

CD99 Antibody [#C12164]

Catalog Number:	C12164
Concentration:	1mg/ml
Swiss-Prot No.:	P14209
Other Names:	CD99 antigen, T-cell surface glycoprotein E2, E2 antigen, Protein MIC2, 12E7, MIC2, MIC2X, MIC2Y
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 CD99.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CD99 antibody detects endogenous levels of total CD99 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:5000
References:	Gelin C., EMBO J. 8:3253-3259(1989). The MGC Project Team, Genome Res. 14:2121-2127(2004). Darling S.M., Cold Spring Harb. Symp. Quant. Biol. 51:205-212(1986).