

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	50 μ g
------------------	------------

Manuals, SDS & CofA

[View Online »](#)

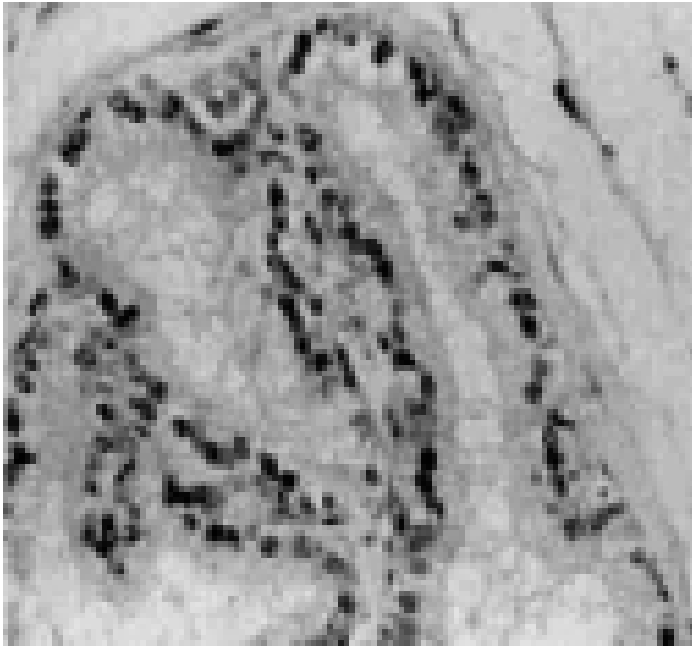


Figure: Immunohistochemistry of ERe in rat prostate using PAb to Estrogen Receptor a (Prod. No. ALX-210-201)

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

UniProt ID P03372

Worry-free Guarantee

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

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)