

## CDYL2 Antibody [#C15112]

<b>Catalog Number:</b>	C15112
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q8N8U2
<b>Other Names:</b>	Chromodomain Y-like protein 2; CDY-like 2
<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 CDYL2.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	CDYL2 antibody detects endogenous levels of total CDYL2 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 ELISA: 1:40000
<b>References:</b>	Ota T., Nat. Genet. 36:40-45(2004).