

CD3EAP Antibody [#C15463]

Catalog Number:	C15463
Concentration:	1mg/ml
Swiss-Prot No.:	O15446
Other Names:	antisense to ERCC-1; CAST; CD3-epsilon-associated protein; CD3e-associated protein; CD3EAP; DNA-directed RNA polymerase I RPA34; RNA polymerase I-associated factor PAF49; RPA34
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 CD3EAP.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CD3EAP antibody detects endogenous levels of total CD3EAP protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Whitehead C.M., Chromosoma 106:493-502(1997). Yamazaki T., J. Biol. Chem. 274:18173-18180(1999). The MGC Project Team; Genome Res. 14:2121-2127(2004).