



# eBlot™ L2 Fast Protein Transfer Device

## Quick Guide



GenScript USA Inc.

Email: [product@genscript.com](mailto:product@genscript.com)

Address: 860 Centennial Ave, Piscataway NJ, 08854 USA

Tel: 1-732-885-9188 | 1-877-436-7274 (Toll-Free)

Fax: 1-732-518-5150



### Step 1.

#### Reagents and Materials Preparation

- Pre-run PAGE gel
- eBlot L2 Transfer Buffer, 1X (300 mL per standard transfer)
- eBlot L2 Equilibrium Buffer, 1X (60 mL per membrane in the plastic container provided)
- Distilled water (250 mL per transfer in the silver tray provided)
- PVDF or Nitrocellulose membrane
- Dry transfer sponges (2 pieces per transfer)

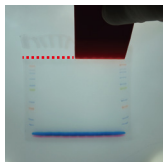
#### Note:

- Regularly empty the waste container.

### Step 3.

Wet PVDF/NC membrane with eBlot L2 Equilibrium buffer for 1 minute (for PVDF membrane: pre-wet the membrane with methanol, ethanol, or isopropanol before equilibrium).

After gel electrophoresis, rinse the gel cassette with water (5 seconds) to remove any remaining running buffer. Open the gel cassette and cut off the gel well fingers. Place the gel in the eBlot L2 Equilibrium buffer for 1 minute\*.



#### Note:

- Cut off any uneven part of the gel
- \* The gel equilibrium time should NOT exceed 5 minutes

### Step 2.

Place the transfer cassette on the bench.



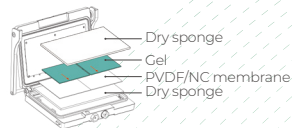
- Release the cassette latch by rotating it upward.
- Lift the latch and fully open the cassette cover to approximately >90°.

#### Note:

- The anode side of the transfer cassette is marked with a red sealing strip.

### Step 4.

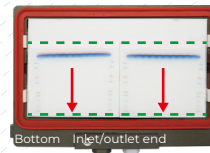
On the red anode side, assemble the transfer sandwich in the following order: **dry sponge, equilibrated membrane, equilibrated gel**. Use the roller to roll out bubbles between the membrane and the gel, then place another piece of **dry sponge** on top of the gel.



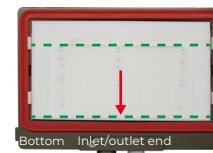
#### Note:

- The sponge doesn't need to be pre-wet
- Note the correct size of the components: Sponge ≥ Membrane ≥ Gel
- Place the gel and membrane close to the bottom, with large molecular proteins close to the inlet/outlet end.
- Avoid overlapping between two mini gels or two membranes.

Transfer 2 mini gels in one cassette



Transfer 1 midi gel in one cassette





# eBlot™ L2 Fast Protein Transfer Device

## Quick Guide



GenScript USA Inc.

Email: [product@genscript.com](mailto:product@genscript.com)

Address: 860 Centennial Ave, Piscataway NJ, 08854 USA

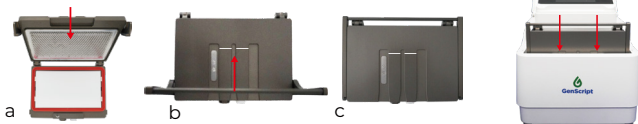
Tel: 1-732-885-9188 | 1-877-436-7274 (Toll-Free)

Fax: 1-732-518-5150



### Step 5.


Close the transfer cassette, then insert the assembled cassette into the selected channel.

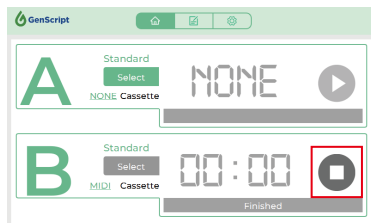


Note the proper orientation and loading of the cassette:


- The latch and scale lines are facing toward the user when inserting the cassette, as shown in Fig. c.
- The transfer cassette is fully inserted into the bottom of the instrument, with the scale lines of the cassette completely covered.

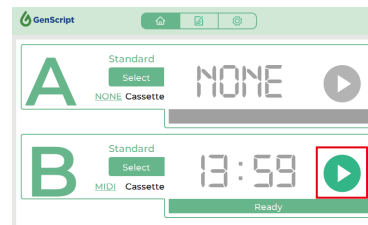
### Step 7.

The device will beep when the countdown reaches 0. Press the stop button  to return to the initial interface.



### Step 6.

Press  for the corresponding channel to initiate the transfer program. The screen will display a countdown timer.



### Step 8.

Remove and open the transfer cassette. Immediately place the transfer sandwich into a tray containing distilled water. Disassemble the transfer sandwich and take out the membrane for the next steps. After each transfer, rinse the cassette with water for 30-60 seconds and set it dry.

Note:

If the PVDF membrane is dry after transfer, it must be reactivated in applicable alcohol before proceeding to the subsequent experiments.

