

## CIITA Antibody [#C10541]

<b>Catalog Number:</b>	C10541
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P33076
<b>Other Names:</b>	MHC class II transactivator, CIITA, CIITA, MHC2TA
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), 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 CIITA.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	CIITA antibody detects endogenous levels of total CIITA protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000   ELISA: 1:10000
<b>References:</b>	Steimle V., Cell 75:135-146(1993). Al-Kandari W., J. Mol. Biol. 369:1175-1187(2007). Al-Kandari W., Mol. Immunol. 44:311-321(2007). Croager E.J., Biochim. Biophys. Acta 1353:231-235(1997).