



Product Information Sheet

Product Name:	Rabbit Anti-chicken IgY (H+L), HiLyte Fluor™ 680-labeled
Catalog Number:	29709-1-H680
Size:	1 mg
Concentration:	1mg/mL
Degree of Substitution: (DOS)	Dye:protein molar ratio is specified on the vial.
Fluorescence:	Excitation/Emission wavelength= 678 nm/699 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™ 680-IgY conjugate has been optimized in fluorophore/protein labeling ratio to ensure high fluorescent signal and uncompromised IgY function. The fluorescence of HiLyte Fluor™ 680 is only slightly red-shifted compared to those of Cy5.5 dyes, resulting in an optimal match to filters designed for Cy5.5 dyes. Fluorescence emission of HiLyte Fluor™ 680 is well separated from that of other commonly used red fluorophores, such as TAMRA, R-phycoerythrin and HiLyte Fluor™ 647 dyes, making it ideal for three- and four-color labeling.</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