

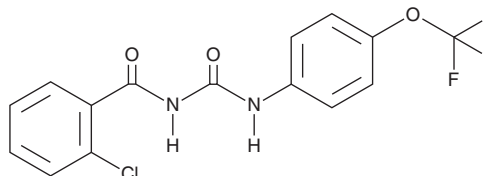
# PRODUCT INFORMATION



## Triflumuron

Item No. 30858

**CAS Registry No.:** 64628-44-0  
**Formal Name:** 2-chloro-N-[[[4-(trifluoromethoxy)phenyl]amino]carbonyl]-benzamide  
**Synonym:** OMS-2015  
**MF:** C<sub>15</sub>H<sub>10</sub>ClF<sub>3</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 358.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 250 nm  
**Supplied as:** A crystalline 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

Triflumuron is supplied as a crystalline solid. A stock solution may be made by dissolving the triflumuron in the solvent of choice, which should be purged with an inert gas. Triflumuron is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of triflumuron in these solvents is approximately 10 and 30 mg/ml respectively.

### Description

Triflumuron is an insecticide and a chitin synthesis inhibitor.<sup>1</sup> It reduces cuticle thickness in *T. molitor* pupae when administered at doses of 1 and 2 µg/insect.<sup>2</sup> Topical application of triflumuron (1 µg/insect) to female *M. domestica* adults reduces egg hatching by greater than 95%.<sup>1</sup> It induces mortality in *A. aegypti*, *C. quinquefasciatus*, and *A. stephensi* third instar larvae with EC<sub>50</sub> values of 0.2, 0.3, and 0.1 ppb, respectively.<sup>3</sup>

### References

- Howard, J. and Wall, R. The effects of triflumuron, a chitin synthesis inhibitor, on the housefly, *Musca domestica* (Diptera: Muscidae). *B. Entomol. Res.* **85(1)**, 71-77 (1995).
- Soltani, N., Soltani-Mazouni, N., and Delachambre, J. Evaluation of triflumuron, a benzoylphenylurea derivative, on *Tenebrio molitor* pupae (Col., Tenebrionidae): Effects on cuticle. *J. Appl. Entomol.* **120(1-5)**, 627-629 (1996).
- Batra, C.P., Mittal, P.K., Adak, T., et al. Efficacy of IGR compound starycide 480 SC (triflumuron) against mosquito larvae in clear and polluted water. *J. Vector Borne Dis.* **42(3)**, 109-116 (2005).

#### 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, 07/25/2023

#### 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](http://WWW.CAYMANCHEM.COM)