

00-8877 Mini PAP PEN Lot No.

Mini PAP Pen

APPLICATION:

The Mini PAP Pen is most useful for immunohistochemical and fluorescent staining methods. It prevents the waste of valuable reagents by creating a water repellent circle around the section.

RECOMMENDED USAGE:

1. FOR PARAFFIN-EMBEDDED SECTIONS:

- Deparaffinize and rehydrate tissue sections.
- Rinse tissue sections with PBS and wipe off excess liquid around the tissue.
- Draw a circle on the glass slide around the section with Mini PAP pen.
- Dry the circle for 1-2 minutes at room temperature.
- Continue the normal immunohistostaining process.

2. FOR FROZEN SECTIONS:

- Preclean glass slide.
- Coat slide with Histogrip[™] (Cat no. 00-8050) or other adhesive media (Optional).
- Cut and mount tissue sections on the slides.
- Rinse tissue sections with PBS and wipe off excess liquid around the tissue.
- Draw a circle on the glass slide around the section with Mini PAP pen.
- Dry the circle for 1-2 minutes at room temperature.
- Continue the normal immunohistostaining process.

Note: The water repellent circle is removable in xylene or xylene substitutes. Since it is not influenced by water soluble media, it will not interfere with microscopic study.

STORAGE:

Store at room temperature (15-30°C). Store with cap on tight and tip pointed down. Press down on tip when pen starts to dry out.

Product of Japan

For research use only. CAUTION: Not intended for human or animal therapeutic or diagnostic use.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 E-mail: techsupport@invitrogen.com

PI008877 (Rev 10/09) DCC-09-14

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com. By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.



00-8877 Mini PAP PEN Lot No.

Mini PAP Pen

APPLICATION:

The Mini PAP Pen is most useful for immunohistochemical and fluorescent staining methods. It prevents the waste of valuable reagents by creating a water repellent circle around the section.

RECOMMENDED USAGE:

3. FOR PARAFFIN-EMBEDDED SECTIONS:

- Deparaffinize and rehydrate tissue sections.
- Rinse tissue sections with PBS and wipe off excess liquid around the tissue.
- Draw a circle on the glass slide around the section with Mini PAP pen.
- Dry the circle for 1-2 minutes at room temperature.
- Continue the normal immunohistostaining process.

4. FOR FROZEN SECTIONS:

- Preclean glass slide.
- Coat slide with Histogrip[™] (Cat no. 00-8050) or other adhesive media (Optional).
- Cut and mount tissue sections on the slides.
- Rinse tissue sections with PBS and wipe off excess liquid around the tissue.
- Draw a circle on the glass slide around the section with Mini PAP pen.
- Dry the circle for 1-2 minutes at room temperature.
- Continue the normal immunohistostaining process.

Note: The water repellent circle is removable in xylene or xylene substitutes. Since it is not influenced by water soluble media, it will not interfere with microscopic study.

STORAGE:

Store at room temperature (15-30°C). Store with cap on tight and tip pointed down. Press down on tip when pen starts to dry out.

Product of Japan

For research use only. CAUTION: Not intended for human or animal therapeutic or diagnostic use.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 E-mail: techsupport@invitrogen.com

08877 (Rev 10/09) DCC-09-1480

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.