

# Estrogen receptor $\alpha$ polyclonal antibody

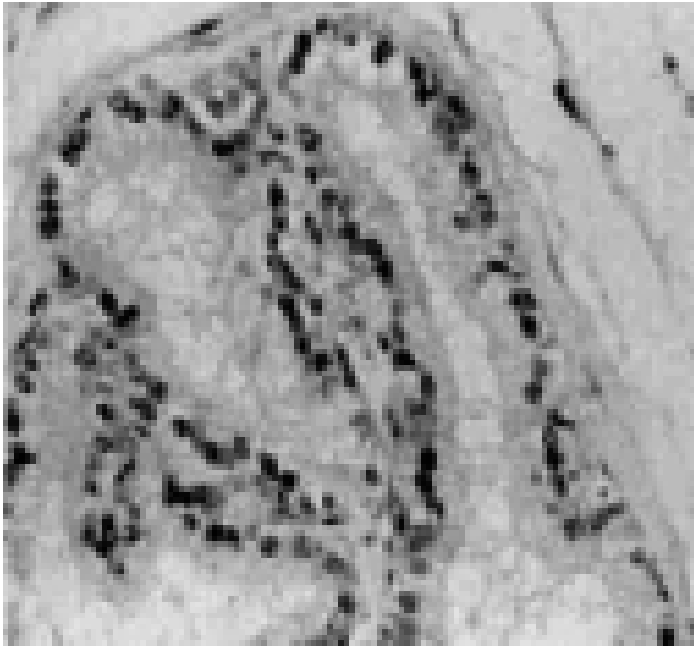
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 4

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



**Figure:** Immunohistochemistry of ER $\alpha$  in rat prostate using PAb to Estrogen Receptor  $\alpha$  (Prod. No. ALX-210-201)

## 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** ER $\alpha$

**Application** IHC (FS), WB

**Application Notes** Detects a band of ~64kDa by Western blot.

**Crossreactivity** Does not cross-react with ER $\beta$ .

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

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to aa 1-21 (M1TMTLHTKASGMALLHQIQGN21) of N-terminal human ER $\alpha$  (estrogen receptor  $\alpha$ ). This sequence is completely conserved in rat and mouse ER $\alpha$  and is highly conserved in sheep, pig, dog, bovine and chicken ER $\alpha$ .

**Purity Detail** Epitope-affinity purified IgG.

**Recommendation Dilutions/Conditions** Immunohistochemistry (frozen sections, 1-5 $\mu$ g/ml)Western Blot (1 $\mu$ g/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Human, Rat

**Specificity** Recognizes endogenous ER $\alpha$ .

**UniProt ID** P03372

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)