

# Nitrotyrosine monoclonal antibody (7A5)

The enzymatic conversion of arginine to citrulline by nitric oxide synthase (NO Synthase) leads to the formation of nitric oxide (NO). Peroxynitrite is formed following reaction of NO with superoxide, producing a potent oxidant that reacts with a variety of biological molecules, including tyrosine residues in proteins. Tyrosine nitration can inactivate enzymes and receptors that depend on tyrosine residues for their activity. Nitration prevents phosphorylation of tyrosine residues important for signal transduction.

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

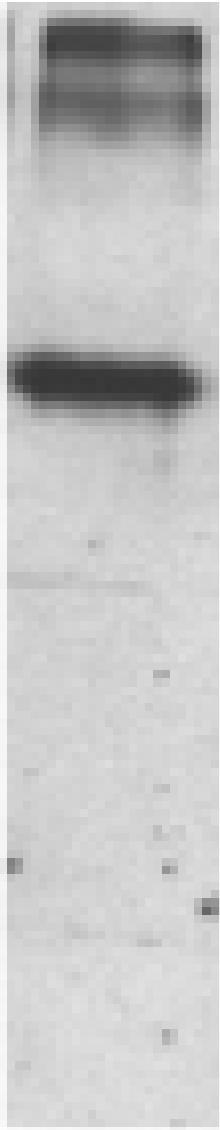
Citations: 3

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

67 kDa -



Western blot analysis of Nitrotyrosine-modified BSA probed with Nitrotyrosine mAb (7A5) at 1 $\mu$ g/ml.

## Handling & Storage

<b>Use/Stability</b>	Stable at -80°C up to 1 year, at 4°C up to 3 months.
<b>Handling</b>	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Application</b>	ELISA, WB
<b>Clone</b>	7A5
<b>Formulation</b>	Lyophilized from 1ml of 2x PBS containing 0.09% sodium azide, PEG, and sucrose.
<b>Host</b>	Mouse
<b>Immunogen</b>	KLH modified with Nitrotyrosine.
<b>Isotype</b>	IgG1
<b>Purity Detail</b>	Thiophilic adsorption and size exclusion chromatography purified.
<b>Recommendation Dilutions/Conditions</b>	ELISA (0.05µg/ml)Western Blot (0.5µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Reconstitution</b>	Reconstitute with 1ml water (15 minutes at room temperature).
<b>Species Reactivity</b>	Dog, Human, Mouse, Rat
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

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)