

See More With Less

The New Immun-Star HRP Detection Kit Brings You Stronger Signal with Less Sample

Chemiluminescent western blot detection is a highly sensitive alternative to isotopic detection. Instead of radioactively labeled antibodies, enzyme-conjugated antibodies convert a chemiluminescent substrate to produce a light signal. This light signal can be captured on film or with dedicated imaging equipment.

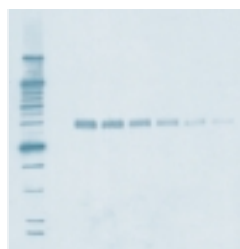
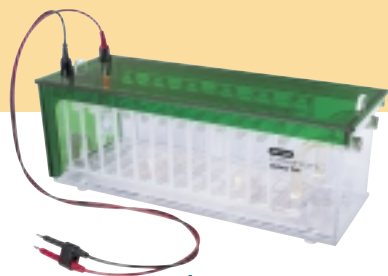
Bio-Rad has expanded its Immun-Star line of western blot products to include the Immun-Star HRP kit for HRP-conjugated secondary antibodies. The chemical reaction involves a peroxidase-catalyzed oxidation of luminol, resulting in the production of light. The addition of an enhancer, included in the kit, intensifies the light signal and increases its duration. The Immun-Star HRP kit is therefore a highly sensitive, fast-acting alternative to isotopic detection.

Bio-Rad's Immun-Star™ HRP
Chemiluminescent Detection Kit

BIO-RAD



See More with Bio-Rad's Immun-Star HRP Chemiluminescent Detection Kit



Ordering Information

Catalog #	Description	Substrate	Antibody	TBS	Tween 20	Blocker
170-5040	Substrate only, 500 ml	•				
170-5041	Substrate only, 100 ml	•				
170-5042	GAR detection reagents*, 500 ml substrate	•	•			
170-5043	GAM detection reagents*, 500 ml substrate	•	•			
170-5044	GAM detection kit**, 500 ml substrate	•	•	•	•	•
170-5045	GAR detection kit**, 500 ml substrate	•	•	•	•	•
170-5046	GAR HRP conjugate, 2 ml		•			
170-5047	GAM HRP conjugate, 2 ml		•			

Conjugates listed above are specially formulated to be compatible with the Immun-Star HRP kit. Tween is a trademark of ICI Americas, Inc.

* Detection reagents also include 2 ml of conjugate.

** Detection kits also include 10x TBS, 1 L, nonfat dry milk, 75 g, and Tween 20, 15 ml.

Supporting Products

Catalog #	Description	Catalog #	Description
165-6001	Criterion™ Cell		Nitrocellulose Membrane (0.2 µm)
170-4070	Criterion Blotter with Plate Electrodes	162-0112	Roll, 33 cm x 3 m, 3
170-4071	Criterion Blotter with Wire Electrodes	162-0146	Sheets, 7 x 8.4 cm, 10
		162-0147	Sheets, 13.5 x 16.5 cm, 10
		162-0150	Sheets, 20 x 20 cm, 5
			Nitrocellulose/Filter Paper Sandwich (0.2 µm)
		162-0232	Sheets, 8.5 x 13.5 cm, 20
		162-0233	Sheets, 8.5 x 13.5 cm, 50
		162-0212	Sheets, 7 x 8.5 cm, 20
		162-0213	Sheets, 7 x 8.5 cm, 50
			Supported Nitrocellulose Membrane (0.45 µm)
		162-0090	Sheets, 7 x 8.4 cm, 10
		162-0091	Sheets, 10 x 15 cm, 10
		162-0092	Sheets, 15 x 15 cm, 10
		162-0093	Sheets, 20 x 20 cm, 10
		162-0094	Roll, 30 cm x 3 m, 1
			Supported Nitrocellulose Membrane (0.2 µm)
		162-0095	Sheets, 7 x 8.4 cm, 10
		162-0096	Sheets, 15 x 15 cm, 10
		162-0097	Roll, 30 cm x 3 m, 1
			Nitrocellulose Membrane (0.45 µm)
		162-0113	Sheets, 20 x 20 cm, 5
		162-0114	Sheets, 15 x 9.2 cm, 10
		162-0115	Roll, 33 cm x 3 m, 1
		162-0116	Sheets, 15 x 15 cm, 10
		162-0117	Sheets, 9 x 12 cm, 10
		162-0145	Sheets, 7 x 8.4 cm, 10
		162-0148	Sheets, 11.5 x 16 cm, 10
			Nitrocellulose/Filter Paper Sandwich (0.45 µm)
		162-0234	Sheets, 8.5 x 13.5 cm, 20
		162-0235	Sheets, 8.5 x 13.5 cm, 50
		162-0214	Sheets, 7 x 8.5 cm, 20
		162-0215	Sheets, 7 x 8.5 cm, 50

Features of Immun-Star HRP Detection Kit:

- Low picogram sensitivity levels
- Stability of working solutions for 24 hours at room temperature
- Signal duration of 6–8 hours
- Compatibility with both PVDF and nitrocellulose membranes
- Reusability of blots (stripping and reprobing)
- Short exposure times (usually ~30 sec)

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 507 6191
Canada (905) 712-2771 China (86-21) 63052255 Denmark 45 44 52-1000 Finland 358 (0)9 804 2200 France 01 47 95 69 65 Germany 089 318 84-177
Hong Kong 852-2789-3300 India (91-124) 6398112/113/114, 6450092/93 Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270
Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 52 5 534 2552 to 54 The Netherlands 0318-540666 New Zealand 64-9-4152280
Norway 47-23-38-41-30 Portugal 351-21-472-7700 Russia 7 095 721 1404 Singapore 65-2729877 South Africa 00 27 11 4428508 Spain 590 5200
Sweden 46 (0)8-55 51 27 00 Switzerland 061 717-9555 United Kingdom 0800-181134