

ERGI3 Antibody [#C11927]

Catalog Number:	C11927
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y282
Other Names:	ERGIC and golgi 3, ERGIC3, Erv46, NY-BR-84, Serologically defined breast cancer antigen NY-BR-84
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 ERGI3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ERGI3 antibody detects endogenous levels of total ERGI3 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
References:	Scanlan M.J., Cancer Immun. 1:4-4(2001). Lai C.-H., Genome Res. 10:703-713(2000). Ota T., Nat. Genet. 36:40-45(2004).