

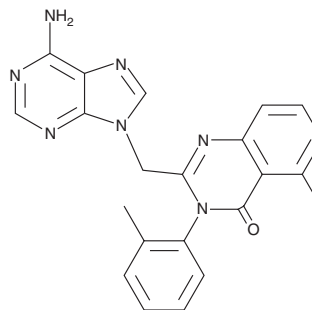
PRODUCT INFORMATION



IC-87114

Item No. 11589

CAS Registry No.: 371242-69-2
Formal Name: 2-[(6-amino-9H-purin-9-yl)methyl]-5-methyl-3-(2-methylphenyl)-4(3H)-quinazolinone
MF: C₂₂H₁₉N₇O
FW: 397.4
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 262, 310 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

IC-87114 is supplied as a crystalline solid. A stock solution may be made by dissolving the IC-87114 in the solvent of choice, which should be purged with an inert gas. IC-87114 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of IC-87114 in these solvents is approximately 30 mg/ml.

IC-87114 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, IC-87114 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. IC-87114 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

IC-87114 is a cell-permeable selective inhibitor of the PI3K catalytic subunit p110δ (IC₅₀ = 0.5 μM).¹ It less effectively inhibits p110γ and p110β (IC₅₀ = 29 and 75 μM, respectively) and has no significant effect on p110α and several other kinases.^{1,2} This product is used to elucidate the role of p110δ in cells, including neutrophils, natural killer cells, and other types of leukocytes.¹⁻⁵ It has also been used in mice.⁶

References

1. Sadhu, C., Masinovsky, B., Dick, K., *et al.* Essential role of phosphoinositide 3-kinase δ in neutrophil directional movement. *J. Immunol.* **170**(5), 2647-2654 (2003).
2. Sadhu, C., Dick, K., Tino, W.T., *et al.* Selective role of PI3Kδ in neutrophil inflammatory responses. *Biochem. Biophys. Res. Commun.* **308**(4), 764-769 (2003).
3. Saudemont, A., Garçon, F., Yadi, H., *et al.* p110γ and p110δ isoforms of phosphoinositide 3-kinase differentially regulate natural killer cell migration in health and disease. *Proc. Natl. Acad. Sci. USA* **106**(14), 5795-5800 (2009).
4. Niedermeier, M., Hennessy, B.T., Knight, Z.A., *et al.* Isoform-selective phosphoinositide 3'-kinase inhibitors inhibit CXCR4 signaling and overcome stromal cell-mediated drug resistance in chronic lymphocytic leukemia: A novel therapeutic approach. *Blood* **113**(22), 5549-5557 (2009).
5. Sujobert, P., Bardet, V., Cornillet-Lefebvre, P., *et al.* Essential role for the p110δ isoform in phosphoinositide 3-kinase activation and cell proliferation in acute myeloid leukemia. *Blood* **106**(3), 1063-1066 (2005).
6. Lee, K.S., Lee, H.K., Hayflick, J.S., *et al.* Inhibition of phosphoinositide 3-kinase δ attenuates allergic airway inflammation and hyperresponsiveness in murine asthma model. *FASEB J.* **20**(3), 455-465 (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, 10/11/2022

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