

NCOA5 Antibody [#C17115]

Catalog Number:	C17115
Concentration:	1mg/ml
Swiss-Prot No.:	Q9HCD5
Other Names:	Nuclear receptor coactivator 5, NCoA-5; Coactivator independent of AF-2; CIA
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 NCOA5.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NCOA5 antibody detects endogenous levels of total NCOA5 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:10000
References:	Sauve F., Mol. Cell. Biol. 21:343-353(2001). Jiang C., J. Biol. Chem. 279:27781-27789(2004). Deloukas P., Nature 414:865-871(2001).