

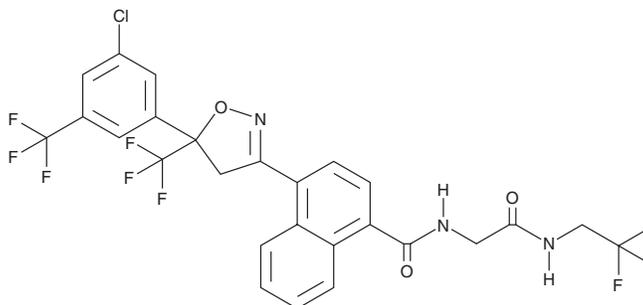
PRODUCT INFORMATION



Afoxolaner

Item No. 33317

CAS Registry No.: 1093861-60-9
Formal Name: 4-[5-[3-chloro-5-(trifluoromethyl)phenyl]-4,5-dihydro-5-(trifluoromethyl)-3-isoxazolyl]-N-[2-oxo-2-[(2,2,2-trifluoroethyl)amino]ethyl]-1-naphthalenecarboxamide
MF: C₂₆H₁₇ClF₉N₃O₃
FW: 625.9
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 221 nm
Supplied as: A crystalline 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

Afoxolaner is supplied as a crystalline solid. A stock solution may be made by dissolving the afoxolaner in the solvent of choice, which should be purged with an inert gas. Afoxolaner is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of afoxolaner in these solvents is approximately 30 mg/ml.

Description

Afoxolaner is an isoxazoline insecticide and acaricide.¹ It is a GABA receptor antagonist that inhibits GABA-induced currents in *P. americana* thoracic neurons (IC₅₀ = 10.8 nM). It induces mortality in 100% of *C. felis* after 24 and 48 hours when used at concentrations of 0.32 and 0.16 µg/ml in an artificial membrane blood feeding device. Afoxolaner (2.5 mg/kg) eradicates *C. felis* infestation in dogs for up to 42 days and *D. variabilis* infestation for up to 30 days.

Reference

1. Shoop, W.L., Hartline, E.J., Gould, B.R., *et al.* Discovery and mode of action of afoxolaner, a new isoxazoline parasiticide for dogs. *Vet. Parasitol.* **201(3-4)**, 179-189 (2014).

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, 07/25/2023

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