

IHC PAP (mini) pen

The IHC PAP (mini) pen provides a hydrophobic barrier when a circle is drawn around a specimen on a glass slide.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-950-232-0001	1ea
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Always keep the pen tightly capped and store horizontally. Press down on the tip when the pen starts to dry out.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Technical Info / Product Notes

The IHC PAP (mini) pen provides a thin, film like water repellant barrier when a circle is drawn around the specimen on the glass slide. This barrier creates the proper surface tension to hold antibody solution within the target area on the slide. This avoids the repeated wipeout around the specimen during the staining procedure. Since the reagent is guarded by the hydrophobic circle around the tissue, it minimizes the undesirable dilution of the reagents with the wash buffer. This pen can be used for PAP, ABC, immunofluorescence method, frozen sections, and in situ hybridization procedures. The IHC PAP (mini) pen is especially useful when working with multiple sections on a single slide because its tip is smaller than IHC PAP pen.

Interpretation of the results will be the sole responsibility of the user.

The pen is applicable to microwave staining and in-situ hybridization. Circles made by the pen are heat resistant up to 120°C, ready to be used directly on wet slides, insoluble in alcohol and acetone, and withstand rehydration and dehydration steps in alcohol. If desired, circles can be removed by xylene after the staining procedure is complete.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)