

PRODUCT INFORMATION



Niguldipine (hydrochloride)

Item No. 19534

CAS Registry No.: 119934-51-9
Formal Name: 1,4-dihydro-2,6-dimethyl-4-(3-nitrophenyl)-3,5-pyridinedicarboxylic acid, 3-[3-(4,4-diphenyl-1-piperidinyl)propyl] 5-methyl ester, monohydrochloride

Synonyms: B 844-39, NSC 617553

MF: $C_{36}H_{39}N_3O_6 \cdot HCl$

FW: 646.2

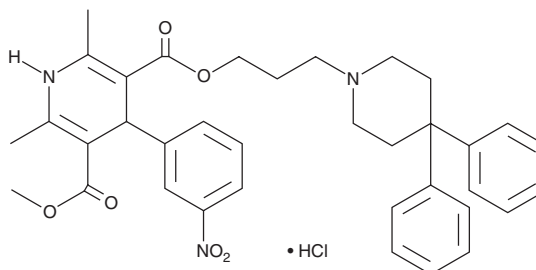
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 354 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Niguldipine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the niguldipine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Niguldipine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of niguldipine (hydrochloride) in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Niguldipine (hydrochloride) is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Niguldipine is a dihydropyridine that acts as a potent, selective α_{1A} -adrenoceptor antagonist ($K_i = 0.16$ nM).¹⁻³ It less potently blocks L- and T-type Ca^{2+} channels (IC_{50} = 0.4 and 0.9 μ M, respectively).^{4,5}

References

1. Li, M.-Y., Tsai, K.-C., and Xia, L. Pharmacophore identification of α_{1A} -adrenoceptor antagonists. *Bioorg. Med. Chem. Lett.* **15**(3), 657-664 (2005).
2. Bylund, D.B. Subtypes of α_1 - and α_2 -adrenergic receptors. *FASEB J.* **6**(3), 832-839 (1992).
3. van Rhee, A.M., Jiang, J.L., Melman, N., et al. Interaction of 1,4-dihydropyridine and pyridine derivatives with adenosine receptors: Selectivity for A_3 receptors. *J. Med. Chem.* **39**(15), 2980-2989 (1996).
4. Klöckner, U. and Isenberg, G. The dihydropyridine niguldipine modulates calcium and potassium currents in vascular smooth muscle cells. *Br. J. Pharmacol.* **97**(3), 957-967 (1989).
5. Perez-Reyes, E., Van Deusen, A.L., and Vitko, I. Molecular pharmacology of human $Ca_v3.2$ T-type Ca^{2+} channels: Block by antihypertensives, antiarrhythmics, and their analogs. *J. Pharmacol. Exp. Ther.* **328**(2), 621-627 (2009).

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, 10/05/2022

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