

## General Detection Buffer

### Tris-buffered saline (TBS), 2 L

20 mM Tris-HCl, 500 mM NaCl (pH 7.5)  
(catalog #170-6435, 1 L, 10x)

|                    |         |
|--------------------|---------|
| Tris base          | 4.84 g  |
| NaCl               | 58.48 g |
| diH <sub>2</sub> O | 1.5 L   |

Adjust pH to 7.5 with HCl.

Adjust volume to 2 L with diH<sub>2</sub>O.

### TTBS wash solution, 1 L

20 mM Tris-HCl, 500 mM NaCl, 0.05% Tween 20 (pH 7.5)  
0.5 ml Tween 20  
1 L TBS

### Citrate-buffered saline (CBS)

20 mM citrate, 500 mM NaCl (pH 5.5)  
Included in Immun-Blot® protein G kits.

### TCBS wash solution, 1 L

20 mM citrate, 500 mM NaCl, 0.05% Tween 20 (pH 5.5)  
0.5 ml Tween 20  
1 L CBS

### Blocking solution, 100 ml

3% gelatin-TBS  
Add 3.0 g gelatin to 100 ml TBS.  
Heat to 50°C; stir to dissolve.

or

3% BSA-TBS  
Add 1.0 g BSA to 100 ml TBS; stir to dissolve.

or

5% nonfat milk-TBS  
Add 5.0 g nonfat dry milk to 100 ml TBS; stir to dissolve.

**Note:** Gelatin can clog membranes and cut off the vacuum flow of microfiltration units; use an alternative blocking solution with the Bio-Dot® or Bio-Dot SF apparatus.

**Note:** Nonfat milk is not recommended for avidin/biotin systems as milk contains endogenous biotin and may cross-react with avidin-containing components in the detection system.

### Antibody dilution buffer, 200 ml

1% gelatin-TTBS  
Add 2.0 g gelatin to 200 ml TTBS.  
Heat to 50°C; stir to dissolve.

or

3% BSA-TTBS  
Add 6.0 g BSA to 200 ml TTBS; stir to dissolve.

or

5% nonfat milk-TTBS  
Add 10.0 g nonfat dry milk to 200 ml TTBS; stir to dissolve.

**Note:** Gelatin can clog membranes and cut off the vacuum flow of microfiltration units; use an alternative blocking solution with the Bio-Dot® or Bio-Dot SF apparatus.

**Note:** Nonfat milk is not recommended for avidin/biotin systems as milk contains endogenous biotin and may cross-react with avidin-containing components in the detection system.

### Antibody buffer (for chemiluminescence, ImmunStar™ AP only)

0.2% nonfat milk-TTBS  
Add 0.4 g nonfat milk to 200 ml TTBS; stir to dissolve.

### Antibody buffer for protein G-HRP, 100 ml

1% gelatin-TCBS  
Add 1.0 g gelatin to 100 ml TCBS.  
Heat to 50°C; stir to dissolve.

### Protein G-HRP conjugate solution, 100 ml

Mix 33 µl protein G conjugate solution in 100 ml 1% gelatin in TCBS.

### Streptavidin-biotinylated AP complex, 100 ml

33 µl streptavidin  
100 ml TTBS  
33 µl biotinylated AP

Incubate the complex 1–3 hr at room temperature before use.

## Total Protein Staining Buffers and Solutions

### Amido black staining solution, 1 L

| For nitrocellulose: |        |
|---------------------|--------|
| Amido black         | 5 g    |
| Methanol            | 400 ml |

Adjust volume to 1 L with diH<sub>2</sub>O.

or

|             |        |
|-------------|--------|
| Amido black | 5 g    |
| Isopropanol | 250 ml |
| Acetic acid | 100 ml |

Adjust volume to 1 L with diH<sub>2</sub>O.

| For PVDF:   |        |
|-------------|--------|
| Amido black | 1 g    |
| Methanol    | 400 ml |
| Acetic acid | 100 ml |

Adjust volume to 1 L with diH<sub>2</sub>O.

### Amido black destain solution, 1 L

| For nitrocellulose: |        |
|---------------------|--------|
| Isopropanol         | 250 ml |
| Acetic acid         | 100 ml |

Adjust volume to 1 L with diH<sub>2</sub>O.

| For PVDF:   |        |
|-------------|--------|
| Methanol    | 400 ml |
| Acetic acid | 100 ml |

Adjust volume to 1 L with diH<sub>2</sub>O.

### Coomassie Blue R-250 staining solution, 1 L

|                      |        |
|----------------------|--------|
| Coomassie Blue R-250 | 1 g    |
| Methanol             | 400 ml |
| Acetic acid          | 100 ml |

Adjust volume to 1 L with diH<sub>2</sub>O.

### Coomassie Blue R-250 destaining solution, 1 L

|             |        |
|-------------|--------|
| Methanol    | 400 ml |
| Acetic acid | 100 ml |

Adjust volume to 1 L with diH<sub>2</sub>O.

### Ponceau S staining solution

|                            |       |
|----------------------------|-------|
| Ponceau S                  | 2 g   |
| Trichloroacetic acid (TCA) | 30 g  |
| Sulfosalicylic acid        | 30 g  |
| diH <sub>2</sub> O         | 80 ml |

### Ponceau S destaining solution

1% acetic acid or PBS

### SYPRO Ruby blot pretreatment solution

|                    |        |
|--------------------|--------|
| Acetic acid        | 70 ml  |
| Methanol           | 100 ml |
| diH <sub>2</sub> O | 830 ml |

### Colloidal gold blot staining solution

Use TTBS wash solution.

## Substrate Buffers and Solutions

### HRP Substrate Buffers

|                             |                                  |        |
|-----------------------------|----------------------------------|--------|
| 4-(chloro-1-naphthol)       | 4CN                              | 60 mg  |
|                             | Methanol                         | 20 ml  |
| Protect mixture from light. |                                  |        |
|                             | 3% H <sub>2</sub> O <sub>2</sub> | 600 µl |
|                             | Substrate solution               | 100 ml |

Mix the two solutions together.

Use immediately. Alternatively, use HRP conjugate substrate solution in kit format.

### HRP conjugate solution

Dissolve contents of premixed substrate color development buffer in diH<sub>2</sub>O to 1 L

|                     |        |
|---------------------|--------|
| Color reagent B     | 600 µl |
| Development buffer  | 100 ml |
| HRP color reagent A | 20 ml  |

Use immediately.

### Diaminobenzidine (DAB)

|                                  |        |
|----------------------------------|--------|
| DAB                              | 50 mg  |
| TBS                              | 100 ml |
| 3% H <sub>2</sub> O <sub>2</sub> | 100 µl |

Use immediately.

### AP Substrate Buffers

|                             |                    |         |
|-----------------------------|--------------------|---------|
| AP color development buffer | MgCl <sub>2</sub>  | 0.233 g |
|                             | Tris base          | 12.1 g  |
|                             | diH <sub>2</sub> O | 800 ml  |

Adjust pH to 9.5 with HCl; adjust volume to 1 L with diH<sub>2</sub>O.

### 5-bromo-4-chloroindolyl phosphate/nitroblue tetrazolium (BCIP/NBT)

|                    |        |
|--------------------|--------|
| Dimethylformamide  | 0.7 ml |
| diH <sub>2</sub> O | 0.3 ml |
| NBT                | 30 mg  |
| Dimethylformamide  | 1 ml   |
| BCIP               | 15 mg  |

Add both solutions to 100 ml AP color development buffer. Use immediately. Alternatively, use AP conjugate substrate solution in kit format.

### Immun-Star™ AP substrate solution (kit format)

Use 5 ml chemiluminescent substrate per 100 cm<sup>2</sup>.

For nitrocellulose membrane blots: Add 500 µl enhancer reagent to 10 ml Immun-Star chemiluminescent substrate. Store at 4°C for up to 1 week.

For PVDF membrane blots: Immun-Star AP generates a very fast light signal on PVDF membrane; therefore, the use of an enhancer is not necessary. The substrate is provided ready to use.

### Immun-Star HRP substrate solution (kit format)

For nitrocellulose and PVDF membrane blots: A 1:1 mixture of luminol/ enhancer to peroxide buffer is recommended. Use 10 ml per 100 cm<sup>2</sup> of membrane (12 ml for one 8.5 × 13.5 cm Criterion™ blot).

## Stripping Buffer

### Acidic glycine stripping buffer

|   |        |
|---|--------|
| Glycine   | 7.5 g  |
| Mg(CH <sub>3</sub> COO) <sub>2</sub> ·4H <sub>2</sub> O | 4.3 g  |
| KCl   | 3.7 g  |
| diH <sub>2</sub> O                                      | 800 ml |
| Adjust pH to 2.2 with HCl.                              |        |
| diH <sub>2</sub> O                                      | to 1 L |

This is an excerpt from Bio-Rad's comprehensive Protein Blotting Guide (Bulletin 2895).



**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 11 5044 5699  
**Canada** 905 364 3435 **China** 86 21 6169 8500 **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 9532 220 **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** 64 9 415 2280 **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 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000

---