

SEPT1 Antibody [#C11420]

Catalog Number:	C11420
Concentration:	1mg/ml
Swiss-Prot No.:	Q8WYJ6
Other Names:	Septin-1, LARP, Serologically defined breast cancer antigen NY-BR-24, Peanut-like protein 3, SEPT1, DIFF6, PNUTL3
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 SEPT1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SEPT1 antibody detects endogenous levels of total SEPT1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Tu Q., Submitted (AUG-1998) to the EMBL/GenBank/DDBJ databases. The MGC Project Team, Genome Res. 14:2121-2127(2004). Scanlan M.J., Cancer Immun. 1:4-4(2001).