Estrogen receptor α monoclonal antibody (33)

Steroid receptors are a large superfamily of ligand-dependent intracellular proteins that stimulate transcription of specific genes by binding to specific DNA sequences following activation by the appropriate hormone. Estrogen receptor (ER) has been classified into two distinct isoforms, α and β . ER α is known to interact in a ligand dependent and independent manner with several known coactivators including Receptor-Interacting Protein 140 (RIP140/ERAP140), ERAP160, and Steroid Receptor Coactivator-1 (SRC-1) to enhance transcriptional activity of target genes.

This antibody is covered by our Worry-Free Guarantee.

Citations: 5

View Online »

Ordering Information

Order Online »

ALX-803-004-C050

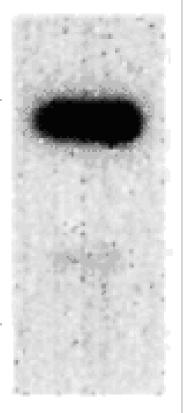
50µg

Manuals, SDS & CofA

View Online »

Fig. 1

67 kDa



56 kDa

Figure: Western blot of recombinant ERa using MAb to Estrogen Receptor α (33) (Prod. No. <u>ALX-803-004</u>).""",Estrogen receptor α monoclonal antibody (33) Western blot,Estrogen receptor α monoclonal antibody (33) Western blot,yes

ALX-803-005,803-

005_color.gif,http://static.enzolifesciences.com/fileadmin/files/image/803-

005_color.gif, **Figure:** Immunohistochemical staining of PRL receptor in rat Islets of Langerhans using MAb to Prolactin Receptor (U5) (Prod. No. ALX-803-005)."""""

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 ERa

Application Flow Cytometry, ICC, IF, IHC (FS), IHC (PS), IP, WB

Application Notes Immunohistochemistry: Staining of ERα in mouse Leydig

cells results in both nuclear and cytoplasmic staining.

Detects a band of ~67kDa by Western blot.

Does not inhibit binding of ER to nuclei and DNA.

Clone 33

Crossreactivity Does not cross-react with androgen receptor, PR

(progesterone receptor), GR (glucocorticoid receptor) or

ERβ.

Formulation Liquid. In Dulbecco's PBS, containing 0.05% sodium

azide, 1mg/mL BSA and 30% glycerol.

Host Mouse

Immunogen Synthetic peptide corresponding to aa 247-261

(E247VGMMKGGIRKDRRG261) of the DNA binding domain of ER α (estrogen receptor α). This sequence is completely conserved in many species including mouse,

rat, pig, bovine, sheep, chicken and frog.

lsotype lgG1

Recommendation Dilutions/Conditions Immunofluorescence (10µg/mL)Immunohistochemistry

(1:20-1:100)Western Blot (4 µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined

individually for each application.

Species Reactivity

Bovine, Human, Monkey, Mouse, Rat

Specificity

Recognizes both the steroid occupied and unoccupied

forms of ERα.

UniProt ID

P03372

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



ENZO LIFE SCIENCES,

Phone: 800.942.0430

info-