

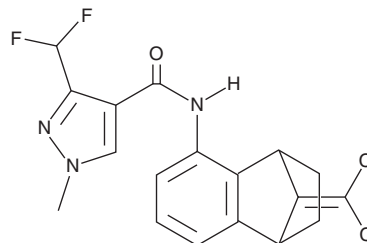
PRODUCT INFORMATION



Benzovindiflupyr

Item No. 34595

CAS Registry No.: 1072957-71-1
Formal Name: N-[9-(dichloromethylene)-1,2,3,4-tetrahydro-1,4-methanonaphthalen-5-yl]-3-(difluoromethyl)-1-methyl-1H-pyrazole-4-carboxamide
MF: C₁₈H₁₅Cl₂F₂N₃O
FW: 398.2
Purity: ≥90%
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

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

Description

Benzovindiflupyr is a fungicide.¹ It inhibits mitochondrial complex II, also known as succinate dehydrogenase (SDH; IC₅₀ = 5.2 nM), and mycelial growth of *S. sclerotiorum* (EC₅₀ = 0.011 µg/ml). Benzovindiflupyr (60 µg/ml) completely protects eggplant leaves from *S. sclerotiorum* infection. Formulations containing benzovindiflupyr have been used to control various fungal diseases in agriculture.

Reference

1. Gao, Y., He, L., Zhu, J., *et al.* The relationship between features enabling SDHI fungicide binding to the Sc-Sdh complex and its inhibitory activity against *Sclerotinia sclerotiorum*. *Pest Manag. Sci.* **76(8)**, 2799-2808 (2020).

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, 10/18/2022

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