

# POLYVIEW<sup>®</sup> PLUS AP (anti-mouse) reagent

High sensitivity, low background  
nanopolymer detection reagent for use with  
HIGHDEF<sup>®</sup> chromogens in ISH and IHC  
applications

The POLYVIEW<sup>®</sup> PLUS AP (anti-mouse) reagent is a ready-to-use, biotin-free, one-step detection reagent suitable for immunohistochemical detection of antigens in formalin-fixed paraffin-embedded tissues and frozen sections. The detection reagent may also be used with blood smears, cytospins, and cell preparations. In addition, the POLYVIEW<sup>®</sup> PLUS AP (anti-mouse) reagent can be adapted for *in situ* hybridization detection.

## Ordering Information

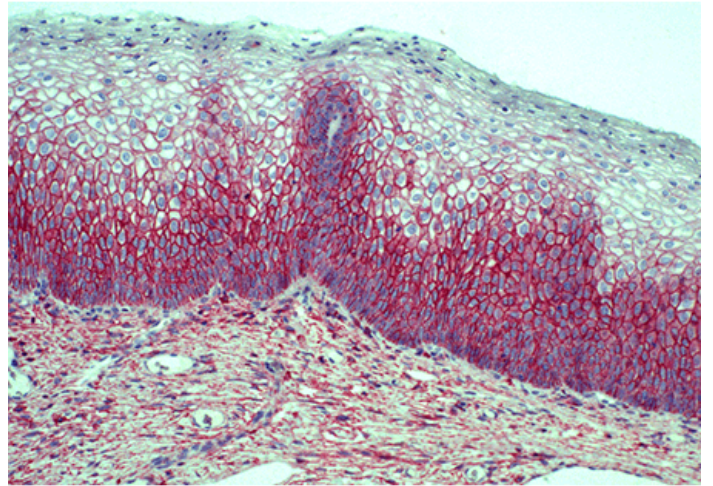
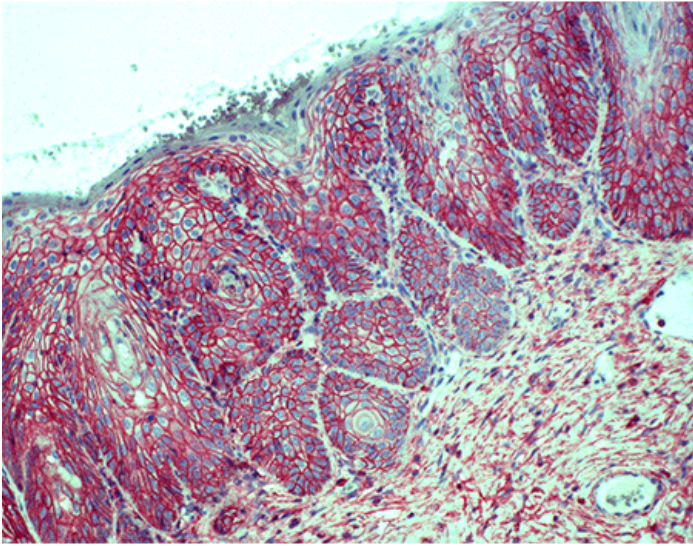
[Order Online »](#)

ENZ-ACC114-0150	150 tests
-----------------	-----------

## Manuals, SDS & CofA

[View Online »](#)

- Ready-to-use reagent
- Biotin-free nanopolymer detection circumvents endogenous biotin background
- High intensity color development delivers sharp, crisp staining
- Compatible with both automated and manual platforms
- The Antibody Blocker/Diluent further reduces background (sold separately)
- Customize with HIGHDEF<sup>®</sup> chromogens of your choice (sold separately)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Do not freeze.
Long Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

**Application** IHC, ISH (in situ hybridization)

**Technical Info / Product Notes** The POLYVIEW<sup>®</sup> PLUS AP (anti-mouse) reagent has been developed by directly labeling anti-mouse immunoglobulins with a nanopolymer of alkaline phosphatase (AP) using proprietary technology. This ensures consistent and reproducible immunodetection of mouse primary antibodies against nuclear, cytoplasmic and membrane antigens in different types of tissues. The single step methodology enables faster staining procedures than traditional two-step methods using biotin and avidin/streptavidin conjugates, with significantly lower background.

The POLYVIEW<sup>®</sup> PLUS AP (anti-mouse) reagent is suitable for use with mouse IgG antibodies. The reagent can be used for manual staining or with automated staining instruments and are well suited for multiplex immunohistochemical staining assays.

The Antibody Blocker/Diluent (ENZ-ACC108) further reduces background and is suitable for diluting the POLYVIEW<sup>®</sup> PLUS AP (anti-mouse) reagent.



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)