

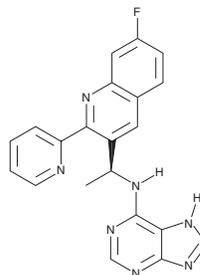
PRODUCT INFORMATION



AMG 319

Item No. 17383

CAS Registry No.: 1608125-21-8
Formal Name: 7-fluoro- α S-methyl-N-9H-purin-6-yl-2-(2-pyridinyl)-3-quinolinemethanamine
MF: C₂₁H₁₆FN₇
FW: 385.4
Purity: \geq 98%
UV/Vis.: λ_{max} : 267 nm
Supplied as: A crystalline solid
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

AMG 319 is supplied as a crystalline solid. A stock solution may be made by dissolving the AMG 319 in the solvent of choice, which should be purged with an inert gas. AMG 319 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AMG 319 in DMSO is approximately 5 mg/ml and approximately 3 mg/ml in ethanol and DMF.

AMG 319 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AMG 319 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AMG 319 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AMG 319 is a potent inhibitor of phosphatidylinositol 3-kinase (PI3K) isoform δ (PI3K δ ; IC₅₀ = 18 nM) that less effectively targets the α , β , and γ isoforms (IC₅₀s = 33, 2.7, and 0.85 μ M, respectively).¹ It blocks anti-IgM/CD40L-induced proliferation (IC₅₀ = 8.6 nM) and reduces Akt phosphorylation (IC₅₀ = 1.5 nM) in B cells.¹ AMG 319 is orally bioavailable and reduces inflammation in two animal models.¹

Reference

1. Cushing, T.D., Hao, X., Shin, Y., *et al.* Discovery and in vivo evaluation of (S)-N-(1-(7-fluoro-2-(pyridin-2-yl)quinolin-3-yl)ethyl)-9H-purin-6-amine (AMG319) and related PI3K δ inhibitors for inflammation and autoimmune disease. *J. Med. Chem.* **58**(1), 480-511 (2015).

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, 11/10/2022

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