

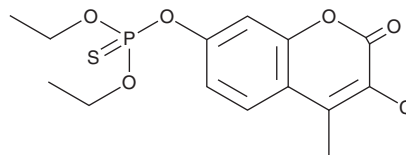
# PRODUCT INFORMATION



## Coumaphos

Item No. 24230

**CAS Registry No.:** 56-72-4  
**Formal Name:** O-(3-chloro-4-methyl-2-oxo-2H-1-benzopyran-7-yl), phosphorothioic acid, O,O-diethyl ester  
**Synonym:** NSC 8944  
**MF:** C<sub>14</sub>H<sub>16</sub>ClO<sub>5</sub>PS  
**FW:** 362.8  
**Purity:** ≥98%  
**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

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

### Description

Coumaphos is an organophosphate pesticide.<sup>1</sup> It is converted into the corresponding oxon metabolite *in vivo*, similar to other organophosphate pesticides, that inhibits acetylcholinesterase (AChE). It is active against adult, but not arrested stage, *O. ostertagi* helminths.<sup>2</sup> Coumaphos is toxic to *A. stephensi* and *A. aegypti* mosquitoes when applied topically, with median lethal doses of 0.002 and 0.012 µg per female mosquito, respectively, but not when used as a contact insecticide.<sup>3</sup> It is lethal to rats (LD<sub>50</sub>s = 41 and 16 mg/kg for male and female rats, respectively).<sup>4</sup> Formulations containing coumaphos have been used to control pests in livestock.

### References

1. Guo, J.-X., Wu, J.J.-Q., Wight, J.B., *et al.* Mechanistic insight into acetylcholinesterase inhibition and acute toxicity of organophosphorus compounds: A molecular modeling study. *Chem. Res. Toxicol.* **19(2)**, 209-216 (2006).
2. Prichard, R.K. Anthelmintics and control. *Vet. Parasitol.* **27(1-2)**, 97-109 (1988).
3. Hadaway, A.B. and Barlow, F. The toxicity of some organophosphorus compounds to adult *Anopheles stephensi*. *Bull. World Health Organ.* **28(1)**, 55-61 (1963).
4. Gaines, T.B. Acute toxicity of pesticides. *Toxicol. Appl. Pharmacol.* **14(3)**, 515-534 (1969).

#### 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, 09/07/2023

#### 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