

A26C2/3 Antibody [#C17976]

Catalog Number:	C17976
Concentration:	1mg/ml
Swiss-Prot No.:	Q6S5H5/Q6S545
Other Names:	Prostate, ovary, testis-expressed protein on chromosome 14; POTE-14
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 A26C2/3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	A26C2/3 antibody detects endogenous levels of total A26C2/3 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Bera T.K., Gene 337:45-53(2004).