

p53 (Ab-20) Antibody [#B0529]

Catalog Number:	B0529
Concentration:	1mg/ml
Swiss-Prot No.:	P04637
Other Names:	Antigen NY-CO-13, Cellular tumor antigen p53, Phosphoprotein p53, TP53, Tumor suppressor p53
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 p53 around the phosphorylation site of serine 20 (T-F-S ^P -D-L).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	p53 (Ab-20) antibody detects endogenous levels of total p53 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Lloyd M. Alderson, Cancer Res., Mar 1995; 55: 999 - 1001. WE Mercer, PNAS, Aug 1990; 87: 6166. Timothy J. Underwood, AACR Meeting Abstracts, Apr 2006; 2006: 1167. Leandro Cerchietti, Blood (ASH Annual Meeting Abstracts), Nov 2006; 108: 1413.