

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.