

AN30A Antibody [#C11638]

Catalog Number:	C11638
Concentration:	1mg/ml
Swiss-Prot No.:	Q9BXX3
Other Names:	Serologically defined breast cancer antigen NY-BR-1, Ankyrin repeat domain-containing protein 30A, ANKRD30A
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 AN30A.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	AN30A antibody detects endogenous levels of total AN30A protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Jaeger D., Cancer Res. 61:2055-2061(2001). Deloukas P., Nature 429:375-381(2004). Hahn Y., Gene 366:238-245(2006).