

ARMX3 Antibody [#C11839]

Catalog Number:	C11839
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UH62
Other Names:	ALEX3, DKFZp781N1954, KIAA0443, MGC12199, dJ545K15.2
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 ARMX3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ARMX3 antibody detects endogenous levels of total ARMX3 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	Kurochkin I.V., Biochem. Biophys. Res. Commun. 280:340-347(2001). Zhang Q.-H., Genome Res. 10:1546-1560(2000). Clark H.F., Genome Res. 13:2265-2270(2003).