

# AMPIVIEW<sup>®</sup> Dig (anti-mouse) Detection AP-Red Kit

## Complete staining kit with highest sensitivity, lowest background

AMPIVIEW<sup>®</sup> Dig (Anti-Mouse) Detection AP-Red Kit is intended for AMPIVIEW<sup>®</sup> RNA probes ISH staining. The kit includes DIGX<sup>®</sup> mouse anti-digoxigenin, POLYVIEW<sup>®</sup> PLUS AP IHC reagent, HIGHDEF<sup>®</sup> 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

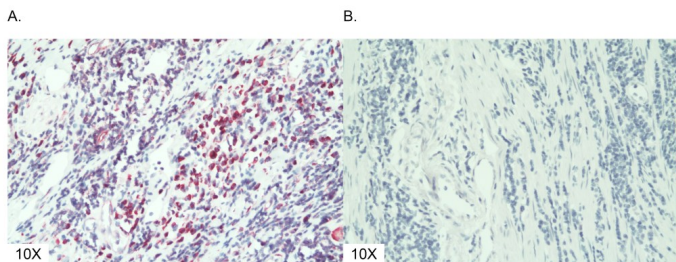
[Order Online »](#)

ENZ-ACC172-0020	20 tests
-----------------	----------

### Manuals, SDS & CofA

[View Online »](#)

- Highest sensitivity, complete staining kit for digoxigenin-labeled AMPIVIEW<sup>®</sup> RNA Probes
- POLYVIEW<sup>®</sup> PLUS reagents' biotin free nanopolymer detection circumvents endogenous background while providing specific antigen localization
- HIGHDEF<sup>®</sup> chromogens provides high intensity color development to deliver sharp, crisp staining
- Compatible with both manual and automated platforms



Ki67 (red) detected in ovarian cancer tissue specimens with A. AMPIVIEW® Ki67 (AS) Dig RNA probes and B. AMPIVIEW® NSP Dig RNA probes. Tissue samples were pre-treated with antigen retrieval reagent, pH 9 (ENZ-ACC113), Proteinase K (ENZ-33801), and AMPIVIEW® Ki67 (AS) Dig RNA probes (ENZ-GEN271) were amplified with DIGX® mouse anti-digoxigenin linker (RTU) (ENZ-ACC304) and detected with POLYVIEW® PLUS (anti-mouse) AP (ENZ-ACC114), combined with HIGHDEF® Red AP chromogen kit (ENZ-ACC158) and hematoxylin (ENZ-ACC106).

## 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<sup>®</sup> PLUS AP (Anti-Mouse) Reagent (Ready-To-Use) (ENZ-ACC114), HIGHDEF<sup>®</sup> Red AP Chromogen/Substrate, HIGHDEF<sup>®</sup> Hematoxylin (ENZ-ACC106), Antigen Retrieval Reagent, pH 9 (10X) (ENZ-ACC113), DIGX<sup>®</sup> Mouse anti-digoxigenin linker (ENZ-ABS304), Proteinase K (ENZ-33801), AMPIVIEW<sup>®</sup> Wash Buffer 1, AMPIVIEW<sup>®</sup> Wash Buffer 2 (ENZ-ACC161)

**Technical Info / Product Notes** AMPIVIEW<sup>®</sup> Dig (anti-mouse) Detection AP-Red Kit is a convenient detection kit for AMPIVIEW<sup>®</sup> RNA probes with reagents from antigen retrieval to counterstain. AMPIVIEW<sup>®</sup> RNA Probes are uniquely designed with the precision of targeted, sequence specific RNA, powered by Enzo's LoopRNA<sup>™</sup> ISH technology to deliver superior sensitivity. Sample specimens are pre-treated with antigen retrieval and Proteinase K treatment before digoxigenin-labeled AMPIVIEW<sup>®</sup> RNA probes are amplified with DIGX<sup>®</sup> anti-digoxigenin linker and detected with POLYVIEW<sup>®</sup> PLUS detection reagents combined with HIGHDEF<sup>®</sup> chromogens and counterstain. Spatial biology of target genes can be visualized within its morphological context under the light microscope.



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)