PRODUCT INFORMATION



Fluazuron

Item No. 22283

CAS Registry No.:	86811-58-7	
Formal Name:	N-[[[4-chloro-3-[[3-chloro-5-	
	(trifluoromethyl)-2-pyridinyl]oxy]phenyl] amino]carbonyl]-2,6-difluoro-benzamide	
Synonym:	CGA 157419	
MF:	$C_{20}H_{10}CI_{2}F_{5}N_{3}O_{3}$	
FW:	506.2	
Purity:	≥98%	F ₃ C CI
UV/Vis.:	λ _{max} : 257 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

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

Fluazuron is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, fluazuron should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Fluazuron has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Fluazuron is an acaricide widely-used to control tick populations in cattle, sheep, and pigs.^{1,2} Fluazuron inhibits chitin formation limiting viable egg production and prevents immature ticks from moulting.¹ It is effective against populations of R. microplus, R. sanguineus, S. scabei, and Y. pestis at host doses ranging from 1.5 to 100 mg/kg.^{1,3-6} Fluazuron does not effect populations of F. candida or O. gazella, alleviating concerns about non-specific insecticide activity.^{2,7} Resistance to fluazuron in certain R. microplus strains may be emerging.8

References

- 1. Bull, M.S., Swindale, S., Overend, D., et al. Aust. Vet. J. 74(6), 468-470 (1996).
- 2. Zortéa, T., Segat, J.C., Maccari, A.P., et al. Chemosphere 173(1), 460-465 (2017).
- 3. Gomes, L.V.C., Lopes, W.D.Z., Cruz, B.C., et al. Exp. Parasitol. 153(1), 22-28 (2015).
- 4. de Oliveira, P.R., Calligaris, I.B., Roma, G.C., et al. Exp. Parasitol. 131(1), 35-39 (2012).
- 5. Pasay, C., Rothwell, J., Mounsey, K., et al. Parasit. Vectors. 5:40, (2012).
- 6. Davis, R.M., Cleugh, E., Smith, R.T., et al. J. Vector Ecol. 33(2), 278-284 (2008).
- 7. Kryger, U., Deschodt, C., Davis, A.L.V., et al. Vet. Parasitol. 143(3-4), 380-384 (2007).
- 8. Reck, J., Klafke, G.M., Webster, A., et al. Vet. Parasitol. 201(1-2), 128-163 (2014).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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, 11/14/2022

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