

HMGB1 Antibody [#C10316]

Catalog Number:	C10316
Concentration:	1mg/ml
Swiss-Prot No.:	P09429
Other Names:	HMG-1, HMG1, High mobility group protein 1, SBP-1, amphoterin, high-mobility group box 1, sulfoglucuronyl carbohydrate binding protein high-mobility group (nonhistone chromosomal) protein 1
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 HMGB1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	HMGB1 antibody detects endogenous levels of total HMGB1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Wen L., Nucleic Acids Res. 17:1197-1214(1989). Ferrari S., Genomics 35:367-371(1996). Rasmussen R.K., Electrophoresis 18:588-598(1997).