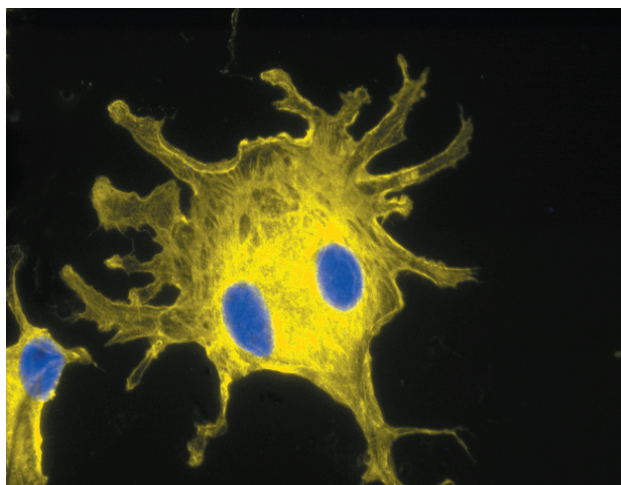


Streptavidin & Conjugates

Brilliant streptavidin conjugates for your research



Actins in bovine pulmonary artery endothelial cells were detected with biotin-conjugated phalloidin, visualized with R-phycoerythrin-conjugated streptavidin (Cat#60669). The nuclei were stained with Hoechst-33342 (Cat# 83218)

- Guaranteed quality and performance
- High purity Streptavidin (Cat#60659)
- Conjugated with AnaSpec's proprietary HiLyte Fluor™ dyes, enzymes and fluorescent proteins
- Superior fluorescence intensity and photostability compared with FITC and Cy dyes
- Application: Flow cytometry, Immunofluorescence, *in situ* Hybridization and Microarrays
- Competitive pricing

HiLyte Fluor™ Labeled Streptavidin

Product	HiLyte Fluor 488	HiLyte Fluor 555	HiLyte Fluor TR	HiLyte Fluor 594	HiLyte Fluor 647	HiLyte Fluor 680	HiLyte Fluor 750
Streptavidin	60665	60666	60671	60672-H594	60667	60659-H680	60659-H750

Other Streptavidin Conjugates

Product	FITC	5-FAM	5-TAMRA	B-PE	R-PE	Crosslinked Allophycocyanin	AP	HRP
Streptavidin	60659-FITC	60664	60670	60662	60669	60663	60660	60668

HiLyte Fluor™ Labeled Streptavidin Sampler Kits

These sampler kits contain 3 different HiLyte Fluor™ labeled streptavidin - HiLyte Fluor™ 488, 555, and 647.

Size	HiLyte Fluor Labeled Streptavidin Sampler Kits
3x20µg	72003-20
3x200µg	72003-200

Custom Service Available
Antibody Labeling
Assay Development

