

## CPE Antibody [#C14961]

<b>Catalog Number:</b>	C14961
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P16870
<b>Other Names:</b>	carboxypeptidase e; carboxypeptidase e preproprotein; cbpe; cpe
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human CPE.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	CPE antibody detects endogenous levels of total CPE protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%) Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Manser E., Biochem. J. 267:517-525(1990). Utsunomiya N., Diabetologia 41:701-705(1998). Sjoebloom T., Science 314:268-274(2006).