

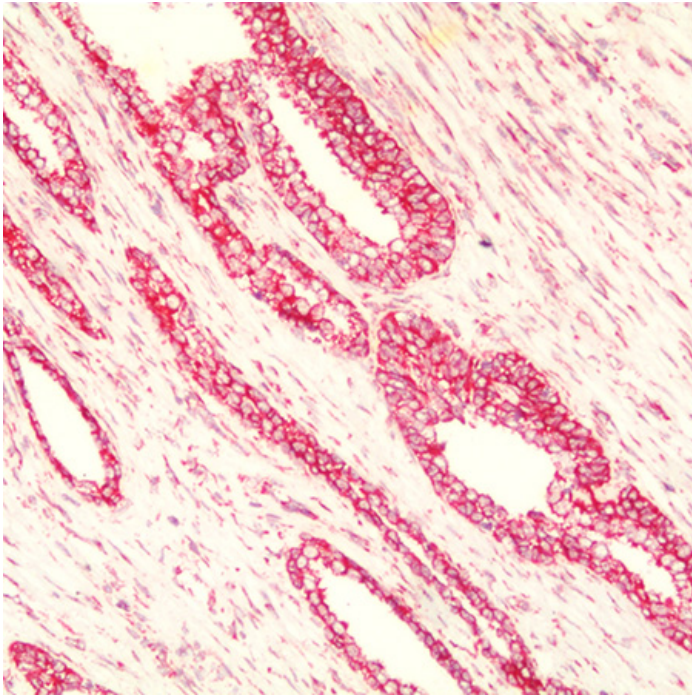
AMPIVIEW® GAPDH and NSP Dig Control Probes Kit

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

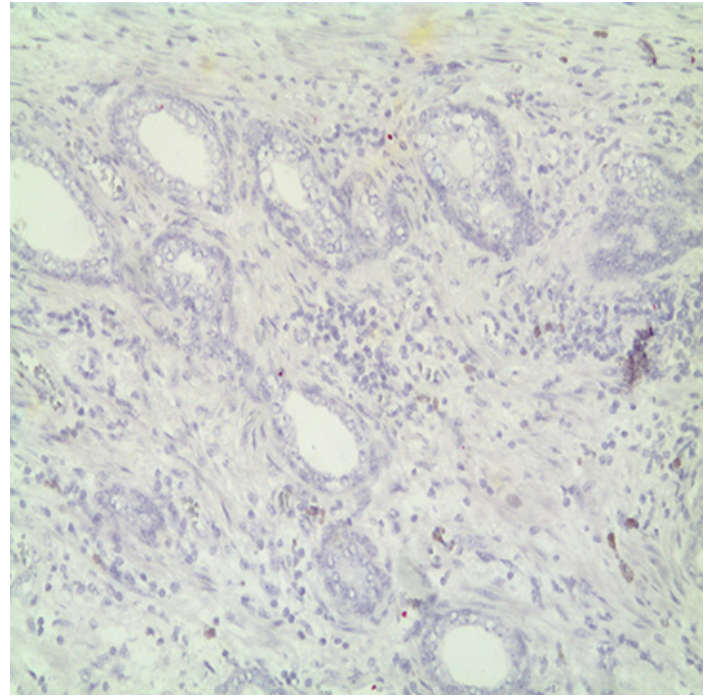
AMPIVIEW® GAPDH (AS) Dig RNA Probes are digoxigenin-labeled RNA probes targeting GAPDH nucleic acid (DNA/RNA or RNA) for *in situ* hybridization in tissues. GAPDH has been used as a housekeeping gene due to its ubiquitous nature. AMPIVIEW® NSP Dig RNA probes have been designed with a non-specific sequence to use as a negative control. The probes are formulated in a buffered formamide solution with hybridization enhancers.

Manuals, SDS & CofA

[View Online »](#)



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



No signal detected in prostate tissue with AMPIVIEW[®],_φ 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.

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[®] GAPDH and NSP Dig Control Probes Kit contains AMPIVIEW[®] GAPDH (AS) Dig RNA probes (positive control) and NSP Dig RNA probes (negative control) for *in situ* hybridization in tissues and cells. AMPIVIEW[®] GAPDH (AS) Dig RNA probes have been designed to target human glyceraldehyde-3-phosphate dehydrogenase (GAPDH) in tissues and cells. AMPIVIEW[®] NSP Dig RNA probes have been designed with a non-specific sequence to use as a negative control. Both probes have a digoxigenin label and have been optimized to produce clear results with Enzo's DIGX[®] anti-digoxigenin linker and nanopolymer-based detection systems such as 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[®] GAPDH (AS) Dig RNA Probes (2µg/mL)
2 x 1 mL AMPIVIEW[®] NSP Dig RNA Probes (RTU)
1 x 2 mL AMPIVIEW[®] Hybridization Buffer (1X)

Glyceraldehyde-3-phosphate dehydrogenase (GAPDH) is a glycolytic and key regulatory enzyme and most commonly used as a housekeeping gene.

AMPIVIEW[®] GAPDH (AS) Dig RNA Probes are optimized for detection of GAPDH 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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com