

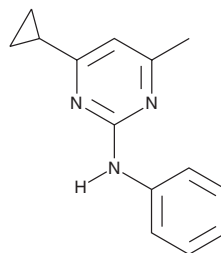
PRODUCT INFORMATION



Cyprodinil

Item No. 24233

CAS Registry No.: 121552-61-2
Formal Name: 4-cyclopropyl-6-methyl-N-phenyl-2-pyrimidinamine
MF: C₁₄H₁₅N₃
FW: 225.3
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

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

Description

Cyprodinil is an anilino-pyrimidine broad-spectrum fungicide that inhibits the biosynthesis of methionine in phytopathogenic fungi.¹ It inhibits mycelial cell growth of *B. cinerea*, *P. herpotrichoides*, and *H. oryzae* on amino acid-free media (IC₅₀s = 0.44, 4.8, and 0.03 μM, respectively), an effect that is reversed by addition of methionine or homocysteine. In an MDA-kb2 assay, cyprodinil acts as an androgen receptor (AR) agonist (EC₂₀ = 1.91 μM) in the absence of the AR agonist DHT and inhibits the androgenic effect of DHT (IC₂₀ = 15.1 μM).² It is cytotoxic in a yeast antiandrogen screen (YAS; EC₂₀ = 27.8 μM) but not in an MDA-kb2 assay (EC₂₀ = >50 μM). Cyprodinil increases proliferation of estrogen receptor-expressing BG-1 ovarian cancer cells when used at low micromolar concentrations in combination with 17β-estradiol.³ It also increases tumor mass in a BG-1 ovariectomized mouse xenograft model after 70 days when administered at a dose of 3 mg/kg every three days. Formulations containing cyprodinil have been used in the control of fungi in agriculture.

References

1. Masner, P., Muster, P., and Schmid, J. Possible methionine biosynthesis inhibition by pyrimidinamine fungicides. *Pestic. Sci.* **42**(3), 163-166 (1994).
2. Orton, F., Rosivatz, E., Scholze, M., et al. Widely used pesticides with previously unknown endocrine activity revealed as *in vitro* antiandrogens. *Environ. Health Perspect.* **119**(6), 794-800 (2011).
3. Go, R.-E., Kim, C.-W., and Choi, K.-C. Effect of fenhexamid and cyprodinil on the expression of cell cycle- and metastasis-related genes via an estrogen receptor-dependent pathway in cellular and xenografted ovarian cancer models. *Toxicol. Appl. Pharmacol.* **289**(1), 48-57 (2015).

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, 11/15/2022

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