

AMPIVIEW[®] 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[®] 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 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 »](#)

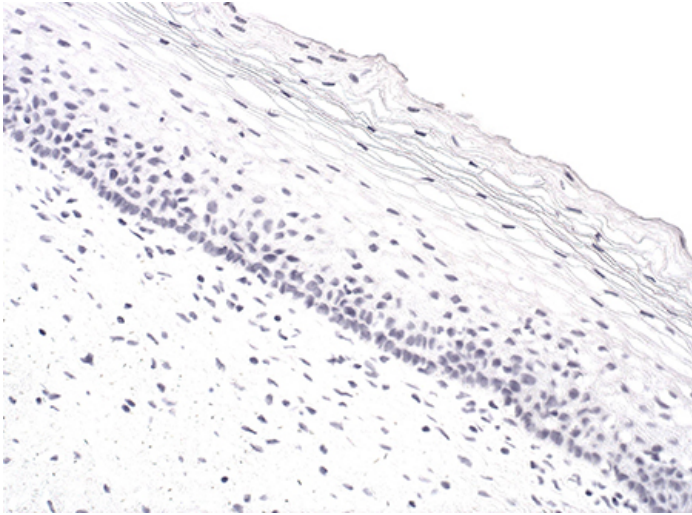
Ordering Information

[Order Online »](#)

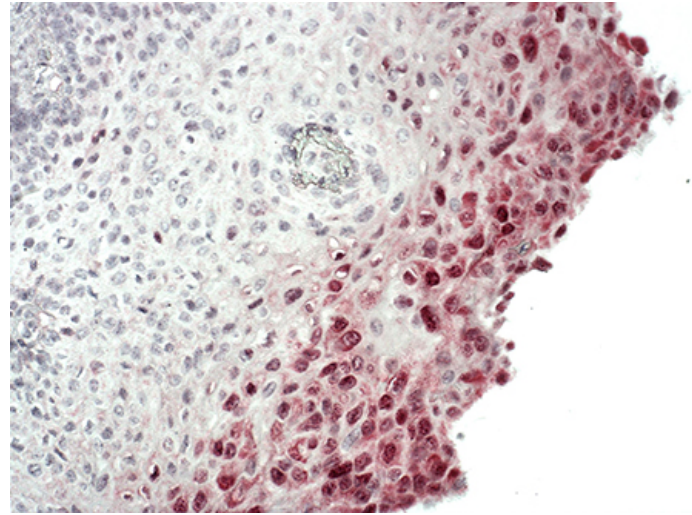
ENZ-GEN148-2000	2ml
-----------------	-----

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	<p>AMPIVIEW[®] 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.</p> <p>AMPIVIEW[®] 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[®] Plus, DIGX[®] anti-digoxigenin linker and , high-quality chromogens and counterstains, HIGHDEF[®] to produce clear results that can be visualized with a light microscope.</p>
Contents	<p>AMPIVIEW[®] HPV High-Risk RNA Probes</p> <p>AMPIVIEW[®] Hybridization Buffer (ENZ-ACC152)</p>
Technical Info / Product Notes	<p>AMPIVIEW[®] HPV High-Risk 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, included in this set.</p>



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)