

IRS-1 Antibody [#B0668]

Catalog Number:	B0668
Concentration:	1mg/ml
Swiss-Prot No.:	P35568
Other Names:	IRS-1, IRS1, Insulin receptor substrate-1
Storage/Stability:	Store at -20°C/1 year
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.
Immunogen:	The antiserum was produced against synthesized peptide derived from Internal of human IRS-1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	IRS-1 antibody detects endogenous levels of total IRS-1protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:5000
References:	Joëlle Dupont, Am J Physiol Endocrinol Metab, Feb 1998; 274: E309 - E316. DG Baskin, Endocrinology, Apr 1994; 134: 1952. Zhefu Ma, Mol. Cell. Biol., Dec 2006; 26: 9338 - 9351. Alexandros Tzatsos, J. Biol. Chem., Jun 2007; 282: 18069 - 18082.