

## TSPAN8 Antibody [#C10423]

<b>Catalog Number:</b>	C10423
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P19075
<b>Other Names:</b>	Tetraspanin-8, Tspan-8, Transmembrane 4 superfamily member 3, Tumor-associated antigen CO-029, TSPAN8, TM4SF3
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human TSPAN8.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	TSPAN8 antibody detects endogenous levels of total TSPAN8 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Szala S., Proc. Natl. Acad. Sci. U.S.A. 87:6833-6837(1990). The MGC Project Team, Genome Res. 14:2121-2127(2004).