

ELAC2 Antibody [#C11594]

Catalog Number:	C11594
Concentration:	1mg/ml
Swiss-Prot No.:	Q9BQ52
Other Names:	Zinc phosphodiesterase ELAC protein 2, EC 3.1.26.11, Ribonuclease Z 2, RNase Z 2, tRNase Z 2, tRNA 3 endonuclease 2, ElaC homolog protein 2, Heredity prostate cancer protein 2, ELAC2, HPC2
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 ELAC2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ELAC2 antibody detects endogenous levels of total ELAC2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Tavtigian S.V., Nat. Genet. 27:172-180(2001). Ota T., Nat. Genet. 36:40-45(2004). Takaku H., Nucleic Acids Res. 31:2272-2278(2003).