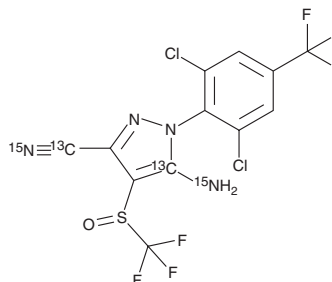


# PRODUCT INFORMATION



**Fipronil-<sup>13</sup>C<sub>2</sub>, <sup>15</sup>N<sub>2</sub>**  
Item No. 39587

**Formal Name:** 5-(amino-<sup>15</sup>N)-1-(2,6-dichloro-4-(trifluoromethyl)phenyl)-4-((trifluoromethyl)sulfinyl)-1H-pyrazole-5-<sup>13</sup>C-3-carbonitrile-<sup>13</sup>C-<sup>15</sup>N  
**MF:** C<sub>10</sub>[<sup>13</sup>C]<sub>2</sub>H<sub>4</sub>Cl<sub>2</sub>F<sub>6</sub>N<sub>2</sub>[<sup>15</sup>N]<sub>2</sub>OS  
**FW:** 441.1  
**Purity:** ≥95%  
**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

Fipronil-<sup>13</sup>C<sub>2</sub>, <sup>15</sup>N<sub>2</sub> is supplied as a solid. A stock solution may be made by dissolving the fipronil-<sup>13</sup>C<sub>2</sub>, <sup>15</sup>N<sub>2</sub> in the solvent of choice, which should be purged with an inert gas. Fipronil-<sup>13</sup>C<sub>2</sub>, <sup>15</sup>N<sub>2</sub> is slightly soluble in methanol and sparingly soluble in DMSO.

## Description

Fipronil-<sup>13</sup>C<sub>2</sub>, <sup>15</sup>N<sub>2</sub> is intended for use as an internal standard for the quantification of fipronil (Item No. 24128) by GC- or LC-MS. Fipronil is an insecticide that acts as a selective antagonist of insect GABA receptors (IC<sub>50</sub>s = 28 and 1,660 nM for cockroach and rat receptors, respectively).<sup>1</sup> It also inhibits desensitizing and non-desensitizing glutamate-induced chloride currents in cockroach neurons (IC<sub>50</sub>s = 800 and 10 nM, respectively). Fipronil induces activity of the cytochrome P450 (CYP) isoforms CYP1A1/2, CYP2B1/2, and CYP3A1/2 in isolated rat liver microsomes.<sup>2</sup> Formulations containing fipronil have been used as insecticides in veterinary care.

## References

1. Narahashi, T., Zhao, X., Ikeda, T., *et al.* Glutamate-activated chloride channels: Unique fipronil targets present in insects but not in mammals. *Pestic. Biochem. Physiol.* **97**(2), 149-152 (2010).
2. Caballero, M.V., Ares, I., Martínez, M., *et al.* Fipronil induces CYP isoforms in rats. *Food Chem. Toxicol.* **83**, 215-221 (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, 09/06/2023

## 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