

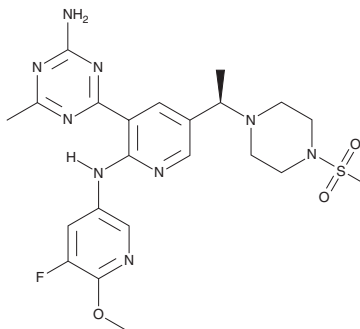
PRODUCT INFORMATION



AMG 511

Item No. 41853

CAS Registry No.: 1253573-53-3
Formal Name: 4-[2-[(5-fluoro-6-methoxy-3-pyridinyl)amino]-5-[(1R)-1-[4-(methylsulfonyl)-1-piperazinyl]ethyl]-3-pyridinyl]-6-methyl-1,3,5-triazin-2-amine
MF: C₂₂H₂₈FN₉O₃S
FW: 517.6
Purity: ≥98%
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

AMG 511 is supplied as a solid. A stock solution may be made by dissolving the AMG 511 in the solvent of choice, which should be purged with an inert gas. AMG 511 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

An inhibitor of class I PI3Ks (K_i s = 4, 6, 2, and 1 nM for PI3K α , PI3K β , PI3K δ , and PI3K γ , respectively).¹ It is selective for class I PI3Ks over 372 other protein kinases at 1 μ M. AMG 511 inhibits phosphorylation of Akt in U87MG glioblastoma cells (IC_{50} = 4 nM) and reduces tumor growth in a U87MG mouse xenograft model when administered at doses of 1, 3, and 10 mg/kg per day.

Reference

1. Norman, M.H., Andrews, K.L., Bo, Y.Y., *et al.* Selective class I phosphoinositide 3-kinase inhibitors: Optimization of a series of pyridyltriazines leading to the identification of a clinical candidate, AMG 511. *J. Med. Chem.* **55**(17), 7796-7816 (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, 11/01/2024

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