

Product Information

APC Conjugated Antibodies

Catalog No.	Product Description
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

Unit Size

100 uL, 500 uL

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

Form: Blue solution

Spectral Property

$\lambda_{\text{abs}}/\lambda_{\text{em}}$ (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

The products are prepared by labeling IgG (H+L) 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.

For most fluorescent secondary antibody applications, a concentration of 1 - 10 µg/mL is sufficient; however, optimal conditions should be determined empirically.

Other APC products

Cat.#	Product Name	Unit Size
29048-200uL	Streptavidin, APC Conjugate	200 uL
29048-1mL	Streptavidin, APC Conjugate	1 mL
92036	Mix-n-Stain APC Antibody Labeling Kit	25-50ug labeling
92037	Mix-n-Stain APC Antibody Labeling Kit	50-100ug labeling

Other Related Products

Cat.#	Product Name	Unit Size
40061-T	RedDot™2 Far Red Nuclear Counterstain, 200X in DMSO, Trial Size (15-20 tests)	25 uL
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 ^{HT} Pap Pen 2.5 mm tip, ~400 uses	1 pen
22006	Super ^{HT} Pap Pen 4 mm tip, ~800 uses	1 pen
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 secondary antibodies and other conjugates featuring bright and photostable fluorescent CF™ dyes, plus hundreds of other fluorescent reagents and assay kits for life science research.

For research use only.

TWEEN is a registered trademark of Uniqema Americas LLC.