

Estrogen Receptor β antibody (mAb)

Catalog Nos: 39767, 39768

RRID: AB_2793335

Clone: CO1531

Isotype: IgG2a

Application(s): IHC, IP, WB

Reactivity: Human

Quantities: 100 μ g, 10 μ g

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 μ g/ μ l

Molecular Weight: 59 kDa

Background: ER β (estrogen receptor β) is a ligand-activated transcription factor composed of several domains important for hormone binding, DNA binding and activation of transcription. ER β localizes to the nucleus where it may form a homodimer or a heterodimer with ER alpha. 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 Estrogen Receptor beta antibody was raised against a recombinant protein corresponding to amino acids 212-477 of human Estrogen Receptor beta.

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:

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.