

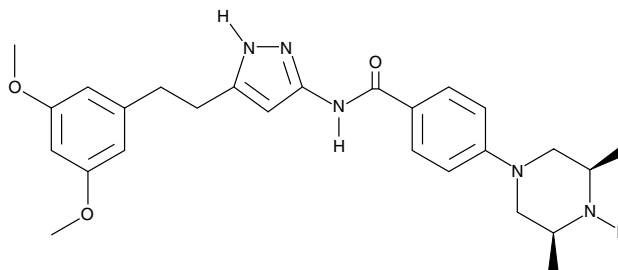
Product Information



AZD 4547

Item No. 13111

CAS Registry No.: 1035270-39-3
Formal Name: *rel*-N-[5-[2-(3,5-dimethoxyphenyl)ethyl]-1H-pyrazol-3-yl]-4-[(3R,5S)-3,5-dimethyl-1-piperazinyl]-benzamide
MF: C₂₆H₃₃N₅O₃
FW: 463.6
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 308 nm



Laboratory Procedures

For long term storage, we suggest that AZD 4547 be stored as supplied at -20°C. It should be stable for at least two years.

AZD 4547 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 4547 in the solvent of choice. AZD 4547 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of AZD 4547 in ethanol is approximately 1.6 mg/ml and approximately 25 mg/ml in DMSO and DMF.

AZD 4547 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AZD 4547 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AZD 4547 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.

AZD 4547 is a selective inhibitor of the fibroblast growth factor receptor (FGFR) tyrosine kinase with IC₅₀ values of 0.2, 2.5, and 1.8 nM for FGFR1, 2, and 3, respectively.¹ It displays weaker activity against FGFR4 (IC₅₀ = 165 nM), VEGFR2 (IC₅₀ = 24 nM), and a panel of other tyrosine and serine/threonine kinases such as insulin-like growth factor 1 receptor (IC₅₀ = 581 nM), cyclin-dependent kinase 2 (IC₅₀ = >10 μM), and p38 (IC₅₀ = >100 μM).¹ AZD 4547 has demonstrated antiproliferative activity in tumor cell lines expressing deregulated FGFRs (IC₅₀s = 18-281 nM).¹ In an FGFR-driven human tumor xenograft mouse model, oral administration of AZD 4547 at concentrations up to 12.5 mg/kg/day resulted in potent dose-dependent antitumor activity.¹

Reference

1. Gavine, P.R., Mooney, L., Kilgour, E., *et al.* AZD4547: An orally bioavailable, potent, and selective inhibitor of the fibroblast growth factor receptor tyrosine kinase family. *Cancer Res.* **72**(8), 2045-2056 (2012).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13111

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 11/10/2014

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com