

Product Information

Streptavidin Conjugates

Cat. No.	Product Name	Unit Size
L155A	Andy Fluor 350 Streptavidin	100 µg
L155B	Andy Fluor 350 Streptavidin	500 µg
L156A	Andy Fluor 488 Streptavidin	100 µg
L156B	Andy Fluor 488 Streptavidin	500 µg
L157A	Andy Fluor 555 Streptavidin	100 µg
L157B	Andy Fluor 555 Streptavidin	500 µg
L158A	Andy Fluor 594 Streptavidin	100 µg
L158B	Andy Fluor 594 Streptavidin	500 µg
L159A	Andy Fluor 647 Streptavidin	100 µg
L159B	Andy Fluor 647 Streptavidin	500 µg
L160A	Cy3 Streptavidin	100 µg
L160B	Cy3 Streptavidin	500 µg
L161A	Cy5 Streptavidin	100 µg
L161B	Cy5 Streptavidin	500 µg

Spectral Properties:

Product Name	Ex nm	Em nm
Andy Fluor 350 Streptavidin	345	440
Andy Fluor 488 Streptavidin	495	520
Andy Fluor 555 Streptavidin	553	565
Andy Fluor 594 Streptavidin	590	615
Andy Fluor 647 Streptavidin	650	665
Cy3 Streptavidin	550	565
Cy5 Streptavidin	650	667

Storage upon receipt:

- -20 °C
- Protect from light
- Avoid freeze-thaw cycles

Product Description

Applied BioProbes offers a variety of streptavidin conjugates with our outstanding series of Andy Fluor™ dyes. Andy Fluor™ dyes are superior to other fluorescent dyes for protein labeling by having advantages in brightness, photostability, specificity and novel features ideal for in vivo imaging.

Streptavidin conjugates are commonly used as secondary reagent to localize antigens in cells and tissues, and to detect biomolecules in immunoassays and DNA hybridization techniques.

Product Specification

Physical State: Lyophilized power

Buffer: PBS, pH 7.4

Stabilizer: 0.1% BSA

Preservative: 0.02% Sodium Azide

Reconstitution and Storage: Store lyophilized power at 2-8°C. When ready to use, rehydrate with dH₂O (50 µL for 100 µg antibody or 250 µL for 500 µg antibody) to make 2 mg/mL solution and centrifuge if not clear. Product is stable for about 6 months at 2-8°C as an undiluted liquid. Prepare working dilution fresh each day. For extended storage after rehydration, add an equal volume of glycerol for a final concentration of 50%, and store at -20 °C as a liquid.

Guidelines for Use

Streptavidin conjugates are used as secondary detection reagents to detect biotinylated probes in histochemical application, flow cytometry, blot analysis, and immunoassays. The following are commonly used methods for employing streptavidin conjugates as a secondary detection reagent.

A biotinylated primary probe such as an antibody, single-stranded nucleic acid probe, or lectin is bound to tissues, cells, or other surfaces. Excess protein is removed by washing, and detection is mediated by streptavidin conjugate.

Centrifuge the protein conjugate solution briefly in a microcentrifuge before use. Add only the supernatant to the experiment. This step eliminates any protein aggregates that may have formed during storage, thereby reducing nonspecific background staining.

Because staining protocols vary with application, determine appropriate dilutions of streptavidin conjugates empirically. A final concentration of 1–10 µg/mL should be satisfactory for most histochemical applications.