

Revised: July 20, 2020



CF® Dye Streptavidin Conjugates

Glowing Products for Science"

Catalog no.	Conjugate	Ex/Em (nm)
29031	CF®350	347/448
29032	CF®405S	404/431
29033	CF®405M	408/452
29056	CF®405L	395/545
29065	CF®430	426/498
29066	CF®440	490/515
29034	CF®488A	490/515
29081	CF®514	516/548
29030	CF®532	527/558
29043	CF®543	541/560
29038	CF®555	555/565
29035	CF®568	562/583
29086	CF®583R	586/609
29036	CF®594	593/614
29037	CF®633	630/651
29041	CF®640R	642/662
29039	CF®647	650/665
29040	CF®660R	663/682
29072	CF®680R	680/701

Unit Size: 1 mg

Concentration: 2 mg/mL with 0.01% sodium azide upon addition of 0.5 mL PBS.

Form: Lyophilized powder

Storage and Handling

Product is stable for at least 2 years at -20°C with desiccant. Upon reconstitution of the lyophilized powder in 0.5 mL PBS, store at 4°C and protect from light.

Product Description

Biotium offers a variety of streptavidin products conjugated to our outstanding series of CF® dyes and other labels. CF® dyes are superior to other fluorescent dyes for protein labeling by having combined advantages in brightness, photostability, specificity and novel features ideal for *in vivo* imaging.

Streptavidin 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. For most fluorescent streptavidin applications, a concentration of 1 - 2 ug/mL is sufficient; however, optimal conditions should be determined empirically.

A full selection of CF® dye labeled products including secondary antibodies, antibody labeling kits, and other bioconjugates such as phalloidins, annexin V and α -bungarotoxin are also available for many CF® dyes.

Related Products

Catalan		
Catalog number	Product	
29049	HRP Streptavidin, 1 mg	
29071	Alkaline Phosphatase Streptavidin	
29048	APC Streptavidin	
29044	R-PE Streptavidin	
40061	RedDot™2 Far Red Nuclear Counterstain, 200X in DMSO	
23001	EverBrite™ Mounting Medium, 10 mL	
23002	EverBrite™ Mounting Medium with DAPI, 10 mL	
23003	EverBrite™ Hardset Mounting Medium, 10 mL	
23004	EverBrite™ Hardset Mounting Medium with DAPI, 10 mL	
23005	CoverGrip™ Coverslip Sealant, 15 mL	
22005	Mini Super ^н Pap Pen 2.5 mm tip, ~400 uses	
22006	Super ^{н⊤} Pap Pen 4 mm tip, ~800 uses	
22015	Fixation Buffer, 100 mL	
22016	Permeabilization Buffer, 100 mL	
22017	Permeabilization and Blocking Buffer, 100 mL	
22010	10% Fish Gelatin Blocking Buffer, 100 mL	
22011	Fish Gelatin Powder, 2 X 50 g	
22014	30% Bovine Serum Albumin Solution, 100 mL	
22002	Tween®-20, 50 mL	

Please visit www.biotium.com to view our full selection of CF® dye and R-PE conjugates, including secondary antibodies, phalloidin, Annexin V, and α-bungarotoxin, as well as Mix-n-Stain™ antibody labeling kits. Biotium also offers a variety of apoptosis and cell viability assays for flow cytometry analysis, including mitochondrial membrane potential dyes and NucView® 488 Caspase-3 Substrate for live cells.

CF dye technology is covered by pending US and international patents.