

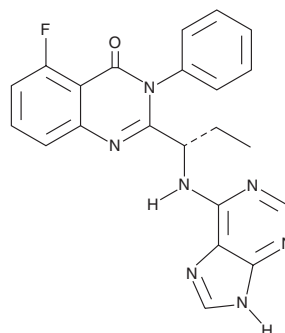
PRODUCT INFORMATION



CAL-101

Item No. 15279

CAS Registry No.: 870281-82-6
Formal Name: 5-fluoro-3-phenyl-2-[(1S)-1-(9H-purin-6-ylamino)propyl]-4(3H)-quinazolinone
Synonyms: GS-1101, Idelalisib
MF: C₂₂H₁₈N₇O
FW: 415.4
Purity: ≥98%
UV/Vis.: λ_{max}: 269, 311 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

CAL-101 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAL-101 in the solvent of choice, which should be purged with an inert gas. CAL-101 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAL-101 in these solvents is approximately 15, 25, and 30 mg/ml, respectively.

CAL-101 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAL-101 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAL-101 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

The phosphatidylinositol 3-kinase (PI3K) signaling pathway has central roles in cell growth, development, and survival.^{1,2} CAL-101 is a cell permeable inhibitor of the PI3K catalytic subunit p110δ (IC₅₀ = 2.5 nM) that demonstrates 40- to 300-fold selectivity against other PI3K class I enzymes (IC₅₀s = 820, 565, and 89 nM for p110α, β, and γ, respectively).³ At 1 μM it can block constitutive PI3K signaling in malignant B-cell lines and primary patient tumor cells, resulting in decreased phosphorylation of Akt and other downstream effectors, an increase in poly(ADP-ribose) polymerase, and apoptosis.³ It has also been shown to inhibit the chemotaxis of chronic lymphocytic leukemia cells and to downregulate the secretion of chemokines triggered by B-cell receptor signaling.⁴

References

1. Hennessy, B.T., Smith, D.L., Ram, P.T., *et al.* Exploiting the PI3K/AKT pathway for cancer drug discovery. *Nat. Rev. Drug Discov.* **4**(12), 988-1004 (2005).
2. Hirsch, E., Ciruolo, E., Ghigo, A., *et al.* Taming the PI3K team to hold inflammation and cancer at bay. *Pharmacol. Ther.* **118**(2), 192-205 (2008).
3. Lannutti, B.H., Meadows, S.A., Hermann, S.E.M., *et al.* CAL-101, a p110δ selective phosphatidylinositol-3-kinase inhibitor for the treatment of B-cell malignancies, inhibits PI3K signaling and cellular viability. *Blood* **117**(2), 591-594 (2011).
4. Hoellenriegel, J., Meadows, S.A., Sivina, M., *et al.* The phosphoinositide 3'-kinase delta inhibitor, CAL-101, inhibits B-cell receptor signaling and chemokine networks in chronic lymphocytic leukemia. *Blood* **118**(13), 3603-3612 (2011).

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, 11/07/2022

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