

PRODUCT INFORMATION

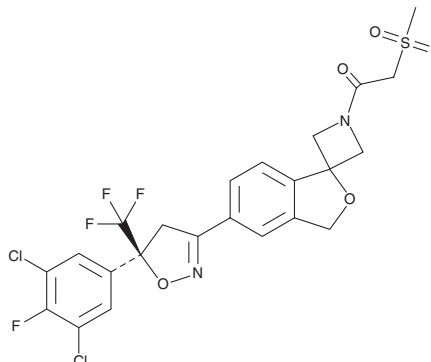


Sarolaner

Item No. 37321

CAS Registry No.: 1398609-39-6
Formal Name: 1-[5'-[(5S)-5-(3,5-dichloro-4-fluorophenyl)-4,5-dihydro-5-(trifluoromethyl)-3-isoxazolyl]spiro[azetidine-3,1'(3'H)-isobenzofuran]-1-yl]-2-(methylsulfonyl)-ethanone

Synonym: PF-6450567
MF: C₂₃H₁₈Cl₂F₄N₂O₅S
FW: 581.4
Purity: ≥98%
UV/Vis.: λ_{max}: 261 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

Sarolaner is supplied as a solid. A stock solution may be made by dissolving the sarolaner in the solvent of choice, which should be purged with an inert gas. Sarolaner is slightly soluble in ethanol, DMSO, and dimethyl formamide.

Description

Sarolaner is an isoxazoline ectoparasiticide.¹ It inhibits GABA-induced chloride currents in CHO-K1 cells expressing either the gene for susceptible *C. felis* flea resistance-to-dieldrin subunits containing alanine at residue 285 (CfRDL-A285; IC₅₀ = 135 nM) or the gene for CfRDL subunits containing the alanine-to-serine resistance mutation (CfRDL-S285; IC₅₀ = 136 nM). It induces mortality in *C. felis* fleas (LC₈₀ = 0.3 µg/ml) and *O. turicata* soft ticks (LC₁₀₀ = 0.003 µg/ml). Sarolaner (2.5 mg/kg) eliminates live *C. felis* fleas, as well as *R. sanguineus*, *I. ricinus*, and *D. reticulatus* ticks, on dogs after 48 hours and reduces the number of fleas and ticks for at least 35 days. Formulations containing sarolaner have been used in the treatment and control of flea and tick infestations in dogs.

Reference

1. McTier, T.L., Chubb, N., Curtis, M.P., *et al.* Discovery of sarolaner: A novel, orally administered, broad-spectrum, isoxazoline ectoparasiticide for dogs. *Vet. Parasitol.* **222**, 3-11 (2016).

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, 12/21/2022

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