

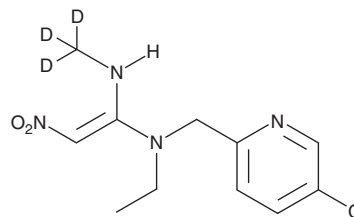
# PRODUCT INFORMATION



## Nitenpyram-d<sub>3</sub>

Item No. 26783

**Formal Name:** (1E)-N-[(6-chloro-3-pyridinyl)methyl]-N-ethyl-N'-methyl-d<sub>3</sub>-2-nitro-1,1-ethenediamine  
**MF:** C<sub>11</sub>H<sub>12</sub>ClD<sub>3</sub>N<sub>4</sub>O<sub>2</sub>  
**FW:** 273.7  
**Chemical Purity:** ≥98% (Nitenpyram)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>3</sub>); ≤1% d<sub>0</sub>  
**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

Nitenpyram-d<sub>3</sub> is intended for use as an internal standard for the quantification of nitenpyram (Item No. 21271) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Nitenpyram-d<sub>3</sub> is supplied as a solid. A stock solution may be made by dissolving the nitenpyram-d<sub>3</sub> in the solvent of choice, which should be purged with an inert gas. Nitenpyram-d<sub>3</sub> is soluble in organic solvents such as methanol and DMSO.

### Description

Nitenpyram is a neonicotinoid insecticide.<sup>1</sup> It is an antagonist of housefly neuronal-type nicotinic acetylcholine receptors (nAChRs; IC<sub>50</sub> = 2 nM for housefly head homogenates).<sup>2</sup> Nitenpyram induces mortality in brown planthopper (*N. lugens*) larvae when used at a concentration of 0.5 ppm.<sup>3</sup> Orally administered nitenpyram (11.4, 11.4, and 57 mg/animal, respectively) eliminates live adult fleas on cats, small dogs, and large dogs after six hours of treatment.<sup>4</sup> It also reduces the number of aphids, ladybugs, ladybug larvae, lacewings, and syrphid larvae on adult cotton plants when applied to cotton seeds at a concentration of 4 g/kg of seeds.<sup>5</sup> Formulations containing nitenpyram have been used in the control of fleas and ticks on companion animals.

### References

1. Vo, D.T., Hsu, W.H., Abu-Basha, E.A., *et al.* Insect nicotinic acetylcholine receptor agonists as flea adulticides in small animals. *J. Vet. Pharmacol. Ther.* **33**(4), 315-322 (2010).
2. Nauen, R., Ebbinghaus, U., and Tietjen, K. Ligands of the nicotinic acetylcholine receptor as insecticides. *Pest. Manag. Sci.* **55**(5), 608-610 (1999).
3. Minamida, I., Iwanaga, K., Tabuchi, T., *et al.* Synthesis and insecticidal activity of acyclic nitroethene compounds containing a heteroarylmethylamino group. *J. Pesticide Sci.* **18**, 41-48 (1993).
4. Dobson, P., Tinembart, O., Fisch, R.D., *et al.* Efficacy of nitenpyram as a systemic flea adulticide in dogs and cats. *Vet. Rec.* **147**(25), 709-713 (2000).
5. Zhang, Z., Zhang, X., Wang, Y., *et al.* Nitenpyram, dinotefuran, and thiamethoxam used as seed treatments act as efficient controls against *Aphis gossypii* via high residues in cotton leaves. *J. Agric. Food Chem.* **64**(49), 9276-9285 (2016).

#### 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, 08/04/2023

#### 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)