

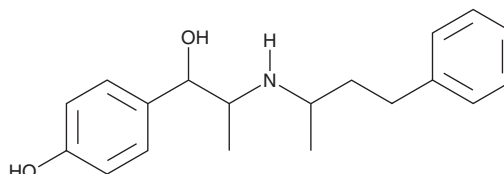
# PRODUCT INFORMATION



## Nylidrin

Item No. 33232

**CAS Registry No.:** 447-41-6  
**Formal Name:** 4-hydroxy- $\alpha$ -[1-[(1-methyl-3-phenylpropyl)amino]ethyl]-benzenemethanol  
**Synonym:** Buphenine  
**MF:** C<sub>19</sub>H<sub>25</sub>NO<sub>2</sub>  
**FW:** 299.4  
**Purity:**  $\geq$ 85%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 226 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

Nylidrin is supplied as a crystalline solid. A stock solution may be made by dissolving the nylidrin in the solvent of choice, which should be purged with an inert gas. Nylidrin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of nylidrin in these solvents is approximately 30 mg/ml.

Nylidrin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, nylidrin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Nylidrin has a solubility of approximately 0.12 mg/ml in a 1:7 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Nylidrin is an agonist of  $\beta$ -adrenergic receptors and antagonist of NR1A/2B subunit-containing NMDA receptors.<sup>1,2</sup> It binds to  $\beta$ -adrenergic receptors ( $K_d = 0.37 \mu\text{M}$ ) and activates adenylate cyclase with a  $K_a$  value of  $1.3 \mu\text{M}$ .<sup>1</sup> Nylidrin is selective for NR1A/2B over NR1A/2A and NR1A/2C subunit-containing NMDA receptors ( $\text{IC}_{50}\text{s} = 0.18, 32, \text{ and } 42 \mu\text{M}$ , respectively, for the recombinant receptors expressed in *Xenopus* oocytes).<sup>2</sup> It inhibits NMDA-induced currents in primary rat cortical neurons ( $\text{IC}_{50} = 0.22 \mu\text{M}$ ). It also decreases blood pressure and increases heart rate in spontaneously hypertensive rats (SHRs) with a minimum effective dose (MED) of 0.5 mg/kg.<sup>3</sup> Nylidrin (20 and 100  $\mu\text{M}$ ) inhibits influenza hemagglutinin 2-mediated membrane fusion in Vero E6 cells.<sup>4</sup> It inhibits infection of MDCK cells by H1N1 and H3N2 influenza isolates *in vitro* ( $\text{EC}_{50}\text{s} = 7.2$  and  $12.1 \mu\text{M}$ , respectively) and prevents infection in a mouse model of mouse-adapted H1N1 infection.

### References

1. Bilezikian, J.P., Dornfeld, A.M., and Gammon, D.E. Structure-binding-activity analysis of beta-adrenergic amines—I. Binding to the beta receptor and activation of adenylate cyclase. *Biochem. Pharmacol.* **27**(10), 1445-1454 (1978).
2. Whittemore, E.R., Ilyin, V.I., Konkoy, C.S., *et al.* Subtype-selective antagonism of NMDA receptors by nylidrin. *Eur. J. Pharmacol.* **337**(2-3), 197-208 (1997).
3. Yen, T.T. and Pearson, D.V. Nylidrin: A potent antihypertensive agent in hypertensive rats. *Res. Commun. Chem. Pathol. Pharmacol.* **23**(1), 11-28 (1979).
4. Jang, Y., Shin, J.S., Lee, J.-Y., *et al.* In vitro and in vivo antiviral activity of nylidrin by targeting the hemagglutinin 2-mediated membrane fusion of influenza A virus. *Viruses* **12**(5), 581 (2020).

#### 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, 12/08/2022

#### 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