

DLX3 Antibody [#C10136]

Catalog Number:	C10136
Concentration:	1mg/ml
Swiss-Prot No.:	O60479
Other Names:	Homeobox protein DLX-3, DLX3
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 DLX3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	DLX3 antibody detects endogenous levels of total DLX3 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
References:	Price J.A., Hum. Mol. Genet. 7:563-569(1998). Sumiyama K., Proc. Natl. Acad. Sci. U.S.A. 99:780-785(2002). Dong J., Am. J. Med. Genet. A 133:138-141(2005).