

Nitric oxide synthase (endothelial) polyclonal antibody

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

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-509-R100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

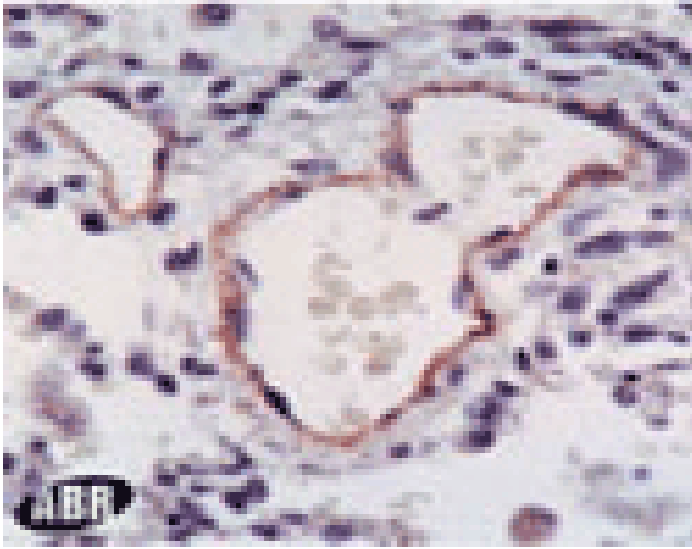


Figure: Immunolocalization of endothelial nitric oxide synthase (eNOS; NOS III) in human lung tissue using PAb to eNOS (Prod. No. ALX-210-509).

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 eNOS, NOS III

Application IHC (FS), IHC (PS), WB

Application Notes Detects a band of ~140kDa by Western blot.

Crossreactivity Does not cross-react with human iNOS (NOS II) or rat nNOS (NOS I).

Formulation Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 1179-1194 (S1179LQERQLRGAVPWAFD1194) of human endothelial nitric oxide synthase (eNOS; NOS III).

Purity Detail Epitope-affinity purified IgG.

Recommendation Dilutions/Conditions Immunohistochemistry (frozen sections, 1:100)Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Bovine, Dog, Human, Mouse, Porcine, Rat

UniProt ID P29474

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)