

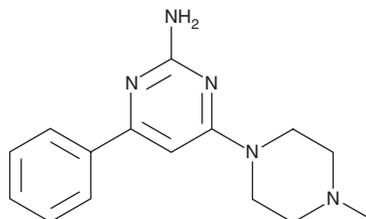
# PRODUCT INFORMATION



**VUF10460**

Item No. 14665

**CAS Registry No.:** 1028327-66-3  
**Formal Name:** 4-(4-methyl-1-piperazinyl)-6-phenyl-2-pyrimidinamine  
**MF:** C<sub>15</sub>H<sub>19</sub>N<sub>5</sub>  
**FW:** 269.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 245, 316 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

VUF10460 is supplied as a crystalline solid. A stock solution may be made by dissolving the VUF10460 in the solvent of choice, which should be purged with an inert gas. VUF10460 is soluble in the organic solvent DMSO at a concentration of approximately 100 mg/ml.

## Description

VUF10460 is a histamine H<sub>4</sub> receptor agonist (K<sub>i</sub> = 5.01 for the human receptor).<sup>1,2</sup> It is selective for the rat histamine H<sub>4</sub> receptor over the rat histamine H<sub>3</sub> receptor (K<sub>i</sub>s = 34.67 and 1,778.28 nM, respectively).<sup>2</sup> VUF10460 (10 mg/kg) potentiates the formation of HCl-induced ulcers in the rat gastric mucosa, an effect that is not blocked by the histamine H<sub>4</sub> receptor antagonist JNJ-7777120 (Item No. 10011925).

## References

1. Cowart, M.D., Altenbach, R.J., Liu, H., *et al.* Rotationally constrained 2,4-diamino-5,6-disubstituted pyrimidines: A new class of histamine H<sub>4</sub> receptor antagonists with improved druglikeness and in vivo efficacy in pain and inflammation models. *J. Med. Chem.* **51(20)**, 6547-6557 (2008).
2. Coruzzi, G., Adami, M., Pozzoli, C., *et al.* Selective histamine H<sub>3</sub> and H<sub>4</sub> receptor agonists exert opposite effects against the gastric lesions induced by HCl in the rat stomach *Eur J. Pharmacol.* **669(1-3)**, 121-127 (2011).

### 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, 12/13/2022

## 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](http://WWW.CAYMANCHEM.COM)