

## Enos (Ab-494) Antibody [#B7072]

<b>Catalog Number:</b>	B7072
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P29474
<b>Other Names:</b>	cNOS; Constitutive NOS; EC 1.14.13.39; EC-NOS; ECNOS; Endothelial NOS; Nitric-oxide synthase, endothelial; NOS, type III; NOS3; NOSIII
<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 non-phosphopeptide derived from human Enos around the phosphorylation site of threonine 494 (K-K-T <sup>P</sup> -F-K).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Enos (Ab-494) antibody detects endogenous levels of total Enos 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>	Janssens S.P., J. Biol. Chem. 267:14519-14522(1992). Marsden P.A., FEBS Lett. 307:287-293(1992). Marsden P.A., J. Biol. Chem. 268:17478-17488(1993).