

Nicardipine . hydrochloride

Ca²⁺ channel inhibitor

Ca²⁺ channel antagonist. Antihypertensive.

Citations: 2

[View Online »](#)

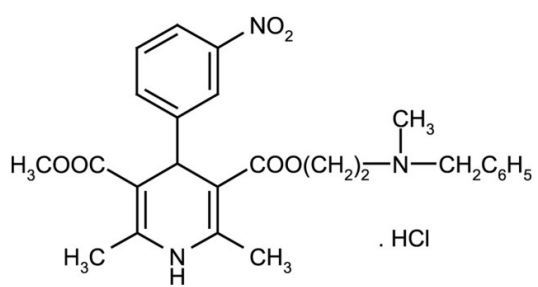
Ordering Information

[Order Online »](#)

ALX-550-273-G001	1g
ALX-550-273-G005	5g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	YC-93 . HCl, 1,4-Dihydro-2,6-dimethyl-4-(3-nitrophenyl)-3,5-pyridinedicarboxylic acid 2-[methyl(phenylmethyl)amino]ethyl ester . HCl
Appearance	Light yellow solid.
CAS	54527-84-3
Couple Target	Calcium channel
Couple Type	Inhibitor
Formula	$C_{26}H_{29}N_3O_6 \cdot HCl$
Identity	Determined by IR.
MI	14: 6495
MW	479.5 . 36.5
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (1mg/ml), ethanol:water 25:75-70:30, propylene glycol, or methanol; slightly soluble in acetone, 100% ethanol, chloroform and water; insoluble in 0.1M NaOH.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

Last modified: May 29, 2024