

AMPIVIEW® HPV 31/33/51 (AS) DNP RNA Probes Set

High-risk HPV 31/33/51 RNA in situ hybridization probe

The AMPIVIEW® HPV 31/33/51 (AS) DNP RNA Probes set is a ready-to-use in situ hybridization (ISH) probe, powered by Enzo's LoopRNA ISH™ technology, for chromogenic detection of Human Papillomavirus types 31, 33, and 51 (HPV 31/33/51), high-risk oncogenic strains associated with HPV-related cervical, oropharyngeal, and anogenital conditions. This probe set extends high-risk HPV coverage for comprehensive ISH-based genotyping in research specimens. in formalin-fixed paraffin-embedded (FFPE) tissue sections and cells. The probe is DNP-labeled and detected using an unconjugated anti-DNP linker followed by a polymer-based POLYVIEW® PLUS detection system combined with HIGHDEF® chromogens and counterstain, visualized under a standard light microscope.

Supplied as 2 x 1.0 mL ready-to-use probe (ENZ-GEN316) with 1 x 2.0 mL AMPIVIEW® Hybridization Buffer (Fast Hyb Solution) (ENZ-ACC187).

AMPIVIEW® RNA probes are designed to detect the RNA or DNA/RNA of Human Papillomavirus types 31, 33, and 51 (HPV 31/33/51), high-risk oncogenic strains associated with HPV-related cervical, oropharyngeal, and anogenital conditions. This probe set extends high-risk HPV coverage for comprehensive ISH-based genotyping in research specimens. In situ hybridization (ISH) is a molecular technique that uses labeled nucleic acid probes to detect and localize specific RNA or DNA sequences directly within tissue sections or cells. Powered by Enzo's LoopRNA ISH™ technology, AMPIVIEW® RNA probes deliver superior sensitivity for chromogenic detection of RNA, DNA, or RNA/DNA targets on formalin-fixed paraffin-embedded (FFPE) tissue and cells, and are compatible with Enzo's complete immunohistochemistry detection workflow, from antigen retrieval to counterstain.

Ordering Information

[Order Online »](#)

ENZ-GEN316-2000	
-----------------	--

Manuals, SDS & CofA

[View Online »](#)

- Powered by Enzo's LoopRNA ISH™ technology for superior sensitivity
- Ready-to-use, DNP-labeled probe compatible with AMPIVIEW® Rabbit Anti-DNP Linker, POLYVIEW® PLUS detection and HIGHDEF® chromogens
- 1-hour hybridization with AMPIVIEW® Hybridization Buffer (Fast Hyb Solution) (ENZ-ACC187) for fast, same-day results
- Compatible with manual staining and automated platforms (LEICA BOND, Roche DISCOVERY/ULTRA) with prior optimization

Handling & Storage

Use/Stability	Stable under recommended storage conditions up to the expiration date indicated on the vial label. Avoid freeze-thaw cycles.
Handling	Protect from light and humidity. Avoid freeze-thaw cycles. Avoid RNases; wear gloves and work in a clean environment.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ISH
Contents	2 x 1.0 mL AMPIVIEW® HPV 31/33/51 (AS) DNP RNA Probes set (ENZ-GEN316); 1 x 2.0 mL AMPIVIEW® Hybridization Buffer (ENZ-ACC187)
Quantity	2 x 1.0 mL probe + 1 x 2.0 mL hybridization buffer
Recommendation Dilutions/Conditions	Ready to use solution for manual use; Optimization may be required for automated staining systems.
Specificity	HPV 31, HPV 33, HPV 51 (antisense)
Technical Info / Product Notes	Ready-to-use probe for manual staining. Prior optimization required for automated staining systems. AMPIVIEW® RNA probes have been optimized with Enzo's AMPIVIEW® and POLYVIEW® Detection Systems combined with HIGHDEF® chromogens. Use of other linkers or polymer-based reagents will require prior optimization. Autostainers such as LEICA BOND or Roche DISCOVERY/ULTRA can be used but may require prior optimization



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)