

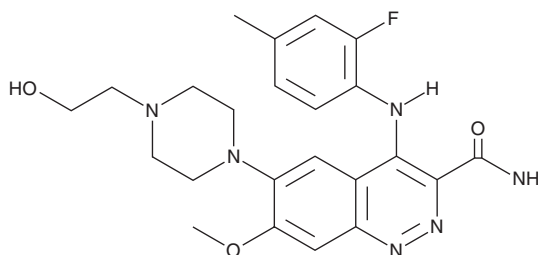
PRODUCT INFORMATION



AZD 7507

Item No. 37485

CAS Registry No.: 1041852-85-0
Formal Name: 4-[(2-fluoro-4-methylphenyl)amino]-6-[4-(2-hydroxyethyl)-1-piperazinyl]-7-methoxy-3-cinnolinecarboxamide
MF: C₂₃H₂₇FN₆O₃
FW: 454.5
Purity: ≥98%
UV/Vis.: λ_{max}: 292, 365 nm
Supplied as: A 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 7507 is supplied as a solid. A stock solution may be made by dissolving the AZD 7507 in the solvent of choice, which should be purged with an inert gas. AZD 7507 is soluble in the organic solvent DMSO.

Description

AZD 7507 is an inhibitor of CSF-1 receptor tyrosine kinase (FMS; IC₅₀ = 3 nM).¹ It is selective for FMS over c-KIT, aurora kinase 5 (ARK-5), FMS-related tyrosine kinase 4 (FLT4), p90 ribosomal S6 kinase (RSK1), and 77 other kinases at 1 μM. AZD 7507 (100 nM) induces apoptosis in isolated Fms⁺ mouse bone marrow cells but not Fms⁻ mouse bone marrow cells. *In vivo*, AZD 7507 (100 mg/kg per day) reduces the levels of tumor-associated Fms⁺ macrophages, increases the number of tumor-associated CD8⁺ T cells, decreases tumor weight, and increases survival in a UN-KPC-960 murine model of pancreatic ductal adenocarcinoma (PDAC). It decreases α-smooth muscle actin (α-Sma) levels and increases E-cadherin levels on isolated UN-KPC-960 tumor tissue, as well as reduces the levels of Ifn-γ, Il-10, Il-17a, chemokine (C-X-C motif) ligand 2 (Cxcl2), chemokine (C-C motif) ligand 2 (Ccl2), Ccl7, and Ccl12 in protein lysates from the same tumors.

Reference

1. Candido, J.B., Morton, J.P., Bailey, P., *et al.* CSF1R⁺ macrophages sustain pancreatic tumor growth through T cell suppression and maintenance of key gene programs that define the squamous subtype. *Cell Rep.* 23(5), 1448-1460 (2018).

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, 10/12/2023

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