

CD88/C5aR (Ab-338) Antibody [#B0853]

Catalog Number:	B0853
Concentration:	1mg/ml
Swiss-Prot No.:	P21730
Other Names:	C5AR, C5AR1, C5R1, C5a anaphylatoxin chemotactic receptor, C5a-R, CD88 antigen
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 CD88/C5aR around the phosphorylation site of serine 338 (T-R-S ^P -T-V).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CD88/C5aR (Ab-338) antibody detects endogenous levels of total CD88/C5aR protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	W Fureder, J. Immunol., Sep 1995; 155: 3152. T Werfel, J. Immunol., Aug 1996; 157: 1729. Anthony A. Floreani, J. Immunol., May 1998; 160: 5073. M Burg, J. Immunol., Dec 1996; 157: 5574.