

POLYVIEW[®] PLUS

HRP (anti-rabbit)

reagent

High sensitivity, low background nanopolymer detection reagent for use with

HIGHDEF[®] chromogens in IHC and IHC applications

The POLYVIEW[®] PLUS HRP (anti-rabbit) 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[®] PLUS HRP (anti-rabbit) reagent can be adapted for *in situ* hybridization detection.

Application Note

[Tumor Phenotyping and Immune Cell Regulation using Multiplex IHC](#)

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

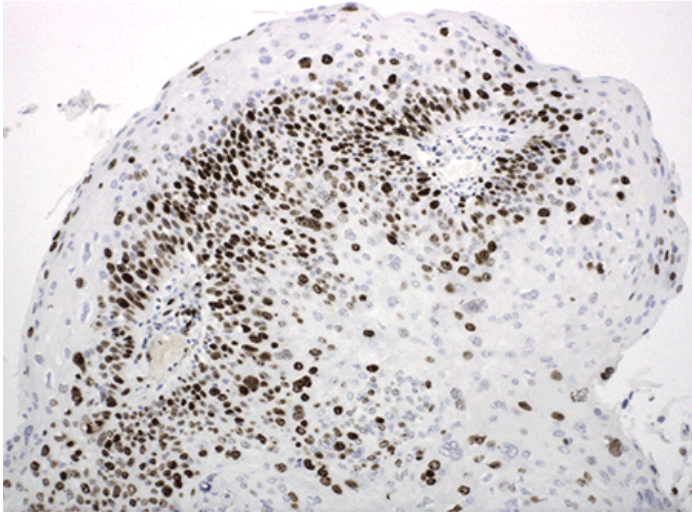
ENZ-ACC103-0150

150 tests

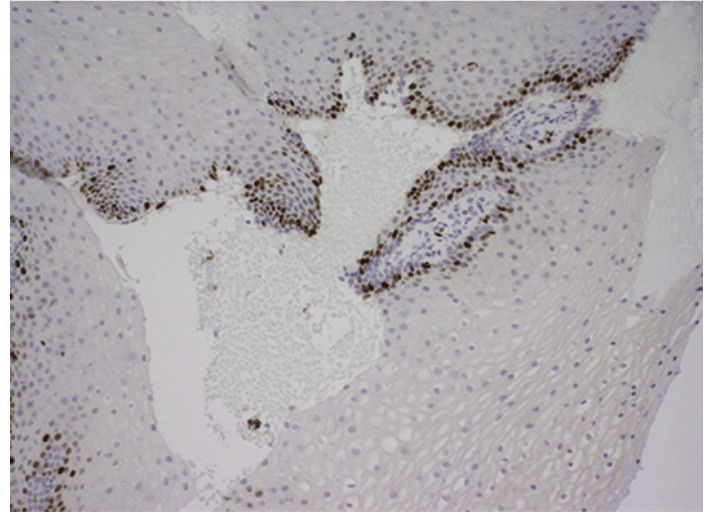
Manuals, SDS & CofA

[View Online »](#)

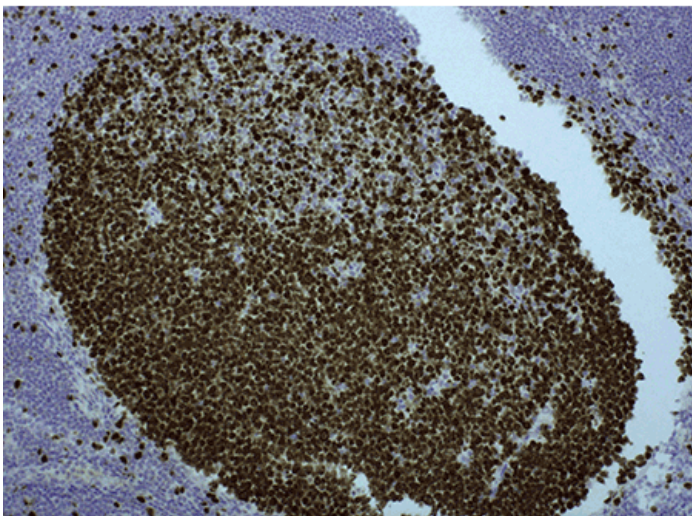
- Ready-to-use reagent
- Biotin-free nanopolymer detection circumvents endogenous biotin background
- High intensity color development delivers sharp, crisp staining
- Compatible with both automated and manual platforms
- The Antibody Blocker/Diluent further reduces background (sold separately)
- Customize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



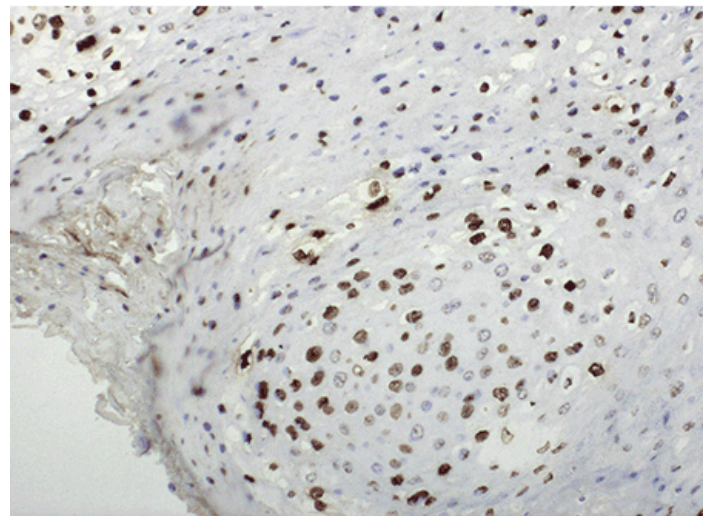
Cervical tissue stained with anti-Ki-67 antibody (rabbit), followed by POLYVIEW® PLUS HRP (anti-rabbit)



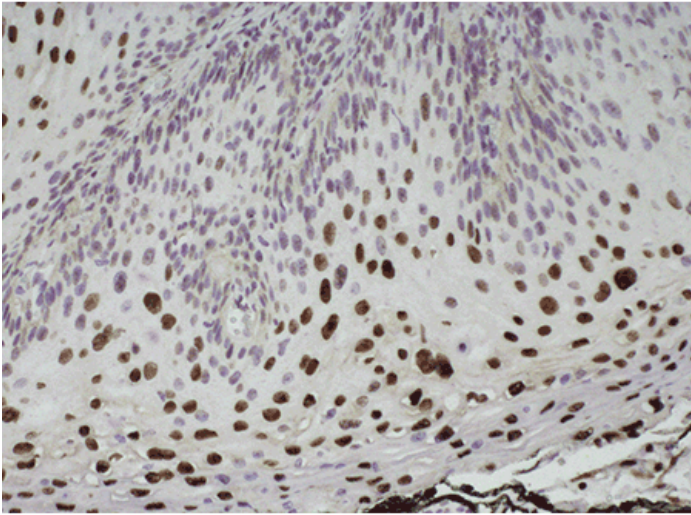
Cervical tissue stained with anti-Ki-67 antibody (rabbit), followed by POLYVIEW® PLUS HRP (anti-rabbit)



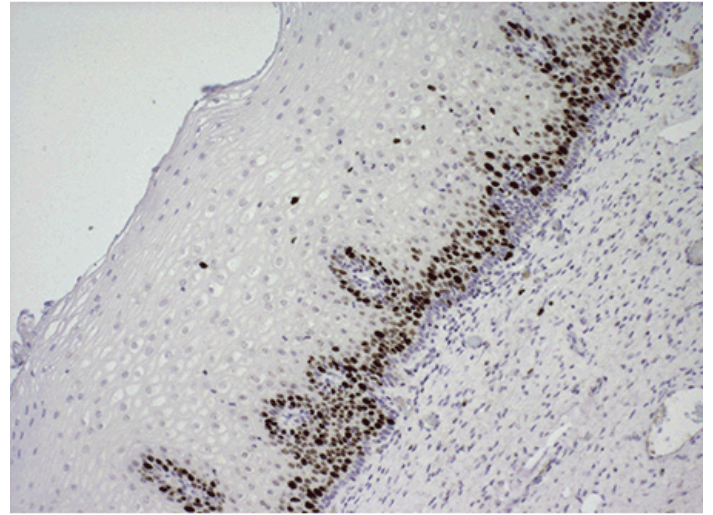
Tonsil tissue stained with anti-Ki-67 antibody (rabbit), followed by POLYVIEW® PLUS HRP (anti-rabbit) reagent



Cervical tissue screened with PATHO-GENE® HPV Type 16/18/31/33/51 (ENZ-32882), then rabbit-anti-biotin linker (ENZ-32892), followed by POLYVIEW® PLUS HRP (anti-rabbit)



Cervical tissue screened with PATHO-GENE[®] HPV Type 16/18/31/33/51 (ENZ-32882), then rabbit-anti-biotin linker (ENZ-32892), followed by POLYVIEW[®] PLUS HRP (anti-rabbit)



Cervical tissue stained with anti-Ki-67 antibody (rabbit), followed by POLYVIEW[®] PLUS HRP (anti-rabbit)

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 HRP (anti-rabbit) reagent has been developed by directly labeling anti-rabbit immunoglobulins with a nanopolymer of horseradish peroxidase (HRP) using proprietary technology. This ensures consistent and reproducible immunodetection of rabbit 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[®] PLUS HRP (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 HRP (anti-rabbit) 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)