

POLYVIEW® PLUS

HRP-DAB (Anti-Rabbit) Kit

Complete staining kit with highest sensitivity, lowest background

The POLYVIEW® PLUS HRP-DAB (Anti-Rabbit) 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 may also be used with blood smears, cytosmeears, and cell preparations.

Citations: 7

[View Online »](#)

Ordering Information

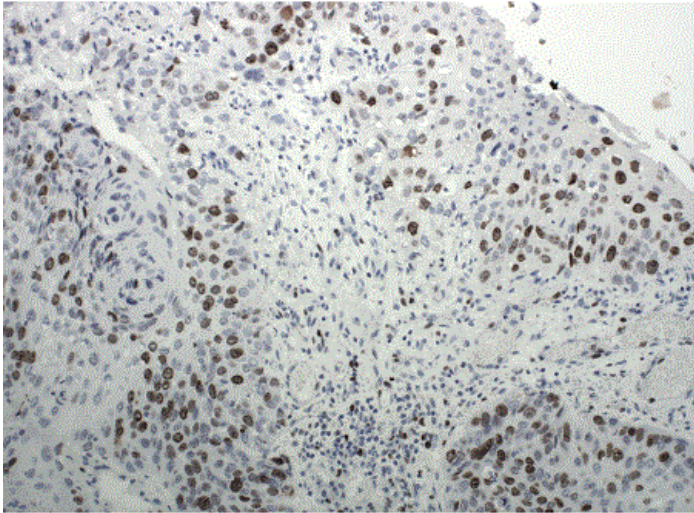
[Order Online »](#)

ENZ-KIT159-0150	150 tests
-----------------	-----------

Manuals, SDS & CofA

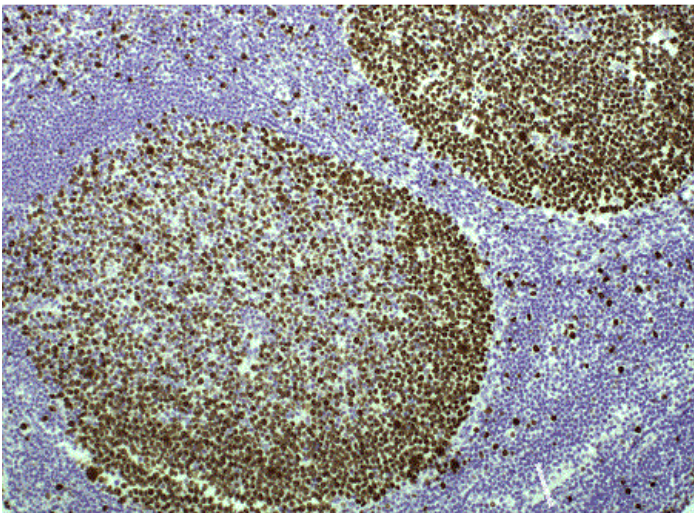
[View Online »](#)

- 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



POLYVIEW® PLUS HRP-DAB (Anti-Rabbit) Kit used to detect Ki-67 on cervical tissue.

.....



POLYVIEW® PLUS HRP-DAB (Anti-Rabbit) Kit used to stain tonsil tissue.

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-Rabbit) 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-Rabbit) Kit contains POLYVIEW[®] PLUS HRP (Anti-Rabbit) Reagent that has been developed by directly labeling anti-rabbit immunoglobulins with enzymes using a proprietary tandem hyper-labelling technology in ready-to-use format. This ensures consistent and reproducible immunodetection of rabbit antibodies against nuclear, cytoplasmic and membrane antigens in different types of tissues. The single step POLYVIEW[®] PLUS HRP (Anti-Rabbit) Reagent enables faster staining procedures 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 blockers included in the kit to give minimal background.

The POLYVIEW[®] PLUS HRP-DAB (Anti-Rabbit) Kit is suitable for use with all rabbit antibodies, both monoclonal and polyclonal. The reagent from this kit can be used for manual staining or with automated staining instruments and is well suited for multiplex immunohistochemical staining assays.



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)