



Product Information Sheet

Product Name:	Rabbit Anti-Mouse IgG (H+L), AMCA-labeled
Catalog Number:	28164-1-AMCA
Size:	1 mg
Concentration:	1 mg/mL
Degree of Substitution: (DOS)	Dye/protein mole ratio is labeled on each vial.
Fluorescence:	Blue fluorescence. Excitation/Emission wavelength= 344 nm/ 446 nm
Storage buffer:	10 mM phosphate, 150 mM NaCl, pH 7.2, with BSA and 2 mM sodium azide as preservatives.
Storage:	The dye-protein conjugate 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.
Instructions:	The recommended concentration for most immunofluorescent staining is 1-10 µg/mL. If the non-specific binding background is high, you may decrease the concentration to minimize the background. You may also centrifuge the conjugate briefly and use the supernatant only for staining. This Ig G has minimal cross-reaction to human serum proteins.

For *in vitro* research use only