



Product Information Sheet

Product Name:	Rabbit Polyclonal Anti-His Tag IgG, HiLyte Fluor™ 555-labeled
Catalog Number:	29673-H555
Size:	0.1 mg
Concentration:	1mg/ml
Degree of Substitution: (DOS)	Dye/protein mole ratio is labeled on each vial.
Fluorescence:	Orange fluorescence. Excitation/Emission wavelength= 553 nm/568 nm
Storage buffer:	10 mM phosphate, 150 mM NaCl, pH 7.2, with BSA and 2 mM sodium azide.
Storage:	The dye-IgG 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 rabbit polyclonal anti-His tag IgG is raised against 6x His synthetic peptide. It recognizes 6x His tag attached to both amino and carboxy terminus of the proteins being expressed in the cells. The recommended concentration is 0.5-5 µg/mL for immunostaining, immunoprecipitation, and other applications.

For *in vitro* research use only