



Product Information Sheet

Product Name:	Rabbit Anti-mouse IgG (H+L), Hilyte Fluor™ 750-labeled
Catalog Number:	28164-1-H750
Size:	1 mg
Concentration:	1mg/mL
Degree of Substitution: (DOS)	Dye:protein molar ratio is specified on the vial.
Fluorescence:	Excitation/Emission wavelength = 754 nm/778 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:	<p>HiLyte Fluor™ 750-IgG conjugate has been optimized in fluorophore/protein labeling ratio to ensure high fluorescent signal and uncompromised IgG function. Spectrally similar to Cy7 dye, HiLyte Fluor™ 750 is the longest-wavelength HiLyte Fluor™ dye currently available. Its fluorescence emission maximum at 778 nm is well separated from commonly used far-red fluorophores such as HiLyte Fluor™ 647, HiLyte Fluor™ 680 or allophycocyanin (APC), facilitating multicolor analysis. With a peak excitation at ~754 nm, conjugates of HiLyte Fluor™ 750 dyes are well excited by a xenon-arc lamp or dye-pumped lasers operating in the 720–750 nm range.</p> <p>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.</p>

For *in vitro* research use only