

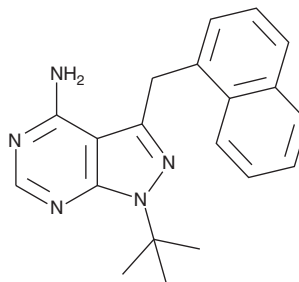
PRODUCT INFORMATION



1-NM-PP1

Item No. 13330

CAS Registry No.: 221244-14-0
Formal Name: 1-(1,1-dimethylethyl)-3-(1-naphthalenylmethyl)-1H-pyrazolo[3,4-d]pyrimidin-4-amine
Synonym: PP1 Analog II
MF: C₂₀H₂₁N₅
FW: 331.4
Purity: ≥95%
UV/Vis.: λ_{max}: 224, 282 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

1-NM-PP1 is supplied as a crystalline solid. A stock solution may be made by dissolving the 1-NM-PP1 in the solvent of choice, which should be purged with an inert gas. 1-NM-PP1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 1-NM-PP1 in these solvents is approximately 2, 20, and 30 mg/ml, respectively.

1-NM-PP1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1-NM-PP1 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 1-NM-PP1 has a solubility of approximately 0.15 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

1-NM-PP1 is a cell permeable inhibitor of kinases that have been mutated, by a single base substitution, to become 'analog sensitive' (as), as compared to the wild-type kinase. 1-NM-PP1 was first developed to optimally inhibit v-Src-as1, with an I338G substitution, preferentially over v-Src (IC₅₀ = 4.2 nM versus 28 μM, respectively).¹ The homologous mutation in other kinases generated similar analog sensitivity (e.g., IC₅₀ = 3.2 nM for c-Fyn-as1 versus 1.0 μM for c-Fyn; 5.0 nM for Cdk2-as1 versus 29 μM for Cdk2; 8.0 nM for CAMKII-as1 versus 24 μM for CAMKII).² This approach has been used to elucidate functions of several kinases in both mammalian and yeast systems.²⁻⁶

References

1. Bishop, A.C., Kung, C.y., Shah, K., et al. *J. Am. Chem. Soc.* **121**(4), 627-631 (1999).
2. Bishop, A.C., Ubersax, J.A., Petsch, D.T., et al. *Nature* **407**(6802), 395-401 (2000).
3. Weiss, E.L., Bishop, A.C., Shokat, K.M., et al. *Nat. Cell Biol.* **2**(10), 677-685 (2000).
4. Ira, G., Pelliccioli, A., Balijja, A., et al. *Nature* **431**(7011), 1011-1017 (2004).
5. Larochelle, S., Merrick, K.A., Terret, M.E., et al. *Mol. Cell* **25**(6), 839-850 (2007).
6. Van den Broeke, C., Radu, M., Deruelle, M., et al. *Proc. Natl. Acad. Sci. USA* **106**(21), 8707-8712 (2009).

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