

# **Product Information**

## **HRP Conjugated Secondary Antibodies**

Product	Cat. No.	Unit Size
Goat Anti-Mouse IgG (H+L)	20400-100uL	100 uL
	20400-1mL	1 mL
	20400-1mg	1 mg
Goat Anti-Mouse IgG (H+L), Highly Cross-	20401-100uL	100 uL
Adsorbed	20401-1mL	1 mL
Min X Bv, Hs, Hu, Rb, Sw	20401-1mg	1 mg
	20402-100uL	100 uL
Goat Anti-Rabbit IgG (H+L)	20402-1mL	1 mL
	20402-1mg	1 mg
Coat Anti Dakkit In C (Hall) 1 Patt 1 Coas	20403-100uL	100 uL
Goat Anti-Rabbit IgG (H+L), Highly Cross- Adsorbed	20403-1mL	1 mL
Min X Hu, Ms, Rt	20403-1mg	1 mg
Dankov Anti Mayan InC (IIII ) Highly Cross	20404-100uL	100 uL
Donkey Anti-Mouse IgG (H+L), Highly Cross- Adsorbed	20404-1mL	1 mL
Min X Bv, Ch, Gt, GP, Hs, Hu, Rb, Sh, SHm	20404-1mg	1 mg
Dankov Anti Dahhit InC (II.I.) Highly Cross	20405-100uL	100 uL
Donkey Anti-Rabbit IgG (H+L), Highly Cross- Adsorbed	20405-1mL	1 mL
Min X Bv, Ch, Gt, GP, Hs, Hu, Ms, Rt, Sh, SHm	20405-1mg	1 mg
	20406-100uL	100 uL
Goat Anti-Rat IgG (H+L), Highly Cross-Adsorbed Min X Hu, Bv, Hs, Rb	20406-1mL	1 mL
WIII A FIU, DV, FIS, ND	20406-1mg	1 mg
Goat Anti-Human IgG (H+L), Highly Cross-	20470-100uL	100 uL
Adsorbed	20470-1mL	1 mL
Min X Bv, Hs, Ms	20470-1mg	1 mg
Coot Anti Chiokon lay (U.I.) Highly Cross	20474-100uL	100 uL
Goat Anti-Chicken IgY (H+L), Highly Cross- Adsorbed	20474-1mL	1 mL
Min X Bv, Gt, Gp, SHm, Hs, Hu, Ms, Rb, Rt, Sh	20474-1mg	1 mg
	20475-100uL	100 uL
Goat Anti-Llama IgG (H+L)	20475-1mL	1 mL
	20475-1mg	1 mg
Chicken Anti-Goat IgG (H+L)	20839-100uL	100 uL
		1 ml
Chicken Anti-Goat IgG (H+L)	20839-1mL	1 mL
Chicken Anti-Goat IgG (H+L)	20839-1mL 20839-1mg	1 mg
Chicken Anti-Goat IgG (H+L)  Goat Anti-Digoxigenin (DIG)		

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**

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

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

## Storage and Handling

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

Reconstitution (lyophilized format only): add 1 mL pH ~7.4 1X PBS and mix gently to dissolve. Reconstituted HRP conjugate can be stored in aliquots at -20°C for at least 6 months. Do not add azide to HRP conjugates.

### **Product Description**

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

## **Recommended Dilution Range**

For most secondary antibody applications, a concentration of 1-10 ug/mL is sufficient; however, appropriate dilutions should be determined empirically. For more detailed protocols for immunofluorescence staining for fluorescence microscopy, flow cytometry, or fluorescence-based western detection, please visit our website.

#### **Related Products**

Cat. No.	Product
10061	10-Acetyl-3,7-dihydroxyphenoxazine (Amplex® Red)
30015	DAB Substrate Kit
92300-92302, 92437	Mix-n-Stain™ HRP Antibody Labeling Kit
40061	RedDot™2 Far Red Nuclear Counterstain, 200X in DMSO
23001	EverBrite™ Mounting Medium
23002	EverBrite™ Mounting Medium with DAPI
23003	EverBrite™ Hardset Mounting Medium
23004	EverBrite™ Hardset Mounting Medium with DAPI
23005	CoverGrip™ Coverslip Sealant
22005	Mini Super <sup>HT</sup> Pap Pen 2.5 mm tip, ∼400 uses
22006	Super HT Pap Pen 4 mm tip, ~800 uses
22015	Fixation Buffer
22016	Permeabilization Buffer
22017	Permeabilization and Blocking Buffer
22010	10% Fish Gelatin Blocking Buffer
22011	Fish Gelatin Powder
22014	30% Bovine Serum Albumin Solution
22002	Tween®-20

Please visit www.biotium.com to view our full selection of products including a large selection of primary and secondary antibodies labeled with bright and photostable fluorescent CF® Dyes. We also carry Mix-n-Stain™ antibody labeling kits to label antibodies with HRP, CF® Dyes, fluorescent proteins, and other labels.

Tween is a registered trademark of Uniqema Americas LLC, Amplex Red is a registered trademark of Thermo Fisher Scientific.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.