

ZNF287 Antibody [#C11719]

Catalog Number:	C11719
Concentration:	1mg/ml
Swiss-Prot No.:	Q9HBT7
Other Names:	Zinc finger protein with KRAB and SCAN domains 13, Zinc finger protein 287, ZNF287, ZKSCAN13
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 ZNF287.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ZNF287 antibody detects endogenous levels of total ZNF287 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
References:	Romero-Pastrana F., Submitted (DEC-1999) to the EMBL/GenBank/DDBJ databases. Ebert L., Submitted (JUN-2004) to the EMBL/GenBank/DDBJ databases. The MGC Project Team, Genome Res. 14:2121-2127(2004).