

PRODUCT INFORMATION



Pyribencarb

Item No. 42362

CAS Registry No.: 799247-52-2

Formal Name: N-[[2-chloro-5-[(1E)-1-[[[6-methyl-2-pyridinyl)methoxy]imino]ethyl]phenyl)methyl]-carbamic acid, methyl ester

Synonyms: KIF 7767, KUF-1204

MF: $C_{18}H_{20}ClN_3O_3$

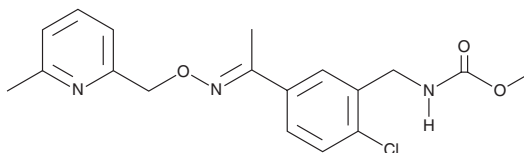
FW: 361.8

Purity: $\geq 95\%$

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

Pyribencarb is supplied as a solid. A stock solution may be made by dissolving the pyribencarb in the solvent of choice, which should be purged with an inert gas. Pyribencarb is slightly soluble (0.1-1 mg/ml) in ethanol and DMSO.

Description

Pyribencarb is a carbamate fungicide.¹ It inhibits the production of aflatoxins in *A. parasiticus* ($IC_{50} = 0.43 \mu\text{M}$).² It inhibits wild-type *B. cinerea* and strobilurin-resistant *B. cinerea* mitochondrial complex II, also known as succinate-cytochrome c reductase (IC_{50} s = 0.079 and $5.1 \mu\text{M}$, respectively) and wild-type *B. cinerea* mitochondrial complex III, also known as decylubiquinol-cytochrome c reductase ($IC_{50} = 0.023 \mu\text{M}$).³ Pyribencarb ($3 \mu\text{g/ml}$) prevents and reduces *B. cinerea* lesion development on cucumber seedlings and inhibits *B. cinerea* mycelial growth ($EC_{50} = 0.006 \mu\text{g/ml}$).¹

References

1. Takagaki, M., Kataoka, S., Kida, K., et al. Disease-controlling effect of a novel fungicide pyribencarb against *Botrytis cinerea*. *J. Pestic. Sci.* **35**(1), 10-14 (2010).
2. Sakuda, S., Prabowo, D.F., Takagi, K., et al. Inhibitory effects of respiration inhibitors on aflatoxin production. *Toxins (Basel)* **6**(4), 1193-1200 (2014).
3. Kataoka, S., Takagaki, M., Kaku, K., et al. Mechanism of action and selectivity of a novel fungicide, pyribencarb. *J. Pestic. Sci.* **35**(2), 99-106 (2010).

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, 12/06/2024

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