

CDH18 Antibody [#C12093]

Catalog Number:	C12093
Concentration:	1mg/ml
Swiss-Prot No.:	Q13634
Other Names:	Cadherin-18 [Precursor], Cadherin-14, CDH18, CDH14
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 CDH18.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CDH18 antibody detects endogenous levels of total CDH18 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Shibata T., J. Biol. Chem. 272:5236-5240(1997).