

ZNF232 Antibody [#C11899]

Catalog Number:	C11899
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UNY5
Other Names:	Zinc finger protein 232, Zinc finger and SCAN domain-containing protein 11, ZNF232, ZSCAN11
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from internal of human ZNF232.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ZNF232 antibody detects endogenous levels of total ZNF232 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:5000
References:	Mavrogiannis L.A., Biochim. Biophys. Acta 1518:300-305(2001). Sjoebloom T., Science 314:268-274(2006). The MGC Project Team; Genome Res. 14:2121-2127(2004).