

# PRODUCT INFORMATION



## Deltamethrin-d<sub>5</sub>

Item No. 33710

**Formal Name:** (1R,3R)-3-(2,2-dibromoethenyl)-2,2-dimethyl-cyclopropanecarboxylic acid, cyano (3-phenoxy-d<sub>5</sub>)phenyl methyl ester

**Synonym:** Decamethrin-d<sub>5</sub>

**MF:** C<sub>22</sub>H<sub>14</sub>Br<sub>2</sub>D<sub>5</sub>NO<sub>3</sub>

**FW:** 510.2

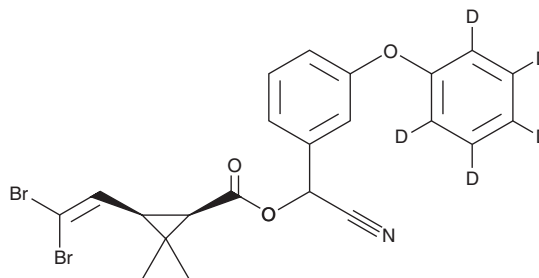
**Chemical Purity:** ≥95% (Deltamethrin; mixture of diastereomers)

**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>5</sub>); ≤1% d<sub>0</sub>

**Supplied as:** A semi-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

Deltamethrin-d<sub>5</sub> is intended for use as an internal standard for the quantification of deltamethrin (Item No. 24172) 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).

Deltamethrin-d<sub>5</sub> is supplied as a semi-solid. A stock solution may be made by dissolving the deltamethrin-d<sub>5</sub> in the solvent of choice, which should be purged with an inert gas. Deltamethrin-d<sub>5</sub> is soluble in acetonitrile, methanol, DMSO, and dimethyl formamide.

### Description

Deltamethrin is a type II pyrethroid insecticide and a modulator of voltage-gated sodium channels (Na<sub>v</sub>).<sup>1</sup> It activates Na<sub>v</sub>1.8 channels and induces a tail current delay in *Xenopus* oocytes expressing the rat channel when used at a concentration of 10 μM. Deltamethrin induces mortality in *A. sinensis* (LC<sub>50</sub> = 4.522 ppm) and *T. urticae* (LC<sub>50</sub> = 0.06 g/L).<sup>2,3</sup> It decreases sperm count, motility, and viability and increases the percentage of morphologically abnormal sperm in rats when administered at a dose of 5 mg/kg.<sup>4</sup> Formulations containing deltamethrin have been used as domestic and agricultural insecticides.

### References

1. Choi, J.S. and Soderlund, D.M. Structure-activity relationships for the action of 11 pyrethroid insecticides on rat Na<sub>v</sub>1.8 sodium channels expressed in *Xenopus* oocytes. *Toxicol. Appl. Pharmacol.* **211**(3), 233-244 (2006).
2. Chang, K.-S., Yoo, D.-H., Shin, E.-H., *et al.* Susceptibility and resistance of field populations of *Anopheles sinensis* (Diptera: Culicidae) collected from Paju to 13 insecticides. *Osong Public Health Res. Perspect.* **4**(2), 76-80 (2013).
3. Penman, D.R., Chapman, R.B., and Bowie, M.H. Direct toxicity and repellent activity of pyrethroids against *Tetranychus urticae* (Acari: Tetranychidae). *J. Econ. Entomol.* **79**(5), 1183-1187 (1986).
4. Abdallah, F.B., Slima, A.B., Dammak, I., *et al.* Comparative effects of dimethoate and deltamethrin on reproductive system in male mice. *Andrologia* **42**(3), 182-186 (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, 03/22/2023

#### 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