

eNOS (Ab-1176) Antibody [#B7071]

- Catalog Number:** B7071-1, B7071-2, B7071
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P29474
- All Names:** Constitutive NOS, EC 1.14.13.39, EC-NOS, ECNOS, Endothelial NOS, NOS, type III, NOS3, NOSIII, Nitric-oxide synthase, endothelial, cNOS
- 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.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human eNOS around the phosphorylation site of serine 1176(T-Q-S^P-F-S).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** eNOS (Ab-1176) antibody detects endogenous levels of eNOS protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 ELISA: 1:40000
- References:** Yong Chool Boo, Am J Physiol Heart Circ Physiol, Nov 2002; 283: 1819.
Christian Pott, Am J Physiol Cell Physiol, May 2006; 290: C1437 - C1445.
Monica Montagnani, J. Biol. Chem., Aug 2001; 276: 30392 - 30398.
Peter F. Mount, Am J Physiol Renal Physiol, Nov 2005; 289: F1103 - F1115.