

Ordering Information

Precision Protein Conjugates and Standards

Catalog #	Description
161-0380	Precision Protein StrepTactin-HRP Conjugate, 0.3 ml
161-0382	Precision Protein StrepTactin-AP Conjugate, 0.3 ml
161-0363	Precision Plus Protein Unstained Standards, 1.0 ml
161-0381	Precision Protein StrepTactin-HRP Conjugate, 0.125 ml
161-0376	Precision Plus Protein WesternC Standards, 0.25 ml

Blotting Reagents

170-6435	Premixed Tris Buffered Saline, 10x, 1 L
161-0782	1x TBS/1% Casein, 1 L
170-6537	Gelatin, blotting grade, 200 ml
170-6531	Tween 20, blotting grade, 100 ml
162-0115	Nitrocellulose Membrane (0.45 μ m), 33 cm x 3 m
162-0112	Nitrocellulose Membrane (0.2 μ m), 33 cm x 3 m
170-5040	Immun-Star HRP Substrate, 500 ml
170-5041	Immun-Star HRP Substrate, 100 ml
170-5070	Immun-Star WesternC Kit, 100 ml

Strep-tag technology for western blot detection is covered by US patent 5,506,121 and UK patent 2,272,698. StrepTactin is covered by German patent application P 19641876.3. StrepTactin and *Strep*-tag are trademarks of Institut für Bioanalytik GmbH. Bio-Rad Laboratories is licensed by Institut für Bioanalytik GmbH to sell these products for research use only. Tween is a trademark of ICI Americas, Inc.

Bio-Rad Laboratories, Inc.

2000 Alfred Nobel Drive, Hercules, CA 94547 USA
(510) 741-1000 4110000 Rev D

Instruction Manual

Precision Protein™ StrepTactin-HRP Conjugate

Catalog #161-0380

Catalog #161-0381

Precision Protein™ StrepTactin-AP Conjugate

Catalog #161-0382



Introduction

Precision Protein StrepTactin conjugates are specifically intended for use with Precision Plus Protein™ unstained standards (catalog #161-0363) and Precision Plus Protein™ WesternC™ standards (catalog #161-0376). These standards contain an integral *Strep*-tag amino acid sequence that is recognized by StrepTactin conjugates. StrepTactin is a modified form of streptavidin that binds with greater affinity and specificity to the *Strep*-tag sequence.

Precision Protein StrepTactin conjugates are labeled with either horseradish peroxidase (HRP) (catalog #161-0380, 161-0381) or with alkaline phosphatase (AP) (catalog #161-0382). The conjugates are compatible with chemiluminescent detection methods (Bio-Rad's Immun-Star™ system) and with colorimetric detection systems (Bio-Rad's Immun-Blot® system).

Specifications

	HRP Conjugate (161-0380, 161-0381)	AP Conjugate (161-0382)
Contents	0.3 ml/0.125 ml	0.3 ml
Number of applications		150 blots/60 blots
Storage	4°C	4°C
Shelf life	1 year at 4°C	1 year at 4°C
Load volume of	1–6 µl colorimetric;	1–4 µl colorimetric;
Precision Plus Protein	5–10 µl chemiluminescent*	5–10 µl chemiluminescent**
unstained standards		
Load volume of	5–10 µl chemiluminescent	5–10 µl chemiluminescent
Precision Plus Protein		
WesternC standards		
Recommended	Colorimetric Kits:	Colorimetric Kits:
working dilution	1:5000	1:5000
	Immun-Star HRP Kit:	Immun-Star AP Kit:
	1:10,000–1:15,000	1:5,000–1:10,000
	Immun-Star WesternC Kit:	
	1:10,000–1:50,000	
Buffers	1x PBS, 5 mg/ml BSA, 0.01%, thimerosal, 3.5 mg/ml, StrepTactin-HRP	50 mM Tris, 0.1 M NaCl, 1 mM MgCl ₂ , 0.1% NaN ₃ , 10 mg/ml BSA, 0.5 mg/ml StrepTactin-AP

* 1:30 – 1:60 dilution of standards

** 1:15 – 1:30 dilution of standards

Abbreviated Procedure

For complete instructions, order any of the Immun-Blot assay kits or the Immun-Star chemiluminescent detection kits. Reading the entire instruction manual is advised for optimal results and avoidance of common problems inherent to western blotting procedures.

NOTE: All steps are to be done at room temperature. Wear gloves and protective clothing when preparing and working with the solutions in the assay.

1. Prepare the membrane blot, i.e., electro-phoretic blotting, passive blotting, or filter lifting. If the antigen of interest is present at very low levels, it is best to separate the Precision Plus Protein standards lane with the rest of the blot and perform the detection in parallel, or perform multiple exposures during chemiluminescent detection. This will avoid overdevelopment of the standards.
2. Block the membrane with blocking solutions of choice for 1 hr with gentle agitation (we recommend 3% BSA, 3% gelatin, or 1% casein; milk solutions have a high biotin content that tends to decrease the signal). To minimize detection of endogenous biotin carrier proteins, add 50 µg/ml avidin to the blocking solution.
3. Decant the blocking solution and wash the membrane with TTBS (0.05% Tween 20 in TBS) for 5 min with gentle agitation.
4. Decant the wash solution and incubate with primary antibody (diluted in TTBS) for 1 hr with gentle agitation.
5. Decant primary antibody solution and wash the membrane in TTBS for 5 min. Repeat twice for a total wash time of 15 min.
6. Decant the wash solution and add diluted StrepTactin-conjugate (dilute according to the recommended working dilution for the kit you are using in TTBS) and the appropriate conjugated second-

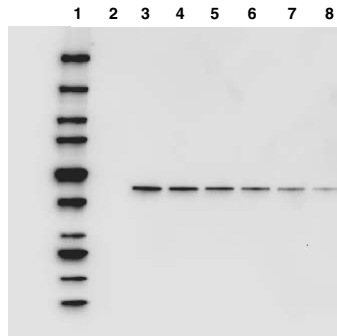
ary antibody, diluted in the same buffer.

Incubate for 1 hr with gentle agitation.

7. Decant the conjugate solution and wash in TTBS for 5 min. Repeat.
8. Perform a final 5 min wash in TBS to remove residual detergent.
9. Proceed to the development step (either the color development step if using Bio-Rad's Immun-Blot assay kits, or the blot development step if using Bio-Rad's Immun-Star kits).

NOTE: When using chemiluminescent detection kits, exposure times may vary. To obtain optimal results, it may be necessary to separate the markers from the rest of the blot and expose them separately, or perform multiple exposures of the standards and unknown lanes.

This product is sold under license from Life Technologies Corporation, Carlsbad, CA, for use only by the buyer of the product. The buyer is not authorized to sell or resell this product or its components.



- Lane 1:** Precision Plus Protein WesternC standards
- Lane 2:** Blank
- Lanes 3–8:** Actin, serial dilution, 200, 150, 100, 75, 50, 25 ng; detection used, Immun-Star WesternC chemiluminescent kit (catalog #170-5070)