

iNOS (human), (recombinant) (His- tag)

Nitric oxide synthases are a family of enzymes catalyzing the production of nitric oxide (NO) from L-arginine. NO is an important cellular signaling molecule. It helps modulate vascular tone, insulin secretion, airway tone, and peristalsis, and is involved in angiogenesis and neural development. It may function as a retrograde neurotransmitter. Nitric oxide is mediated in mammals by the calcium-calmodulin controlled isoenzymes eNOS (endothelial NOS) and nNOS (neuronal NOS). The inducible isoform, iNOS, is involved in immune response, binds calmodulin at physiologically relevant concentrations, and produces NO as an immune defense mechanism, as NO is a free radical with an unpaired electron. It is the proximate cause of septic shock and may function in autoimmune disease.

Ordering Information

[Order Online »](#)

ALX-201-852-0010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -80°C. Keep stock vial on ice during experiments.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	NOS II, Nitric oxide synthase (inducible)
Formulation	Liquid. In 50 mM TRIS, pH 7.4, containing 0.1 mM EDTA, 10% glycerol, 100 mM NaCl, 0.1 mM DTT, 10 µM tetrahydrobiopterin.
MW	~130kDa/subunit; homodimer.
Purity	≥90% (SDS-PAGE)
Source	Produced in <i>E. coli</i> . Human iNOS is fused to an N-terminal His-tag.
Specific Activity	≥100 U/mg. 1 U is defined as the amount of enzyme required to produce 1 nmol of NO (nitric oxide) per minute at 37°C. NO synthesis is measured by the conversion of oxyhemoglobin to methemoglobin.
UniProt ID	P35228

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-
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)