



Glowing products for science

Streptavidin Conjugates

A high quality biotin-binding protein conjugated to Biotium's signature bright and photostable fluorescent CF® Dyes, and a selection of other labels.



Product Description

CF® Dyes conjugates of Streptavidin are high quality biotin-binding proteins labeled with the superior CF® Dyes or a selection of other labels. These conjugates are typically used as secondary reagents to detect biotinylated probes such as primary antibodies for flow cytometry, western blotting, immunofluorescence staining, and other applications.

- Conjugates with wide choice of CF® Dye colors, plus HRP, R-PE, APC, PerCP, and Alkaline Phosphatase
- Superior CF® Dyes are bright, photostable, and water-soluble
- Ideal for flow cytometry, western blotting, and immunofluorescence staining

Blue dyes have lower fluorescence and can give higher non-specific background than other dye colors. Conjugates of blue fluorescent dyes like CF®350, CF®405S, and CF®405M are not recommended for detecting low abundance targets.

Superior CF® Dyes

Biotium's next-generation CF® Dyes were designed to be highly water-soluble with advantages in brightness and photostability compared to Alexa Fluor®, DyLight®, and other fluorescent dyes. Learn more about [CF® Dyes](#).

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Colors	Blue, Green, Orange, Red, Far-red, Near-infrared
Detection method/readout	Fluorescence microscopy, Live cell imaging, Flow cytometry, Near-IR imager, In vivo near-IR imaging
Antibody reactivity (target)	Biotin
Antibody/conjugate formulation	AP conjugates: 1 mg/mL in 50 mM Tris buffer (pH 8.0) with 1 mM MgCl ₂ , 50% glycerol, and 0.05% sodium azide, APC Conjugate: 0.5 mg/mL in PBS, 50% glycerol, 2 mg/mL BSA (IgG-free and protease-free), and 0.05% sodium azide, CF® Dye Conjugates: 2 mg/mL in 1X PBS with 0.01% sodium azide after reconstitution, HRP conjugate: 1 mg/mL PBS/10 mg/mL trehalose after reconstitution, PerCP Conjugate: 0.5 mg/mL in PBS, 2 mg/mL BSA (IgG-free and protease-free), and 0.05% sodium azide, R-PE Conjugate: 0.5 mg/mL in PBS, 2 mg/mL BSA (IgG-free and protease-free), and 0.05% sodium azide
Product origin	Alkaline Phosphatase: Recombinant, E. coli, HRP: Horseradish (Armoracia rusticana), Streptavidin, recombinant from E. coli, APC from algae, BSA from bovine serum (Bos taurus) or recombinant BSA produced in Chinese hamster ovary cells., PerCP from algae

Streptavidin Conjugates

Product	Conjugation	Ex/Em	Size	Catalog number
CF®350 Streptavidin	CF®350	347/448 nm	1 mg	29031
CF®405S Streptavidin	CF®405S	404/431 nm	1 mg	29032
CF®405M Streptavidin	CF®405M	408/452 nm	1 mg	29033
CF®405L Streptavidin	CF®405L	395/545 nm	1 mg	29056
CF®430 Streptavidin	CF®430	426/498 nm	1 mg	29065
CF®440 Streptavidin	CF®440	440/515 nm	1 mg	29066
CF®488A Streptavidin	CF®488A	490/515 nm	1 mg	29034
CF®514 Streptavidin	CF®514	516/548 nm	1 mg	29081
CF®532 Streptavidin	CF®532	527/558 nm	1 mg	29030
CF®543 Streptavidin	CF®543	541/560 nm	1 mg	29043
CF®555 Streptavidin	CF®555	555/565 nm	1 mg	29038
CF®568 Streptavidin	CF®568	562/583 nm	1 mg	29035
CF®583R Streptavidin	CF®583R	586/609 nm	1 mg	29086
CF®594 Streptavidin	CF®594	593/614 nm	1 mg	29036
CF®633 Streptavidin	CF®633	630/650 nm	1 mg	29037
CF®640R Streptavidin	CF®640R	642/662 nm	1 mg	29041
CF®647 Streptavidin	CF®647	650/665 nm	1 mg	29039
CF®660R Streptavidin	CF®660R	663/682 nm	1 mg	29040
CF®680R Streptavidin	CF®680R	680/701 nm	1 mg	29072
CF®740 Streptavidin	CF®740	742/767 nm	1 mg	29129
Alkaline Phosphatase Streptavidin (1 mg/mL)	Alkaline Phosphatase	N/A	100 uL	29071-100uL
1 mL	29071-1mL			
HRP Streptavidin	HRP	N/A	1 mg	29049
R-PE Streptavidin (0.5 mg/mL)	R-PE	496, 546, 565/578 nm	200 uL	29044-200uL
1 mL	29044-1mL			
APC Streptavidin (0.5 mg/mL)	APC	650/660 nm	200 uL	29048-200uL
1 mL	29048-1mL			
PerCP Streptavidin (0.5 mg/mL)	PerCP	482/675 nm	200 uL	29140-200uL
1 mL	29140-1mL			

Storage and Handling

Liquid formats:

Streptavidin, R-PE Conjugate and PerCP Conjugate: Store at 4°C, protected from light. Storage at -20°C is not recommended for R-PE or PerCP conjugates. Product is stable for at least 6 months from date of receipt when stored as recommended.

Streptavidin, APC Conjugate: Store at -20°C, protected from light. Product is stable for at least 6 months from date of receipt when stored as recommended.

Streptavidin, Alkaline Phosphatase Conjugate: Store at -20°C. Product is stable for at least 6 months from date of receipt when stored as recommended.

Lyophilized format:

CF® Dye Streptavidin Conjugates: Product is stable for at least 2 years at -20°C with desiccant. Add 0.5 mL of PBS to 1 mg lyophilized CF® Dye Streptavidin Conjugate and mix gently to dissolve.

Streptavidin, HRP Conjugate: add 1 mL pH ~7.4 1X PBS and mix gently to dissolve. Reconstituted HRP conjugate can be stored in aliquots at -20°C for at least 6 months. Do not add azide to HRP conjugates.

CF is a registered trademark of Biotium

References

Download a list of [CF® dye references](#).

This datasheet was generated on February 3, 2026 at 07:23:32 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/streptavidin-conjugates/>