

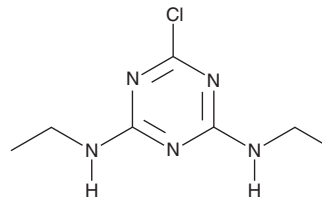
PRODUCT INFORMATION



Simazine

Item No. 36491

CAS Registry No.: 122-34-9
Formal Name: 6-chloro-N²,N⁴-diethyl-1,3,5-triazine-2,4-diamine
Synonym: NSC 25999
MF: C₇H₁₂ClN₅
FW: 201.7
Purity: ≥95%
UV/Vis.: λ_{max}: 222 nm
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

Simazine is supplied as a solid. A stock solution may be made by dissolving the simazine in the solvent of choice, which should be purged with an inert gas. Simazine is slightly soluble in DMSO and dimethyl formamide.

Description

Simazine is a triazine herbicide.¹⁻³ Pre-emergent application of simazine (1 kg/ha) controls annual and perennial monocot and dicot weeds and increases crop yield in sorghum crops.¹ It also controls weed emergence in kharif maize crops when applied at a concentration of 2 kg/ha.² Simazine (200 and 400 mg/kg) induces spleen cell apoptosis and reduces the proliferation of B and T cells in mice.³ Formulations containing simazine have been used in the control of broadleaf weeds and annual grasses in agriculture.

References

1. Patro, G.K., Tosh, G.C., and Nayak, B.C. Weed control in sorghum through cultural and chemical methods. *Indian J. Agric. Sci.* **42(12)**, 1128-1131 (1972).
2. Jadhav, S.N. and Khuspe, V.S. Relative efficacy of varying levels and times of application of simazine and atrazine on the control of weeds and yield of kharif maize (*Zea mays* Linn). *J. Maharashtra Agric. Univ.* **6(2)**, 169-171 (1981).
3. Ren, R., Sun, D.-J., Yan, H., *et al.* Oral exposure to the herbicide simazine induces mouse spleen immunotoxicity and immune cell apoptosis. *Toxicol. Pathol.* **41(1)**, 63-72 (2013).

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, 11/30/2022

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