

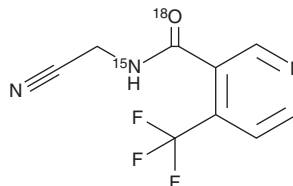
PRODUCT INFORMATION



Flonicamid-¹⁵N,¹⁸O

Item No. 39578

CAS Registry No.: 2483830-60-8
Formal Name: N-(cyanomethyl)-4-(trifluoromethyl)-3-Pyridinecarboxamide-¹⁵N-¹⁸O
MF: C₉H₆F₃N₂[¹⁵N][¹⁸O]
FW: 232.2
Purity: ≥98%
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Flonicamid-¹⁵N,¹⁸O is supplied as a solution in methanol. A stock solution may be made by dissolving the flonicamid-¹⁵N,¹⁸O in the solvent of choice, which should be purged with an inert gas. Flonicamid-¹⁵N,¹⁸O is soluble in methanol.

Description

Flonicamid-¹⁵N,¹⁸O is intended for use as an internal standard for the quantification of flonicamid (Item No. 25608) by GC- or LC-MS. Flonicamid is a pyridinecarboxamide insecticide.¹ It inhibits the brown planthopper (*N. lugens*) inward-rectifying potassium channel K_{ir}1 (IC₅₀ = 0.64 μM). Flonicamid (100 μg/ml) inhibits aphid (*M. persicae*) salivation and sap ingestion during feeding on Japanese radish leaves.² It induces toxicity in the aphid species *A. gossypii*, *R. padi*, *S. graminum*, and *L. erysimi* (LC₅₀s = 0.64, 0.74, 0.78, and 2.01 μg/ml, respectively) and rats (LD₅₀ = ≥844 mg/kg).^{2,3} Formulations containing flonicamid have been used for the control of pest insects in agriculture.

References

1. Ren, M., Niu, J., Hu, B., *et al.* Block of Kir channels by flonicamid disrupts salivary and renal excretion of insect pests. *Insect Biochem. Mol. Biol.* **99**, 17-26 (2018).
2. Morita, M., Ueda, T., Yoneda, T., *et al.* Flonicamid, a novel insecticide with a rapid inhibitory effect on aphid feeding. *Pest. Manag. Sci.* **63**(10), 969-973 (2007).
3. Morita, M., Yoneda, T., and Akiyoshi, N. Research and development of a novel insecticide, flonicamid. *J. Pestic. Sci.* **39**(3), 179-180 (2014).

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, 05/08/2026

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