

NCoR1 Antibody [#C0358]

Catalog Number:	C0358
Concentration:	1mg/ml
Swiss-Prot No.:	O75376
Other Names:	KIAA1047, N-CoR1
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 NCoR1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	NCoR1 antibody detects endogenous levels of total NCoR1 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
References:	Igor Girault, Clin. Cancer Res., Apr 2003; 9: 1259. Xinjian Peng, Carcinogenesis, May 2007; 28: 1000 - 1007. Satoshi Mikawa, Genome Res., May 2007; 17: 586 - 593. Claire M. Banwell, Clin. Cancer Res., Apr 2006; 12: 2004 - 2013.