

AMPIVIEW® Dig (Anti-Rabbit) Detection AP-Red Kit

Complete staining kit with highest sensitivity, lowest background

AMPIVIEW® Dig (anti-rabbit) Detection AP-Red Kit is intended for AMPIVIEW® RNA probes ISH staining. The kit includes DIGX® rabbit anti-digoxigenin, POLYVIEW® PLUS AP IHC reagent, HIGHDEF® chromogen and counterstain, and it is suitable visualization of the spatial localization of target genes in formalin-fixed paraffin-embedded (FFPE) tissue specimens without disrupting morphology.

Ordering Information

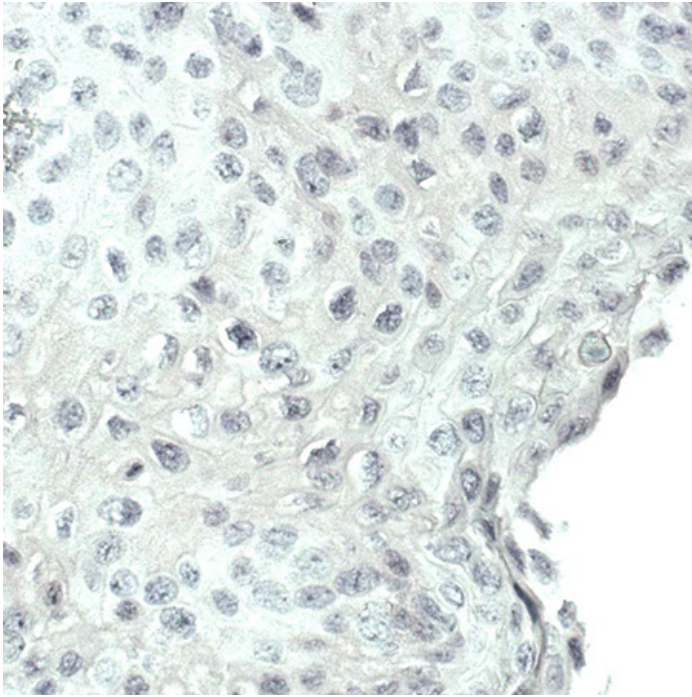
[Order Online »](#)

ENZ-ACC160-0020	20 assays
-----------------	-----------

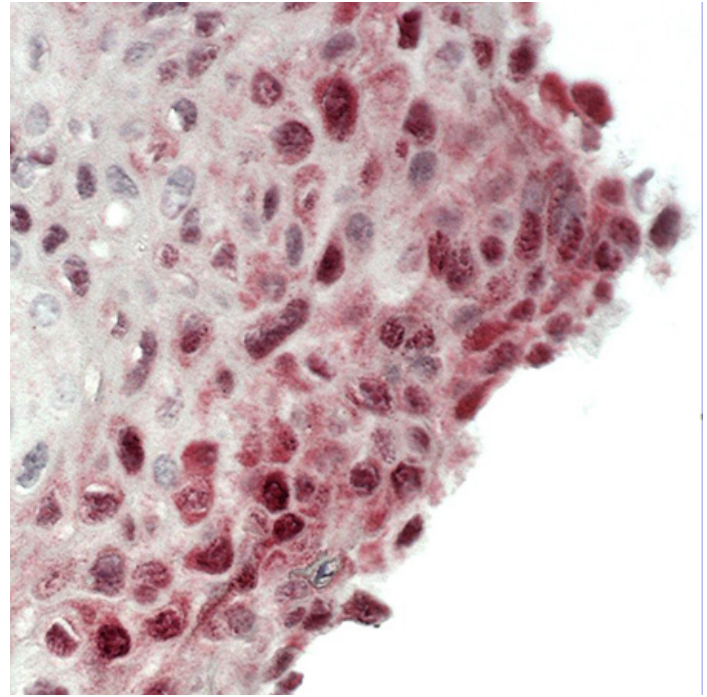
Manuals, SDS & CofA

[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® chromogens provides high intensity color development to deliver sharp, crisp staining
- Compatible with both manual and automated platforms



High-risk HPV not detected with AMPIVIEW NSP-DIG RNA probes (negative control probes) developed with POLYVIEW PLUS AP (anti-rabbit) and HIGHDEF Red AP chromogen/substrate.



High-risk HPV detected (red) with AMPIVIEW HPV High-Risk RNA probes developed with POLYVIEW PLUS AP (anti-rabbit) and HIGHDEF Red AP chromogen/substrate.

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 AP (Anti-Rabbit) Reagent (Ready-To-Use) (ENZ-ACC110), HIGHDEF[®] Red Chromogen/Substrate, HIGHDEF[®] Hematoxylin (ENZ-ACC106), Antigen Retrieval Reagent, pH 9 (10X) (ENZ-ACC113), DIGX[®] Rabbit anti-digoxigenin linker (ENZ-ABS303), Proteinase K (ENZ-33801), AMPIVIEW[®] Wash Buffer 1, AMPIVIEW[®] Wash Buffer 2

Technical Info / Product Notes AMPIVIEW[®] Dig (anti-rabbit) Detection AP-Red Kit contains POLYVIEW[®] PLUS AP (anti-r) Reagent that has been developed by directly labeling anti-rabbit immunoglobulins with enzymes using a proprietary tandem hyper-labeling 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. POLYVIEW[®] PLUS AP(Anti-rabbit) 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 reagent is combined with HIGHDEF[®] chromogens and counterstains.



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)