

Zinc Stain and Destain Kit

Catalog Number 161-0440

Instructions:

As with any chemical product, wear safety glasses and gloves while using Zinc Stain.

1. Prepare the working solution by diluting Zinc Stain Solution A (Imidazole) and Solution B (Zinc Sulfate) 1:10 with distilled, deionized (DDI) water in separate containers.
2. Place the SDS-PAGE gel in a container with diluted Imidazole, Solution A. Place the container on an orbital mixing platform set to a low mix speed for 10 minutes.
3. Transfer the gel to the diluted Zinc Sulfate, Solution B. Completely immerse the gel for even staining. Allow 30–60 seconds for the gel to develop.
4. To stop the reaction, transfer the gel to a container filled with DDI water and rinse for 3 minutes. Discard this water and replace with fresh DDI water.
5. To visualize the bands, place the gel against the black background on the reverse side of this card. The protein bands will be visible as dark bands against the opaque white gel.
6. Zinc Destaining Protocol: consult Table 1 in the Zinc Stain Kit manual for destain dilutions and wash times dependent on gel thickness.

For complete instructions, consult the instruction manual.

Bio-Rad Laboratories, 2000 Alfred Nobel Drive, Hercules, CA 94547

