

# PRODUCT DATA SHEET



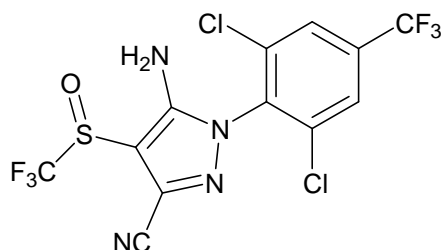
PRODUCT: Fipronil  
5-Amino-1-[2,6-dichloro-4-(trifluoromethyl)phenyl]-4-[(trifluoromethyl)sulfinyl]-1H-pyrazole-3-carbonitrile

CAS NO. 120068-37-3

CATALOG NO.: CM-102

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid (m.p. 200-201°C)

MOLECULAR FORMULA:  $C_{12}H_4Cl_2F_6N_4OS$  MOLECULAR WEIGHT: 437.2

PURITY: 98%(TLC: 100% Methylene chloride;  $R_f = 0.55$ )

SOLUBILITY: Soluble in DMSO (25 mg/ml) and ethanol (25 mg/ml)

STORAGE: Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.

APPLICATION NOTES: GABA gated chloride channel blocker. The product is not sterile.

REFERENCES: 1. A.M.Hosie *et al.* *Br.J.Pharmacol.* 1995 **115** 909

**The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.**