



Product Information Sheet

Product Name:	Anti-Actin (Smooth Muscle Specific), HiLyte Fluor™ 488-labeled
Catalog Number:	29553-H488
Size:	0.05 mg
Concentration:	1mg/ml
Degree of Substitution: (DOS)	Dye/protein mole ratio is labeled on each vial.
Fluorescence:	Green fluorescence. Excitation/Emission wavelength= 499 nm/523 nm. The fluorescence of HiLyte Fluor™ 488 is brighter and more resistant to photobleaching compared to other green fluorophores, such as FITC.
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 anti-actin polyclonal antibody was raised against a synthetic peptide corresponding to N-terminus of human α -smooth muscle actin. This peptide sequence is identical in rat, mouse, chicken, bovine, etc. The other actin isoforms, such as α - and γ -cytoplasmic, α -skeletal, α -cardiac actins, are not stained by this antibody. The recommended concentration is 1-10 μ g/mL for immunostaining.

For *in vitro* research use only