

Product Information

Streptavidin, APC Conjugates

Cat. No.	Product Description	Unit Size
29048-1mL	Streptavidin, Allophycocyanin crosslinked conjugate	1 mL
29048-200uL	Streptavidin, Allophycocyanin crosslinked conjugate	200 uL

Concentration: 0.5 mg/mL in pH ~7.4 PBS containing 50% glycerol, 2 mg/ml bovine serum albumin (IgG-free and protease-free) and 0.05% sodium azide.

Form: Blue solution

Spectral Property

$\lambda_{abs}/\lambda_{em}$ (in pH 7.4 PBS buffer): 633, 640 nm/660 nm

Storage and Handling

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

Product Description

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 - 10 µg/mL is sufficient; however, optimal conditions should be determined empirically.

The products are prepared by labeling streptavidin with allophycocyanine (APC). APC belongs to the family of phycobiliproteins, which are derived from cyanobacteria and eukaryotic algae. APC has absorption maxima at 650 and emission at 660 nm. Because of the high fluorescence quantum yield (~0.68), APC is by far the brightest dye for the 650 nm wavelength region.

Other APC products

Cat. No.	Product Name
20411	Goat Anti-Mouse IgG (H+L)
20412	Goat Anti-Rabbit IgG (H+L)
20413	Goat Anti-Rat IgG (H+L), highly cross-adsorbed Min X Hu, Bv, Hs and Rb
20414	F(ab') ₂ fragment of Goat anti-Mouse IgG (H+L)
20415	F(ab') ₂ fragment Goat Anti-Rabbit IgG (H+L)

Bv: bovine; Ch: chicken; Gt: goat; GP: Guinea pig; Hs: horse; Hu: human; Ms: mouse; Rb: rabbit; Sh: sheep; SHm: Syrian hamster; Sw: swine; Rt: rat

Related Products

CF® dye streptavidin conjugates

Cat. No.	Product Description	Unit Size
29031	Streptavidin, CF®350 Conjugate	1 mg
29032	Streptavidin, CF®405S Conjugate	1 mg
29033	Streptavidin, CF®405M Conjugate	1 mg
29034	Streptavidin, CF®488A Conjugate	1 mg
29038	Streptavidin, CF®555 Conjugate	1 mg
29035	Streptavidin, CF®568 Conjugate	1 mg
29036	Streptavidin, CF®594 Conjugate	1 mg
29037	Streptavidin, CF®633 Conjugate	1 mg
29039	Streptavidin, CF®647 Conjugate	1 mg
29040	Streptavidin, CF®660R Conjugate	1 mg

CF®dye conjugated mouse monoclonal anti-biotin antibody

Cat. No.	Product Description	Unit Size
20203	CF®405S mouse monoclonal anti-biotin, 2 mg/mL	0.25 mL
20203-1	CF®405S mouse monoclonal anti-biotin, 2 mg/mL	50 uL
20204	CF®488A mouse monoclonal anti-biotin, 2 mg/mL	0.25 mL
20204-1	CF®488A mouse monoclonal anti-biotin, 2 mg/mL	50 uL
20205	CF®594 mouse monoclonal anti-biotin, 2 mg/mL	0.25 mL
20205-1	CF®594 mouse monoclonal anti-biotin, 2 mg/mL	50 uL
20206	CF®633 mouse monoclonal anti-biotin, 2 mg/mL	0.25 mL
20206-1	CF®633 mouse monoclonal anti-biotin, 2 mg/mL	50 uL
20207	CF®640R mouse monoclonal anti-biotin, 2 mg/mL	0.25 mL
20207-1	CF®640R mouse monoclonal anti-biotin, 2 mg/mL	50 uL

Biotium offers a wide selection of fluorescent dye labeled products including secondary antibodies, antibody labeling kits, and other bioconjugates, featuring our proprietary line of exceptionally bright and photostable CF® dyes. Please visit our website at www.biotium.com for details.