

## MED24 Antibody [#C10176]

<b>Catalog Number:</b>	C10176
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O75448
<b>Other Names:</b>	100 kDa thyroid hormone receptor associated protein, ARC100, Activator-recruited cofactor 100 kDa component, DRIP100, KIAA0130, THRAP4, TR100, Thyroid hormone receptor-associated protein complex 100 kDa component, Vitamin D3 receptor-interacting protein complex component DRIP100
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human MED24.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	MED24 antibody detects endogenous levels of total MED24 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Yuan C.-X., Proc. Natl. Acad. Sci. U.S.A. 95:7939-7944(1998). Rachez C., Genes Dev. 12:1787-1800(1998). Nagase T., DNA Res. 2:167-174(1995).