

PRODUCT INFORMATION

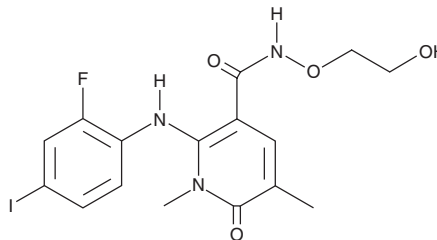


AZD 8330

Item No. 11588

CAS Registry No.: 869357-68-6
Formal Name: 2-[(2-fluoro-4-iodophenyl)amino]-1,6-dihydro-N-(2-hydroxyethoxy)-1,5-dimethyl-6-oxo-3-pyridinecarboxamide

Synonym: ARRY-424704
MF: C₁₆H₁₇FIN₃O₄
FW: 461.2
Purity: ≥98%
UV/Vis.: λ_{max}: 245, 321 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

AZD 8330 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 8330 in the solvent of choice, which should be purged with an inert gas. AZD 8330 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AZD 8330 in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

AZD 8330 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AZD 8330 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AZD 8330 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AZD 8330 is a selective allosteric inhibitor of MEK1/2 (IC₅₀ = 7 nM). It blocks phosphorylation of ERK1/2 and inhibits proliferation of certain cancer cells.¹ AZD 8330 is orally bioavailable, inhibits ERK phosphorylation *in vivo*, and reduces tumor growth in a Calu-6 nude rat xenograft model.²

References

1. Baranski, Z., Booij, T.H., Kuijjer, M.L., *et al.* MEK inhibition induces apoptosis in osteosarcoma cells with constitutive ERK1/2 phosphorylation. *Genes Cancer* **6**(11-12), 503-512 (2015).
2. Cohen, R.B., Aamdal, S., Nyakas, M., *et al.* A phase I dose-finding, safety and tolerability study of AZD8330 in patients with advanced malignancies. *Eur. J. Cancer* **49**(7), 1521-1529 (2013).

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, 10/31/2022

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