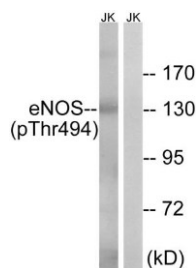


eNOS(Phospho-Thr494)Antibody [#8A7072]

Catalog Number:	8A7072
Concentration:	1mg/ml
Molecular Weight:	133 Kd
Origin of Antibody:	Rabbit
Swiss-Prot No.:	P29474
NCBI Gene Symbol:	NOS3
Other Names:	cNOS; Constitutive NOS; EC-NOS; ECNOS; Endothelial NOS; Nitric-oxide synthase, endothelial; NOS, type III; NOS3; NOSIII
All Sites:	H:T495 M:T494 R:T494
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 human eNOS around the phosphorylation site of Thr494.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
Specificity:	eNOS (Phospho-Thr494) Antibody detects endogenous levels of eNOS only when phosphorylated at Thr494.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Janssens S.P., J. Biol. Chem. 267:14519-14522(1992). Janssens S.P., J. Biol. Chem. 267:22694-22694(1992). Marsden P.A., FEBS Lett. 307:287-293(1992).

Application Images:

Figure1 (WB)



Western blot analysis of lysates from Jurkat cells, using eNOS (Phospho-Thr494) Antibody. The lane on the right is blocked with the phospho peptide