## NMDAR1 (Ab-890) Antibody [#B0692]

Catalog Number:B0692Concentration:1mg/mlSwiss-Prot No.:Q05586

Other Names: LURZ1, GRIN1, Glutamate [NMDA] receptor subunit zeta 1 precursor, N-methyl-D-aspartate

receptor, NMD-R1, NMDZ1, NMZ1, NR1

Storage/Stability: Store at -20°C/1 year

Form of Antibody: 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.

**Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human

NMDAR1 around the phosphorylation site of serine 890 (A-S-S<sup>P</sup>-F-K).

**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-

specific immunogen.

**Specificity:** NMDAR1 (Ab-890) antibody detects endogenous levels of total NMDAR1 protein.

Reactivity: Human, Mouse, Rat

**Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000

**References:** Guang Bai, J. Biol. Chem., Jan 1998; 273: 1086.

Guang Bai, J. Biol. Chem., Mar 1995; 270: 7737. Mai T. Dang, PNAS, Oct 2006; 103: 15254 - 15259.

Technical Support: <a href="mailto:support@aatbio.com">support@aatbio.com</a>; 408-733-1055