

ANKRD26 Antibody [#C14497]

Catalog Number:	C14497
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UPS8
Other Names:	ankyrin repeat domain 26; ANR26
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 ANKRD26.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ANKRD26 antibody detects endogenous levels of total ANKRD26 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Kikuno R., DNA Res. 6:197-205(1999). Deloukas P., Nature 429:375-381(2004). Ota T., Nat. Genet. 36:40-45(2004).