

# PRODUCT INFORMATION

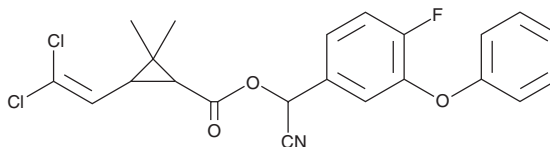


## Cyfluthrin

Item No. 24231

**CAS Registry No.:** 68359-37-5  
**Formal Name:** 3-(2,2-dichloroethenyl)-2,2-dimethyl-cyclopropanecarboxylic acid, cyano(4-fluoro-3-phenoxyphenyl)methyl ester

**Synonym:**  $\beta$ -Cyfluthrin  
**MF:** C<sub>22</sub>H<sub>18</sub>Cl<sub>2</sub>FNO<sub>3</sub>  
**FW:** 434.3  
**Purity:**  $\geq$ 98%  
**Supplied as:** A 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

Cyfluthrin is supplied as a solid. A stock solution may be made by dissolving the cyfluthrin in the solvent of choice, which should be purged with an inert gas. Cyfluthrin is soluble in methanol (heated) and is slightly soluble in chloroform.

### Description

Cyfluthrin is a pyrethroid insecticide and a modulator of voltage-gated sodium channels (Na<sub>v</sub>).<sup>1,2</sup> It slowly activates rat recombinant Na<sub>v</sub>1.8 channels, delays inactivation by longer than 40 ms, and induces persistent tail currents in channels expressed in *X. laevis* oocytes.<sup>1</sup> It also decreases the mean burst rate in rat primary neurons (IC<sub>50</sub> = 0.99  $\mu$ M, respectively).<sup>2</sup> Cyfluthrin is toxic to various insects, including *A. melinus*, *G. ashmeadi*, *E. eremicus*, and *E. formosa* (LC<sub>50</sub>s = 7, 67, 96, and 63 ng/ml, respectively) and the *A. sinensis* mosquito (LC<sub>50</sub> = 0.446 ppm).<sup>3,4</sup> It is also toxic to *A. mellifera* honeybees (LD<sub>50</sub> = 0.22 ppm), affecting locomotor activity and wing fanning behavior with an increase in the mean bout duration of time spent upside down, indicating disruption of the righting reflex, and a decrease in wing fanning behavior when administered at a dose of 10 ng/bee.<sup>5</sup> Formulations containing cyfluthrin have been used for the control of insects in agriculture and for non-commercial purposes.

### 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. Baskar, M.K. and Murthy, P.B. *In vitro* evaluation of pyrethroid-mediated changes on neuronal burst parameters using microelectrode arrays. *Neurotoxicology* **57**, 270-281 (2016).
3. Prabhaker, N., Morse, J.G., Castle, S.J., *et al.* Toxicity of seven foliar insecticides to four insect parasitoids attacking citrus and cotton pests. *J. Econ. Entomol.* **100(4)**, 1053-1061 (2007).
4. 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).
5. Oliver, C.J., Softley, S., Williamson, S.M., *et al.* Pyrethroids and nectar toxins have subtle effects on the motor function, grooming and wing fanning behaviour of honeybees (*Apis mellifera*). *PLoS One* **10(8):e0133733**, (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, 11/11/2022

#### 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