

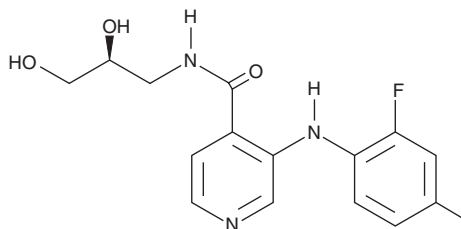
PRODUCT INFORMATION



AS-703026

Item No. 11226

CAS Registry No.: 1236699-92-5
Formal Name: N-[(2S)-2,3-dihydroxypropyl]-3-[(2-fluoro-4-iodophenyl)amino]-4-pyridinecarboxamide
Synonyms: MSC1936369B, Pimasertib
MF: C₁₅H₁₅FIN₃O₃
FW: 431.2
Purity: ≥95%
UV/Vis.: λ_{max}: 204, 284, 353 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

AS-703026 is supplied as a crystalline solid. A stock solution may be made by dissolving the AS-703026 in the solvent of choice, which should be purged with an inert gas. AS-703026 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AS-703026 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

AS-703026 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AS-703026 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AS-703026 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

The mitogen-activated protein kinases, MEK1 and 2, are dual-specificity threonine/tyrosine kinases that play key roles in the activation of the Ras-Raf-MEK-ERK pathway and are often upregulated in a variety of cancer cell types.¹ AS-703026 is an orally bioavailable small molecule that selectively binds to and inhibits MEK1/2, preventing the activation of downstream effector proteins and transcription factors. It potently inhibits growth and survival of human INA-6 multiple myeloma cells and cytokine-induced osteoclast differentiation with IC₅₀ values of 10 and 18.2 nM, respectively.² In mice bearing H929 MM xenograft tumors, 30 mg/kg AS-703026 reduced tumor growth significantly, which correlated with downregulated ERK1/2 activity, induced PARP cleavage, and decreased microvessels *in vivo*.² At 10 μM, AS-703026 suppressed proliferation and transformation of K-Ras mutated colorectal cancer cells resistant to EGFR antibody therapy.³

References

1. VanScyoc, W.S., Holdgate, G.A., Sullivan, J.E., *et al.* Enzyme kinetics and binding studies on inhibitors of MEK protein kinase. *Biochemistry* (2008).
2. Kim, K., Kong, S.-Y., Fulciniti, M., and *et al.* Blockade of the MEK/ERK signalling cascade by AS703026, a novel selective MEK1/2 inhibitor, induces pleiotropic anti-myeloma activity *in vitro* and *in vivo*. *Br. J. Haematol.* **149**(4), 537-549 (2010).
3. Yoon, J., Koo, K.-H., and Choi, K.-Y. MEK1/2 inhibitors AS703026 and AZD6244 may be potential therapies for KRAS mutated colorectal cancer that is resistant to EGFR monoclonal antibody therapy. *Cancer Res.* **71** (445), (2011).

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, 10/17/2022

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