

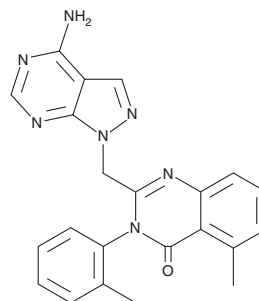
PRODUCT INFORMATION



PIK-293

Item No. 21283

CAS Registry No.: 900185-01-5
Formal Name: 2-[[4-amino-1H-pyrazolo[3,4-d]pyrimidin-1-yl)methyl]-5-methyl-3-(2-methylphenyl)-4(3H)-quinazolinone
MF: C₂₂H₁₉N₇O
FW: 397.4
Purity: ≥98%
UV/Vis.: λ_{max}: 227, 271, 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

PIK-293 is supplied as a crystalline solid. A stock solution may be made by dissolving the PIK-293 in the solvent of choice, which should be purged with an inert gas. PIK-293 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PIK-293 in ethanol is approximately 0.25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

PIK-293 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PIK-293 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. PIK-293 has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

PIK-293 is a cell-permeable, selective inhibitor of the phosphatidylinositol 3-kinase (PI3K) catalytic subunit p110δ (IC₅₀s = 100, 25, 10, and 0.24 μM for p110 subunit isoforms α, β, γ, and δ, respectively).¹

Reference

1. Knight, Z.A., Gonzalez, B., Feldman, M.E., *et al.* A pharmacological map of the PI3-K family defines a role for p110α in insulin signaling. *Cell* **125**, 733-747 (2006).

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, 11/18/2022

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