

DMC1 Antibody [#C13043]

Catalog Number:	C13043
Concentration:	1mg/ml
Swiss-Prot No.:	Q14565
Other Names:	Meiotic recombination protein DMC1/LIM15 homolog; DMC1; DMC1H; LIM15
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 DMC1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	DMC1 antibody detects endogenous levels of total DMC1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 85%, Positives = 100%);
Applications:	IHC: 1:50~1:100 ELISA: 1:10000
References:	Sato S., DNA Res. 2:183-185(1995). Habu T., Nucleic Acids Res. 24:470-477(1996). Collins J.E., Genome Biol. 5:RESEARCH84.1-RESEARCH84.11(2004).