

TACD1 Antibody [#C10390]

Catalog Number:	C10390
Concentration:	1mg/ml
Swiss-Prot No.:	P16422
Other Names:	Tumor-associated calcium signal transducer 1 [Precursor], Major gastrointestinal tumor-associated protein GA733-2, Epithelial cell surface antigen, Epithelial glycoprotein, EGP, Adenocarcinoma-associated antigen, KSA, KS 1/4 antigen, Cell surface glycoprotein Trop-1, CD326 antigen, TACSTD1, GA733-2, M1S2, M4S1, MIC18, TROP1
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 TACD1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TACD1 antibody detects endogenous levels of total TACD1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Strnad J., Cancer Res. 49:314-317(1989). Simon B., Proc. Natl. Acad. Sci. U.S.A. 87:2755-2759(1990). Perez M.S., J. Immunol. 142:3662-3667(1989).