

Product Information

HRP Conjugated Antibodies

Cat. No.	Product Description
20400	Goat Anti-Mouse IgG (H+L)
20401	Goat Anti-Mouse IgG (H+L), highly cross-adsorbed Min X Bv, Hs, Hu, Rb, Sw
20402	Goat Anti-Rabbit IgG (H+L)
20403	Goat Anti-Rabbit IgG (H+L), highly cross-adsorbed Min X Hu, Ms, Rt
20404	Donkey Anti-Mouse IgG (H+L), highly cross-adsorbed Min X Bv, Ch, Gt, GP, Hs, Hu, Rb, Sh, SHm
20405	Donkey Anti-Rabbit IgG (H+L), highly cross-adsorbed Min X Bv, Ch, Gt, GP, Hs, Hu, Ms, Rt, Sh, SHm
20406	Goat Anti-Rat IgG (H+L), highly cross-adsorbed Min X Hu, Bv, Hs, Rb
20470	Goat Anti-Human IgG (H+L), highly cross-adsorbed Min X Bv, Hs, Ms
20474	Goat Anti-Chicken IgY (H+L), highly cross-adsorbed Min X Bv, Gt, Gp, SHm, Hs, Hu, Ms, Rb, Rt, Sh
20475	Goat Anti-Llama IgG (H+L)
20839	Chicken Anti-Goat 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

HRP Streptavidin

Cat. No.	Product Name	Unit Size
29049	HRP Streptavidin	1 mg

Unit Size

Whole IgG / Streptavidin: 100uL or 1 mL (liquid format), or 1 mg (lyophilized)

Concentration

Liquid format: 1 mg/mL in pH ~7.4 PBS containing 50% glycerol and 15 mg/ml bovine serum albumin (IgG-free and protease-free).

Lyophilized format (after reconstitution): 1 mg/mL in pH ~7.4 PBS containing 10 mg/ml trehalose.

Color and Form

Liquid format: light brown solution

Lyophilized format: light brown 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.

Reconstitution (lyophilized format only): add 1mL pH ~7.4 1X PBS and mix gently to dissolve.

Product Description

This product is prepared by labeling IgG (H+L) or recombinant streptavidin with horseradish peroxidase. HRP conjugates could be used for Western Blot, ELISA, immunohistochemistry and other standard immunoassay applications.

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

HRP antibody conjugates

PSF006

Related Products

Cat. No.	Product Name	Unit Size
10061	10-Acetyl-3,7-dihydroxyphenoxazine (Amplex Red)	5 mg
30015	DAB substrate kit	1 kit
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.