



Product Information Sheet

Product Name:	Goat Anti-Mouse IgG (H+L), Alkaline Phosphatase Conjugated
Catalog Number:	28174
Size:	0.5 mg
Concentration:	1mg/ml
Storage buffer:	10 mM Tris-HCl, 250 mM NaCl, pH 8.0, with 2 mM sodium azide as preservative and BSA as stabilizer.
Storage:	The product is stable for six months at 4°C.
Specification:	The antibody was immunoaffinity purified from the serum of goat immunized with purified mouse IgG, and then conjugated with alkaline phosphatase. The recommended dilution range is 1:1000 –1:10,000 for ELISA and Western blot and 1:200 – 1:1000 for immunohistochemistry. The optimal working concentration should be determined experimentally. The antibody conjugate should be diluted in Tris-buffered saline with BSA or equivalent as stabilizer. Phosphate-buffered saline should not be used to dilute alkaline phosphatase-antibody conjugate.

For *in vitro* research use only