

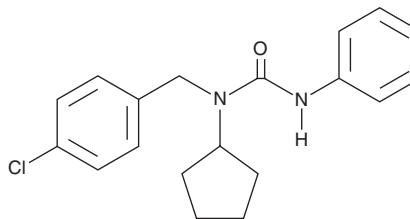
PRODUCT INFORMATION



Pencycuron

Item No. 36127

CAS Registry No.: 66063-05-6
Formal Name: N-[(4-chlorophenyl)methyl]-N-cyclopentyl-N'-phenyl-urea
MF: C₁₉H₂₁ClN₂O
FW: 328.8
Purity: ≥95%
UV/Vis.: λ_{max}: 241 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

Pencycuron is supplied as a solid. A stock solution may be made by dissolving the pencycuron in the solvent of choice, which should be purged with an inert gas. Pencycuron is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of pencycuron in these solvents is approximately 10 and 25 mg/ml, respectively.

Description

Pencycuron is a phenylurea fungicide.¹ It is active against the plant pathogenic fungi *G. graminis*, *R. oryzae*, and *R. solani*, but not *P. intermedium* or *P. irregulare*, when used at concentrations of 1 and 10 µg/ml. In plant studies, pencycuron (125, 250, and 500 ppm) reduces the incidence of *R. solani*-induced sheath blight on rice.²

References

1. Duffy, B. Combination of pencycuron and *Pseudomonas fluorescens* strain 2-79 for integrated control of rhizoctonia root rot and take-all of spring wheat. *Crop Prot.* **19(1)**, 21-25 (2000).
2. Yamada, Y., Saito, J., and Takase, I. Development of a new fungicide, pencycuron. *J. Pest. Sci.* **13(2)**, 375-387 (1988).

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, 07/14/2025

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