

ELF5 Antibody [#C11875]

Catalog Number:	C11875
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UKW6
Other Names:	ETS-related transcription factor Elf-5, E74-like factor 5, Epithelium-specific Ets transcription factor 2, ESE-2, Epithelium-restricted ESE-1-related Ets factor, ELF5, ESE2
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 ELF5.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ELF5 antibody detects endogenous levels of total ELF5 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Taylor T.D., Nature 440:497-500(2006). Ota T., Nat. Genet. 36:40-45(2004). Oettgen P., J. Biol. Chem. 274:29439-29452(1999).