

# AMPIVIEW<sup>®</sup> DNP (anti-rabbit) Detection HRP-DAB Kit

**Complete staining kit with highest sensitivity, lowest background**

AMPIVIEW<sup>®</sup> DNP (Anti-Rabbit) Detection HRP-DAB Kit is intended for AMPIVIEW<sup>®</sup> RNA probes ISH staining. The kit includes AMPIVIEW<sup>®</sup> Rabbit Anti-DNP, Enzo's POLYVIEW<sup>®</sup> PLUS detection system, HIGHDEF<sup>®</sup> 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.

## Ordering Information

[Order Online »](#)

ENZ-ACC177-0020	20 tests
-----------------	----------

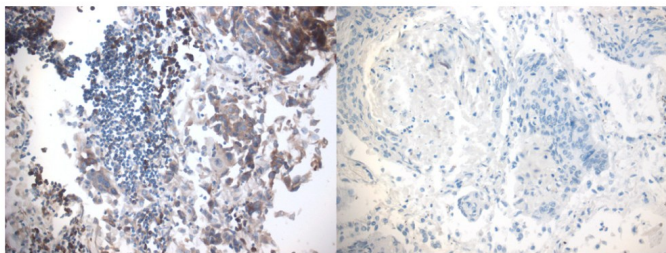
## Manuals, SDS & CofA

[View Online »](#)

- Highest sensitivity, complete staining kit for DNP-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> DAB (Brown) HRP provides high intensity color development to deliver sharp, crisp staining
- Compatible with both manual and automated platforms

A.

B.



PD-L1 (brown) detected in lung carcinoma tissue with A. AMPIVIEW® PD-L1 (AS) DNP RNA Probes and B. AMPIVIEW® NSP DNP RNA Probes. Tissue samples were pre-treated with antigen retrieval reagent, pH 9 (ENZ-ACC113), Proteinase K (ENZ-33801), and AMPIVIEW® PD-L1 (AS) DNP RNA probes (ENZ-GEN291) were amplified with AMPIVIEW® anti-DNP Linker (RTU) (ENZ-ACC928) and detected with POLYVIEW® PLUS (anti-rabbit) HRP (ENZ-ACC103), combined with HIGHDEF® DAB Chromogen/Substrate (ENZ-ACC105) 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 HRP (Anti-Rabbit) Reagent (Ready-To-Use) (ENZ-ACC103), HIGHDEF<sup>®</sup> DAB Chromogen/Substrate (ENZ-ACC105), HIGHDEF<sup>®</sup> Hematoxylin (ENZ-ACC106), Antigen Retrieval Reagent, pH 9 (10X) (ENZ-ACC113), AMPIVIEW<sup>®</sup> Rabbit anti-DNP linker (ENZ-ABS928), 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> DNP (anti-rabbit) Detection HRP-DAB 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 dinitrophenol (DNP)-labeled AMPIVIEW<sup>®</sup> RNA probes are amplified with AMPIVIEW<sup>®</sup> anti-DNP 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)