

Nitric oxide synthase (inducible) polyclonal antibody

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 13

[View Online »](#)

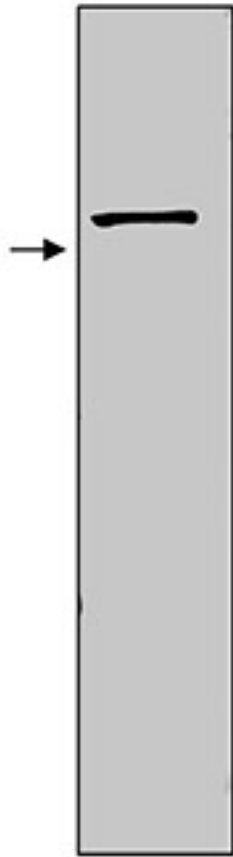
Ordering Information

[Order Online »](#)

BML-SA200-0025	25µl
BML-SA200-0100	100µl

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of lysate from mouse macrophage RAW 264.7 cell line, which was induced with IFN γ and LPS. The nitrocellulose membrane was probed with Anti-iNOS [Prod. no. BML-SA200] at 1/1000 dilution. Blot was developed with "Vectastain ABC Kit" and BCIP/NBT. Arrow points to MW ~130kDa.

Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	iNOS, NOS II
Application	ICC, IHC, WB
Application Notes	Detects a band of ~130kDa by Western blot.
Crossreactivity	Does not cross-react with eNOS or nNOS.
Formulation	Liquid. In PBS containing 0.05% sodium azide.
Gene/Protein Identifier	NM_010927 (RefSeq)
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 1131-1141 of mouse macrophage NOS.
Recommendation Dilutions/Conditions	Immunocytochemistry (1:500)Immunohistochemistry (1:500)Western Blot (1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	P29477
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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)