Endothelin A receptor polyclonal antibody

Endothelins are a family of peptides that are potent vasoconstrictors produced by mammalian vascular endothelial cells. The receptors for the endothelins are found on various tissues including, but not restricted to, lung and intestine.

This antibody is covered by our Worry-Free Guarantee.

Citations: 12

View Online »

Ordering Information

Order Online »

ALX-210-507A-C250

250µg

Manuals, SDS & CofA

View Online »

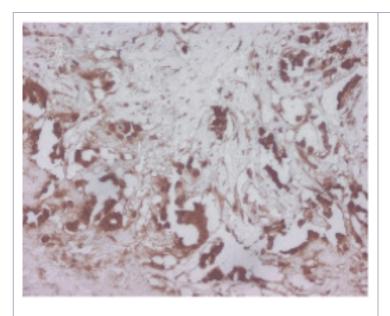


Figure: Immunohistochemical staining of invasive ductal carcinoma (G2) using PAb to Endothelin A Receptor (Prod. No. <u>ALX-210-507A</u>). Expression can be seen in the cytoplasma and in the cell membrane.""",Endothelin A receptor polyclonal antibody Immunohistochemistry,Endothelin A receptor polyclonal antibody Immunohistochemistry,yes ALX-210-509,210-

509_1.gif,http://static.enzolifesciences.com/fileadmin/files/image/210-509_1.gif,**Figure:** Immunolocalization of endothelial nitric oxide synthase (eNOS; NOS III) in human lung tissue using PAb to eNOS (Prod. No. <u>ALX-210-509</u>)."""""

Handling & Storage

Handling Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name ETA-receptor

Application IHC, WB

Application NotesAs this antibody is generated against the cytosolic portion

of endothelin A receptor, permeabilization is required for

immunochemical analysis.

Detects bands of 34-52kDa by Western blot.

Does not block the binding of endothelins to the receptor.

Formulation Liquid. In TRIS-citrate-phosphate, pH 7-8, containing 0.1%

 $so dium\ azide.$

Host Sheep

Immunogen Synthetic peptide corresponding to aa 410-422

(Q410EQNHNTERSSHK422) of C-terminal rat ETA-

receptor (endothelin A receptor).

Purity Detail Epitope-affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections, 4µg/ml)Western

Blot (10-20µg/ml using ECL)Suggested

dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Human, Mouse, Rat

Specificity

Recognizes an epitope on the cytoplasmic side of the $\mathsf{ET}_\mathtt{A}$ -receptor.

UniProt ID

P26684

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-