

## ZNF638 Antibody [#C11052]

<b>Catalog Number:</b>	C11052
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q14966
<b>Other Names:</b>	Zinc finger protein 638, Nuclear protein 220, Zinc finger matrin-like protein, Cutaneous T-cell lymphoma-associated antigen se33-1, CTCL tumor antigen se33-1, ZNF638, NP220, ZFML
<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 ZNF638.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ZNF638 antibody detects endogenous levels of total ZNF638 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
<b>References:</b>	Inagaki H., J. Biol. Chem. 271:12525-12531(1996). Eichmueller S., Proc. Natl. Acad. Sci. U.S.A. 98:629-634(2001). Hillier L.W., Nature 434:724-731(2005).