

# IHC PAP pen

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

Citations: 5

[View Online »](#)

## Ordering Information

[Order Online »](#)

ADI-950-233-0001	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 pen provides a thin, film-like water repellent barrier when a circle is drawn around a specimen on a glass slide. This barrier creates the proper surface tension to hold antibody solution within the target area on the slide. This avoids repeated wipeout around the specimen during staining procedures. 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. Two specimens separated by two circles can be applied to the same slide.

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-  
be@enzolifesciences.com](mailto:info-be@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)