

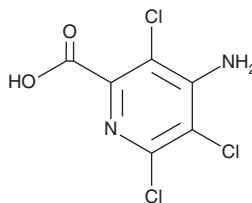
PRODUCT INFORMATION



Picloram

Item No. 45895

CAS Registry No.: 1918-02-1
Formal Name: 4-amino-3,5,6-trichloro-2-pyridinecarboxylic acid
Synonym: NSC 233899
MF: $C_6H_3Cl_3N_2O_2$
FW: 241.5
Purity: $\geq 95\%$
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

Picloram is supplied as a solid. A stock solution may be made by dissolving the picloram in the solvent of choice, which should be purged with an inert gas. Picloram is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Picloram is an herbicide.¹⁻³ It increases membrane lipid peroxidation, reduces catalase (CAT), peroxidase (POD), and superoxide dismutase (SOD) activity, and induces mitochondrial membrane and chloroplast injury in *E. adenophorum* when applied at 120-240 g/hectare.² Picloram inhibits root growth in *A. thaliana* ($IC_{50} = 5.8 \mu\text{M}$).¹ Residual picloram has been detected in soil more than two years after application.³ Formulations containing picloram have been used for weed control in pastures and agricultural areas.

References

1. Shi, R.-C., Liu, Q., Wei, W., *et al.* Study on design, synthesis, and herbicidal activity of 4-amino-3,5-dichloro-6-indolyl-2-picolinic acids. *J. Agric. Food Chem.* **74**(22), 17551-17561 (2026).
2. Liu, X., Qi, C., Wang, Z., *et al.* Effect of picloram herbicide on physiological responses of *Eupatorium adenophorum* Spreng. *Chil. J. Agric. Res.* **74**(4), 438-444 (2014).
3. Braz, G.B.P., Souza, M.F., Almeida, D.P., *et al.* Residue of picloram in new agricultural areas: Impacts, detection, and mitigation measures. *ACS Omega* **11**(19), 27745-27750 (2026).

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, 06/30/2026

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