

Integrin β 1 (Ab-788) Antibody [#B0445]

Catalog Number:	B0445
Concentration:	1mg/ml
Swiss-Prot No.:	P05556
Other Names:	CD29, FNRB, Fibronectin receptor beta subunit, ITB1, Integrin VLA-4 beta subunit, Integrin beta-1 precursor, integrin beta-1
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 Integrin β 1 around the phosphorylation site of threonine 788 (A-V-T ^P -T-V).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Integrin β 1 (Ab-788) antibody detects endogenous levels of total Integrin β 1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Yves Benninger, J. Neurosci., Jul 2006; 26: 7665 - 7673. Deepak K. Sharma, Cancer Res., Sep 2005; 65: 8233 - 8241. Lia Scotti Campos, J. Biol. Chem., Feb 2006; 281: 5300 - 5309.