AMPIVIEW® DNP (anti-rabbit) Detection HRP-DAB Kit

Complete staining kit with highest sensitivity, lowest background

AMPIVIEW[®] DNP (Anti-Rabbit) Detection HRP-DAB Kit is intended for AMPIVIEW[®] RNA probes ISH staining. The kit includes AMPIVIEW[®] Rabbit Anti-DNP, Enzo's POLYVIEW[®] PLUS detection system, HIGHDEF [®] chromogens and countertains, 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 »

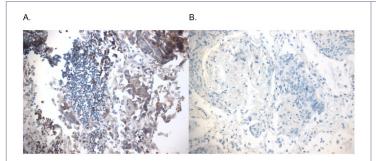
ENZ-ACC177-0020

20 tests

Manuals, SDS & CofA

View Online »

- Highest sensitivity, complete staining kit for DNP-labeled AMPIVIEW[®] 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



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® PLUS HRP (Anti-Rabbit) Reagent (Ready-

To-Use) (ENZ-ACC103), HIGHDEF® DAB

Chromogen/Substrate (ENZ-ACC105), HIGHDEF[®]
Hematoxylin (ENZ-ACC106), Antigen Retrieval Reagent,

pH 9 (10X) (ENZ-ACC113), AMPIVIEW[®] Rabbit anti-DNP linker (ENZ-ABS928), Proteinase K (ENZ-33801).

AMPIVIEW® Wash Buffer 1, AMPIVIEW® Wash Buffer 2

(ENZ-ACC161)

Technical Info / Product Notes

AMPIVIEW[®] DNP (anti-rabbit) Detection HRP-DAB Kit is a convenient detection kit for AMPIVIEW[®] RNA probes with reagents from antigen retrieval to counterstain. AMPIVIEW [®] RNA Probes are uniquely designed with the precision of targeted, sequence specific RNA, powered by Enzo's LoopRNA[™] ISH technology to deliver superior sensitivity. Sample specimens are pre-treated with antigen retrieval and Proteinase K treatment before dinitrophenol (DNP)-labeled AMPIVIEW[®] RNA probes are amplified with AMPIVIEW[®] anti-DNP linker and detected with POLYVIEW[®] PLUS detection reagents combined with HIGHDEF[®] chromogens and counterstain. Spatial biology of target genes can be visualized within its morphological context under the light microscope.

Last modified: January 13, 2025



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg
Phone: +32 3 466 0420
infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com