

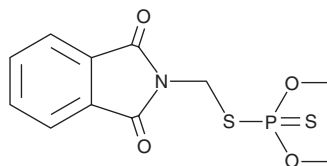
PRODUCT INFORMATION



Phosmet

Item No. 25812

CAS Registry No.: 732-11-6
Formal Name: S-[(1,3-dihydro-1,3-dioxo-2H-isoindol-2-yl)methyl] phosphorodithioic acid, O,O-dimethyl ester
Synonym: Phthalophos
MF: C₁₁H₁₂NO₄PS₂
FW: 317.3
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

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

Description

Phosmet is an organophosphate insecticide and acaricide.¹⁻³ It reduces apple damage by a large variety of insects, including apple maggots, codling moths, and obliquebanded leafrollers when used as either a border or cover spray at a concentration of 1.9 kg AI/hectare.² Phosmet is effective in controlling *S. scabiei* in pigs when applied as a 20% pour-on solution.³ It is toxic to rats *via* oral administration (LC₅₀ = 230 mg/kg).⁴ Formulations containing phosmet have been used in the control of insects and mites in agriculture.

References

1. Abdollahi, M. and Karami-Mohajeri, S. A comprehensive review on experimental and clinical findings in intermediate syndrome caused by organophosphate poisoning. *Toxicol. Appl. Pharmacol.* **258(3)**, 309-314 (2012).
2. Trimble, R.M. and Vickers, P.M. Evaluation of border sprays for managing the codling moth (Tortricidae: Lepidoptera) and the apple maggot (Tephritidae: Diptera) in Ontario apple orchards. *J. Econ. Entomol.* **93(3)**, 777-787 (2000).
3. Hewett, G.R. Phosmet for the systemic control of pig mange in growing pigs. *Vet. Parasitol.* **18(3)**, 265-268 (1985).
4. European Food Safety Authority. Conclusion on the peer review of the pesticide risk assessment of the active substance phosmet. *EFSA J.* **9(5)**, 2162 (2011).

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, 11/22/2022

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