

## Integrin $\alpha$ 4 (Ab-1027) Antibody [#B1071]

<b>Catalog Number:</b>	B1071
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P13612
<b>Other Names:</b>	CD49d, ITA4, Integrin alpha-4 precursor, Integrin alpha-IV, VLA-4
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized non-phosphopeptide derived from human Integrin $\alpha$ 4 around the phosphorylation site of serine 1027 (R-D-S <sup>P</sup> -W-S).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Integrin $\alpha$ 4 (Ab-1027) antibody detects endogenous levels of total Integrin $\alpha$ 4 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Takada Y., EMBO J. 8:1361-1368(1989). Rubio M., Eur. J. Immunol. 22:1099-1102(1992). Szabo M.C., Mol. Immunol. 32:1453-1454(1995).