

7-Nitroindazole

7-Nitroindazole is a selective and BBB-penetrable inhibitor of nitric oxide synthase

7-Nitroindazole is a selective and BBB-penetrable inhibitor of nitric oxide synthase (NOS). 7-Nitroindazole has anti-injurious and neuroprotective properties.

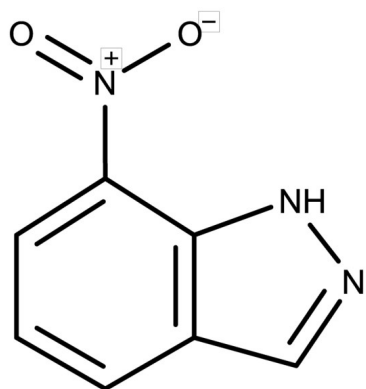
Ordering Information

[Order Online »](#)

ENZ-CHM465-0001	1g
-----------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1H-Indazole, 7-nitro

Appearance Light yellow to yellow solid.

CAS 2942-42-9

Couple Target NOS

Couple Type Inhibitor

Formula $C_7H_5N_3O_5$

MW 163.14

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (100mg/ml).



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