

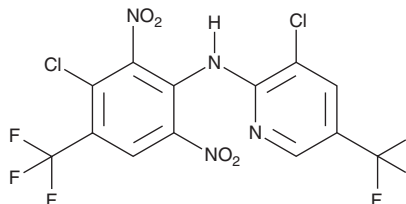
PRODUCT INFORMATION



Fluazinam

Item No. 36221

CAS Registry No.: 79622-59-6
Formal Name: 3-chloro-N-[3-chloro-2,6-dinitro-4-(trifluoromethyl)phenyl]-5-(trifluoromethyl)-2-pyridinamine
Synonyms: ASC 66825, ASC 67178, IKF 1216
MF: C₁₃H₄Cl₂F₆N₄O₄
FW: 465.1
Purity: ≥95%
UV/Vis.: λ_{max}: 238 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

Fluazinam is supplied as a solid. A stock solution may be made by dissolving the fluazinam in the solvent of choice, which should be purged with an inert gas. Fluazinam is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of fluazinam in ethanol is approximately 15 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of fluazinam can be prepared by directly dissolving the solid in aqueous buffers. The solubility of fluazinam in PBS (pH 7.2) is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Fluazinam is a fungicide.¹ It inhibits mycelial growth of 35 plant pathogenic fungi, including *S. sclerotiorum*, *P. infestans*, *C. rolfisii*, *B. cinerea*, and *Fusarium* (EC₅₀s = <0.1-42.1 ppm). Fluazinam (0.56 kg/ha) reduces the incidence of Sclerotinia blight in peanut fields.² It decreases the incidence of leaf drop in lettuce fields experimentally infected with *S. minor* or *S. sclerotiorum* when applied at a concentration of 561 g AI/ha. Formulations containing fluazinam have been used as fungicides in agriculture.

References

1. Komyoji, T., Sugimoto, K., Mitani, S., *et al.* Biological properties of a new fungicide, fluazinam. *J. Pesticide Sci.* **20**(2), 129-135 (1995).
2. Smith, F.D., Phipps, P.M., and Stipes, R.J. Fluazinam: A new fungicide for control of Sclerotinia blight and other soilborne pathogens of peanut. *Peanut Sci.* **19**(2), 115-120 (1992).

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, 03/27/2023

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