

Isradipine

Ca²⁺ channel blocker

Potent and selective L-type calcium channel blocker. A clinically useful antihypertensive agent which has anti-atherosclerotic effects and improves endothelium-mediated NO-dependent vasodilation. Displays neuroprotective activity in a rat model of focal ischemia.

Citations: 2

[View Online »](#)

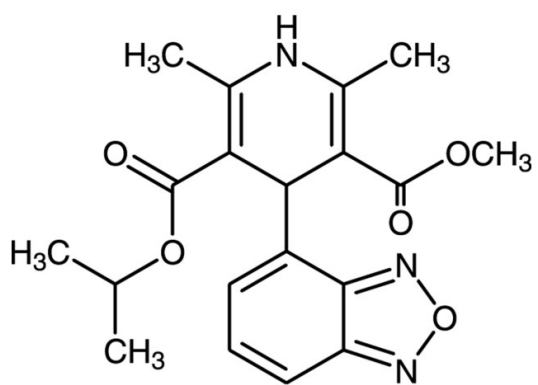
Ordering Information

[Order Online »](#)

BML-CA213-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Yellow solid.
CAS	75695-93-1
Couple Target	Calcium channel
Couple Type	Blocker
Formula	$C_{19}H_{21}N_3O_5$
MW	371.4
Purity	≥98% (HPLC)
RTECS	U57968632
Solubility	Soluble in DMSO (25 mg/ml) or 100% ethanol (15 mg/ml).

Last modified: May 29, 2024



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-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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

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