

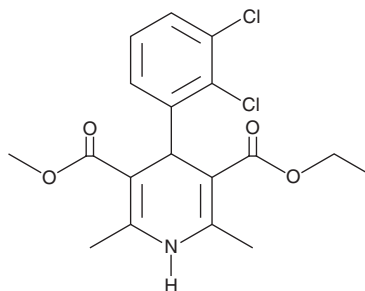
# PRODUCT INFORMATION



## Felodipine

Item No. 17535

**CAS Registry No.:** 72509-76-3  
**Formal Name:** 4-(2,3-dichlorophenyl)-1,4-dihydro-2,6-dimethyl-3,5-pyridinedicarboxylic acid, 3-ethyl 5-methyl ester  
**Synonym:** H 154/82  
**MF:** C<sub>18</sub>H<sub>19</sub>Cl<sub>2</sub>NO<sub>4</sub>  
**FW:** 384.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 236, 360 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

Felodipine is supplied as a crystalline solid. A stock solution may be made by dissolving the felodipine in the solvent of choice, which should be purged with an inert gas. Felodipine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of felodipine in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Felodipine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, felodipine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Felodipine has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Felodipine is an inhibitor of L-type calcium channels.<sup>1</sup> It induces relaxation of precontracted isolated porcine coronary artery segments (EC<sub>50</sub> = 0.15 nM), which highly express L-type calcium channels.<sup>2</sup> Felodipine is selective for L-type calcium channels over N-, R-, and P/Q-type channels at 10 μM, as well as the T-type Ca<sub>v</sub>3.2 channel (IC<sub>50</sub> = 6.8 μM).<sup>1,3</sup> Felodipine preferentially inhibits L-type calcium channels in isolated rat portal vein over rat left ventricle (IC<sub>50</sub>s = 33.9 and 3,981 nM, respectively).<sup>4</sup> It decreases mean arterial blood pressure and total peripheral resistance in a rabbit model of hypertension induced by renal artery ligation when administered intravenously at doses of 30 and 100 nmol/kg.<sup>5</sup>

### References

1. Furukawa, T., Yamakawa, T., Midera, T., et al. Selectivities of dihydropyridine derivatives in blocking Ca<sup>2+</sup> channel subtypes expressed in *Xenopus* oocytes. *J. Pharmacol. Exp. Ther.* **291**(2), 464-473 (1999).
2. Johnson, J.D. and Fugman, D.A. Calcium and calmodulin antagonists binding to calmodulin and relaxation of coronary segments. *J. Pharmacol. Exp. Ther.* **226**(2), 330-334 (1983).
3. Perez-Reyes, E., Van Deusen, A.L., and Vitko, I. Molecular pharmacology of human Ca<sub>v</sub>3.2 T-type Ca<sup>2+</sup> channels: Block by antihypertensives, antiarrhythmics, and their analogs. *J. Pharmacol. Exp. Ther.* **328**(2), 621-627 (2009).
4. Ljung, B. Vascular selectivity of felodipine: Experimental pharmacology. *J. Cardiovasc. Pharmacol.* **15**(Suppl 4), S11-S16 (1990).
5. Bolt, G.R. and Saxena, P.R. Acute systemic and regional hemodynamic effects of felodipine, a new calcium antagonist, in conscious renal hypertensive rabbits. *J. Cardiovasc. Pharmacol.* **6**(4), 707-712 (1984).

#### 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, 11/17/2022

#### 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