

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

Complete staining kit with highest sensitivity, lowest background

AMPIVIEW® Dig (Anti-Rabbit) Detection HRP-DAB Kit is intended for AMPIVIEW® RNA probes ISH staining. The kit includes DIGX® rabbit 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.

Ordering Information

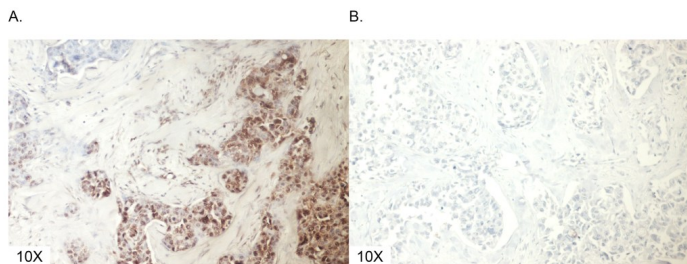
[Order Online »](#)

ENZ-ACC171-0020	20 tests
-----------------	----------

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



NFκB (brown) detected in breast cancer with A. AMPIVIEW® NFκB (AS) RNA Probes and B. AMPIVIEW® NSP Dig RNA probes. Tissue samples are pretreated with antigen-retrieval reagent, pH 9 (ENZ-ACC113), Proteinase K (ENZ-33801), and AMPIVIEW® RNA Probes are amplified with DIGX® rabbit anti-digoxigenin (ENZ-ABS303) and detected with POLYVIEW® PLUS (anti-r) 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), DIGX[®] Rabbit anti-digoxigenin linker (ENZ-ABS303), Proteinase K (ENZ-33801), AMPIVIEW[®] Wash Buffer 1, AMPIVIEW[®] Wash Buffer 2 (ENZ-ACC161)

Technical Info / Product Notes AMPIVIEW[®] Dig (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 digoxigenin-labeled AMPIVIEW[®] RNA probes are amplified with DIGX[®] anti-digoxigenin 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.



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)