



Product Information Sheet

Product Name:	Goat Anti-Mouse IgG (H+L), Biotinylated
Catalog Number:	28175-05
Size:	0.5 mg
Concentration:	1mg/mL
Degree of Substitution: (DOS)	Biotin/protein mole ratio is labeled on each vial.
Buffer:	10 mM phosphate, 150 mM NaCl, pH 7.6, BSA and 2 mM sodium azide.
Storage:	The biotinylated IgG is stable for 2~3 months at 4°C. For long-term storage, divide the solution into aliquots and store at -20°C or add an equal volume of glycerol (ACS grade or higher) and store the solution at -20°C without aliquoting. Avoid multiple thaw-freeze cycles. The product is stable for 1 year at -20°C.
Instruction:	The recommended concentration for Western blot is 1 µg/mL, and for immunoprecipitation and immunostaining is 5-10 µg/mL.

For *in vitro* research use only