

AMPIVIEW[®] Ubiquitin (AS) Dig RNA Probes Set

A digoxigenin-labeled RNA probe for the *in situ* hybridization (ISH) detection ubiquitin

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. The probes are formulated in a buffered formamide solution with hybridization enhancers.

Ordering Information

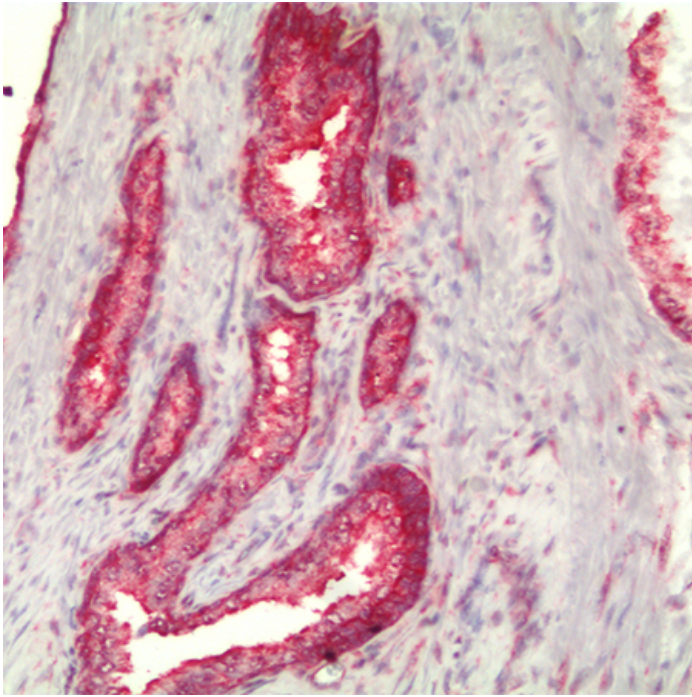
[Order Online »](#)

ENZ-GEN125-2000	2ml
-----------------	-----

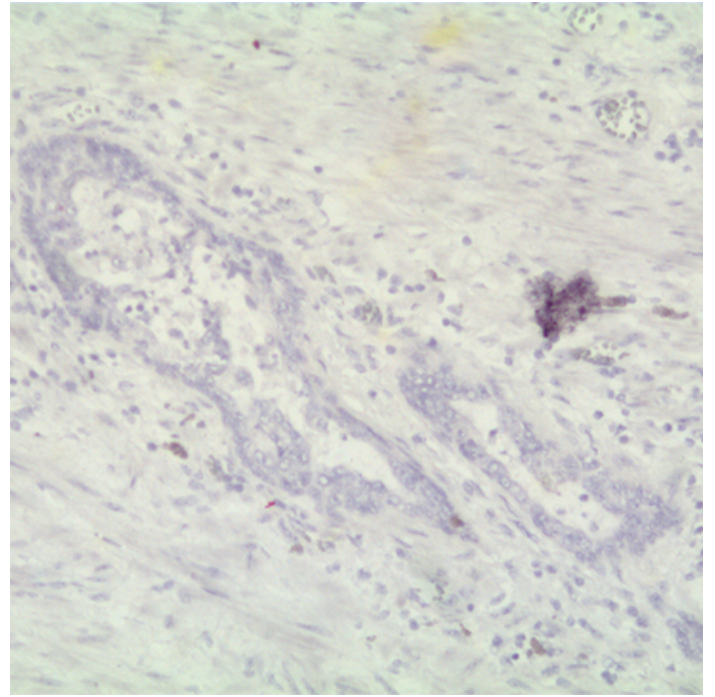
Manuals, SDS & CofA

[View Online »](#)

- Validated on the [Roche DISCOVERY ULTRA Platform](#)



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.



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.

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	<p>AMPIVIEW® Ubiquitin (AS) Dig RNA probes have been designed to target ubiquitin in tissues and cells.</p> <p>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.</p>
Contents	<p>2 x 1 mL AMPIVIEW® Ubiquitin (AS) Dig RNA Probes (2µg/mL)</p> <p>1 x 2 mL AMPIVIEW® Hybridization Buffer (1X)</p>

Ubiquitin is a regulatory protein found in nearly all eukaryotic organism. 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 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 set.

Protocol for the Roche DISCOVERY ULTRA Platform:

[Download the protocol](#)

Last modified: October 17, 2024



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-
be@enzolifesciences.com](mailto:info-be@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)