

Mucin-14 Antibody [#C11879]

Catalog Number:	C11879
Concentration:	1mg/ml
Swiss-Prot No.:	Q9ULC0
Other Names:	Endomucin [Precursor], Endomucin-2, Mucin-14, Gastric cancer antigen Ga34, EMCN, EMCN2, MUC14
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 Mucin-14.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Mucin-14 antibody detects endogenous levels of total Mucin-14 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:5000
References:	Kinoshita M., FEBS Lett. 499:121-126(2001). Kuhn A., J. Invest. Dermatol. 119:1388-1393(2002). The MGC Project Team, Genome Res. 14:2121-2127(2004).