

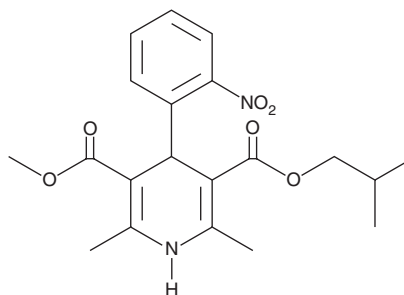
# PRODUCT INFORMATION



## Nisoldipine

Item No. 20998

**CAS Registry No.:** 63675-72-9  
**Formal Name:** 1,4-dihydro-2,6-dimethyl-4-(2-nitrophenyl)-3,5-pyridinedicarboxylic acid, 3-methyl 5-(2-methylpropyl) ester  
**Synonyms:** (±)-BAY-K-5552, (±)-Nisoldipine  
**MF:** C<sub>20</sub>H<sub>24</sub>N<sub>2</sub>O<sub>6</sub>  
**FW:** 388.4  
**Purity:** ≥97%  
**UV/Vis.:** λ<sub>max</sub>: 235, 330 nm  
**Supplied as:** A crystalline solid  
**Storage:** Room temperature  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Nisoldipine is supplied as a crystalline solid. A stock solution may be made by dissolving the nisoldipine in the solvent of choice. Nisoldipine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of nisoldipine in ethanol is approximately 3 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Nisoldipine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, nisoldipine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Nisoldipine has a solubility of approximately 0.1 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Nisoldipine is a calcium channel inhibitor.<sup>1</sup> It binds to calcium channels in isolated rat ventricular membranes (K<sub>d</sub> = 0.04 nM) and inhibits calcium uptake by smooth muscle cells.<sup>1,2</sup> Nisoldipine inhibits acetylcholine-induced contraction of isolated rabbit coronary arteries (IC<sub>50</sub> = 0.03 nM).<sup>1</sup> *In vivo*, nisoldipine (3 mg/kg) reduces ventricular tachycardia and fibrillation and increases survival in a rat model of ventricular arrhythmias induced by myocardial ischemia.<sup>3</sup> Dietary administration of nisoldipine (50-100 mg/kg) reduces systolic blood pressure in spontaneously hypertensive rats.<sup>4</sup> Formulations containing nisoldipine have been used in the treatment of hypertension.

### References

1. Knorr, A. The pharmacology of nisoldipine. *Cardiovasc. Drugs Ther.* **1(4)**, 393-402 (1987).
2. Janis, R.A., Shrikhande, A.V., Greguski, R., et al. Review of nisoldipine binding studies. *Nisoldipine 1987*. Hugenholtz, P.G., and Meyer, J., editors, 1st edition, Springer-Verlag (1987).
3. Fagbemi, O. and Parratt, J.R. Suppression by orally-administered nifedipine, nisoldipine and niludipine of early, life-threatening ventricular arrhythmias resulting from acute myocardial ischaemia. *Br. J. Pharmacol.* **74(1)**, 12-14 (1981).
4. Stasch, J.-P., Kazda, S., Hirth, C., et al. Role of nisoldipine on blood pressure, cardiac hypertrophy, and atrial natriuretic peptides in spontaneously hypertensive rats. *Hypertension* **10(3)**, 303-307 (1987).

#### 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, 10/11/2022

#### 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