

# AMPIVIEW® p53 (AS) Dig RNA Probes Set

A digoxigenin-labeled RNA probe for the detection of human p53 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® p53 (AS) Dig RNA Probes are digoxigenin-labeled probes targeting Human p53 for *in situ* hybridization. The probes are formulated in a buffered formamide solution with hybridization enhancers.

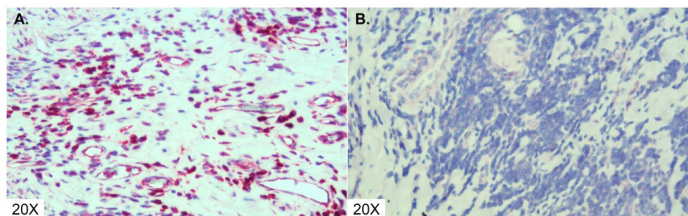
## Ordering Information

[Order Online »](#)

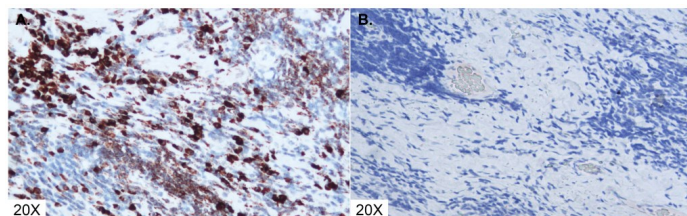
ENZ-GEN279-2000	2ml
-----------------	-----

## Manuals, SDS & CofA

[View Online »](#)



p53 (brown) detected in Mantle Lymphoma tissue with A. AMPIVIEW® p53 (AS) Dig RNA probes (Human) and B. AMPIVIEW® NSP Dig RNA Probes amplified with DIGX® rabbit anti-digoxigenin linker (ENZ-ABS303) and detected with POLYVIEW® PLUS (anti-r) AP reagent (ENZ-ACC110) combined with HIGHDEF® Red Chromogen Kit (ENZ-ACC158) and counterstained with HIGHDEF® hematoxylin (ENZ-ACC106).



p53 (brown) detected in Mantle Lymphoma tissue with A. AMPIVIEW® p53 (AS) Dig RNA probes (Human) and B. AMPIVIEW® NSP Dig RNA Probes amplified with DIGX® mouse anti-Digoxigenin linker (ENZ-ABS304) and detected with POLYVIEW® PLUS (anti-m) HRP reagent (ENZ-ACC104) combined with HIGHDEF® DAB chromogen/substrate (ENZ-ACC105) and counterstained with HIGHDEF® hematoxylin (ENZ-ACC106).

# Handling & Storage

Use/Stability	Aliquot and store at -20°C. Long term storage recommended at -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® p53 (AS) Dig RNA probes have been designed to target the nucleic acid of Tumor Protein 53 (TP53), which encodes the p53 protein, for its detection in tissues and cells. p53 protein plays a crucial role in maintaining genomic stability and prevenitng cancer by detecting DNA damage, allowing DNA repair, triggering apoptosis if damage if irreparable and preventing the proliferation of mutated cells. AMPIVIEW® p53 (AS) Dig RNA probes are conjugated with digoxigenin and have been optimized to be amplified with DIGX® anti-Digoxigenin 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® p53 (AS) Dig RNA Probes (ENZ-GEN278) AMPIVIEW® Hybridization Buffer (ENZ-ACC152)
Technical Info / Product Notes	AMPIVIEW® p53 (AS) Dig RNA Probes are optimized for amplification with DIGX® anti-digoxigenin 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-  
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)