the solvent of choice. BMS 794833 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of BMS 794833 in these solvents is approximately 17 mg/ml.

Description

BMS 794833 is an inhibitor of c-Met and VEGF receptor 2 (VEGFR2) with IC₅₀ values of 1.7 and 15 nM, respectively, that also inhibits Axl, Flt3, and Ron.^{1,2} It inhibits proliferation of GTL-16 gastric carcinoma cells (IC₅₀ = 39 nM) and induces tumor stasis in GTL-16 and U87 mouse xenograft models when administered at a dose of 25 mg/kg.¹

References

1. Borzilleri, R.M. and Cai, Z.-W. 4-Pyridinone compounds and their use for cancer. *Bristol-Myers Squibb Company*. **WO2009094417A1** (2009).
2. Moran-Jones, K. The therapeutic potential of targeting the HGF/cMET axis in ovarian cancer. *Mol. Diagn. Ther.* **20(3)**, 199-212 (2016).

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, 12/01/2022

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