

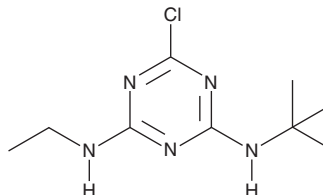
PRODUCT INFORMATION



Terbuthylazine

Item No. 34192

CAS Registry No.: 5915-41-3
Formal Name: 6-chloro-N²-(1,1-dimethylethyl)-N⁴-ethyl-1,3,5-triazine-2,4-diamine
Synonyms: GS-13529, Terbutazine
MF: C₉H₁₆ClN₅
FW: 229.7
Purity: ≥98%
UV/Vis.: λ_{max}: 223 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

Terbuthylazine is supplied as a solid. A stock solution may be made by dissolving the terbuthylazine in the solvent of choice, which should be purged with an inert gas. Terbuthylazine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of terbuthylazine in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

Terbuthylazine is a triazine herbicide.¹ Terbuthylazine is active against a variety of broadleaf and grass weeds, including *A. retroflexus*, *D. stramonium*, *S. nigrum*, and *S. halepense*, when used in combination with (S)-metolachlor.² It has been found in surface groundwater and soil samples.^{3,4} Formulations containing terbuthylazine have been used for weed control in agriculture and as algicides in water coolant systems.

References

1. Štěpánová, S., Plhalová, L., Doleželová, P., et al. The effects of subchronic exposure to terbuthylazine on early developmental stages of common carp. *Sci. World J.* 615920 (2012).
2. Meseldžija, M. and Dudić, M. Terbuthylazine application with herbicides of different mode of action in maize crop. *IX International Scientific Agriculture Symposium* (2018).
3. Guzzella, L., Pozzoni, F., and Giuliano, G. Herbicide contamination of surficial groundwater in Northern Italy. *Environ. Pollut.* **142**(2), 344-353 (2006).
4. Sahid, I.B. and Teoh, S.S. Persistence of terbuthylazine in soils. *Bull. Environ. Contam. Toxicol.* **52**(2), 226-230 (1994).

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, 11/18/2022

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