

AMPIVIEW™ RNA Probes

Powered by Enzo's LoopRNA ISH™ Technology

AMPIVIEW™ RNA probes are uniquely designed with the precision of targeted, sequence-specific RNA probes, powered by Enzo's LoopRNA ISH™ technology to deliver superior sensitivity. When AMPIVIEW™ RNA probes hybridize to nucleic acid targets, loops form exposing biotin or digoxigenin labels, making these probes not only the most sensitive means of detection, but also adaptable with existing immunohistochemistry (IHC) detection solutions and chromogens. Additionally, sample morphology is preserved and the signal of the targeted biomarker can be seen under the light microscope.

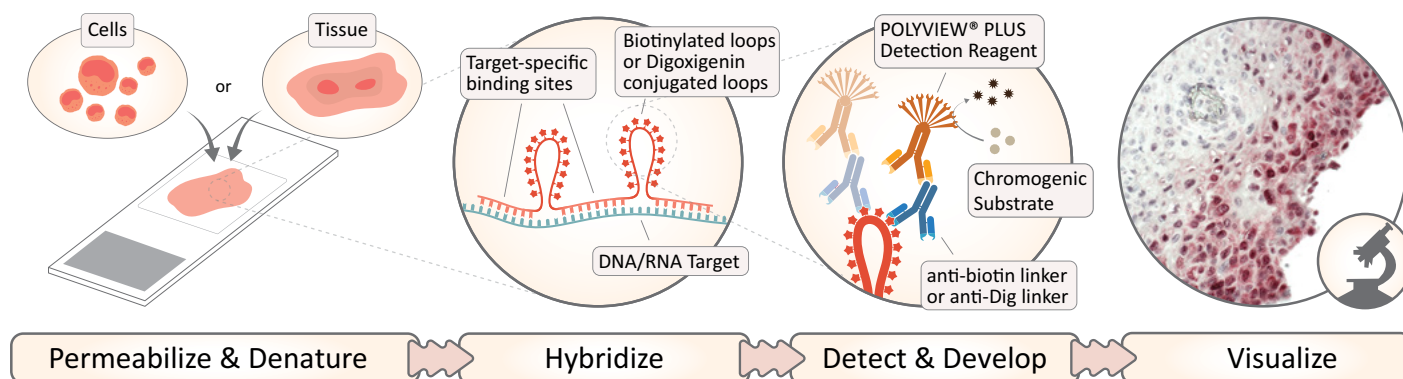
Main Features

- Detection of unique nucleic acid targets (DNA/RNA) down to a single cell level
- High sensitivity and reliability
- Protocols for formalin-fixed, paraffin-embedded (FFPE) tissues and cells
- Mild protocols to preserve tissue morphology
- End-user can use existing ISH set ups and protocols
- Optimized with Enzo's cost-effective IHC/ISH detection portfolio
- Flexible products ready for any workflow (Manual or Automated)

APPLICATIONS

- Basic Research
- Biomarkers
- Cancer
- Developmental Biology
- Drug Discovery and Development
- Infectious Diseases
- Neuroscience
- Pathogen Identification
- And more...

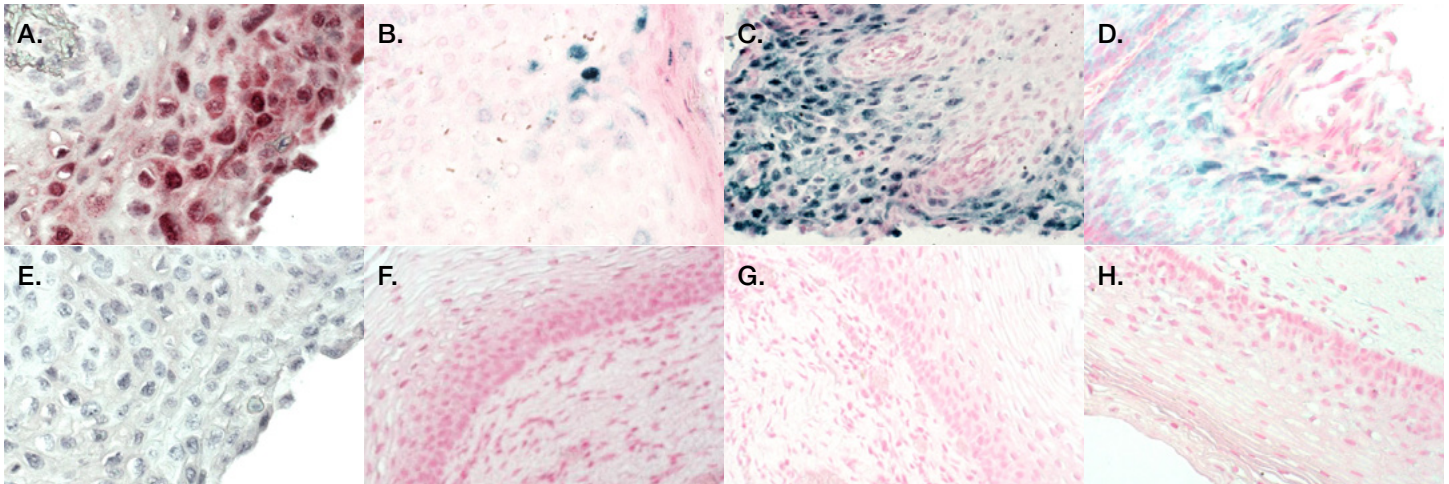
Visualize the Spatial Biology of Nucleic Acids in Tissues and Cells with a Light Microscope



AMPIVIEW™ RNA Probes



Highly Specific and Sensitive RNA Probes Powered by Enzo's LoopRNA ISH™ Technology



Strong signal can be viewed with AMPIVIEW™ HPV RNA probes on HPV infected cervical tissue (A-D). A. AMPIVIEW™ HPV High-Risk RNA probes signal (red) B. AMPIVIEW™ HPV 6/11 RNA probes signal (blue) C. AMPIVIEW™ HPV 16/18 RNA probes signal (blue) D. AMPIVIEW™ HPV 31/33/51 RNA probes signal (green). No specific signal can be viewed when AMPIVIEW™ HPV RNA probes were tested on non-infected cervical tissue (E-H). E. AMPIVIEW™ HPV High-Risk RNA probes signal F. AMPIVIEW™ HPV 6/11 RNA probes signal G. AMPIVIEW™ HPV 16/18 RNA probes signal H. AMPIVIEW™ HPV 31/33/51 RNA probes signal.

Ordering Information

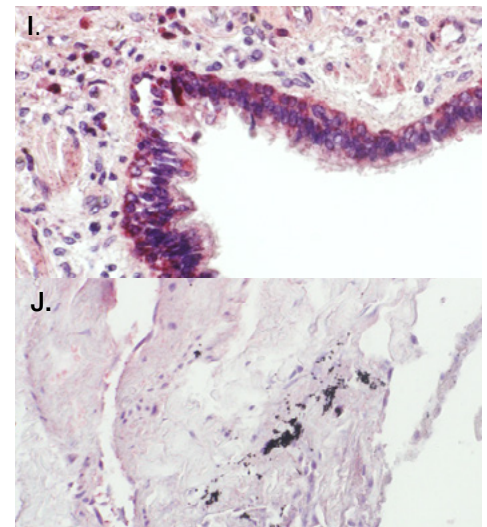
AMPIVIEW™ RNA Probes Sets

Product Name	Product #
AMPIVIEW™ HPV 6/11 RNA Probes Set	ENZ-GEN145
AMPIVIEW™ HPV 16/18 RNA Probes Set	ENZ-GEN146
AMPIVIEW™ HPV 31/33/51 RNA Probes Set	ENZ-GEN147
AMPIVIEW™ HPV High-Risk RNA Probes Set	ENZ-GEN148
AMPIVIEW™ SARS-CoV-2 RNA Probes Set	ENZ-GEN149
AMPIVIEW™ HER2 (AS) RNA Probes Set	ENZ-GEN161

AMPIVIEW™ RNA Probes sets come with targeted probes and AMPIVIEW™ Hybridization Buffer.

For more product and detection systems information, visit enzolifesciences.com/AMPIVIEW

For Research Use Only. Not for Use in Diagnostic Procedures.



I. Strong signal (red) is evident in SARS-CoV-2 infected tissue with AMPIVIEW™ SARS-CoV-2 RNA probes combined with Enzo's IHC detection system. J. No specific labeling is evident in normal lung tissue incubated with AMPIVIEW™ SARS-CoV-2 RNA probes combined with Enzo's IHC detection system.

ASK US ABOUT CUSTOM PROBES

Enzo can design probes for practically ANY gene in ANY genome to be used in ANY tissue or cells.



Global Headquarters
ENZO LIFE SCIENCES, INC.
 Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES (ELS) AG
 Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg
 Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
 Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
 Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
 Phone (UK customers):
 0845 601 1488
 Phone: +44 1392 825900
info-uk@enzolifesciences.com

For our global distributor network and detailed product information visit us online: www.enzolifesciences.com
 All product names, logos, and brands are the property of their respective owners.