

ZNF148 Antibody [#C11914]

Catalog Number:	C11914
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UQR1
Other Names:	Beta enolase repressor factor 1, G-rich box-binding protein, Transcription factor BFCOL1, Transcription factor ZBP-89, ZBP89, ZFP148, ZN148, ZNF148, Zinc finger DNA binding protein 89, Zinc finger protein 148
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 ZNF148.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ZNF148 antibody detects endogenous levels of total ZNF148 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
References:	Beausoleil S.A., Proc. Natl. Acad. Sci. U.S.A. 101:12130-12135(2004). Boopathi E., J. Biol. Chem. 279:35242-35254(2004). Beausoleil S.A., Nat. Biotechnol. 24:1285-1292(2006).