

Estrogen Receptor α antibody (mAb)

Catalog No: 61035

RRID: AB_2793482

Clone: F3

Isotype: IgG1

Application(s): ChIP, IP, WB

Reactivity: Human

Quantity: 100 μ g

Purification: Protein G Chromatography

Host: Mouse

Concentration: 0.61 μ g/ μ l

Molecular Weight: 66 kDa

Background: ER- α (estrogen receptor α) is a ligand-activated transcription factor composed of several domains important for hormone binding, DNA binding and activation of transcription. ER- α localizes to the nucleus where it may form a homodimer or a heterodimer with estrogen receptor 2. Estrogen and its receptors are essential for sexual development and reproductive function, but also play a role in other tissues such as bone. Estrogen receptors are also involved in pathological processes including breast cancer, endometrial cancer and osteoporosis.

Immunogen: This ER-alpha antibody was raised against a peptide within the C-terminal region of human Estrogen Receptor alpha.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

WB: 1:500 - 1:1,000 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.