

## 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 attributes

<b>Colors</b>	Blue, Green, Orange, Red, Far-red, Near-infrared
<b>Detection method/readout</b>	Fluorescence microscopy, Live cell imaging, Flow cytometry, Near-IR imager, In vivo near-IR imaging
<b>Antibody reactivity (target)</b>	Biotin
<b>Product origin</b>	Streptavidin, recombinant from E. coli

## 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, 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](#).

## Streptavidin Conjugates

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

CF is a registered trademark of Biotium, Inc. Alexa Fluor, Texas Red, and DyLight are registered trademarks of Thermo Fisher Scientific.

## References

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