

## ZP1 Antibody [#C19635]

<b>Catalog Number:</b>	C19635
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P60852
<b>Other Names:</b>	mgc87693; zona pellucida glycoprotein 1; zona pellucida glycoprotein 1 (sperm receptor); zp1; zp1 protein
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), 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 ZP1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ZP1 antibody detects endogenous levels of total ZP1 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000   ELISA: 1:10000
<b>References:</b>	Hughes D.C., Biochim. Biophys. Acta 1447:303-306(1999). Taylor T.D., Nature 440:497-500(2006).