

## IRS-1 (Ab-636/639) Antibody [#B7122]

- Catalog Number:** B7122-1, B7122-2, B7122
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P35568
- All Names:** IRS-1, IRS1, Insulin receptor substrate-1
- 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.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human IRS-1 around the phosphorylation site of serine 636 or serine 639 (M-S-P-K-S).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** IRS-1 (Ab-636/639) antibody detects endogenous levels of total IRS-1 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** IHC: 1:50~1:100 ELISA: 1:20000
- References:** Ozes ON, et al. (2001) Proc Natl Acad Sci USA; 98(8): 4640-4645.  
Tzatsos A, et al. (2006) Mol Cell Biol; 26(1): 63-76.  
Kadowaki T, et al. (2000) J Clin Invest; 106(4): 459-465.  
Ozes ON, et al. (2001) Proc Natl Acad Sci USA; 98(8): 4640-4645.  
Szanto I, et al. (2000) Proc Natl Acad Sci USA; 97(5): 2355-2360.