

IR Antibody [#B0492]

Catalog Number:	B0492
Concentration:	1mg/ml
Swiss-Prot No.:	P06213
Other Names:	CD220 antigen, EC 2.7.10.1, IR, insulin receptor, kinase InsR
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 IR.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	IR antibody detects endogenous levels of total IR protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:5000
References:	Ebina Y., Cell 40:747-758(1985). Ullrich A., Nature 313:756-761(1985). Seino S., Diabetes 39:123-128(1990).