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

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

AMPINEW® 33NA probest are uniquely designed with/the procession 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

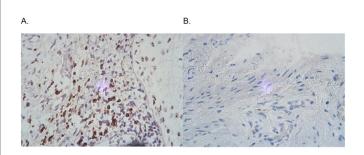
Order Online »

ENZ-GEN320-2000

2ml

Manuals, SDS & CofA

View Online »



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 NotesAMPIVIEW® 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)

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 inlcuded). Dilutions

and concentration optimization can be done with AMPIVIEW® hybridization buffer, included in this set.



Technical Info / Product Notes

ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoBelgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com