

AMPIVIEW® EGFR (AS) Dig RNA Probes Set

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

AMPIVIEW® EGFR (AS) Dig RNA Probes are digoxigenin-labeled RNA probes targeting EGFR 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

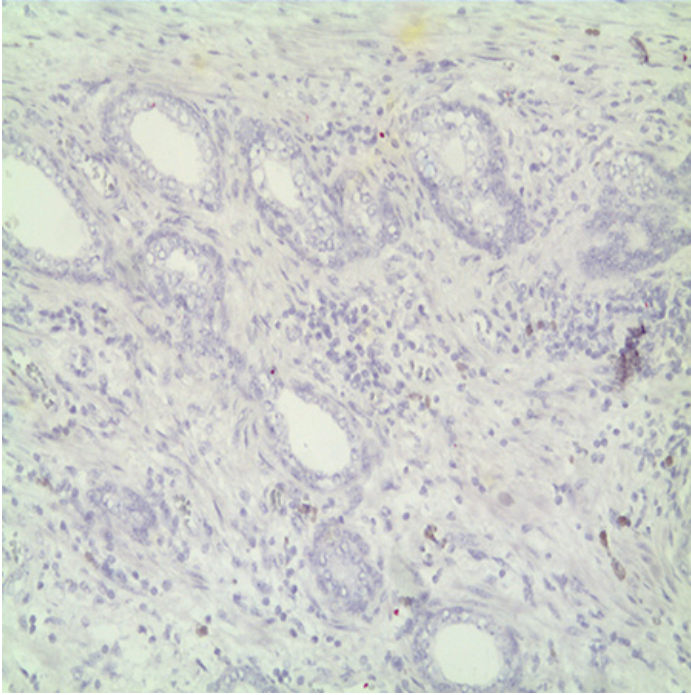
[Order Online »](#)

ENZ-GEN129-2000	2ml
-----------------	-----

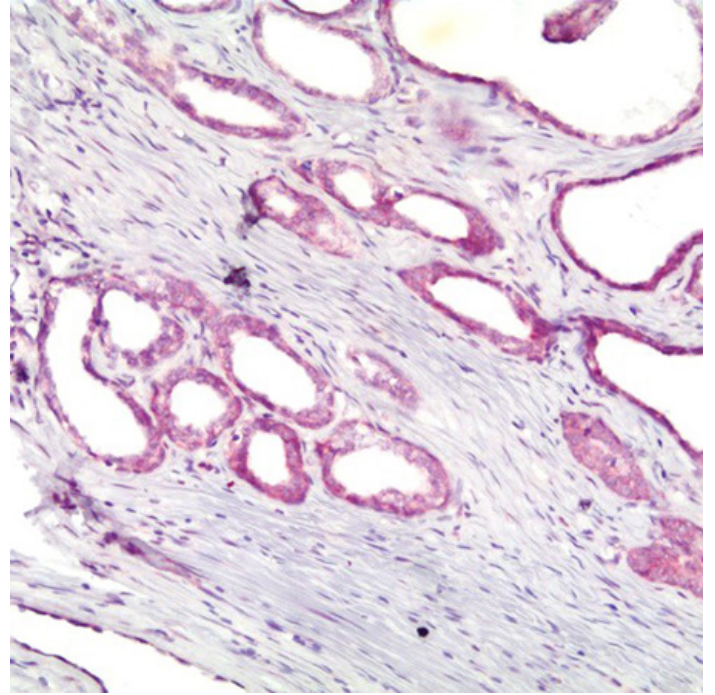
Manuals, SDS & CofA

[View Online »](#)

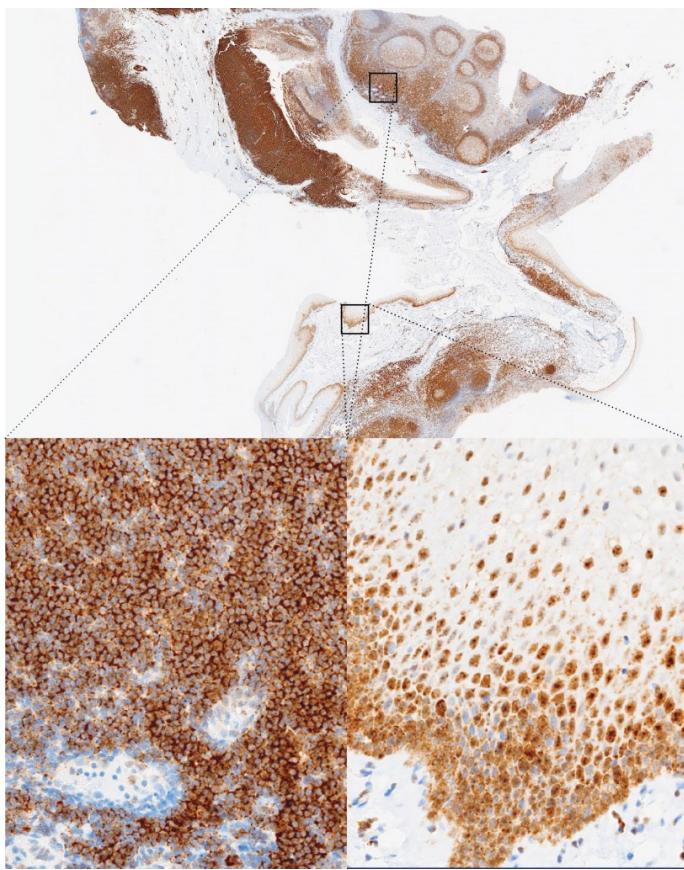
- Validated on the [Roche DISCOVERY ULTRA Platform](#)



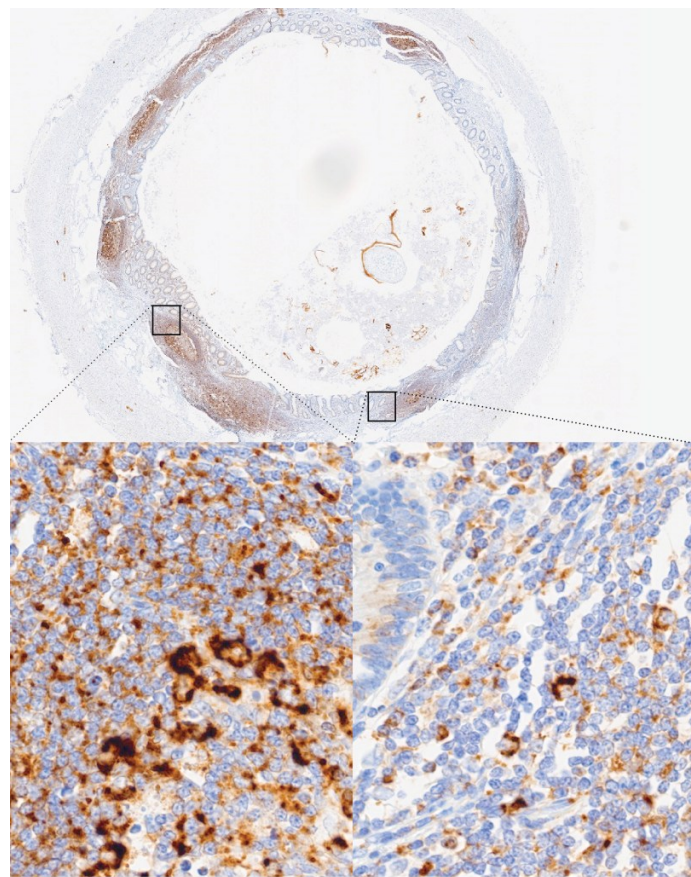
AMPIVIEW® NSP Dig RNA probes (no signal) detected in prostate hyperplasia tissue specimen with DIGX® rabbit anti-digoxigenin linker (ENZ-ABS303), POLYVIEW® PLUS AP (anti-rabbit) (ENZ-ACC110), combined with HIGHDEF® Red Chromogen/Substrate and Hematoxylin counterstain (ENZ-ACC106). AMPIVIEW® NSP Dig RNA probes are negative control probes.



EGFR (red) detected in prostate hyperplasia tissue specimen with AMPIVIEW® EGFR Dig RNA probes, DIGX® rabbit anti-digoxigenin linker (ENZ-ABS303), POLYVIEW® PLUS AP (anti-rabbit) (ENZ-ACC110) combined with HIGHDEF® Red Chromogen/Substrate and Hematoxylin counterstain (ENZ-ACC106).



EGFR (brown) was detected in tonsil tissue with AMPIVIEW® EGFR (AS) Dig RNA Probes (ENZ-GEN129) using Ventana detection reagents on the Ventana Discovery Ultra.



EGFR (brown) was detected in appendix tissue with AMPIVIEW® EGFR (AS) Dig RNA Probes (ENZ-GEN129) using Ventana detection reagents on the Ventana Discovery Ultra.

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 [®] EGFR (AS) Dig RNA probes have been designed to target the EGFR or HER-1 DNA and RNA in tissues and cells. AMPIVIEW [®] EGFR (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 [®] EGFR (AS) Dig RNA Probes (2µg/mL) 1 x 2 mL AMPIVIEW [®] Hybridization Buffer (1X)
Technical Info / Product Notes	AMPIVIEW [®] EGFR (AS) Dig RNA Probes are optimized for detection of EGFR 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)