

Estrogen receptor α polyclonal antibody

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

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-201-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

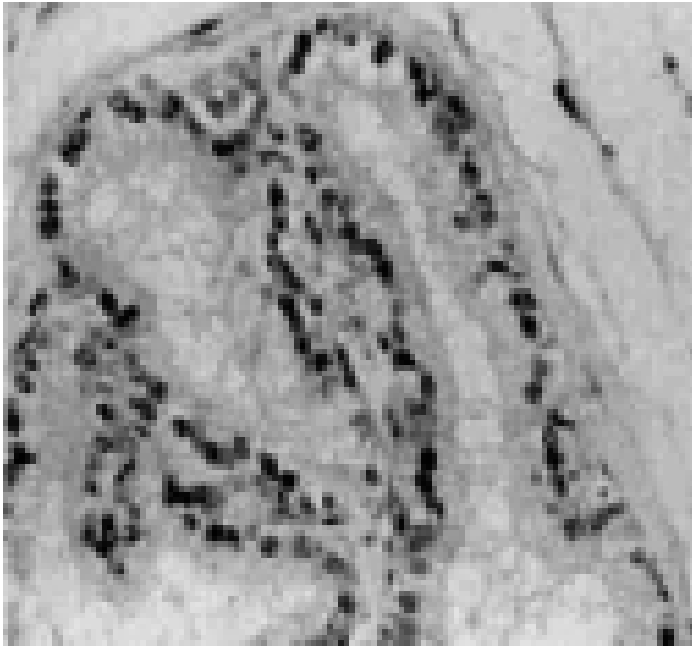


Figure: Immunohistochemistry of ERα in rat prostate using PAb to Estrogen Receptor α (Prod. No. ALX-210-201)

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 ERα

Application IHC (FS), WB

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

Crossreactivity Does not cross-react with ERβ.

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

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 1-21 (M1TMTLHTKASGMALLHQQGN21) of N-terminal human ERα (estrogen receptor α). This sequence is completely conserved in rat and mouse ERα and is highly conserved in sheep, pig, dog, bovine and chicken ERα.

Purity Detail Epitope-affinity purified IgG.

Recommendation Immunohistochemistry (frozen sections, 1-5µg/ml)Western Blot (1µg/ml)Suggested
Dilutions/Conditions 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 ERα.

UniProt ID P03372



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AS
Phone: +41 61 926 8989
info-
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-
be@enzolifesciences.com

France
Phone: +33 472 440 655
info-
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-
de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-
uk@enzolifesciences.com

This antibody is covered by our **Worry-Free Guarantee**.