

ERR alpha Antibody - ChIP Grade

Cat. No. C15410230-100

Type: Polyclonal	Specificity: Human, mouse, rat
Size: 100 µl	Isotype: NA
Concentration: 0.7 µg/µl	Host: Rabbit
Lot No.: 39764	Purity: Affinity purified
Storage buffer: NA	Storage conditions: NA
Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.	

Description

Polyclonal antibody raised in rabbit against **ERR alpha** (estrogen-related receptor alpha), using a KLH-conjugated synthetic peptide.

Applications

Applications	Suggested dilution	References
ChIP assay	5 µg/ChIP	Fig 3
Immunoprecipitation	1:100 - 1:500	Fig 1
Western Blotting	1:500 - 1:3,000	Fig 2, 4, 5

* Optimal dilutions/concentrations should be determined by the researcher.

Validation Data

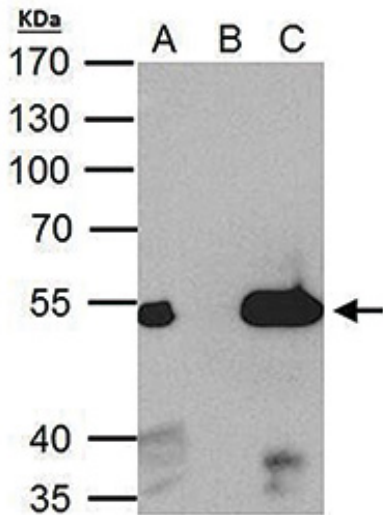


Figure 1. IP

ERR alpha antibody immunoprecipitates ERR alpha protein in IP experiments. IP Sample: 293T whole cell extract A. 40 µg 293T whole cell extract B. Control with 2 µg of preimmune rabbit IgG C. Immunoprecipitation of ERR alpha protein by 2 µg of ERR alpha antibody. The immunoprecipitated ERR alpha protein was detected by western blot with the ERR alpha antibody (Cat. No. C15410230) diluted 1:1,000.

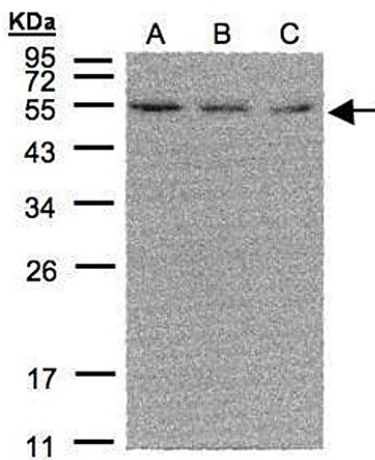


Figure 2. Western blot

Sample: 30 µg of whole cell lysate A: 293T B: A431 C: H1299 12% SDS PAGE C15410230 diluted 1:500

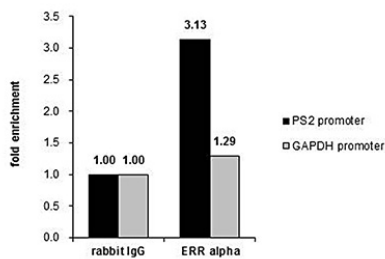


Figure 3. ERR alpha antibody ChIP results

ChIP was performed with MCF-7 chromatin extract and 5 µg of either control rabbit IgG or ERR alpha antibody. The precipitated DNA was detected by PCR with primer set, targeting to PS2 promoter, used as a positive control or GAPDH promoter, used as a negative control.

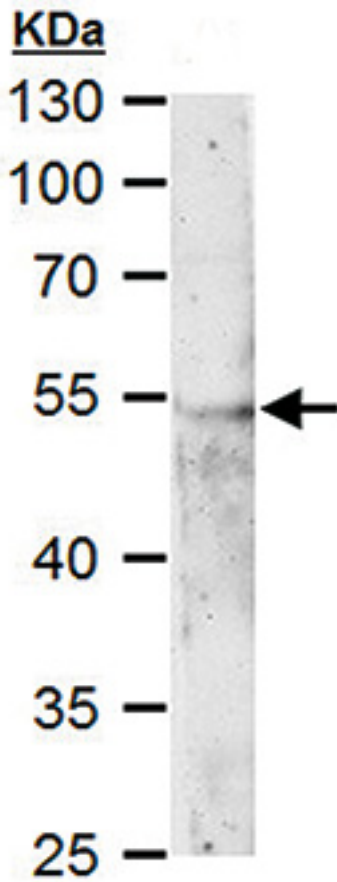


Figure 4. ERR alpha antibody WB results

ERR alpha antibody detects ERR alpha protein by western blot analysis. 50 μ g mouse kidney extract 10 % SDS-PAGE ERR alpha antibody (cat. No. C15410230) dilution 1:500

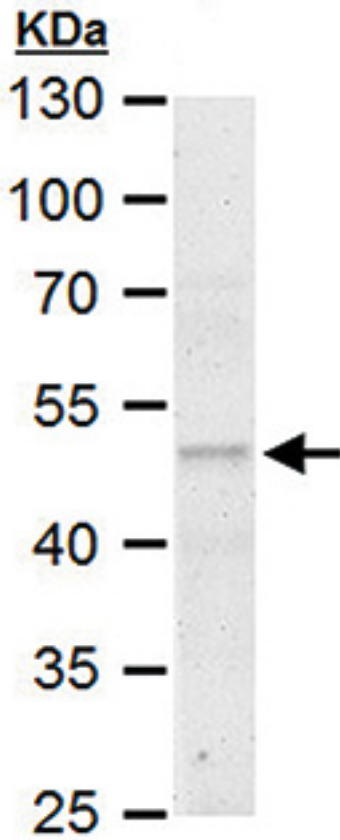


Figure 5. ERR alpha antibody WB results

ERR alpha antibody detects ERR alpha protein by western blot analysis. 50 μ g rat kidney extract 10 % SDS-PAGE ERR alpha antibody (Cat. No. C15410230) dilution 1:500