

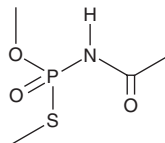
# PRODUCT INFORMATION



## Acephate

Item No. 24038

**CAS Registry No.:** 30560-19-1  
**Formal Name:** N-acetyl-phosphoramidothioic acid, O,S-dimethyl ester  
**Synonym:** RE 12420  
**MF:** C<sub>4</sub>H<sub>10</sub>NO<sub>3</sub>PS  
**FW:** 183.2  
**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

Acephate is supplied as a solid. A stock solution may be made by dissolving the acephate in the solvent of choice. Acephate is soluble in organic solvents such as chloroform and DMSO, which should be purged with an inert gas. Acephate is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Acephate is an organophosphate insecticide that is active against a variety of insects.<sup>1</sup> It inhibits acetylcholinesterase in the house fly with a topical 24-hour LD<sub>50</sub> value of 1.8 µg/g but has millimolar IC<sub>50</sub> values for vertebrate cholinesterases.<sup>2,3</sup> In rats, acephate (140 mg/kg) induces reversible hyperglycemia after two and six hours and increases plasma corticosterone, liver glucose-6-phosphate, and liver tyrosine aminotransferase.<sup>1</sup> It also increases liver glycogen levels by 3.5-fold after six hours. Formulations containing acephate have been used in the control of insects in agriculture and horticulture.

### References

1. Joshi, A.K.R., and Rajini, P.S. Reversible hyperglycemia in rats following acute exposure to acephate, an organophosphorus insecticide: Role of gluconeogenesis. *Toxicology* **257(1-2)**, 40-45 (2009).
2. Rojakovick, A.S., and March, R.B. *In vitro* and *in vivo* inhibition of house fly acetylcholinesterase by the N-acetyl phosphoramidothioate orthene. *J. Econ. Entomol.* **65(5)**, 1505-1507 (1972).
3. Hussain, M.A., Mohamad, R.B., and Oloffs, P.C. Studies on the toxicity, metabolism, and anticholinesterase properties of acephate and methamidophos. *J. Environ. Sci. Health B.* **20(1)**, 129-147 (1985).

#### 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, 10/04/2022

#### 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