

PRODUCT INFORMATION

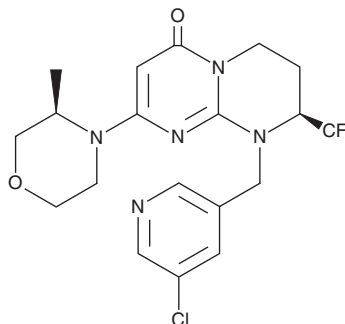


SAR405

Item No. 16979

CAS Registry No.: 1523406-39-4
Formal Name: (8S)-9-[(5-chloro-3-pyridinyl)methyl]-6,7,8,9-tetrahydro-2-[(3R)-3-methyl-4-morpholinyl]-8-(trifluoromethyl)-4H-pyrimido[1,2-a]pyrimidin-4-one

MF: C₁₉H₂₁ClF₃N₅O₂
FW: 443.9
Purity: ≥98%
UV/Vis.: λ_{max}: 235, 278 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

SAR405 is supplied as a crystalline solid. A stock solution may be made by dissolving the SAR405 in the solvent of choice, which should be purged with an inert gas. SAR405 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SAR405 in these solvents is approximately 5, 10, and 15 mg/ml, respectively.

SAR405 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SAR405 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. SAR405 has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SAR405 is an inhibitor of the phosphoinositide 3-kinase (PI3K) class III isoform Vps34 (IC₅₀ = 1.2 nM).¹ It is without effect (IC₅₀ > 10,000 nM) on other PI3K isoforms and mTOR. SAR405 alters vesicle trafficking and inhibits autophagy by blocking autophagosome formation.¹ SAR405 also synergizes with the mTOR inhibitor everolimus (Item No. 11597) in preventing the proliferation of renal tumor cell lines.¹

Reference

1. Ronan, B., Flamand, O., Vescovi, L., *et al.* A highly potent and selective Vps34 inhibitor alters vesicle trafficking and autophagy. *Nat. Chem. Biol.* **10**(12), 1013-1019 (2014).

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, 12/06/2022

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