

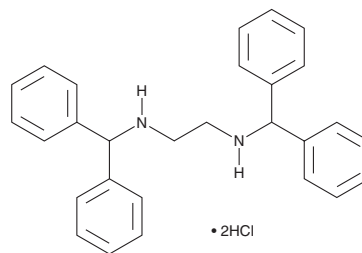
# PRODUCT INFORMATION



## AMN082

Item No. 28779

**CAS Registry No.:** 97075-46-2  
**Formal Name:** N,N'-bis(diphenylmethyl)-1,2-ethanediamine, dihydrochloride  
**MF:** C<sub>28</sub>H<sub>28</sub>N<sub>2</sub> • 2HCl  
**FW:** 465.5  
**Purity:** ≥98%  
**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

AMN082 is supplied as a crystalline solid. A stock solution may be made by dissolving the AMN082 in the solvent of choice, which should be purged with an inert gas. AMN082 is soluble in organic solvents such as ethanol and DMSO. The solubility of AMN082 in these solvents is approximately 1 and 100 mM, respectively. It is also soluble in water. The solubility of AMN082 in water is approximately 5 mM. We do not recommend storing the aqueous solution for more than one day.

### Description

AMN082 is an orally bioavailable allosteric agonist of metabotropic glutamate receptor 7 (mGluR7).<sup>1</sup> It inhibits cAMP accumulation induced by forskolin in CHO cells expressing human mGluR7b (EC<sub>50</sub> = 64 nM). AMN082 (10 μM) is selective for mGluR7a and mGluR7b over mGluR1b, mGluR4, and mGluR8a with 140, 90, 15, 18, and 20% activation, respectively, as well as over mGluR2, mGluR3, mGluR5a, mGluR6, GluR3, and NMDA receptors containing NR1a, NR2A, or NR2B subunits. AMN082 also binds to the norepinephrine transporter (NET) and α<sub>1</sub>-adrenergic receptor (α<sub>1</sub>-AR).<sup>2</sup> It increases the proliferation of neural progenitor cells (NPCs) and induces their differentiation into neurons when used at a concentration of 1 μM.<sup>3</sup> AMN082 (2.5 mg/kg) inhibits apomorphine-induced circling in a rat model of Parkinson's disease induced by 6-hydroxy dopamine (6-OHDA; Item No. 25330).<sup>4</sup>

### References

1. Mitsukawa, K., Yamamoto, R., Ofner, S., *et al.* A selective metabotropic glutamate receptor 7 agonist: Activation of receptor signaling via an allosteric site modulates stress parameters *in vivo*. *Proc. Natl. Acad. Sci. U.S.A.* **102(51)**, 18712-18717 (2005).
2. Sukoff Rizzo, S.J., Leonard, S.K., Gilbert, A., *et al.* The metabotropic glutamate receptor 7 allosteric modulator AMN082: A monoaminergic agent in disguise? *J. Pharmacol. Exp. Ther.* **338(1)**, 345-352 (2011).
3. Tian, Y., Liu, Y., Chen, X., *et al.* AMN082 promotes the proliferation and differentiation of neural progenitor cells with influence on phosphorylation of MAPK signaling pathways. *Neurochem. Int.* **57(1)**, 8-15 (2010).
4. Greco, B., Lopez, S., van der Putten, H., *et al.* Metabotropic glutamate 7 receptor subtype modulates motor symptoms in rodent models of Parkinson's disease. *J. Pharmacol. Exp. Ther.* **332(3)**, 1064-1071 (2010).

#### 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, 11/30/2022

#### 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