

Niguldipine . HCl

Ca²⁺ channel blocker

A classic L-type Ca²⁺ channel blocker (IC₅₀=75nM) which can also inhibit T-type Ca²⁺ channels (IC₅₀=244nM). It is a novel inhibitor of drug transport by p-glycoprotein (both stereoisomers are equipotent at inhibition of p-glycoprotein).

Citations: 4

[View Online »](#)

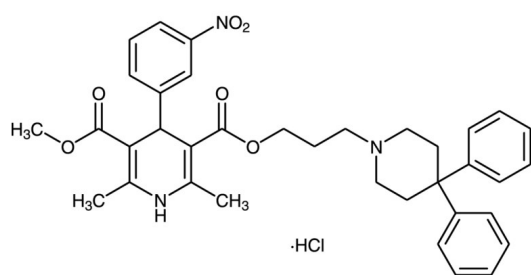
Ordering Information

[Order Online »](#)

BML-CA216-0050	50mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(±)-Niguldipine . hydrochloride, (±)-1,4-Dihydro-2,6-dimethyl-4-(3'-nitrophenyl)-3,5-pyridinedicarboxylic acid 3-(4',4'-diphenyl-1'-piperidiny)propylmethyl ester . HCl
Appearance	Yellow crystalline solid.
CAS	119934-51-9
Couple Target	Calcium channel
Couple Type	Blocker
Formula	$C_{36}H_{39}N_3O_6 \cdot HCl$
MW	646.2
Purity	≥95% (TLC)
Solubility	Soluble in DMSO or methanol.



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-eu@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)