

ZP4 Antibody [#C19638]

Catalog Number:	C19638
Concentration:	1mg/ml
Swiss-Prot No.:	Q12836
Other Names:	ZBP; zona pellucida sperm-binding 4; zona pellucida sperm-binding B; ZP1; ZPB
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 peptide derived from internal of human ZP4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ZP4 antibody detects endogenous levels of total ZP4 protein.
Reactivity:	Human, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Harris J.D., DNA Seq. 4:361-393(1994). Gregory S.G., Nature 441:315-321(2006). The MGC Project Team; Genome Res. 14:2121-2127(2004).