

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

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

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-029-R200	200µl
------------------	-------

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** Immunohistochemistry (frozen sections, 1:100)Western Blot (1:50-1:500)Suggested  
**Dilutions/Conditions** 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



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)  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France

Phone: +33 1 72 440 1335  
[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)