

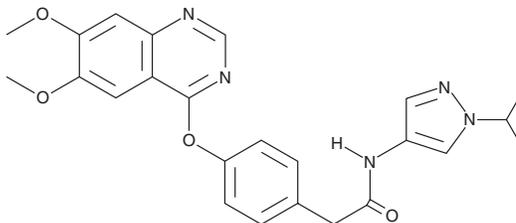
PRODUCT INFORMATION



AZD 2932

Item No. 17741

CAS Registry No.: 883986-34-3
Formal Name: 4-[[6,7-dimethoxy-4-quinazolinyloxy]-N-[1-(1-methylethyl)-1H-pyrazol-4-yl]-benzeneacetamide
MF: C₂₄H₂₅N₅O₄
FW: 447.5
Purity: ≥98%
UV/Vis.: λ_{max}: 239, 320 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 2932 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 2932 in the solvent of choice, which should be purged with an inert gas. AZD 2932 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AZD 2932 in these solvents is approximately 0.25 mg/ml.

Description

AZD 2932 is an inhibitor of VEGFR2, PDGFRβ, CSF1R, c-KIT, and FLT3 receptor tyrosine kinases (IC₅₀s = 4, 8, 100, 9, and 7 nM, respectively).¹ It also inhibits Pdgfra phosphorylation (IC₅₀ = 2 nM) in lysates from C6 rat glial tumors grown in athymic mice. AZD 2932 (12.5-50 mg/kg) reduces tumor growth and inhibits PDGFRβ and VEGFR2 phosphorylation in a 1:1 ratio in a C6 mouse xenograft model. Xenograft models using non-PDGFR expressing tumor cells were insensitive to AZD 2932 treatment.

Reference

1. Plé, P., Jung, F., Ashton, S., *et al.* Discovery of AZD2932, a new quinazoline ether inhibitor with high affinity for VEGFR-2 and PDGFR tyrosine kinases. *Bioorg. Med. Chem. Lett.* **22(1)**, 262-266 (2012).

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, 12/14/2022

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