

AMPIVIEW[®] HPV HR-15 (AS) DNP RNA Probes Set

A mixture of DNP-labeled RNA probes for the detection of HPV

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 HR-15 (AS) DNP RNA Probes are a mixture of DNP-labeled probes targeting HPV types 16, 18, 31, 33, 35, 39, 45, 51, 52, 56, 58, 59, 66, 68 and 82 for *in situ* hybridization. The probes are formulated in a buffered formamide solution with hybridization enhancers.

Ordering Information

[Order Online »](#)

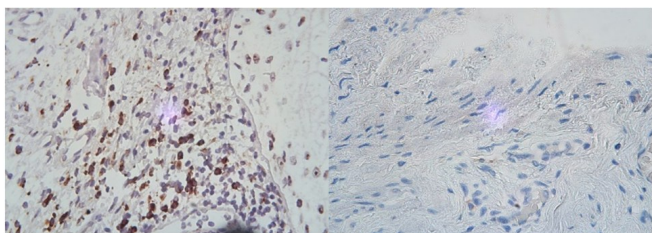
ENZ-GEN320-2000	2ml
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)

A.

B.



HPV 16, 18, 31, 33, 35, 39, 45, 51, 52, 56, 58, 59, 66, 68 and/or 82 (brown) detected with AMPIVIEW® HPV HR-15 (AS) DNP RNA probes in A. HPV infected cervical cancer tissue sample and B. non-infected cervical tissue sample.

Handling & Storage

Use/Stability	Aliquot and store at -20°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	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application ISH (in situ hybridization)

Application Notes AMPIVIEW[®] HPV HR-15 (AS) DNP RNA probes have been designed to target the nucleic acid of E6/E7 for human papillomavirus (HPV) genotype 16, 18, 31, 33, 35, 39, 45, 51, 52, 56, 58, 59, 66, 68 and 82 for its detection in tissues and cells. AMPIVIEW[®] HPV HR-15 (AS) DNP RNA probes are conjugated with dinitrophenol and have been optimized to be amplified with AMPIVIEW[®] Rabbit anti-DNP linker and detected with POLYVIEW[®] PLUS combined with HIGHDEF[®] chromogens and counterstain to produce clear results that can be visualized with a light microscope.

Contents AMPIVIEW[®] HPV HR-15 (AS) DNP RNA Probes (ENZ-GEN319)
AMPIVIEW[®] Hybridization Buffer (ENZ-ACC152)

Technical Info / Product Notes AMPIVIEW[®] HPV HR-15 (AS) DNP RNA Probes are optimized for amplification with AMPIVIEW[®] Rabbit anti-DNP Linker and detection with POLYVIEW[®] PLUS detection reagents combined with 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-
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)