

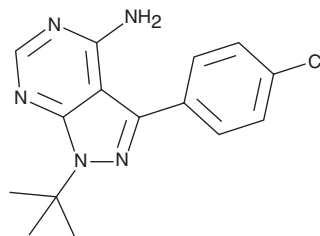
PRODUCT INFORMATION



PP2

Item No. 13198

CAS Registry No.: 172889-27-9
Formal Name: 3-(4-chlorophenyl)-1-(1,1-dimethylethyl)-1H-pyrazolo[3,4-d]pyrimidin-4-amine
Synonym: AGL 1879
MF: C₁₅H₁₆ClN₅
FW: 301.8
Purity: ≥95%
UV/Vis.: λ_{max}: 256, 286 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

PP2 is supplied as a crystalline solid. A stock solution may be made by dissolving the PP2 in the solvent of choice, which should be purged with an inert gas. PP2 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PP2 in these solvents is approximately 0.15, 1.4, and 3 mg/ml, respectively.

PP2 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PP2 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. PP2 has a solubility of approximately 0.1 mg/ml in a 1:9 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 Src family of non-receptor tyrosine kinases regulate cell adhesion, growth, and differentiation through activation of multiple intracellular signaling pathways. Normally inactive, Src kinases are transiently activated during mitosis and constitutively activated by abnormal mutations. PP2 is a potent, reversible, ATP-competitive, and selective inhibitor of the Src family of protein tyrosine kinases. It inhibits p56^{lck} (IC₅₀ = 4 nM), p59^{fynT} (IC₅₀ = 5 nM), Hck (IC₅₀ = 5 nM), and Src (IC₅₀ = 100 nM).¹ PP2 does not significantly affect the activity of EGFR kinase (IC₅₀ = 480 nM), JAK2 (IC₅₀ > 50 μM), or ZAP-70 (IC₅₀ > 100 μM).¹ PP2 inhibits the activation of focal adhesion kinase as well as its phosphorylation at Tyr⁵⁷⁷. PP2 also inhibits anti-CD3-stimulated tyrosine phosphorylation of human T-cells with an IC₅₀ value of 600 nM.¹

Reference

1. Hanke, J.H., Gardner, J.P., Dow, R.L., *et al.* Discovery of a novel, potent, and Src family-selective tyrosine kinase inhibitor. Study of Lck- and FynT-dependent T cell activation. *J. Biol. Chem.* **271**(2), 695-701 (1996).

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, 02/16/2023

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