

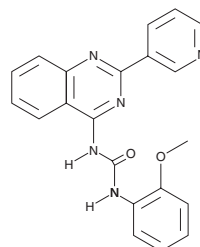
# PRODUCT INFORMATION



## VUF 5574

Item No. 27676

**CAS Registry No.:** 280570-45-8  
**Formal Name:** N-(2-methoxyphenyl)-N'-[2-(3-pyridinyl)-4-quinazoliny]-urea  
**MF:** C<sub>21</sub>H<sub>17</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 371.4  
**Purity:** ≥98%  
**Supplied as:** A 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

VUF 5574 is supplied as a solid. A stock solution may be made by dissolving the VUF 5574 in the solvent of choice, which should be purged with an inert gas. VUF 5574 is soluble in the organic solvent DMSO at a concentration of approximately 4 mM.

### Description

VUF 5574 is an antagonist of the adenosine A<sub>3</sub> receptor (K<sub>i</sub> = 4.03 nM for the recombinant human receptor).<sup>1</sup> It is greater than 2,500-fold selective for adenosine A<sub>3</sub> over A<sub>1</sub> and A<sub>2A</sub> receptors. VUF 5574 (100 nM) decreases phosphorylation of ERK1/2 induced by adenosine (Item No. 21232) in isolated porcine coronary artery smooth muscle cells.<sup>2</sup> It increases oxygen-glucose deprivation-induced reductions in the amplitude of field excitatory postsynaptic potentials (EPSPs) in a rat hippocampal slice model of ischemia when used at a concentration of 100 nM.<sup>3</sup> VUF 5574 (2 µg, intracisternal) reduces sodium nitroprusside-induced heart rate increases in rats.<sup>4</sup>

### References

1. van Muijwijk-Koezen, J.E., Timmerman, H., van der Goot, H., *et al.* Isoquinoline and quinazoline urea analogues as antagonists for the human adenosine A<sub>3</sub> receptor. *J. Med. Chem.* **43(11)**, 2227-2238 (2000).
2. Shen, J., Halenda, S.P., Sturek, M., *et al.* Cell-signaling evidence for adenosine stimulation of coronary smooth muscle proliferation via the A<sub>1</sub> adenosine receptor. *Circ. Res.* **97(6)**, 574-582 (2005).
3. Pugliese, A.M., Coppi, E., Spalluto, G., *et al.* A<sub>3</sub> adenosine receptor antagonists delay irreversible synaptic failure caused by oxygen and glucose deprivation in the rat CA1 hippocampus *in vitro*. *Br. J. Pharmacol.* **147(5)**, 524-532 (2006).
4. El-Mas, M.M., El-Gowilly, S.M., Fouda, M.A., *et al.* Role of adenosine A<sub>2A</sub> receptor signaling in the nicotine-evoked attenuation of reflex cardiac sympathetic control. *Toxicol. Appl. Pharmacol.* **254(3)**, 229-237 (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, 11/01/2022

#### 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