

PRODUCT DATA SHEET



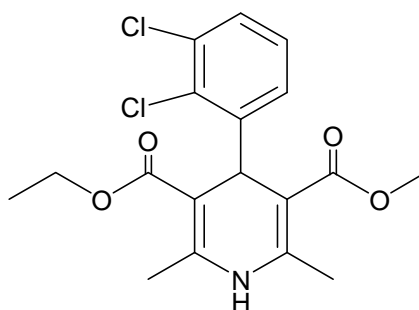
PRODUCT: Felodipine
Ethyl methyl 4-(2,3-dichlorophenyl)-1,4-dihydro-
2,6-dimethyl-3,5-pyridine dicarboxylate

CAS NO.: 72509-76-3

CATALOG NO.: CA-236

LOT NO.:

STRUCTURE:



PHYSICAL APPEARANCE: Pale yellow solid

MOLECULAR FORMULA: $C_{18}H_{19}Cl_2NO_4$

MOLECULAR WEIGHT: 384.3

PURITY: 98% (TLC: 2% Methanol/methylene chloride; $R_f = 0.30$)

SOLUBILITY: Soluble in: DMSO (>25 mg/mL); ethanol (20 mg/mL)

STORAGE: Store as supplied, at room temperature for up to 3 years. Store solutions at $-20^{\circ}C$ for up to 3 months.

APPLICATION NOTES: Felodipine is an L-type Ca^{2+} channel blocker¹. In isolated rat ventricular myocytes, using a whole-cell patch-clamp technique, it displays a K_i of 11 nM². It is a clinically useful antihypertensive agent³ which attenuates the hypertrophic effect of cyclosporine on transplanted hearts⁴.

REFERENCES: 1. T. Furakawa *et al.* *J. Cardiovasc. Pharmacol.* 2005 **45** 241
2. A. Zahradnikova *et al.* *J. Pharmacol. Exp. Ther.* 2007 **322** 638
3. B.A. Manzo *et al.* *Pharmacotherapy* 2003 **23** 1508
4. J Schwitter *et al.* *J. Heart Lung Transplant.* 1999 **18** 1003

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.