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 »

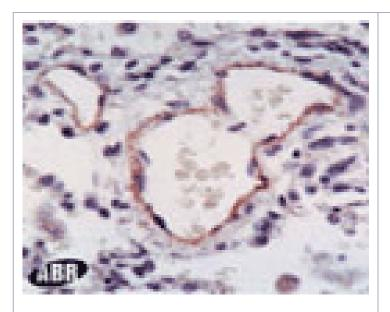


Figure: Immunolocalization of endothelial nitric oxide synthase (eNOS; NOS III) in human lung tissue using PAb to eNOS (Prod. No. ALX-210-509).

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 NotesDetects 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

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com