

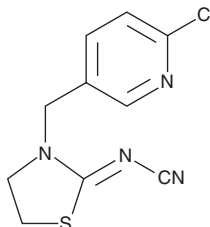
PRODUCT INFORMATION



Thiacloprid

Item No. 25824

CAS Registry No.: 111988-49-9
Formal Name: [N(Z)]-N-[3-[(6-chloro-3-pyridinyl)methyl]-2-thiazolidinylidene]-cyanamide
Synonym: BAY YRC 2894
MF: C₁₀H₉CIN₄S
FW: 252.7
Purity: ≥98%
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

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

Description

Thiacloprid is a neonicotinoid insecticide.¹ It is an antagonist of nicotinic acetylcholine receptors (nAChRs; IC₅₀ = 2.7 nM for the *D. melanogaster* enzyme). Thiacloprid is lethal to *D. melanogaster* (LD₅₀ = 0.03 mg/kg). It eliminates aphids on apple tree leaves in field studies when applied at a concentration of 0.0072% AI/L.² Thiacloprid is toxic to mice (LD₅₀ = 28 mg/kg).¹

References

1. Tomizawa, M., Lee, D.L., and Casida, J.E. Neonicotinoid insecticides: Molecular features conferring selectivity for insect versus mammalian nicotinic receptors. *J. Agric. Food Chem.* **48**(12), 6016-6024 (2000).
2. Elbert, A., Erdelen, C., Kuhnhold, J., et al. Thiacloprid, a novel neonicotinoid insecticide for foliar application. *Proc. Brighton Crop. Prot. Conf. - Pests and Diseases* **1**, 21-26 (2000).

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, 09/13/2023

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