

NKX6.3 Antibody [#C10009]

Catalog Number:	C10009
Concentration:	1mg/ml
Swiss-Prot No.:	A6NJ46
Other Names:	NKX63, Homeobox protein Nkx-6.3, NKX6-3
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 NKX6.3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NKX6.3 antibody detects endogenous levels of total NKX6.3 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
References:	Ensembl, Submitted (SEP-2007) to UniProtKB. Nusbaum C., Nature 439:331-335(2006). Mural R.J., Submitted (SEP-2005) to the EMBL/GenBank/DDBJ databases.