

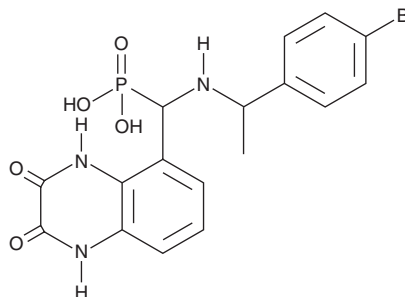
# PRODUCT INFORMATION



## NVP-AAM077

Item No. 30622

**CAS Registry No.:** 2514919-70-9  
**Formal Name:** (((1-(4-bromophenyl)ethyl)amino)(2,3-dioxo-1,2,3,4-tetrahydroquinoxalin-5-yl)methyl)phosphonic acid  
**MF:**  $C_{17}H_{17}BrN_3O_5P$   
**FW:** 454.2  
**Purity:**  $\geq 95\%$  (mixture of isomers)  
**UV/Vis.:**  $\lambda_{max}$ : 216 nm  
**Supplied as:** A crystalline solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

NVP-AAM077 is supplied as a crystalline solid. A stock solution may be made by dissolving the NVP-AAM077 in water. We do not recommend storing the aqueous solution for more than one day.

### Description

NVP-AAM077 is an NMDA receptor antagonist ( $IC_{50} = 0.008 \mu M$ ).<sup>1</sup> It is selective for NMDA receptors containing NR1A/NR2A subunits over NMDA receptors containing NR1A/NR2B subunits ( $IC_{50}s = 0.27$  and  $29.6 \mu M$ , respectively, in *X. laevis* oocytes expressing receptors containing the respective subunits). NVP-AAM077 inhibits seizures induced by maximal electroshock (MES) in mice with an  $ED_{50}$  value of 23 mg/kg. It increases the activity of neuronal nitric oxide synthase (nNOS) in primary hippocampal neurons.<sup>2</sup> NVP-AAM077 (10 mg/kg) inhibits proliferation of neural progenitor cells (NPCs) in the subventricular zone (SVZ) and in the subgranular zone (SGZ) of the hippocampal dentate gyrus of adult mice, an effect not seen in *nNOS*<sup>-/-</sup> mice. It also increases the escape latency and decreases the percentage of time spent in the target quadrant in the Morris water maze.

### References

1. Auberson, Y.P., Allgeier, H., Bischoff, S., *et al.* 5-Phosphonomethylquinoxalinediones as competitive NMDA receptor antagonists with a preference for the human 1A/2A, rather than 1A/2B receptor composition. *Bioorg. Med. Chem. Lett.* **12**(7), 1099-1102 (2002).
2. Hu, M., Sun, Y.-J., Zhou, Q.-G., *et al.* Reduced spatial learning in mice treated with NVP-AAM077 through down-regulating neurogenesis. *Eur. J. Pharmacol.* **622**(1-3), 37-44 (2009).

#### 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, 12/14/2022

#### 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