## Estrogen receptor β polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 8

View Online »

**Ordering Information** 

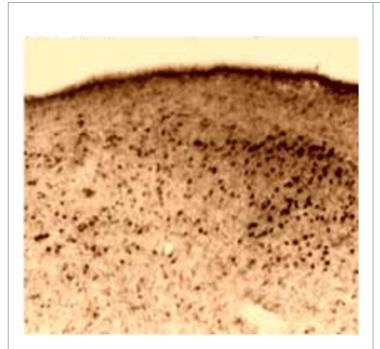
Order Online »

ALX-210-135-C050

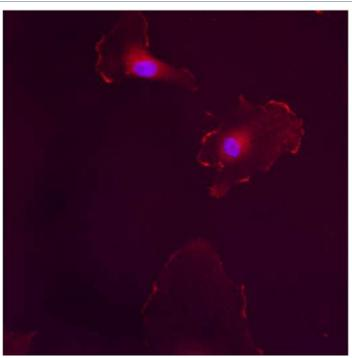
50µg

Manuals, SDS & CofA

**View Online »** 



**Figure 2:** IHC of ER β in primate amygdala using Estrogen receptor β, pAb (Prod. No. ALX-210-135).



**Figure 1**: Estrogen receptor  $\beta$ , pAb (Prod. No. ALX-210-135) used in immunofluorescence.

## **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 ERβ

**Application** Gel Shift, ICC, IHC (FS), IHC (PS), WB

**Application Notes**Detects a band of ~55kDa by Western blot.

Crossreactivity Does not cross-react with ERa.

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

azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to aa 100-115

(A100EPQKSPWCEARSLEH115) of human, mouse and

rat ER $\beta$  (estrogen receptor  $\beta$ ).

Purity Detail Epitope-affinity purified IgG.

Recommendation Dilutions/Conditions Gel Shift Assay (1-5µg/ml)Immunocytochemistry (2-

4µg/ml)Immunohistochemistry (paraffin sections, 1-

2μg/ml)Western Blot (1-2μg/ml)Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Human, Monkey, Mouse, Rat, Sheep

UniProt ID Q92731

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com