

## OVOL1 Antibody [#C10046]

<b>Catalog Number:</b>	C10046
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O14753
<b>Other Names:</b>	Putative transcription factor Ovo-like 1, hOvo1, OVOL1
<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 N-terminal of human OVOL1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	OVOL1 antibody detects endogenous levels of total OVOL1 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	The German cDNA consortium, Submitted (SEP-2004) to the EMBL/GenBank/DDBJ databases. The MGC Project Team, Genome Res. 14:2121-2127(2004). Chidambaram A., Mamm. Genome 8:950-951(1997).