

CD32(Phospho-Tyr292) Antibody [#8A0626]

Catalog Number:	8A0626
Concentration:	1mg/ml
Origin of Antibody:	Rabbit
Swiss-Prot No.:	P31994
Other Names:	CD32, CDW32, FC gamma receptor IIB, FC-gamma RII, FC-gamma RII-B, FC-gamma-RIIB, FCG2, FCG2B, FCGR2, FCGR2B, FCRII, FCgammaRIIB, IGFR2, IGG FC receptor II beta, IGG FC receptor II-B, LY-17, Low affinity immunoglobulin gamma FC region receptor II precursor, Low affinity immunoglobulin gamma FC region receptor II-B precursor.
All Sites:	Human: Tyr292
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 phosphopeptide derived from human CD32 around the phosphorylation site of tyrosine 292 (I-T-Y ^P -S-L).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
Specificity:	CD32 (Phospho-Tyr292) antibody detects endogenous levels of CD32 only when phosphorylated at tyrosine 292.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:1000
References:	Stuart S.G., EMBO J. 8:3657-3666(1989). Brooks D.G., J. Exp. Med. 170:1369-1385(1989). Engelhardt W., Eur. J. Immunol. 20:1367-1377(1990).