

ALX3 Antibody [#C10213]

Catalog Number:	C10213
Concentration:	1mg/ml
Swiss-Prot No.:	O95076
Other Names:	Homeobox protein aristaless-like 3, Proline-rich transcription factor ALX3, ALX3
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 ALX3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ALX3 antibody detects endogenous levels of total ALX3 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
References:	Wimmer K., Genes Chromosomes Cancer 33:285-294(2002).