

eStain™ L1 Protein Staining Device Quick Guide

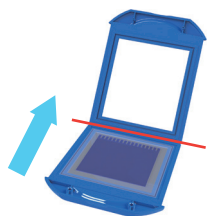
For complete instructions and advanced programming, please refer to the manual



860 Centennial Ave.
Piscataway, NJ 08854 U.S.A.

Tel: 1-732-885-9188
Fax: 1-732-210-0262
Email: product@genscript.com

1. Add distilled water into the tray provided. When gel electrophoresis is complete, carefully remove the gel from the gel cassette and briefly rinse the gel with distilled water for 1 min



2. Open the gel holder and place it on the table with the fabric side down. Carefully place the rinsed gel on the fabric side of the gel holder. Ensure that the upper part of the gel (with comb) point to the axis of the gel holder and is as close to the axis as possible

3. Pre-wet a piece of filter paper (provided) with distilled water and place it on top of the gel



4. Close the gel holder



8. The channel stops flashing and the screen returns to normal. Take out the gel holder and the stained gel. The staining process is complete.



7. The instrument beeps as the program countdown to 0



6. Press the corresponding channel to start the program. The channel timer starts to countdown



5. Choose a Channel and insert the gel holder into the channel.

Note: Please ensure that the fabric side of the gel holder is facing the user before inserting it into the Channel

