

## Estrogen Receptor- $\beta$ (Ab-105) Antibody [#B0922]

<b>Catalog Number:</b>	B0922
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q92731
<b>Other Names:</b>	ESR2; ESTRB; Estrogen receptor beta; NR3A2
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized non-phosphopeptide derived from human Estrogen Receptor- $\beta$ around the phosphorylation site of serine 105 (Q-K-S <sup>P</sup> -P-W).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Estrogen Receptor- $\beta$ (Ab-105) antibody detects endogenous levels of total Estrogen Receptor- $\beta$ protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Ogawa S., Biochem. Biophys. Res. Commun. 243:122-126(1998). Moore J.T., Biochem. Biophys. Res. Commun. 247:75-78(1998). Mosselman S., FEBS Lett. 392:49-53(1996).