

HMG17 Antibody [#C10279]

Catalog Number:	C10279
Concentration:	1mg/ml
Swiss-Prot No.:	P05204
Other Names:	HMG17, high mobility group protein N2, high-mobility group (nonhistone chromosomal) protein 17, high-mobility group nucleosomal binding domain 2, high-mobility group nucleosome binding domain 2, nonhistone chromosomal protein HMG-17
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 HMG17.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	HMG17 antibody detects endogenous levels of total HMG17 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: 1:40000
References:	Landsman D., J. Biol. Chem. 261:7479-7484(1986). Gregory S.G., Nature 441:315-321(2006). Giancotti V., Eur. J. Biochem. 198:211-216(1991).