

SPIC Antibody [#C11346]

Catalog Number:	C11346
Concentration:	1mg/ml
Swiss-Prot No.:	Q8N5J4
Other Names:	Transcription factor Spi-C, SPIC
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 SPIC.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SPIC antibody detects endogenous levels of total SPIC protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Carlsson R., Gene 299:271-278(2002). The MGC Project Team, Genome Res. 14:2121-2127(2004).