

SPZ1 Antibody [#C11631]

Catalog Number:	C11631
Concentration:	1mg/ml
Swiss-Prot No.:	Q9BXG8
Other Names:	Spermatogenic leucine zipper protein 1, Testis-specific protein 1, Testis-specific protein NYD-TSP1, SPZ1, TSP1
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 SPZ1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SPZ1 antibody detects endogenous levels of total SPZ1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Sha J.-H., Asian J. Androl. 5:83-88(2003). Ota T., Nat. Genet. 36:40-45(2004). Schmutz J., Nature 431:268-274(2004).