

POLE4 Antibody [#C15412]

Catalog Number:	C15412
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NR33
Other Names:	DNA polymerase epsilon subunit 4, EC 2.7.7.7; DNA polymerase II subunit 4; DNA polymerase epsilon subunit p12
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 POLE4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	POLE4 antibody detects endogenous levels of total POLE4 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Li Y., J. Biol. Chem. 275:23247-23252(2000). Hillier L.W., Nature 434:724-731(2005). Matsuoka S., Science 316:1160-1166(2007).