

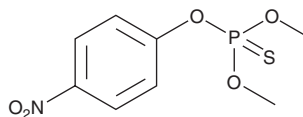
PRODUCT INFORMATION



Methyl Parathion

Item No. 25786

CAS Registry No.: 298-00-0
Formal Name: phosphorothioic acid, O,O-dimethyl
O-(4-nitrophenyl) ester
Synonym: BAY 11405
MF: C₈H₁₀NO₅PS
FW: 263.2
Purity: ≥98%
UV/Vis.: λ_{max}: 272 nm
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

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

Description

Methyl parathion is an organophosphate insecticide.¹ Methyl parathion is lethal to lab strains and field isolates of tobacco budworm larvae (LD₅₀s = 7 and 81.8-128.3 μg/g, respectively).² It reduces the number of stink bugs (*C. sayi*) caught per 100 net sweeps when applied to alfalfa fields at a concentration of 0.4 pounds per acre.³ Methyl parathion increases sister chromatid exchange (SCE) in a concentration-dependent manner and induces cell cycle arrest at the M₁ phase in V79 cells at a concentration of 40 μg/ml.⁴ It is toxic to rats (LD₅₀ = 14 mg/kg).¹

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. Martinez-Carrillo, J.L. and Reynolds, H.T. Dosage-mortality studies with pyrethroids and other insecticides on the tobacco budworm (*Lepidoptera: Noctuidae*) from the Imperial Valley, California. *J. Econ. Entomol.* **76(5)**, 983-986 (1983).
3. Reynolds, H.T., Stern, V.M., Fukoto, T.R., *et al.* Potential use of dylox and other insecticides in a control program for field crop pests in California. *J. Econ. Entomol.* **53(1)**, 72-78 (1960).
4. Chen, H.H., Hsueh, J.L., Sirianni, S.R., *et al.* Induction of sister-chromatid exchanges and cell cycle delay in cultured mammalian cells treated with eight organophosphorus pesticides. *Mutat. Res.* **88(3)**, 307-316 (1981).

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, 07/12/2023

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