

## iNOS (Ab-151) Antibody [#B8233]

<b>Catalog Number:</b>	B8233
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P35228
<b>Other Names:</b>	EC 1.14.13.39; HEP-NOS; hepatocyte NOS; inducible NOS; nitric oxide synthase, inducible; NOS type II; NOS2; NOS2A
<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 iNOS around the phosphorylation site of tyrosine 151 (Q-Y-Y <sup>P</sup> -G-S).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	iNOS (Ab-151) antibody detects endogenous levels of total iNOS protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%); Rat (Identities = 92%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Geller D.A., Proc. Natl. Acad. Sci. U.S.A. 90:3491-3495(1993). Sherman P.A., Biochemistry 32:11600-11605(1993). Charles I.G., Proc. Natl. Acad. Sci. U.S.A. 90:11419-11423(1993).