

ETS1 (Ab-38) Antibody [#B0478]

Catalog Number:	B0478
Concentration:	1mg/ml
Swiss-Prot No.:	P27577
Other Names:	C-ets-1 protein, ETS-1, ETS1, p54
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 non-phosphopeptide derived from human ETS1 around the phosphorylation site of threonine 38 (L-L-T ^P -P-S).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ETS1 (Ab-38) antibody detects endogenous levels of total ETS1 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Tsukasa Higuchi, Mol. Cell. Biol., May 2007; 27: 3353 - 3366. H Suzuki, PNAS, May 1995; 92: 4442. V Fafeur, Cell Growth Differ., Jun 1997; 8: 655.