

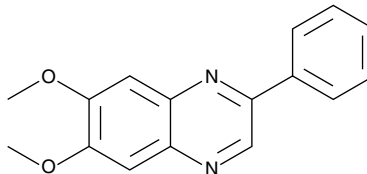
# Product Information



## AG-1296

Item No. 10010592

**CAS Registry No.:** 146535-11-7  
**Formal Name:** 6,7-dimethoxy-2-phenyl-quinoxaline  
**Synonym:** Tyrphostin AG-1296  
**MF:** C<sub>16</sub>H<sub>14</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 266.3  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid  
**UV/Vis.:** λ<sub>max</sub>: 233, 264, 367 nm



### Laboratory Procedures

For long term storage, we suggest that AG-1296 be stored as supplied at -20°C. It should be stable for at least two years.

AG-1296 is supplied as a crystalline solid. A stock solution may be made by dissolving the AG-1296 in an organic solvent purged with an inert gas. AG-1296 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of AG-1296 in these solvents is approximately 3 mg/ml.

AG-1296 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AG-1296 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. AG-1296 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF: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 act as PTK blockers. PTK inhibitors specific for platelet-derived growth factor (PDGF) receptor kinase could help in the treatment of atherosclerosis, restenosis, pulmonary fibrosis, and gliomas.<sup>1</sup> AG-1296 is a potent and selective inhibitor of PDGF receptor kinase with an IC<sub>50</sub> value of about 0.4 μM both *in vitro* and in cells (Swiss 3T3 cells).<sup>2</sup> It inhibits ligand-stimulated DNA synthesis in PDGF receptors and stem cell factor (SCF)/kit receptor transfected cells with an IC<sub>50</sub> values of 1.5 and 1.8 μM, respectively.<sup>2</sup> Treatment of sis oncogene transfected NIH3T3 cells with AG-1296 reverses the transformed phenotype.<sup>2</sup>

### References

- Levitzki, A. and Gazit, A. Tyrosine kinase inhibition: An approach to drug development. *Science* **267**(5205), 1782-1788 (1995).
- Kovalenko, M., Gazit, A., Böhmer, A., *et al.* Selective platelet-derived growth factor receptor kinase blockers reverse sis-transformation. *Cancer Res.* **54**, 6106-6114 (1994).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10010592](http://www.caymanchem.com/catalog/10010592)

### 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)

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

#### MATERIAL 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 Material 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, 01/06/2012