

## MED21 Antibody [#C10986]

<b>Catalog Number:</b>	C10986
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q13503
<b>Other Names:</b>	Mediator of RNA polymerase II transcription subunit 21, Mediator complex subunit 21, RNA polymerase II holoenzyme component SRB7, RNAPII complex component SRB7, hSrb7, MED21, SRB7, SURB7
<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 MED21.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	MED21 antibody detects endogenous levels of total MED21 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 ELISA: 1:20000
<b>References:</b>	Chao D.M., Nature 380:82-85(1996). Maldonado E., Nature 381:86-89(1996). Oelgeschlaeger T., Mol. Cell 1:925-931(1998).