



HiLyte Fluor™ Labeled Streptavidin Sampler Kit

Catalog #	AS-72003-20
Unit Size	1 Kit
Kit Size	3 x 20 µg HiLyte Fluor™ Dye-Streptavidin Conjugates

This kit provides three different HiLyte Fluor™ fluorophore labeled streptavidin conjugates, which can be used in optimizing immunofluorescence staining.

INTRODUCTION

Fluorescent dye conjugated streptavidins have been widely used in immunofluorescence staining, fluorescence activated cell sorting, *in situ* hybridization and other fluorescence-based biological applications.

The HiLyte Fluor™ fluorophores are premium quality dyes that have brighter fluorescence and better photostability compared to FITC, TAMRA and Cy® dyes. Prepared at optimal fluorophore to streptavidin labeling ratio, these HiLyte Fluor™ streptavidin conjugates exhibit high fluorescence signal and uncompromised biotin binding capability. The actual dye/streptavidin ratio or degree of substitution (DOS) is labeled on each vial. Every lot has been validated with immunofluorescence staining to guarantee the best performance.

All fluorescent streptavidin conjugates are also available for purchase individually.

KIT COMPONENTS

Catalog#	Fluorescent Streptavidin Conjugate	Fluorescence	Ex/Em (nm)
60665	Streptavidin HiLyte Fluor™ 488 conjugated	Green	499/523
60666	Streptavidin HiLyte Fluor™ 555 conjugated	Orange	553/568
60667	Streptavidin HiLyte Fluor™ 647 conjugated	Red	653/673

STORAGE AND HANDLING

All fluorescent streptavidin conjugates are supplied as 0.5 mg/mL in 10 mM phosphate, 150 mM NaCl, pH 7.2, with bovine serum albumin and 2 mM sodium azide.

Conjugates are stable for 2~3 months at 4°C. For long-term storage, add an equal volume of glycerol (ACS grade or higher) and store at -20°C.

USE AND INSTRUCTIONS

The recommended concentration for most immunofluorescent staining is 1-10 µg/mL (1:500-1:50 dilution of stock solution). A good starting point is 10 µg/mL. If the non-specific binding background is high, you may decrease the concentration. You may also centrifuge the conjugate briefly and use the supernatant for staining.

AnaSpec, Inc. • 34801 Campus Dr • Fremont, CA 94555

Tel: 1-800-452-5530 • Fax 510-791-9572 • service@anaspec.com • www.anaspec.com