

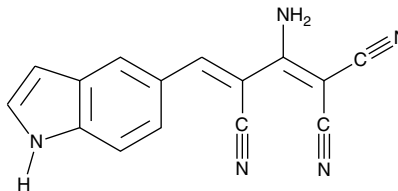
# Product Information



## AG-370

Item No. 10010568

**CAS Registry No.:** 134036-53-6  
**Formal Name:** 3-amino-4-(1H-indol-5-ylmethylene)-2-pentenenitrile  
**Synonym:** NSC 651712  
**MF:** C<sub>15</sub>H<sub>9</sub>N<sub>5</sub>  
**FW:** 259.3  
**Purity:** ≥95% (*cis/trans* mixture)  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid  
**UV/Vis.:** λ<sub>max</sub>: 260, 367 nm



### Laboratory Procedures

For long term storage, we suggest that AG-370 be stored as supplied at -20°C. It should be stable for at least two years. AG-370 is supplied as a crystalline solid. A stock solution may be made by dissolving the AG-370 in an organic solvent purged with an inert gas. AG-370 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AG-370 in these solvents is approximately 30 mg/ml.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. For maximum solubility in aqueous buffers, AG-370 should be dissolved in DMSO and diluted with the aqueous buffer of choice. AG-370 has a solubility of approximately 0.1 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.

Protein tyrosine kinase (PTK) inhibitors are potential antiproliferative agents for diseases caused by the hyperactivity of PTKs. Tyrphostins are a class of antiproliferative compounds which selectively inhibit PTKs of key growth factors such as epidermal growth factor (EGF) or platelet-derived growth factor (PDGF) by blocking the phosphorylation of specific tyrosine residues.<sup>1</sup> AG-370 is a selective inhibitor of PDGF receptor kinase with an IC<sub>50</sub> value of 20 μM in human bone marrow fibroblasts. It displays comparatively weak inhibition of the EGF receptor (IC<sub>50</sub> = 820 μM).<sup>2-4</sup>

### References

1. Gazit, A., Yaish, P., Gilon, C., *et al.* Tyrphostins I: Synthesis and biological activity of protein tyrosine kinase inhibitors. *J. Med. Chem.* **32**(10), 2344-2352 (1989).
2. Bryckaert, M.C., Eldor, A., Fontenay, M., *et al.* Inhibition of platelet-derived growth factor-induced mitogenesis and tyrosine kinase activity in cultured bone marrow fibroblasts by tyrphostins. *Exp. Cell Res.* **199**, 255-261 (1992).
3. Erneux, C. and Takazawa, K. Tyrphostins as molecular tools and potential antiproliferative drugs. *Trends Pharmacol. Sci.* **12**, 171-174 (1991).
4. Gazit, A., App, H., McMahon, G., *et al.* Tyrphostins. 5. Potent inhibitors of platelet-derived growth factor receptor tyrosine kinase: Structure-activity relationships in quinoxalines, quinolines, and indole tyrphostins. *J. Med. Chem.* **39**, 2170-2177 (1996).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10010568](http://www.caymanchem.com/catalog/10010568)

**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, 04/16/2015

### 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

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