

HER3/ErbB3 (Ab-1289) Antibody [#B0945]

Catalog Number:	B0945
Concentration:	1mg/ml
Swiss-Prot No.:	P21860
Other Names:	EC 2.7.10.1, ERB3, ERBB3, HER3, Receptor protein-tyrosine kinase erbB-3 precursor, Tyrosine kinase-type cell surface receptor HER3, c-erbB3, kinase ErbB3
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized non-phosphopeptide derived from human HER3/ErbB3 around the phosphorylation site of tyrosine 1289 (Q-G-Y ^P -E-E).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	HER3/ErbB3 (Ab-1289) antibody detects endogenous levels of total HER3/ErbB3 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	Katoh M., Biochem. Biophys. Res. Commun. 192:1189-1197(1993). Lessor T.J., Mol. Cell. Endocrinol. 175:185-191(2001). Kinugasa Y., Biochem. Biophys. Res. Commun. 321:1045-1049(2004).