



Peroxidase-Conjugated Rabbit anti-Mouse IgG (H+L)

Catalog Number:	28165
Immunogen:	Purified mouse IgG
Host:	Rabbit
Size:	0.5mg
Concentration:	1mg/mL
Degree of Substitution: (DOS)	HRP/protein mole ratio is labeled on each vial.
Specificity:	Affinity purified rabbit anti-mouse IgG (H+L) specific antibody is conjugated with Horseradish Peroxidase (HRP). The antibody reacts with the heavy chains on mouse IgG and with light chains common to most mouse immunoglobulins. It has minimum cross-reaction to human serum proteins.
Application:	<p>This antibody can be used for ELISA, Western blot and IHC. The investigator should determine the working concentrations for specific applications. The results might be affected by several factors, including antigen concentration, sensitivity of detection method, temperature and length of incubations, etc. The following concentration ranges are recommended starting points for this product.</p> <p>ELISA: 1:5,000-1:100,000 Western blot: 1:10,000-1:200,000 with ECL substrates. IHC: 1:500-1:5,000</p>
Storage buffer:	10 mM phosphate, 150 mM NaCl, pH 7.6, with 0.5% Proclin 300 as preservative.
Storage:	The peroxidase conjugate is stable for 1 year at 4°C.

For *in vitro* research use only