

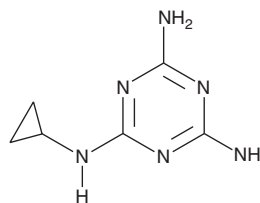
# PRODUCT INFORMATION



## Cyromazine

Item No. 24041

**CAS Registry No.:** 66215-27-8  
**Formal Name:** N<sup>2</sup>-cyclopropyl-1,3,5-triazine-2,4,6-triamine  
**MF:** C<sub>6</sub>H<sub>10</sub>N<sub>6</sub>  
**FW:** 166.2  
**Purity:** ≥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

Cyromazine is supplied as a solid. A stock solution may be made by dissolving the cyromazine in the solvent of choice, which should be purged with an inert gas. Cyromazine is soluble in organic solvents such as methanol and DMSO.

### Description

Cyromazine is a triazine insecticide.<sup>1</sup> It completely inhibits larval development of the Australian sheep blow fly (*L. cuprina*), as measured by adult emergence, when used at a concentration of 0.5 ppm in rearing medium. Cyromazine (1-20 µg/ml) in drinking water reduces egg production by adult blow flies and subsequent larval development. Application of cyromazine (0.025-0.1%) to chicken manure or administration as a chicken feed additive at concentrations of 1.5 and 5 ppm reduces proliferation of the house fly (*M. domestica*).<sup>2</sup> Formulations containing cyromazine have been used to control fly infestations and biting in livestock rearing.

### References

1. Friedel, T. and McDonell, P.A. Cyromazine inhibits reproduction and larval development of the Australian sheep blow fly (diptera: Calliphoridae). *J. Econ. Entomol.* **78(4)**, 868-873 (1985).
2. Williams, R.E. and Berry, J.G. Evaluation of CGA 72662 as a topical spray and feed additive for controlling house flies breeding in chicken manure. *Poult. Sci.* **59(10)**, 2207-2212 (1980).

#### 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, 10/19/2022

#### 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