



## Product Information Sheet

---

Product Name:	Goat Anti-rabbit IgG (H+L), highly cross-adsorbed, HiLyte Fluor™ 680-labeled
Catalog Number:	61056-05-H680
Size:	0.5 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-IgG conjugate has been optimized in fluorophore/protein labeling ratio to ensure high fluorescent signal and uncompromised IgG 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. This IgG is highly adsorbed with minimal cross-reaction to human, mouse, and rat serum protein.</p>

For *in vitro* research use only