

Advanced A-PAP Pen Mini

PAP Pens have been improved! The new formulation of Advanced A-PAP Pens is more easily removed with xylene. Additionally, water repellency and quick-drying are slightly faster than previous liquid blockers. The protocol, volume of the tube, and the diameter of the tip remain the same.

PAP Pen Super-Liquid Blocker Mini is a small-size, ready-to-use hydrophobic barrier pen, designed for immunohistochemistry applications. A minimal amount of antibody is constrained by surface tension within a user-defined border drawn with the pen around a tissue section. The pen is stable up to 129°C in microwave heat.

Ordering Information

[Order Online »](#)

CAC-DAI-APAP-M	1ea
----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Regulatory Status RUO - Research Use Only

Product Details



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)