

## BACE Antibody [#B0052]

<b>Catalog Number:</b>	B0052
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P56817
<b>Other Names:</b>	ASP2, Asp 2, Aspartyl protease 2, BACE1, Beta-secretase precursor, Beta-site APP cleaving enzyme, Beta-site amyloid precursor protein cleaving enzyme, EC 3.4.23.46, Memapsin-2, Membrane-associated aspartic protease 2
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), 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 BACE.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	BACE antibody detects endogenous levels of total BACE protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:100~1:500 ELISA: 1:5000
<b>References:</b>	Vassar R., Science 286:735-741(1999). Sinha S., Nature 402:537-540(1999). Yan R., Nature 402:533-537(1999).