

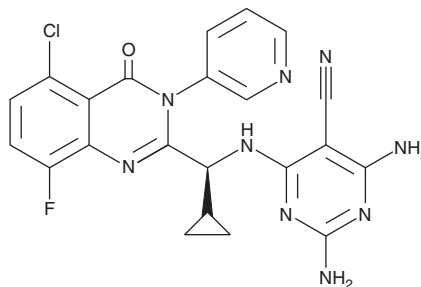
PRODUCT INFORMATION



GS-9901

Item No. 42105

CAS Registry No.: 1640247-87-5
Formal Name: 2,4-diamino-6-[[[(S)-[5-chloro-8-fluoro-3,4-dihydro-4-oxo-3-(3-pyridinyl)-2-quinazoliny]cyclopropylmethyl]amino]-5-pyrimidinicarbonitrile
MF: C₂₂H₁₇ClFN₉O
FW: 477.9
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

GS-9901 is supplied as a solid. A stock solution may be made by dissolving the GS-9901 in the solvent of choice, which should be purged with an inert gas. GS-9901 is sparingly soluble (1-10 mg/ml) in methanol and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

GS-9901 is an inhibitor of PI3K δ (IC₅₀ = 1 nM).¹ It is selective for PI3K δ over PI3K α , PI3K β , and PI3K γ (IC₅₀s = 750, 100, and 190 nM, respectively) and a panel of 456 additional kinases, including 11 lipid and 56 mutant kinases, at 10 μ M. GS-9901 activates basophils in isolated human whole blood (EC₅₀ = 1 nM). *In vivo*, GS-9901 (1 and 3 mg/kg) decreases ankle swelling and disease severity in a rat model of collagen-induced arthritis.

Reference

1. Patel, L., Chandrasekhar, J., Evarts, J., *et al.* Discovery of orally efficacious phosphoinositide 3-kinase δ inhibitors with improved metabolic stability. *J. Med. Chem.* **59(19)**, 9228-9242 (2016).

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, 11/04/2024

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