

ImmEdge® Hydrophobic Barrier PAP Pen

Cat. No.: H-4000

DESCRIPTION

The ImmEdge Pen provides a light blue hydrophobic barrier soluble in all commonly used clearing agents.

INSTRUCTIONS FOR USE:

- 1) Shake well before use.
- 2) Uncap the pen and place the pen on a solid surface with the tip up. Depress the tip with a gloved finger to release air from the pen.
- 3) Place the pen tip on your microscope slide and depress the tip against the slide several times until the tip is saturated. **DO NOT** over-saturate the tip as this will result in excess reagent pooling on the slide.
- 4) Draw a line around each sample on your microscope slide.
- 5) Replace the pen cap and store at room temperature with the tip down.
- 6) Allow the reagent on the slide to dry for 1 – 2 minutes before immersing the slide in an aqueous solution.