

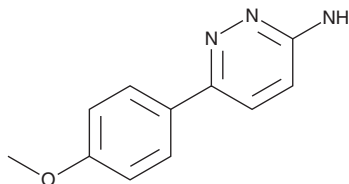
# PRODUCT INFORMATION



## 6-(4-Methoxyphenyl)-3-pyridazinamine

Item No. 17588

**CAS Registry No.:** 4776-87-8  
**Formal Name:** 6-(4-methoxyphenyl)-3-pyridazinamine  
**MF:** C<sub>11</sub>H<sub>11</sub>N<sub>3</sub>O  
**FW:** 201.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 274 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

6-(4-Methoxyphenyl)-3-pyridazinamine is supplied as a crystalline solid. A stock solution may be made by dissolving the 6-(4-methoxyphenyl)-3-pyridazinamine in the solvent of choice, which should be purged with an inert gas. 6-(4-Methoxyphenyl)-3-pyridazinamine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 6-(4-methoxyphenyl)-3-pyridazinamine in ethanol and DMF is approximately 10 mg/ml and approximately 15 mg/ml in DMSO.

6-(4-Methoxyphenyl)-3-pyridazinamine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 6-(4-methoxyphenyl)-3-pyridazinamine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 6-(4-Methoxyphenyl)-3-pyridazinamine has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Ionotropic GABA<sub>A</sub> receptors are ligand-gated ion channels that facilitate the passing of chloride ions across the cell membrane and promote an inhibitory influence on target neurons. These receptors are the major targets for benzodiazepines and related anxiolytic drugs.<sup>1</sup> 6-(4-Methoxyphenyl)-3-pyridazinamine is an aminopyridazine derivative that acts as a GABA<sub>A</sub> receptor antagonist.<sup>2</sup> It can also be used as an intermediate in the synthesis of SR 95531 (Item No. 14585).<sup>2</sup>

### References

1. Karobath, M. and Sperk, G. Stimulation of benzodiazepine receptor binding by  $\gamma$ -aminobutyric acid. *Proc. Natl. Acad. Sci. USA* **76(2)**, 1004-1006 (1979).
2. Wermuth, C.-G., Bourguignon, J.-J., Schlewer, G., et al. Synthesis and structure-activity relationships of a series of aminopyridazine derivatives of  $\gamma$ -aminobutyric acid acting as selective GABA-A antagonists. *J. Med. Chem.* **30(2)**, 239-249 (1987).

#### 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, 02/14/2024

#### 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