

PRODUCT INFORMATION



Bifenthrin-d₅ Item No. 39588

Formal Name: (1R,3R)-rel-3-[(1Z)-2-chloro-3,3,3-trifluoro-1-propen-1-yl]-2,2-dimethylcyclopropanecarboxylic acid, (2-methyl[1,1'-biphenyl-2',3',4',5',6'-d₅]-3-yl)methyl ester

Synonym: (±)-Bifenthrin-d₅

MF: C₂₃H₁₇ClD₅F₃O₂

FW: 427.9

Chemical Purity: ≥95% (Bifenthrin)

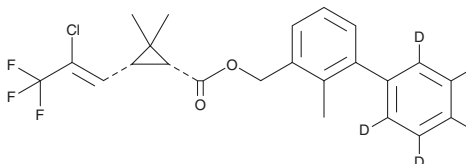
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀

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

Bifenthrin-d₅ is intended for use as an internal standard for the quantification of bifenthrin (Item No. 24057) 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).

Bifenthrin-d₅ is supplied as a solid. A stock solution may be made by dissolving the bifenthrin-d₅ in the solvent of choice, which should be purged with an inert gas. Bifenthrin-d₅ is slightly soluble in chloroform and DMSO.

Description

Bifenthrin is a synthetic pyrethroid insecticide and a slow activator of voltage-gated sodium channel 1.8 (Na_v1.8).¹ It is toxic to *A. gambiae* and *C. quinquefasciatus* mosquitos (LD₅₀s = 0.15 and 0.16 ng/mg, respectively) and reduces the number of insidious flower bugs (*O. insidiosus*) in corn when applied at a concentration of 168 g/ha.^{2,3} Bifenthrin is toxic to *D. magna* (LC₅₀ = 12.4 µg/L) and inhibits *D. magna* reproduction when used at a concentration of 1 µg/L.⁴ It induces fine- and whole-body tremors and decreases motor activity in rats when administered at a dose of 20 mg/kg.⁵ Bifenthrin (8 mg/kg) decreases body mass and locomotor activity, increases memory deficits in the passive avoidance test, as well as decreases the concentration of blood erythrocytes, increases lymphocyte hepatic infiltration, and increases the activity of hepatic alanine transaminase (ALT) in mice.⁶ Formulations containing bifenthrin have been used as insecticides in residential and agricultural settings.

References

1. Choi, J.S. and Soderlund, D.M. *Toxicol. Appl. Pharmacol.* **211**(3), 233-244 (2006).
2. Hougard, J.-M., Duchon, S., Zaim, M., et al. *J. Med. Entomol.* **39**(3), 526-533 (2002).
3. Al-Deeb, M.A., Wilde, G.E., and Zhu, K.Y. *J. Econ. Entomol.* **94**(6), 1353-1360 (2001).
4. Ye, W.H., Wen, Y.Z., Liu, W.P., et al. *J. Environ. Sci. (China)* **16**(5), 843-846 (2004).
5. Wolansky, M.J., McDaniel, K.L., Moser, V.C., et al. *Neurotoxicol. Teratol.* **29**(3), 377-384 (2007).
6. Nieradko-Iwanicka, B., Borzecki, A., and Jodlowska-Jedrych, B. *J. Physiol. Pharmacol.* **66**(1), 129-137 (2015).

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, 09/22/2023

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