

CD147 Antibody [#C10559]

Catalog Number:	C10559
Concentration:	1mg/ml
Swiss-Prot No.:	P35613
Other Names:	CD147, EMMPRIN, M6, OK, TCSF, basigin, basigin (Ok blood group), collagenase stimulatory factor, extracellular matrix metalloproteinase inducer, leukocyte activation M6, tumor cell-derived collagenase stimulatory factor
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 C-terminal of human CD147.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CD147 antibody detects endogenous levels of total CD147 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
References:	Kasinrerk W., J. Immunol. 149:847-854(1992). Miyauchi T., J. Biochem. 110:770-774(1991). Biswas C., Cancer Res. 55:434-439(1995).