

ER- α antibody (mAb)

Catalog No: 61035, 61036

Host: Mouse

Isotype: IgG1

Application: ChIP, EMSA, IF, IP, WB

Reactivity: Human

Purification: Protein G Chromatography

Volume: 100 μ l, 10 μ l

Concentration: 1 μ g/ μ l

Molecular Weight: 66 kDa

Clone: F3

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: Peptide within the region of human Estrogen receptor alpha.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 32 mM glycine, 0.07 mM EDTA, 30% glycerol, and 0.035% sodium azide.

Application Notes:

See references for details.

References:

ChIP: Welboren, W-J. et al (2009) EMBO J. 28:1418-1428.

Storage and Guarantee: Antibodies in solution can be stored at -20°C for 2 years. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of arrival.