

3-Bromo-7-nitroindazole

NOS inhibitor

Potent inhibitor of rat cerebellar nitric oxide synthase (nNOS/NOS I). More potent than 7-nitroindazole.

Citations: 8

[View Online »](#)

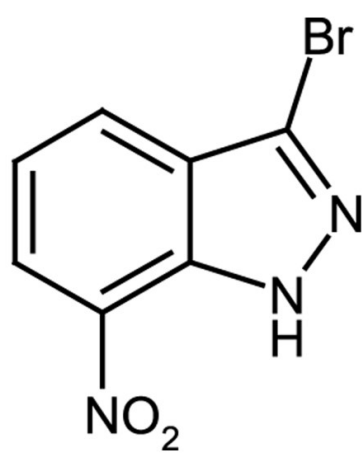
Ordering Information

[Order Online »](#)

ALX-270-009-M005	5mg
------------------	-----

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	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Yellowish solid.
CAS	74209-34-0
Couple Target	NOS
Couple Type	Inhibitor
Formula	$C_7H_4BrN_3O_2$
MW	242.0
Purity	≥98%
Solubility	Soluble in DMSO (50mg/ml), methanol or 100% ethanol; also soluble in water containing 10% DMSO or in a 50:50 mixture of water:DMSO.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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