

EZH1 Antibody [#C11435]

Catalog Number:	C11435
Concentration:	1mg/ml
Swiss-Prot No.:	Q92800
Other Names:	ENX-1, Enhancer of zeste homolog 2
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 EZH1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	EZH1 antibody detects endogenous levels of total EZH1 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
References:	Abel K.J., Genomics 37:161-171(1996). Ogawa M., Biochim. Biophys. Acta 1395:151-158(1998). Nagase T., DNA Res. 4:141-150(1997).