

AMPIVIEW® Dig (Anti-Mouse) Detection HRP-DAB Kit

Complete staining kit with highest sensitivity, lowest background

AMPIVIEW® Dig (Anti-Mouse) Detection HRP-DAB Kit is intended for AMPIVIEW® RNA probes ISH staining. The kit includes DIGX® mouse anti-digoxigenin, Enzo's POLYVIEW® PLUS detection system, HIGHDEF® chromogens and counterstains, and it is suitable for the visualization of the spatial localization of target genes in formalin-fixed paraffin-embedded (FFPE) tissue specimens without disrupting morphology.

Citations: 1

[View Online »](#)

Ordering Information

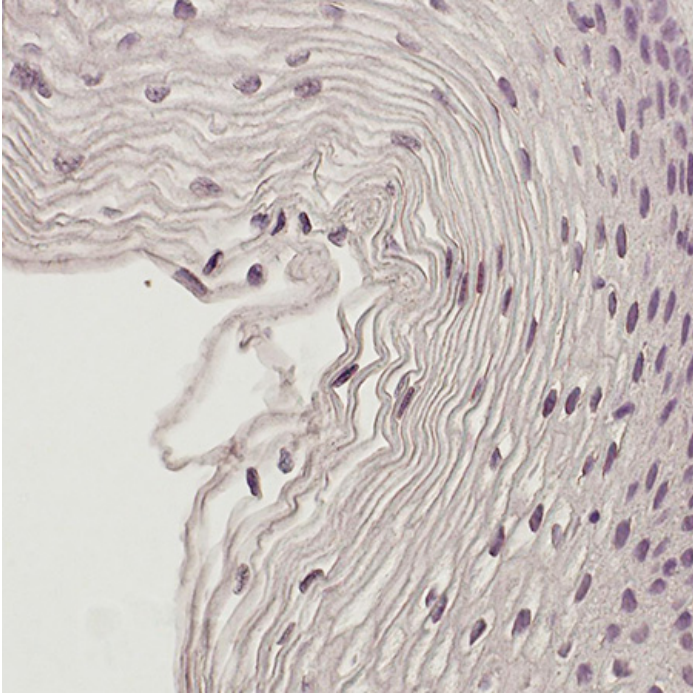
[Order Online »](#)

ENZ-ACC159-0020	20 assays
-----------------	-----------

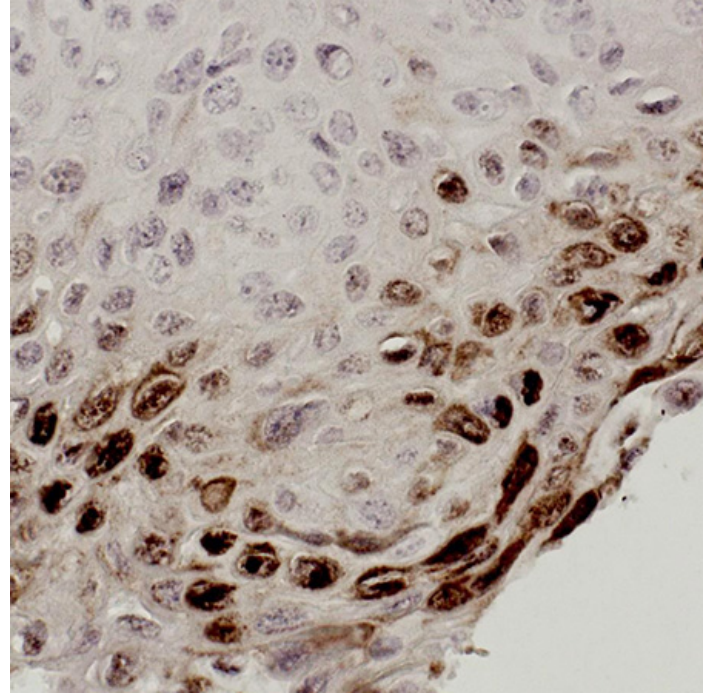
Manuals, SDS & CofA

[View Online »](#)

- Highest sensitivity, complete staining kit for digoxigenin-labeled AMPIVIEW® RNA Probes
- 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



HPV 16/18 not detected in normal cervical tissue with AMPIVIEW 16/18 RNA probes and POLYVIEW (anti-m) HRP/DAB detection kit.



HPV 16/18 detected (brown) in infected cervical tissue with AMPIVIEW HPV 16/18 RNA probes and POLYVIEW (anti-mouse) HRP/DAB detection 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)

Contents POLYVIEW[®] PLUS HRP (Anti-Mouse) Reagent (Ready-To-Use) (ENZ-ACC114), HIGHDEF[®] DAB Chromogen/Substrate (ENZ-ACC105), HIGHDEF[®] Hematoxylin (ENZ-ACC106), Antigen Retrieval Reagent, pH 9 (10X) (ENZ-ACC113), DIGX[®] Mouse anti-digoxigenin linker (ENZ-ABS304), Proteinase K (ENZ-33801), AMPIVIEW[®] Wash Buffer 1, AMPIVIEW[®] Wash Buffer 2

Technical Info / Product Notes

The AMPIVIEW[®] Dig (Anti-Mouse) Detection HRP-DAB 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. 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. Stunning results can be achieved when POLYVIEW[®] PLUS detection reagents are combined with HIGHDEF[®] chromogens and counterstains.

Last modified: October 17, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com