

## DDX5/DEAD-box Protein 5 Antibody [#B0903]

<b>Catalog Number:</b>	B0903
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P17844
<b>Other Names:</b>	DEAD-box protein 5, DEAD-box protein p68, EC 3.6.1.-, HELR, HLR1, Probable RNA-dependent helicase p68
<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 peptide derived from Internal of human DDX5/DEAD-box Protein 5.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	DDX5/DEAD-box Protein 5 antibody detects endogenous levels of total DDX5/DEAD-box Protein 5 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IF: 1:100~1:500 ELISA: 1:20000
<b>References:</b>	Hloch P., Nucleic Acids Res. 18:3045-3045(1990). Roessler O.G., Nucleic Acids Res. 28:932-939(2000). Molina H., Proc. Natl. Acad. Sci. U.S.A. 104:2199-2204(2007).