

AMPIVIEW[®] Ubiquitin and NSP Dig Control Probes Kit

Positive and negative control RNA probes labeled with digoxigenin for *in situ*

AMPIVIEW[®] Ubiquitin (AS) Dig RNA Probes are digoxigenin-labeled RNA probes targeting ubiquitin nucleic acid (DNA/RNA or RNA) for *in situ* hybridization in tissues and cells and can be used as a positive control probe. AMPIVIEW[®] NSP Dig RNA Probes are digoxigenin-labeled probes designed with a non-specific sequence that can be used as a negative control. Both probes are formulated in a buffered formamide solution with hybridization enhancers.

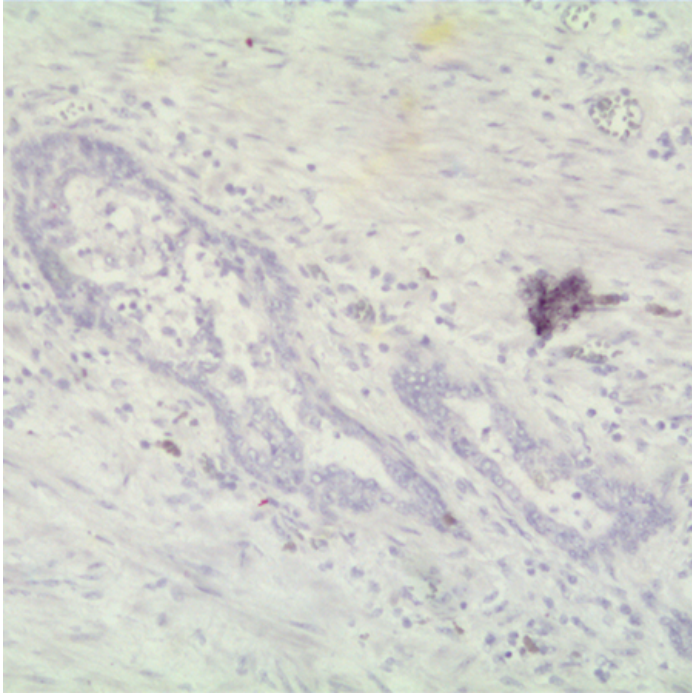
Ordering Information

[Order Online »](#)

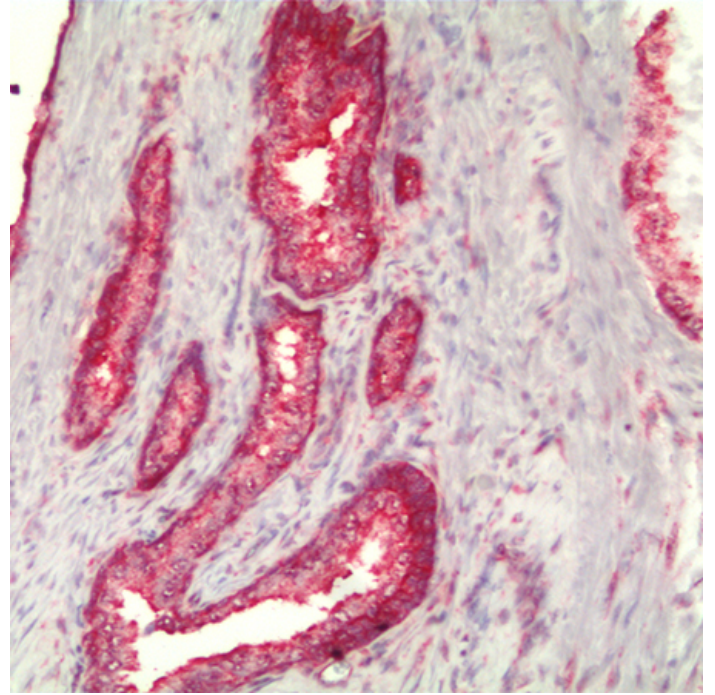
ENZ-KIT223-0040	40 tests
-----------------	----------

Manuals, SDS & CofA

[View Online »](#)



No signal detected in prostate tissue with AMPIVIEW® NSP Dig RNA probes, DIGX® anti-digoxigenin linker, POLYVIEW® PLUS AP Detection reagent combined with HIGHDEF® Red AP IHC chromogen and counterstained with HIGHDEF® Hematoxylin.



Ubiquitin (red) detected in prostate tissue with AMPIVIEW® Ubiquitin (AS) Dig RNA probes, DIGX® anti-digoxigenin linker, POLYVIEW® PLUS AP Detection reagent combined with HIGHDEF® Red AP IHC chromogen and counterstained with HIGHDEF® Hematoxylin.

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 AMPIVIEW[®] Ubiquitin (AS) Dig RNA probes have been designed to target ubiquitin in tissues and cells. AMPIVIEW[®] Ubiquitin (AS) Dig RNA probes are conjugated with digoxigenin and have been optimized to produce clear results with Enzo's DIGX[®] anti-digoxigenin linker and nanopolymer-based detection systems, POLYVIEW[®] PLUS, combined with HIGHDEF[®] chromogens and counterstain to produce clear results that can be visualized with a light microscope.

Contents 2 x 1 mL AMPIVIEW[®] Ubiquitin (AS) Dig RNA Probes (2µg/mL)
2 x 1 mL AMPIVIEW[®] NSP Dig RNA Probes (RTU)
1 x 2 mL AMPIVIEW[®] Hybridization Buffer (1X)

Technical Info / Product Notes Ubiquitin is a regulatory protein found in nearly all eukaryotic organisms. It is highly conserved and virtually identical across all forms of life and ideal to use as a positive control.

AMPIVIEW[®] Ubiquitin (AS) Dig RNA Probes are optimized for detection of ubiquitin RNA or RNA/DNA in FFPE tissues or cells. AMPIVIEW[®] NSP Dig RNA Probes are ready-to-use probes that have been designed with a non-specific sequence. Both probes can be detected with DIGX[®] anti-digoxigenin linker, POLYVIEW[®] PLUS detection reagents combined with HIGHDEF[®] chromogens and counterstains (linker and detection solutions not included). Dilutions and concentration optimization can be done with AMPIVIEW[®] Hybridization Buffer (1X), included in this kit.



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)