

Ku70 (Ab-5) Antibody [#B0449]

Catalog Number:	B0449
Concentration:	1mg/ml
Swiss-Prot No.:	P12956
Other Names:	70 kDa subunit of Ku antigen, ATP-dependent DNA helicase II, 70 kDa subunit, CTC box binding factor 75 kDa subunit, CTC75, CTCBF, G22P1, Lupus Ku autoantigen protein p70, TLAA, Thyroid-lupus autoantigen
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 Ku70 around the phosphorylation site of serine 5 (W-E-S ^P -Y-Y).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Ku70 (Ab-5) antibody detects endogenous levels of total Ku70 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
References:	Nicholas S. Y. Ting, Nucleic Acids Res., Apr 2005; 33: 2090 - 2098. Suparna Mazumder, Mol. Cell. Biol., May 2007; 27: 3511 - 3520. Roopashree Narasimhaiah, Cereb Cortex, Jun 2005; 15: 696 - 707.