

# AMPIVIEW<sup>®</sup> HPV

## High-Risk RNA Probes Set

**A mixture of digoxigenin-labeled RNA probes for the detection of HPV 16/18/31/33/51 for *in situ* hybridization.**

AMPIVIEW<sup>®</sup> RNA probes are uniquely designed with the precision of targeted, sequence-specific RNA probes, powered by Enzo's LoopRNA<sup>™</sup> ISH technology to deliver superior sensitivity. The AMPIVIEW<sup>®</sup> HPV High-Risk RNA Probe is a mixture of digoxigenin-labeled probes targeting HPV type 16, HPV type 18, HPV 31, HPV 33 and HPV 51 for *in situ* hybridization. The probes are formulated in a buffered formamide solution with hybridization enhancers.

Citations: 2

[View Online »](#)

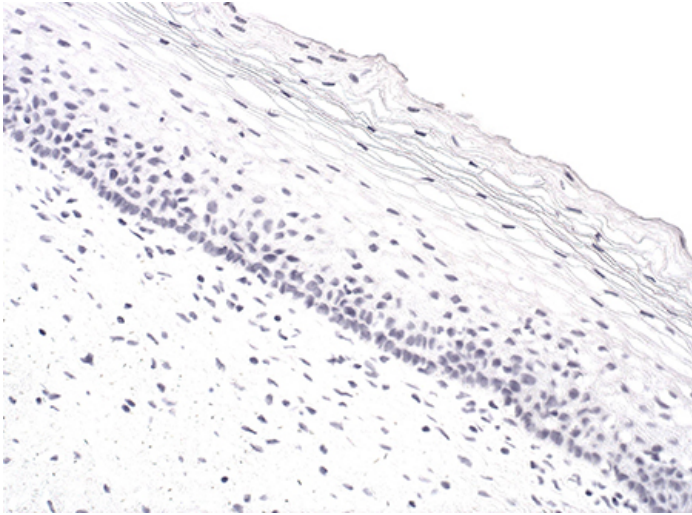
### Ordering Information

[Order Online »](#)

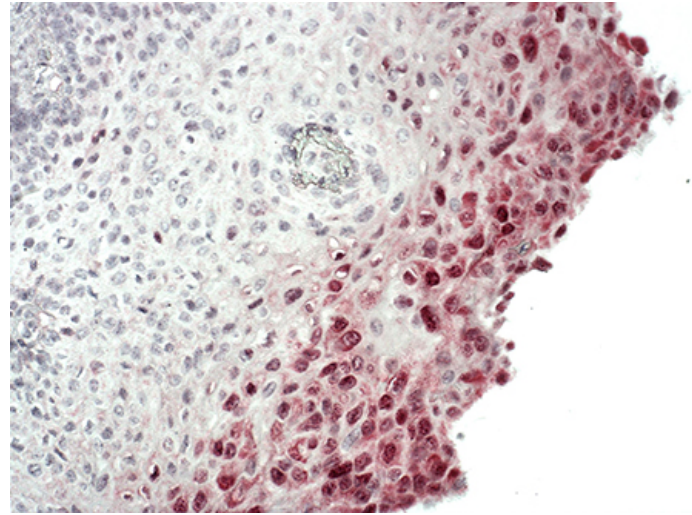
ENZ-GEN148-2000	2ml
-----------------	-----

### Manuals, SDS & CofA

[View Online »](#)



Normal cervical tissue tested negative for HPV 16/18/31/33/51 as no signal is detected with AMPIVIEW™ HPV High Risk RNA Probes (ENZ-GEN148)



HPV infected cervical tissue tested positive for HPV 16/18/31/33/51 (red) detected with AMPIVIEW™ HPV High Risk RNA Probes (ENZ-GEN148)

## Handling & Storage

Use/Stability	Aliquot and store at -20°C or -80°C. Under these conditions, products are stable until its expiration dates.
Handling	Allow contents to warm up to room temperature prior to use.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

**Application** ISH (in situ hybridization)

**Application Notes**

AMPIVIEW<sup>®</sup> HPV High-Risk RNA probes have been designed to target the human papillomavirus (HPV) type 16 and type 18 for its detection in tissues and cells. AMPIVIEW<sup>®</sup> HPV High-Risk RNA probes are conjugated with digoxigenin and have been optimized to produce clear results with Enzo's nanopolymer-based detection system, POLYVIEW<sup>®</sup> Plus, DIGX<sup>®</sup> anti-digoxigenin linker and , high-quality chromogens and counterstains, HIGHDEF<sup>®</sup> to produce clear results that can be visualized with a light microscope.

**Contents**

AMPIVIEW<sup>®</sup> HPV High-Risk RNA Probes  
AMPIVIEW<sup>®</sup> Hybridization Buffer (ENZ-ACC152)

**Technical Info / Product Notes**

AMPIVIEW<sup>®</sup> HPV High-Risk RNA Probes are optimized for detection with either POLYVIEW<sup>®</sup> PLUS detection reagents, HIGHDEF<sup>®</sup> chromogens and counterstains (not included). Dilutions and concentration optimization can be done with AMPIVIEW<sup>®</sup> hybridization buffer, included in this set.

Last modified: October 17, 2024



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-  
fr@enzolifesciences.com](mailto:info-fr@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)