

## eNOS(Phospho-Ser1176)Antibody [#8A7071]

|                            |  |
|----------------------------|--|
| <b>Catalog Number:</b>     | 8A7071   |
| <b>Concentration:</b>      | 1mg/ml   |
| <b>Molecular Weight:</b>   | 133 Kd   |
| <b>Origin of Antibody:</b> | Rabbit   |
| <b>Swiss-Prot No.:</b>     | P29474   |
| <b>NCBI Gene Symbol:</b>   | NOS3   |
| <b>Other Names:</b>        | cNOS; Constitutive NOS; EC-NOS; ECNOS; Endothelial NOS; Nitric-oxide synthase, endothelial; NOS, type III; NOS3; NOSIII  |
| <b>All Sites:</b>          | H:S1177 M:S1176 R:S1176  |
| <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 human eNOS around the phosphorylation site of Ser1176.   |
| <b>Purification:</b>       | 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.                               |
| <b>Specificity:</b>        | eNOS (Phospho-Ser1176) Antibody detects endogenous levels of eNOS only when phosphorylated at Ser1176.   |
| <b>Reactivity:</b>         | Human, Mouse, Rat  |
| <b>Applications:</b>       | WB: 1:500~1:1000 ELISA: 1:20000  |
| <b>References:</b>         | Yong Chool Boo, Am J Physiol Heart Circ Physiol, Nov 2002; 283: 1819.<br>Christian Pott, Am J Physiol Cell Physiol, May 2006; 290: C1437 - C1445.<br>Monica Montagnani, J. Biol. Chem., Aug 2001; 276: 30392 - 30398.<br>Peter F. Mount, Am J Physiol Renal Physiol, Nov 2005; 289: F1103 - F1115. |

### Application Images:

**Figure1 (WB)**

