

## MZF-1 Antibody [#C10502]

<b>Catalog Number:</b>	C10502
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P28698
<b>Other Names:</b>	Myeloid zinc finger 1, MZF-1, Zinc finger protein 42, Zinc finger and SCAN domain-containing protein 6, MZF1, MZF, ZNF42, ZSCAN6
<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 N-terminal of human MZF-1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	MZF-1 antibody detects endogenous levels of total MZF-1 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
<b>References:</b>	Hromas R., J. Biol. Chem. 266:14183-14187(1991). Peterson M.J., Gene 254:105-118(2000). Olsen J.V., Cell 127:635-648(2006).