

# AMPIVIEW® Wnt5a (AS) DNP RNA Probes Set

**A DNP-labeled RNA probe for the detection of Human Wnt5a 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® Wnt5a (AS) DNP RNA Probes are DNP-labeled probes targeting Human Wnt5a for *in situ* hybridization. The probes are formulated in a buffered formamide solution with hybridization enhancers.

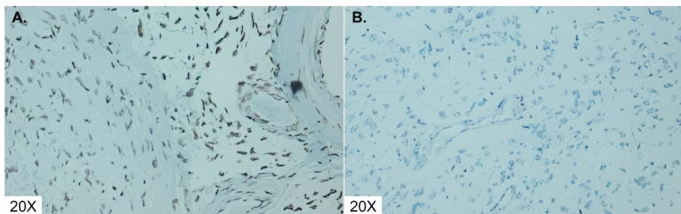
## Ordering Information

[Order Online »](#)

ENZ-GEN306-2000	2ml
-----------------	-----

## Manuals, SDS & CofA

[View Online »](#)



Wnt5a (brown) detected in Glioblastoma tissue with A. AMPIVIEW® Wnt5a (AS) DNP RNA probes (Human) and B. AMPIVIEW® NSP 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

<b>Use/Stability</b>	Aliquot and store at -20°C. Under these conditions, products are stable until its expiration dates.
<b>Handling</b>	Allow contents to warm up to room temperature prior to use.
<b>Short Term Storage</b>	-20°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

**Application** ISH (in situ hybridization)

**Application Notes** AMPIVIEW<sup>®</sup> Wnt5a (AS) DNP RNA probes have been designed to target the nucleic acid for Wnt5a, a protein encoded by the *WNT5A* gene, for its detection in tissues and cells. Wnt5a is a signaling molecule expressed embryonically during development, and has been implicated in many different types of cancer, neurodegeneration, inflammatory diseases and cardiovascular diseases. AMPIVIEW<sup>®</sup> Wnt5a (AS) DNP RNA probes are conjugated with dinitrophenol (DNP) and have been optimized to be amplified with AMPIVIEW<sup>®</sup> Rabbit anti-DNP linker and detected with POLYVIEW<sup>®</sup> PLUS combined with HIGHDEF<sup>®</sup> chromogens and counterstain to produce clear results that can be visualized with a light microscope.

**Contents** AMPIVIEW<sup>®</sup> Wnt5a (AS) DNP RNA Probes (ENZ-GEN305)  
AMPIVIEW<sup>®</sup> Hybridization Buffer (ENZ-ACC152)

**Technical Info / Product Notes** AMPIVIEW<sup>®</sup> Wnt5a (AS) DNP RNA Probes are optimized for amplification with AMPIVIEW<sup>®</sup> Rabbit anti-DNP Linker and detection with POLYVIEW<sup>®</sup> PLUS detection reagents combined with HIGHDEF<sup>®</sup> chromogens and counterstains (not included). Dilutions and concentration optimization can be done with AMPIVIEW<sup>®</sup> 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)