

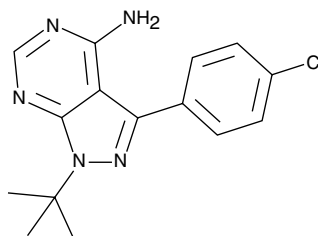
# Product Information



## 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:** C<sub>15</sub>H<sub>16</sub>ClN<sub>5</sub>  
**FW:** 301.8  
**Purity:** ≥95%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that PP2 be stored as supplied at -20°C. It should be stable for at least two years.

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

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<sup>lck</sup> (IC<sub>50</sub> = 4 nM), p59<sup>fyn</sup>T (IC<sub>50</sub> = 5 nM), Hck (IC<sub>50</sub> = 5 nM), and Src (IC<sub>50</sub> = 100 nM).<sup>1</sup> PP2 does not significantly affect the activity of EGFR kinase (IC<sub>50</sub> = 480 nM), JAK2 (IC<sub>50</sub> > 50 μM), or ZAP-70 (IC<sub>50</sub> > 100 μM).<sup>1</sup> PP2 inhibits the activation of focal adhesion kinase as well as its phosphorylation at Tyr<sup>577</sup>. PP2 also inhibits anti-CD3-stimulated tyrosine phosphorylation of human T-cells with an IC<sub>50</sub> value of 600 nM.<sup>1</sup>

### 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).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/13198](http://www.caymanchem.com/catalog/13198)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 02/25/2015

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)