

AMPIVIEW[®] HPV 6/11 RNA Probes Set

A mixture of digoxigenin-labeled HPV 6 and HPV 11-specific RNA probes 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. AMPIVIEW[®] HPV 6/11 RNA probes is a mixture of digoxigenin-labeled probes targeting HPV type 6 and HPV type 11 for *in situ* hybridization. The probes are formulated in a buffered formamide solution with hybridization enhancers.

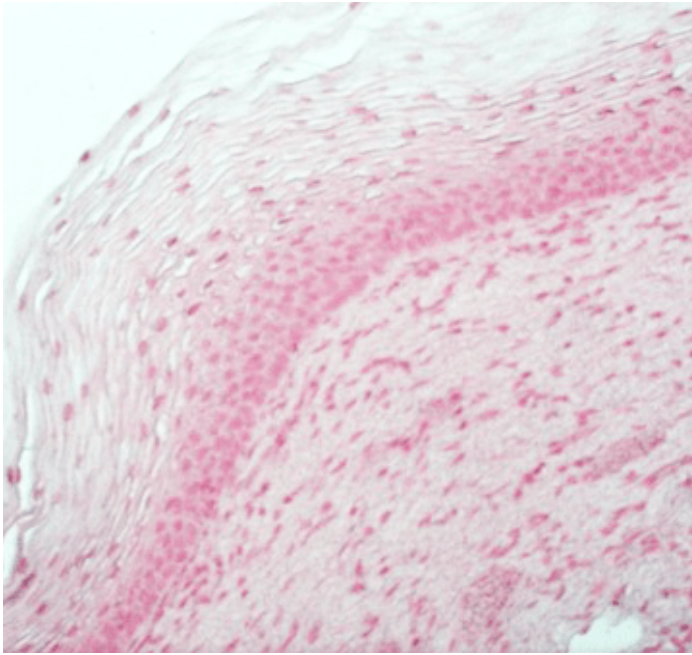
Ordering Information

[Order Online »](#)

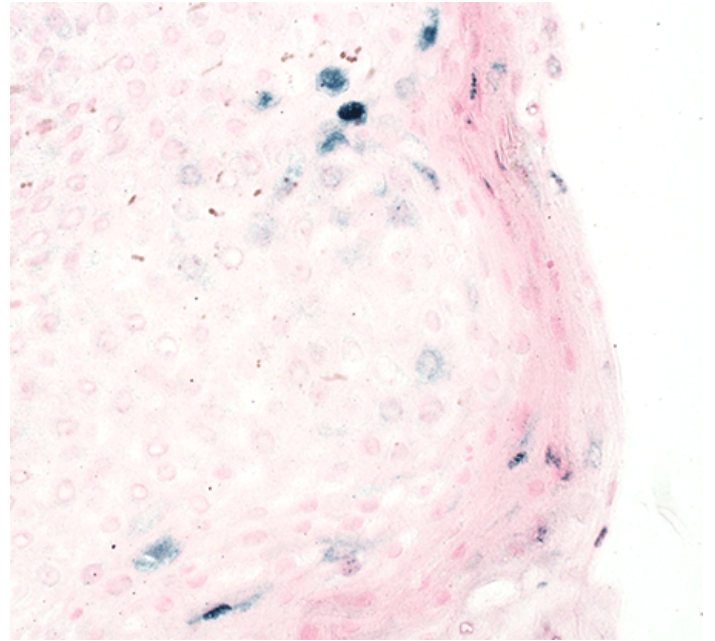
ENZ-GEN145-2000	2ml
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Normal cervical tissue shows no signal when detected with AMPIVIEW™ HPV 6/11 RNA Probes



HPV infected cervical tissue shows infection (blue) with AMPIVIEW™ HPV 6/11 RNA Probes (ENZ-GEN1465)

Handling & Storage

Use/Stability	Upon receipt, aliquot and store at -20°C. For long-term storage, store at -80°C. These products, under these conditions, are stable until the expiration date.
Handling	Allow contents to reach 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 6/11 RNA probes have been designed to target the human papillomavirus (HPV) type 6 and type 11 for its detection in tissues and cells. AMPIVIEW[®] HPV 6/11 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 HIGHDEF[®] chromogens and counterstains to produce clear results that can be visualized with a light microscope.

Contents AMPIVIEW[®] HPV 6/11 RNA Probes
AMPIVIEW[®] Hybridization Buffer (ENZ-ACC152)

Technical Info / Product Notes AMPIVIEW[®] HPV 6/11 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.



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)