



Product Information Sheet

Product Name:	Rabbit Anti-Chicken IgY (H+L), Hilyte Fluor™ 405-labeled
Catalog Number:	29709-05-H405
Size:	0.5 mg
Concentration:	1 mg/mL
Degree of Substitution: (DOS)	Dye/protein molar ratio is written on each vial.
Fluorescence:	Blue fluorescence. Excitation/Emission wavelength= 404 nm/428 nm
Storage buffer:	10 mM phosphate, 150 mM NaCl, pH 7.2, with BSA and 2 mM sodium azide as preservative.
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 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.

For *in vitro* research use only