

Product Information

Biotin Conjugated Antibodies

Catalog No.	Product Description
20189	Donkey Anti-Goat IgG (H+L), highly cross-adsorbed Min X Ch, GP, Hs, Hu, Ms, Rb, Rt, SHm
20190	Donkey Anti-Mouse IgG (H+L), highly cross-adsorbed Min X Bv, Ch, Gt, GP, Hs, Hu, Rb, Sh, SHm
20191	Donkey Anti-Rabbit IgG (H+L), highly cross-adsorbed Min X Bv, Ch, Gt, GP, Hs, Hu, Ms, Rt, Sh, SHm
20183	Goat Anti-Mouse IgG (H+L)
20184	Goat Anti-Mouse IgG (H+L), highly cross-adsorbed Min X Bv, Hs, Hu, Rb, and Sw
20185	Goat Anti-Rabbit IgG (H+L)
20186	Goat Anti-Rabbit IgG (H+L), highly cross-adsorbed Min X Hu, Ms and Rt
20187	F(ab') ₂ fragment of Goat anti-Mouse IgG (H+L)
20188	F(ab') ₂ fragment Goat Anti-Rabbit IgG (H+L)
20471	Biotin Goat Anti-Mouse IgG1 Min X Hu, Bv and Rb
20472	Biotin Goat Anti-Mouse IgG2a Min X Hu, Bv and Rb
20473	Biotin Goat Anti-Mouse IgG2b Min X Hu, Bv and Rb

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

Whole IgG: 50 uL, 0.25mL or 0.5 mL (liquid format), or 1 mg (lyophilized)
F(ab')₂ fragment: 0.25 mL (liquid format)

Concentration

Liquid format: 2 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.

Lyophilized format (after reconstitution): 2 mg/mL in pH ~7.4 PBS containing 15 mg/mL bovine serum albumin (IgG-free and protease-free) and 20 mg/mL trehalose.

Color and Form

Liquid format: colorless solution
Lyophilized format: off-white solid

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. Liquid format antibodies contain 50% glycerol and will not freeze at -20°C.

Reconstitution (lyophilized format only): add 0.5 mL dH₂O and mix gently to dissolve. Store at -20°C. Aliquot to avoid freeze-thaw cycles. Alternatively, glycerol can be added to the antibody so that it will not freeze at -20°C: add 0.25 mL dH₂O to the lyophilized antibody and mix gently to dissolve, then add 0.25 mL glycerol and mix well. Optional: a preservative may be added, such as 0.05% (final concentration) sodium azide.

Note: storage of the antibody for more than a day at final working dilution is not recommended.

Recommended Dilution Range

1-10 µg/mL of the IgG conjugate for most applications (appropriate dilutions of the conjugate should be determined empirically).

Product Description

Biotin secondary antibody conjugates can be used in immunofluorescence or immunohistochemical staining protocols for detection using fluorophore- or enzyme-conjugated streptavidin.

Related Products

A full selection of secondary antibodies, antibody labeling kits, bioconjugates of phalloidin, annexin V and a-bungarotoxin are available for many of our CF™ dyes. Please visit the Biotium website at www.biotium.com for details.

Fluorescent CF™ Dye streptavidin conjugates

Streptavidin conjugated to a selection of Biotium's exceptionally bright and photostable CF Dyes

Cat.#	Product Name	Ex (nm)	Em (nm)	Unit Size
29031	Streptavidin, CF350 Conjugate	347	448	1 mg
29032	Streptavidin, CF405S Conjugate	404	431	1 mg
29033	Streptavidin, CF405M Conjugate	408	452	1 mg
29034	Streptavidin, CF488A Conjugate	490	515	1 mg
29038	Streptavidin, CF555 Conjugate	555	565	1 mg
29035	Streptavidin, CF568 Conjugate	562	583	1 mg
29036	Streptavidin, CF594 Conjugate	593	614	1 mg
29037	Streptavidin, CF633 Conjugate	630	650	1 mg
29039	Streptavidin, CF647 Conjugate	650	665	1 mg
29040	Streptavidin, CF660R Conjugate	663	682	1 mg
29044-200uL	R-PE Streptavidin	496,546	565/578	200uL
29044-1mL				1mL
29048-200uL	APC Streptavidin	633	640/660	200uL
29048-1mL				1mL
29049	HRP Streptavidin	---	---	1mg
29071-100uL	Alkaline Phosphatase Streptavidin	---	---	100uL
29071-1mL				1mL

Biotin Mix-n-Stain™ Antibody Labeling Kits

Conjugate 5-100 ug of your antibody to biotin in 30 minutes

Cat.#	Product Name	Unit Size
92286	Mix-n-Stain Biotin Antibody Labeling Kit	1 x 5-20 ug labeling
92266	Mix-n-Stain Biotin Antibody Labeling Kit	1 x 20-50 ug labeling
92244	Mix-n-Stain Biotin Antibody Labeling Kit	1 x 50-100 ug 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
22014	30% Bovine Serum Albumin Solution	100 mL