

ERF Antibody [#B0477]

Catalog Number:	B0477
Concentration:	1mg/ml
Swiss-Prot No.:	P50548
Other Names:	ETS-domain transcription factor ERF, Ets2 repressor factor
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 ERF.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ERF antibody detects endogenous levels of total ERF protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%);
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Sgouras D.N., EMBO J. 14:4781-4793(1995). Grimwood J., Nature 428:529-535(2004). Olsen J.V., Cell 127:635-648(2006).