

POLYVIEW[®] PLUS AP (anti-rabbit) reagent

High sensitivity, low background
nanopolymer detection reagent for use with
HIGHDEF[®] chromogens in ISH and IHC
applications

The POLYVIEW[®] PLUS AP (anti-rabbit) reagent is a ready-to-use, non-biotin, 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[®] PLUS AP (anti-rabbit) reagent can be adapted for *in situ* hybridization detection.

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

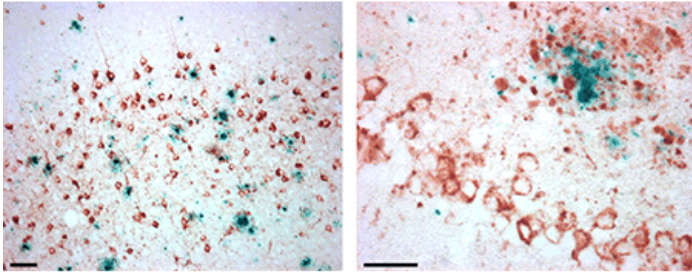
ENZ-ACC110-0150

150 tests

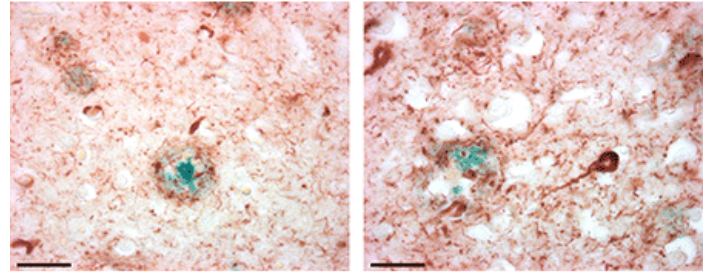
Manuals, SDS & CofA

[View Online »](#)

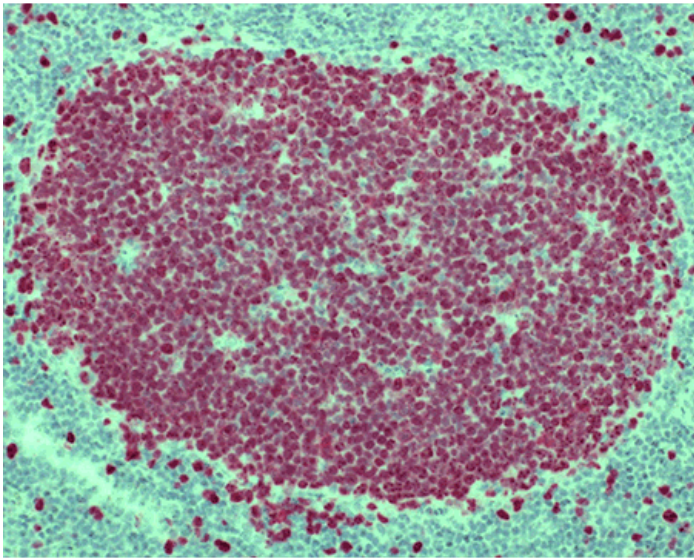
- Ready-to-use reagent
- Biotin-free nanopolymer detection circumvents endogenous biotin background
- High intensity color development delivers sharp, crisp staining
- Compatible with both automated and manual platforms
- The Antibody Blocker/Diluent reduces background further (sold separately)
- Customize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



Dual IHC staining performed on formalin-fixed paraffin-embedded tissue section from human brain with Alzheimer's disease. The tissue sections were stained for alpha-amyloid with POLYVIEW® PLUS AP (anti-rabbit) reagent (Prod. No. ENZ-ACC110) and HIGHDEF® Green AP Chromogen/Substrate (Prod. No. ENZ-ACC130), and phospho-tau with POLYVIEW® PLUS HRP (anti-mouse) reagent (Prod. No. ENZ-ACC104) and HIGHDEF DAB Chromogen/Substrate Set (Prod. No. ENZ-ACC105). Picture courtesy of Dr. Karelle Leroy Laboratoire d'histologie, de neuroanatomie et de neuropathologie, Université Libre de Bruxelles, Bruxelles, Belgium.



Dual IHC staining performed on formalin-fixed paraffin-embedded tissue section from transgenic mouse brain expressing human mutated APP and PS1. The tissue sections were stained for alpha-amyloid with POLYVIEW® PLUS AP (anti-rabbit) reagent (Prod. No. ENZ-ACC110) and HIGHDEF® Green AP Chromogen/Substrate (Prod. No. ENZ-ACC130), and phospho-tau with POLYVIEW® PLUS HRP (anti-mouse) reagent (Prod. No. ENZ-ACC104) and HIGHDEF DAB Chromogen/Substrate Set (Prod. No. ENZ-ACC105). Picture courtesy of Dr. Karelle Leroy Laboratoire d'histologie, de neuroanatomie et de neuropathologie, Université Libre de Bruxelles, Bruxelles, Belgium.



Tonsil tissue stained with anti-Ki-67 antibody (rabbit), followed by POLYVIEW[®] PLUS AP (anti-rabbit) reagent

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[®] PLUS AP (anti-rabbit) reagent has been developed by directly labeling anti-rabbit immunoglobulins with a nanopolymer of alkaline phosphatase (AP) using proprietary technology. This ensures consistent and reproducible immunodetection of rabbit 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[®] PLUS AP (anti-rabbit) reagent is suitable for use with rabbit IgG antibodies, both monoclonal and polyclonal. 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[®] PLUS AP (anti-rabbit) reagent.

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)