

AMPIVIEW® PTEN (AS) DNP RNA Probes Set

**A DNP-labeled RNA probe for the detection
of Human PTEN 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® PTEN (AS) DNP RNA Probes are DNP-labeled probes targeting Human PTEN for *in situ* hybridization. The probes are formulated in a buffered formamide solution with hybridization enhancers.

Ordering Information

[Order Online »](#)

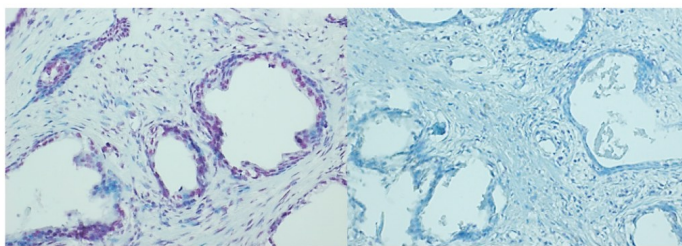
ENZ-GEN131-2000	
-----------------	--

Manuals, SDS & CofA

[View Online »](#)

A.

B.



PTEN (red) detected in prostate cancer tissue with A. AMPIVIEW® PTEN (AS) DNP RNA Probes and B. AMPIVIEW® NSP DNP RNA Probes.

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[®] PTEN (AS) DNP RNA probes have been designed to target the nucleic acid of the phosphatase and tensin homolog (PTEN) that encodes for a phosphatase protein that acts as a tumor suppressor-gene, for its detection in tissues and cells. AMPIVIEW[®] PTEN5 (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[®] PTEN (AS) DNP RNA Probes (ENZ-GEN130)
AMPIVIEW[®] Hybridization Buffer (ENZ-ACC152)

Technical Info / Product Notes AMPIVIEW[®] PTEN (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.



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