

POLYVIEW[®] PLUS

HRP-DAB (Anti-Mouse) Kit

Complete staining kit with highest sensitivity, lowest background

The POLYVIEW[®] PLUS HRP-DAB (Anti-Mouse) Kit is intended for both IHC and ISH staining. The kit is a non-biotin one step detection system suitable for demonstrating antigens in formalin-fixed paraffin-embedded tissues and frozen sections. This kit can be used with blood smears, cytosmeears, and cell preparations.

Citations: 1

[View Online »](#)

Ordering Information

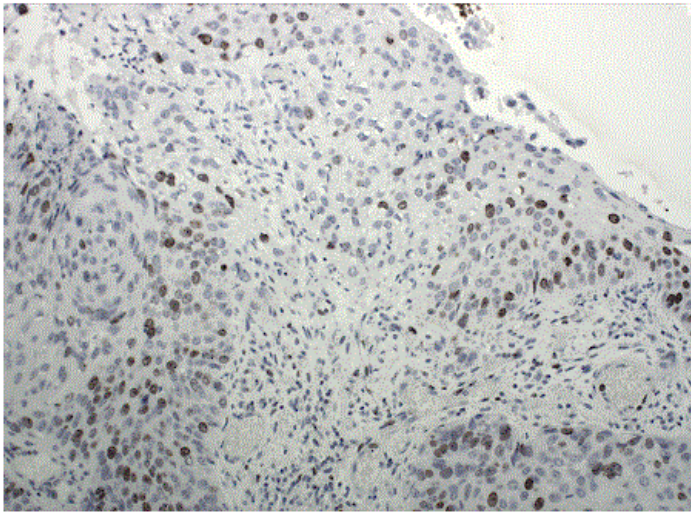
[Order Online »](#)

ENZ-KIT160-0150	150 tests
-----------------	-----------

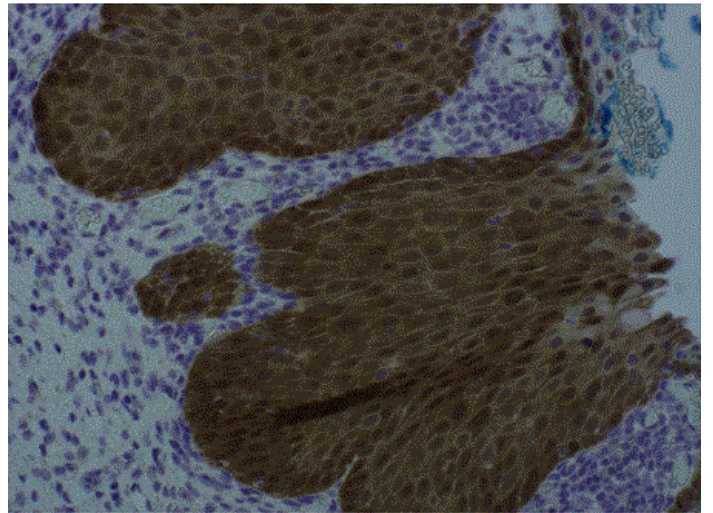
Manuals, SDS & CofA

[View Online »](#)

- Highest sensitivity, complete staining kit for both IHC and ISH
- POLYVIEW[®] PLUS reagents' biotin free nanopolymer detection circumvents endogenous background while providing specific antigen localization
- HIGHDEF[®] DAB (Brown) HRP provides high intensity color development to deliver sharp, crisp staining
- Compatible with both manual and automated platforms



Cervical tissue Ki-67 detected using POLYVIEW® PLUS HRP-DAB (Anti-Mouse) Kit



Cervical tissue p16 detected using POLYVIEW® PLUS HRP-DAB (Anti-Mouse) Kit

Handling & Storage

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)

Assay Time < 2 hours

Contents IHC/ISH Peroxidase Block, Antibody Blocking Buffer, POLYVIEW[®] PLUS HRP (Anti-Mouse) Reagent (Ready-To-Use), HIGHDEF[®] DAB Chromogen, HIGHDEF[®] DAB Substrate Buffer, HIGHDEF[®] Hematoxylin, Antigen Retrieval Reagent, pH 9 (10X)

Technical Info / Product Notes

The POLYVIEW[®] PLUS HRP-DAB (Anti-Mouse) Kit contains POLYVIEW[®] PLUS HRP (Anti-Mouse) Reagent that has been developed by directly labeling anti-mouse immunoglobulins with enzymes using a proprietary tandem hyper-labeling technology in ready-to-use format. This ensures consistent and reproducible immunodetection of mouse antibodies against nuclear, cytoplasmic and membrane antigens in different types of tissues. The single step POLYVIEW[®] PLUS HRP (Anti-Mouse) Reagent allows faster staining procedures as compared to traditional two-step methods using biotin and avidin/streptavidin conjugates, with significantly lower background. Enzo's HIGHDEF[®] DAB (Brown) HRP and HIGHDEF[®] Hematoxylin reagents provide strong signal coupled with the peroxidase and antibody blocking blockers included in the kit to give minimal background.

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)