

AMPIVIEW® Ki67 (AS) DNP RNA Probes Set

A DNP-labeled RNA probe for the detection of human Ki67 for *in situ* hybridization.

AMPIVIEW® RNA probes are uniquely designed with the precision of targeted, sequence-specific RNA probes, powered by Enzo's LoopRNA™ ISH technology to deliver superior sensitivity. AMPIVIEW® Ki67 (AS) DNP RNA Probes are DNP-labeled probes targeting Human Ki67 for *in situ* hybridization. The probes are formulated in a buffered formamide solution with hybridization enhancers.

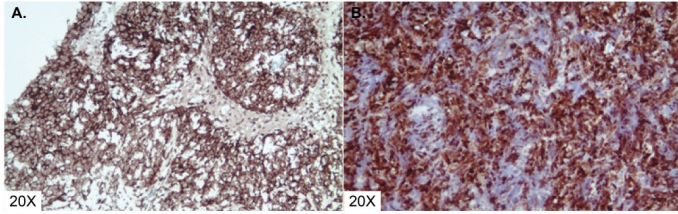
Ordering Information

[Order Online »](#)

ENZ-GEN304-2000	2ml
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Ki-67 (brown) was detected in prostate cancer tissues with A.-B. AMPIVIEW® Ki67 (AS) DNP RNA probes amplified with AMPIVIEW® rabbit anti-DNP linker (ENZ-ABS928) and detected with POLYVIEW® PLUS (anti-r) HRP reagent (ENZ-ACC103) combined with HIGHDEF® DAB chromoge/substrate (ENZ-ACC105) and counterstained with HIGHDEF® hematoxylin (ENZ-ACC106).

Handling & Storage

Use/Stability	Aliquot and store at -20°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	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application ISH (in situ hybridization)

Application Notes

AMPIVIEW® Ki67 (AS) Dig RNA probes have been designed to target the nucleic acid of Marker of Proliferation Ki-67 (MKI67), which encodes the Ki67 protein, for its detection in tissues and cells. Ki67 is a well-known marker for cell proliferation. AMPIVIEW® Ki67 (AS) DNP RNA probes are conjugated with dinitrophenol (DNP) and have been optimized to be amplified with AMPIVIEW® Rabbit anti-DNP linker and detected with POLYVIEW® PLUS combined with HIGHDEF® chromogens and counterstain to produce clear results that can be visualized with a light microscope.

Contents

AMPIVIEW® Ki67 (AS) DNP RNA Probes (ENZ-GEN303)
AMPIVIEW® Hybridization Buffer (ENZ-ACC152)

Technical Info / Product Notes

AMPIVIEW® Ki67 (AS) DNP RNA Probes are optimized for amplification with AMPIVIEW® Rabbit anti-DNP Linker and detection with POLYVIEW® PLUS detection reagents combined with HIGHDEF® chromogens and counterstains (not included). Dilutions and concentration optimization can be done with AMPIVIEW® hybridization buffer, included in this set.



Last modified: July 1, 2025

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

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@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