AMPIVIEW[®] Ubiquitin and NSP Dig Control Probes Kit

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

hypothem at biquities (AS) in the 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

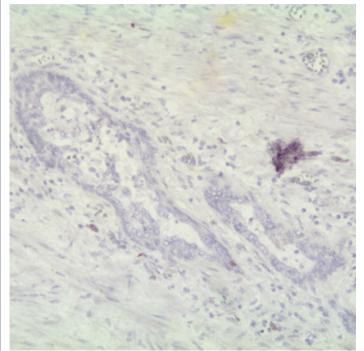
Order Online »

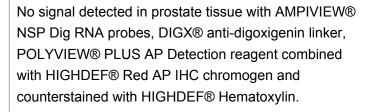
ENZ-KIT223-0040

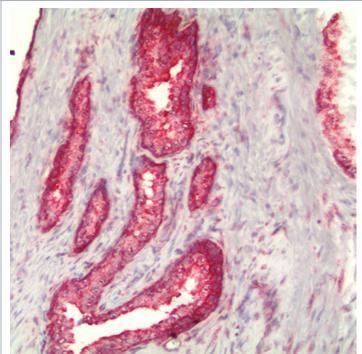
40 tests

Manuals, SDS & CofA

View Online »







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 NotesAMPIVIEW[®] 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 nanopolymner-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

 $(2\mu g/mL)$

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.

Last modified: October 17, 2024

