

53BP1 (Ab-29) Antibody [#B8284]

Catalog Number:	B8284
Concentration:	1mg/ml
Swiss-Prot No.:	Q12888
Other Names:	p53-binding protein 1; P531; TP53B; TP53BP1; TRP53BP1; Tumor suppressor p53-binding 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 peptide derived from Internal of human 53BP1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	53BP1 (Ab-29) antibody detects endogenous levels of total 53BP1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Iwabuchi K., J. Biol. Chem. 273:26061-26068(1998). Nakajima D., Submitted (MAR-2005) to the EMBL/GenBank/DDBJ databases. Bechtel S., BMC Genomics 8:399-399(2007).