

## Materials

---

- Gel running apparatus
- Transfer apparatus
- Power supply
- SDS-PAGE mini gel
- Nitrocellulose membrane cut to size of gel
- Filter Paper (Whatman) cut to size of gel
- Tweezers
- X-ray film
- X-ray processor
- Gel loading pipette tips
- Pipettor, small volumes

## Buffer Formulations

---

### SDS/Running Buffer:

- 25 mM Tris
- 192 mM Glycine
- 0.1% SDS

### Transfer Buffer:

- 20 mM Tris
- 150 mM Glycine
- 20% methanol
- 0.038% SDS

### Blocking Buffer

- 5% non-fat dry milk
- TBS

### Wash Buffer (TBST)

- 125 mM NaCl
- 25 mM Tris pH 8.0
- 0.1% Tween-20

## Procedures

---

### Running Protein Samples onto a Gel

1. Cast a mini SDS-PAGE gel per your labs standard protocols or purchase pre-made gels.
2. Clamp the gel to the apparatus with per manufacturer directions. Add running buffer.
3. Remove the comb gently so as to not disturb the wells.
4. Add 6  $\mu$ L of your select marker to a well.
5. Add 7.5  $\mu$ L of lysate per well (2 mg/mL for 15 $\mu$ g per lane).
6. Apply the anode and cathode wires to the appropriate poles and cover.
7. Run at 130V for 2 hours (or until the dye front is close to the bottom).

### Transfer of Proteins onto Membrane

8. Assemble the “sandwich” for the transfer apparatus per the diagram below.
  - a. Note: Handle gel and membrane with tweezers – do not touch!
9. Pre-wet sponges and filter paper in transfer buffer.

a. Note: Filter papers and membrane should be same size as the gel.

10. Insert the “sandwich” and insert into the transfer apparatus, making sure the gel is on the cathode (-) and the membrane is on the anode (+) side of the apparatus.
11. Transfer the proteins to the membrane at 250mA in transfer buffer for 2 hours
12. Remove the cassette from the apparatus, discard the gel, and place membrane on paper towels to dry. (Wash the membrane in 1xTBST to remove any leftover gel.)

### Blocking

13. Incubate the blot with blocking buffer overnight at 4°C or 2 hours at room temperature with gentle agitation.
14. Remove blot from blocking solution.

### Primary Antibody Incubation

15. Mix antibody with peptide in a 1:1 ratio (e.g. 1µg antibody and 1µg peptide) to ensure a molar excess of peptide and incubate for 30 minutes at 37°C.
16. Incubate the blot with the primary antibody for one hour at room temperature or overnight at 4°C.
17. Wash the blot three (3) times 10 minutes each in washing buffer with gentle agitation.

### Secondary Antibody Incubation

18. Dilute 1 µL anti-rabbit IgG-HRP conjugated secondary (or other appropriate secondary) in 10 mL of blocking buffer to make a 1:10000 dilution
  - a. Note: working dilution of secondary can vary from 1:2000 to 1:10000.
19. Incubate blot with secondary antibody for one (1) hour at room temperature.
20. Wash three (3) times for 10 minutes each in washing buffer with gentle agitation.

### Development

21. Drain wash buffer
22. Add ECL solution (Amersham) per manufacturer directions and develop for 1 minute.
23. Drain the fluid.
24. Cover the blot in plastic wrap.
25. Expose the blot to X-ray film for 1 minute in a dark room.
  - b. If there is no banding, expose the film for 5 minutes, then 30 minutes and up to overnight if the signal is weak.
  - c. If the signal is strong, expose the film for 30 seconds or less.
26. Develop the film in an X-ray processor.

### Notes

27. Optimal dilutions should be determined by each laboratory for each antibody.

#### PromoCell GmbH

Sickingenstr. 63/65  
69126 Heidelberg  
Germany

Email: [info@promocell.com](mailto:info@promocell.com)  
[www.promocell.com](http://www.promocell.com)

#### USA/Canada

Phone: 1 - 866 - 251 - 2860 (toll free)  
Fax: 1 - 866 - 827 - 9219 (toll free)

#### Deutschland

Telefon: 0800 - 776 66 23 (gebührenfrei)  
Fax: 0800 - 100 83 06 (gebührenfrei)

#### France

Téléphone: 0800 90 93 32 (ligne verte)  
Téléfax: 0800 90 27 36 (ligne verte)

#### United Kingdom

Phone: 0800 - 96 03 33 (toll free)  
Fax: 0800 - 169 85 54 (toll free)

#### Other Countries

Phone: +49 6221 - 649 34 0  
Fax: +49 6221 - 649 34 40