

MED14 Antibody [#C10125]

Catalog Number:	C10125
Concentration:	1mg/ml
Swiss-Prot No.:	O60244
Other Names:	ARC150, Activator-recruited cofactor 150 kDa component, CRSP2, CSP2, Cofactor required for Sp1 transcriptional activation subunit 2, DRIP150, EXLM1, Thyroid hormone receptor-associated protein complex 170 kDa component, Transcriptional coactivator CRSP150, Vitamin D3 receptor-interacting protein complex 150 kDa component
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 MED14.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MED14 antibody detects endogenous levels of total MED14 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	Yoshikawa H., Genomics 49:237-246(1998). Ryu S., Nature 397:446-450(1999). Rachez C., Nature 398:824-828(1999).