

AMPIVIEW® HPV 31/33/51 RNA Probes Set

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

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. The AMPIVIEW® HPV 31/33/51 RNA Probe is a mixture of digoxigenin-labeled probes targeting HPV type 31, HPV type 33 and HPV type 51 for *in situ* hybridization. The probes are formulated in a buffered formamide solution with hybridization enhancers.

Citations: 2

[View Online »](#)

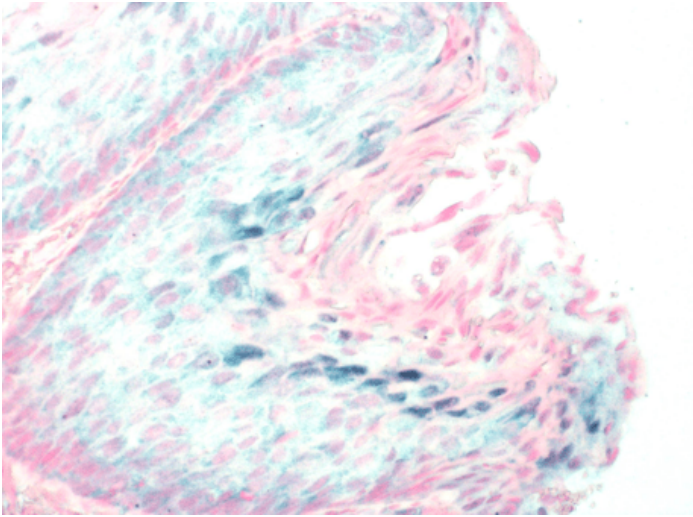
Ordering Information

[Order Online »](#)

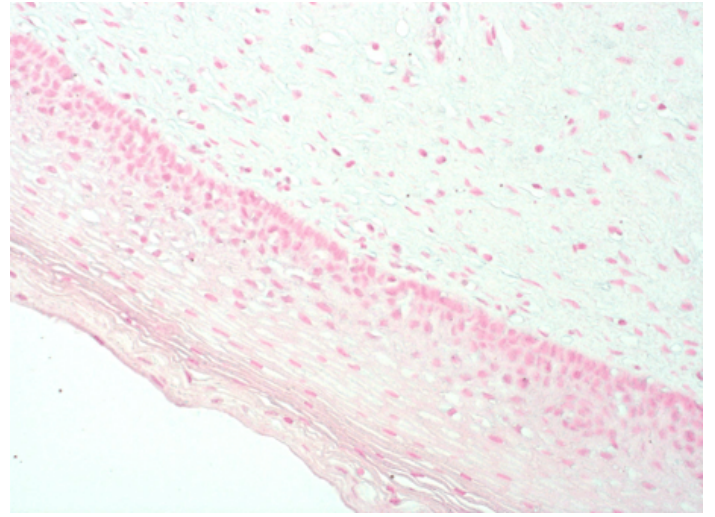
ENZ-GEN147-2000	2ml
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



HPV infected cervical tissue tested positive for HPV 31/33/51 (blue) detected with AMPIVIEW™ HPV 31/33/51 (ENZ-GEN147)



Normal cervical tissue tested negative for HPV 31/33/51 as no signal is detected with AMPIVIEW™ HPV 31/33/51 (ENZ-GEN147)

Handling & Storage

Use/Stability	Stable for at least a year after receipt when stored as recommended. Use probe immediately after dilution with AMPIVIEW® Hybridization Buffer (included).
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® HPV 31/33/51 RNA probes have been designed to target the human papillomavirus (HPV) type 31, type 33 and type 51 for its detection in tissues and cells. AMPIVIEW® HPV 31/33/51 RNA probes are conjugated with digoxigenin and have been optimized to produce clear results with Enzo's nanopolymer-based detection system, POLYVIEW® Plus, DIGX® anti-digoxigenin linker and, high-quality chromogens and counterstains, HIGHDEF® to produce clear results that can be visualized with a light microscope.
Contents	AMPIVIEW® HPV 31/33/51 RNA Probes AMPIVIEW® Hybridization Buffer (ENZ-ACC152)
Technical Info / Product Notes	AMPIVIEW® HPV 31/33/51 RNA Probes are optimized for detection with either POLYVIEW® PLUS detection reagents, HIGHDEF® chromogens and counterstains (not included). Dilutions and concentration optimization can be done with AMPIVIEW® hybridization buffer (ENZ-ACC152), included in this set.



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)