

Smad1 (Ab-465) Antibody [#B0029]

Catalog Number:	B0029
Concentration:	1mg/ml
Swiss-Prot No.:	Q15797
Other Names:	BSP-1; BSP1; Dwarf1-A; Dwf-A; JV4-1; Mad-related protein 1; Mad1; MADH1; MADR1; Mothers against decapentaplegic homolog 1; Mothers against DPP homolog 1; Mothers-against-DPP-related-1; Smad 1; SMAD 1; Transforming growth factor- beta signaling protein-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 Smad1 around the phosphorylation site of serine 465 (I-S-S-V-S ^P).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Smad1 (Ab-465) antibody detects endogenous levels of total Smad1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Riggins G.J., Nat. Genet. 13:347-349(1996). Liu F., Nature 381:620-623(1996). Hoodless P.A., Cell 85:489-500(1996).