

HER2 (Ab-1221/1222) Antibody [#B7103]

- Catalog Number:** B7103-1, B7103-2, B7103
Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl
Form of Antibody: Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Storage/Stability: Store at -20°C/1 year
Immunogen: The antiserum was produced against synthesized non-phosphopeptide derived from human HER2 around the phosphorylation site of tyrosine 1221/1222.
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity: HER2 (Ab-1221/1222) antibody detects endogenous levels of total HER2 protein.
Reactivity: Human
Applications: WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:20000
Swiss-Prot No. : P04626
References: Marone R, et al. (2004) Nat Cell Biol; 6(6): 515-22.
Ren Z, et al. (2002) J Biol Chem; 277(41): 38486-93.
Ouyang X, et al. (1998) Carcinogenesis; 19(11): 2013-9.
Ouyang X, et al. (1998) Exp Cell Res; 241(2): 467-75.
Klapper, L. N. et al. (2000) Cancer Res. 60, 3384-3388.