

Revised: January 21, 2019



# **Product Information**

# Alkaline Phosphatase Conjugated **Antibodies**

Cat. No.	Product Description	
20464-100uL, 1mL	Goat Anti-Mouse IgG (H+L)	
20465-100uL, 1mL	Goat Anti-Rabbit IgG (H+L)	
20466-100uL, 1mL	Donkey Anti-Mouse IgG (H+L), highly cross-adsorbed Min X Bv, Ch, Gt, GP, Hs, Hu, Rb, Sh, SHm	
20467-100uL, 1mL	Donkey Anti-Rabbit IgG (H+L), highly cross-adsorbed Min X Bv, Ch, Gt, GP, Hs, Hu, Ms, Rt, Sh, SHm	

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

# Alkaline Phosphatase Streptavidin

Cat. No.	Product Name	Unit Size
29071-100uL	Alkaline Phosphatase Streptavidin	100 uL
29071-1mL		1mL

#### **Unit Size**

100uL or 1 mL.

#### Concentration

Alkaline phosphatase conjugated antibodies: 1 mg/mL in pH ~7.4 PBS containing 40% glycerol (v/v); 5 mg/ml bovine serum albumin (lgG-free and protease-free); and 0.05% NaN<sub>3</sub> (w/v).

Alkaline phosphatase streptavidin: 1 mg/mL in pH ~7.4 PBS containing 50% glycerol (v/v); 5 mg/ml bovine serum albumin (IgG-free and protease-free); and 0.05% NaN<sub>3</sub> (w/v).

## Color and Form

Colorless, clear solution.

# 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**

This product is prepared by labeling IgG (H+L) or recombinant streptavidin with recombinant alkaline phosphatase. Alkaline phosphatase 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.

Alkaline phosphatase requires phosphate-free buffer with bivalent metal ion (eg. Mg2+) to calatyze the reaction of substrates. After antibody binding, it is recommended to wash the sample with alkaline phosphatase reaction buffer (100 mM Tris-HCl; 100 mM NaCl; 50 mM MgCl<sub>2</sub>; 0.2% Tween-20; pH 9.0) before adding the substrates.

## **Related Products**

Cat No.	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 <sup>HT</sup> Pap Pen 2.5 mm tip, ∼400 uses	1 pen
22006	Super <sup>HT</sup> Pap Pen 4 mm tip, ∼800 uses	1 pen
22020	10X Phosphate-Buffered Saline (PBS)	4 L
22023	Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative	5 x 20 mL
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
10003	BCIP/NBT Kit	1 set
10007	BCIP Pink/NBT Kit	1 set
10005	BCIP Red/NBT Kit	1 set
10009	MUP, free acid	100 mg
10030	FDP (high purity)	5 mg

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 Unigema Americas LLC.