

Nitric oxide synthase (universal) monoclonal antibody (NOS-3F7-B11-B5)

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

Citations: 3

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name NOS (universal)

Application IHC (FS), WB

Application Notes Detects bands of ~130kDa (representing iNOS in samples first induced with interferon (IFN) and lipopolysaccharides (LPS)), ~155kDa (in tissues expressing nNOS) and ~140kDa (in tissues expressing eNOS) by Western blot.

Clone NOS-3F7-B11-B5

Formulation Liquid. Ascites diluted in PBS containing 0.05% sodium azide.

Host Mouse

Immunogen Purified neuronal nitric oxide synthase (nNOS; NOS I) from bovine brain.

Isotype IgM

Recommendation Dilutions/Conditions Immunohistochemistry (frozen sections, 1:100)Western Blot (1:50-1:500)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Bovine, Mouse, Rat

Specificity Recognizes iNOS (NOS II), eNOS (NOS III) and nNOS (NOS I).

UniProt ID E1BK13

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-
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)