

## Trans-Blot® Semi-Dry Quick Reference Guide

Catalog Number 170-3940

Full instructions are provided in the manual and should be read before using the Trans-Blot SD cell.

### I. Steps

1. Prepare transfer buffer and equilibrate acrylamide gel in buffer 20–60 minutes.
2. Cut membrane and filter paper to gel size and wet in buffer 5–10 minutes.
3. Remove safety cover and prepare gel sandwich as follows:

To bottom platinum anode place:

Pre-wet filter paper

Pre-wet membrane

Equilibrated gel

Pre-wet filter paper

**Note: Roll out air bubbles  
between layers.**

4. Secure top stainless steel cathode and safety cover.
5. Run blot.  
Optimized conditions will be needed for different proteins.  
Guide range:\*

Mini-gel      10 V for 30 minutes or 15 V for 15 minutes.

Large gel      25 V for 30 minutes or 15 V for 60 minutes.

**Your optimal conditions:**      Time: \_\_\_\_\_ Volts: \_\_\_\_\_

\*Do not exceed 25 volts and 3 mA/cm<sup>2</sup> current for large gels or 5.5 mA/cm<sup>2</sup> for mini gels.

6. Turn off the power supply, unplug electrodes, and remove blot.

# Exploded View of the Trans-Blot SD with Gel Sandwich

