

# Blotting Starter Kits



Your Complete Blotting Preparation Resource

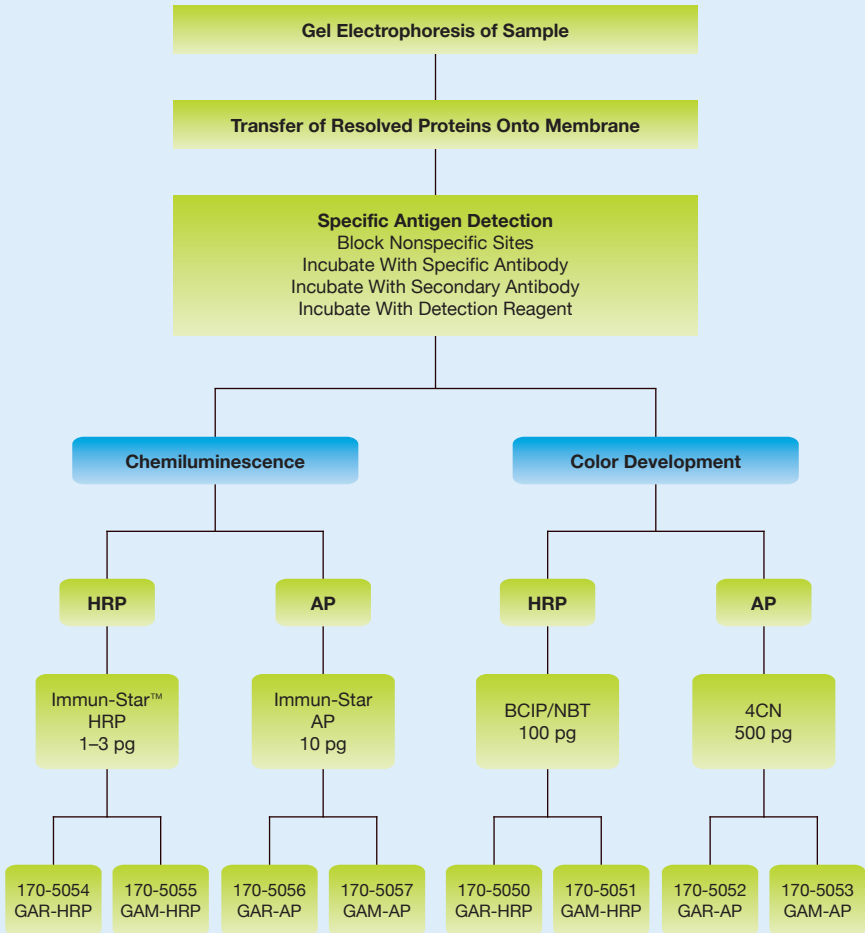
**BIO-RAD**

# Simplify

## Your Blotting Experience

Bio-Rad provides a comprehensive line of reagents and membranes to meet virtually every blotting need.





**Choose an appropriate detection method for your application.** The catalog numbers on the left-hand side of the chart are for chemiluminescent detection, and the ones on the right are for colorimetric detection. Choose your kit based on the sensitivity required and the species of conjugate.

Key: HRP, horseradish peroxidase; AP, alkaline phosphatase; GAR, goat anti-rabbit conjugate; GAM, goat anti-mouse conjugate

# Starter Kit Components

## Ordering Information

Catalog #	Description
170-5050	<b>HRP Color Blot Starter Kit 1</b> , includes 1 L 4CN (HRP substrate), 2 sheets Immun-Blot® PVDF membrane, 2 sheets 0.2 µm nitrocellulose membrane, 50 µl Precision Plus Protein Kaleidoscope™ standards, 1 L 10x TBS, 100 ml 10% Tween 20, 300 g nonfat dry milk, 2 ml GAR-HRP conjugate, 10x Tris/glycine buffer
170-5051	<b>HRP Color Blot Starter Kit 2</b> , same as 170-5050 except with 2 ml GAM-HRP conjugate
170-5052	<b>AP Color Blot Starter Kit 1</b> , includes 1 L BCIP/NBT (AP substrate), 2 sheets Immun-Blot PVDF membrane, 2 sheets 0.2 µm nitrocellulose membrane, 50 µl Precision Plus Protein Kaleidoscope standards, 1 L 10x TBS, 100 ml 10% Tween 20, 300 g nonfat dry milk, 2 ml GAR-AP conjugate, 10x Tris/glycine buffer
170-5053	<b>AP Color Blot Starter Kit 2</b> , same as 170-5052 except with 2 ml GAM-AP conjugate
170-5054	<b>HRP Chemi Blot Starter Kit 1</b> , includes 100 ml Immun-Star HRP substrate, 2 sheets Immun-Blot PVDF membrane, 2 sheets 0.2 µm nitrocellulose membrane, 50 µl Precision Plus Protein unstained standards, 0.3 ml Precision Protein™ StrepTactin-HRP conjugate, 1 L 10x TBS, 100 ml 10% Tween 20, 300 g nonfat dry milk, 2 ml GAR-HRP conjugate, 10x Tris/glycine buffer
170-5055	<b>HRP Chemi Blot Starter Kit 2</b> , same as 170-5054 except with 2 ml GAM-HRP conjugate
170-5056	<b>AP Chemi Blot Starter Kit 1</b> , includes 100 ml Immun-Star AP substrate, 2 sheets Immun-Blot PVDF membrane, 2 sheets 0.2 µm nitrocellulose membrane, 50 µl Precision Plus Protein unstained standards, 0.3 ml Precision Protein StrepTactin-AP conjugate, 1 L 10x TBS, 100 ml 10% Tween 20, 300 g nonfat dry milk, 2 ml GAR-AP conjugate, 10x Tris/glycine buffer
170-5057	<b>AP Chemi Blot Starter Kit 2</b> , same as 170-5056 except with 2 ml GAM-AP conjugate

To order any of the above reagents separately, visit us on the Web at [discover.bio-rad.com](http://discover.bio-rad.com)

StrepTactin is a trademark of Institut für Bioanalytik GmbH. Tween is a trademark of ICI Americas Inc.

### NOTICE TO PURCHASER: LIMITED LICENSE

\*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.



**Bio-Rad**  
**Laboratories, Inc.**

*Life Science  
Group*

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 424 6723 **Australia** 61 2 9914 2800  
**Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 31 3689 6600  
**Canada** 905 364 3435 **China** 86 20 8732 2339  
**Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00  
**Finland** 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 31 884 0  
**Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300  
**Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 03 963 6050  
**Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460  
**Mexico** 52 555 488 7670 **The Netherlands** 0318 540666  
**New Zealand** 0508 805 500 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99  
**Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04  
**Singapore** 65 6415 3188 **South Africa** 27 861 246 723  
**Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55  
**Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000