

PKNOX2 Antibody [#C11510]

Catalog Number:	C11510
Concentration:	1mg/ml
Swiss-Prot No.:	Q96KN3
Other Names:	Homeobox protein PKNOX2, PBX/knotted homeobox 2, Homeobox protein PREP-2, PKNOX2, PREP2
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 PKNOX2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PKNOX2 antibody detects endogenous levels of total PKNOX2 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	The MGC Project Team, Genome Res. 14:2121-2127(2004). Imoto I., Biochem. Biophys. Res. Commun. 287:270-276(2001). Fognani C., Nucleic Acids Res. 30:2043-2051(2002).