

August 8 , 2012

Product Information

Mini Super^{HT} Pap Pen Super^{HT} Pap Pen

Catalog Number:

Mini Super^{HT} Pap Pen: 22005 Super^{HT} Pap Pen: 22006

Unit Size: 1 pen

Size:

Mini Super^{HT} Pap Pen: 2.5 mm tip, ~400 applications Super^{HT} Pap Pen: 4 mm tip, ~800 applications Number of applications varies depending on barrier size

Storage and handling: Store at room temperature. The pens contain organic solvents. Avoid skin and eye contact and keep away from open flame. Cap tightly after each use.

Product Description

Super^{HT} Pap Pens are used to create hydrophobic barriers around tissue sections on glass slides to hold staining solutions in place on the slide. This allows the conservation of antibody staining solution or staining of two sections on the same slide with different solutions. The Mini Super^{HT} Pap Pen has a fine point that is useful for separating multiple sections on the same slide. Super^{HT} Pap Pen barriers are insoluble in aqueous buffers, detergents, alcohol and acetone, but can be removed with xylene. Barriers are stable at temperatures up to 120°C.

Directions

Before first use, fill the tip with barrier fluid. Press the tip down on a clean glass microscope slide to open the tip valve (Figure 1). Keep pressing down until you see the tip fill completely with green liquid. Stop pressing as soon as the tip is completely filled.

Barriers should be created after deparaffinization and rehydration for paraffin sections. For frozen sections, create barriers when slides are dry, before the first buffer incubation step. To create a barrier, dry the area around the section with a cotton swab if necessary. Draw an unbroken circle of Pap Pen fluid around the section (Figure 2). Do not press the valve down while drawing the barrier. Take care to prevent the Pap Pen fluid from touching the tissue section. Let the barrier dry 15-30 seconds before immersing slides or adding buffer. Dry the area immediately outside the barrier if necessary and add just enough buffer to fill the barrier without overflowing. A square of Parafilm® can be placed on top of the section to spread the buffer evenly and prevent evaporation. For overnight incubations, slides should be placed in a humidified chamber to prevent evaporation.



Figure 1. Filling pen tip before first use. A. Pen tip is packaged dry. B. Press tip down against a microscope slide until tip fills completely with green liquid (C).

Parafilm is a registed trademark of Bemis Company. TWEEN is a registered trademark of Uniqema Americas LLC.

Biotium products are high-quality reagents and materials intended for research purposes only. Some products are potentially hazardous chemicals - please read the Material Safety Data Sheet for additional information regarding handling potentially hazardous chemicals. Several of Biotium products and product applications are covered by U.S. and international patents and pending patents. Our products are not available for resale or other commercial uses without a specific agreement from Biotium, Inc. We welcome inquiries about licensing the use of our dyes, trademarks or technologies. Please submit inquiries by e-mail to btinfo@biotium.com.

Related Products

| Cat.# | Product Name | Unit Size |
|---------|--|-----------|
| 23001 | EverBrite™ Mounting Medium | 10 mL |
| 23002 | EverBrite™ Mounting Medium with DAPI | 10 mL |
| 23003 | EverBrite™ Hardset Mounting Medium | 10 mL |
| 23004 | EverBrite™ Hardset Mounting Medium with DAPI | 10 mL |
| 23005 | CoverGrip™ Coverslip Sealant | 15 mL |
| 40061-T | RedDot™2 Far Red Nuclear Counterstain, 200X in DMSO, Trial Size (15-20 tests) | 25 uL |
| 22015 | Fixation Buffer | 100 mL |
| 22016 | Permeabilization Buffer | 100 mL |
| 22017 | Permeabilization and Blocking Buffer (5X) | 100 mL |
| 22010 | 10% Fish Gelatin Blocking Buffer | 100 mL |
| 22011 | Fish Gelatin Powder | 2 x 50 g |
| 22014 | 30% Bovine Serum Albumin Solution | 100 mL |
| 22002 | Tween®-20 | 50 mL |

Please visit www.biotium.com to view our full selection of products featuring bright and photostable fluorescent CF[™] dyes, including secondary antibodies and Mix-n-Stain[™] antibody labeling kits, and R-PE dye conjugates. Biotium also offers a variety of apoptosis and cell viability assays for flow cytometry analysis, including mitochondrial membrane potential dyes, fluorescent Annexin V conjugates, and NucView[™]488 Caspase-3 Substrate for live cells.



Figure 2. Creating barriers. A. Draw an unbroken circle around sections with Pap Pen. B. Pipette buffer into the barrier. Add just enough buffer to fill the barrier without overflowing (C).