

# AMPIVIEW<sup>®</sup> Wnt5a (AS) Dig RNA Probes (Rat) Set

A digoxigenin-labeled RNA probe for the *in situ* hybridization (ISH) detection Wnt5a in

rat cells and tissues

AMPIVIEW<sup>®</sup> Wnt5a (AS) Dig RNA Probes (rat) are digoxigenin-labeled RNA probes targeting Wnt5a for *in situ* hybridization in tissues and cells. The probes are formulated in a buffered formamide solution with hybridization enhancers.

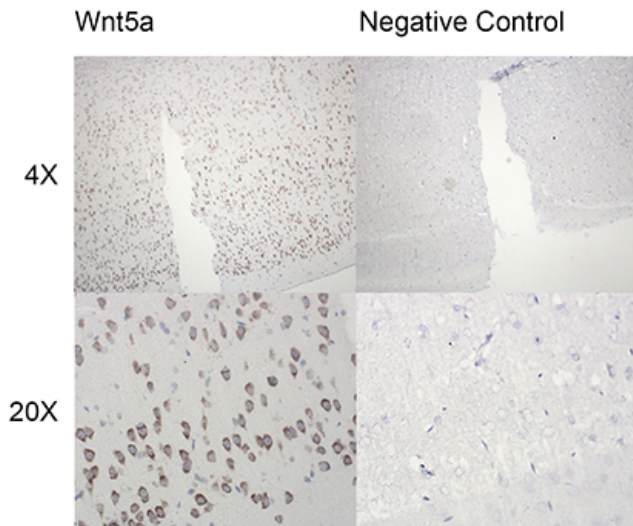
## Ordering Information

[Order Online »](#)

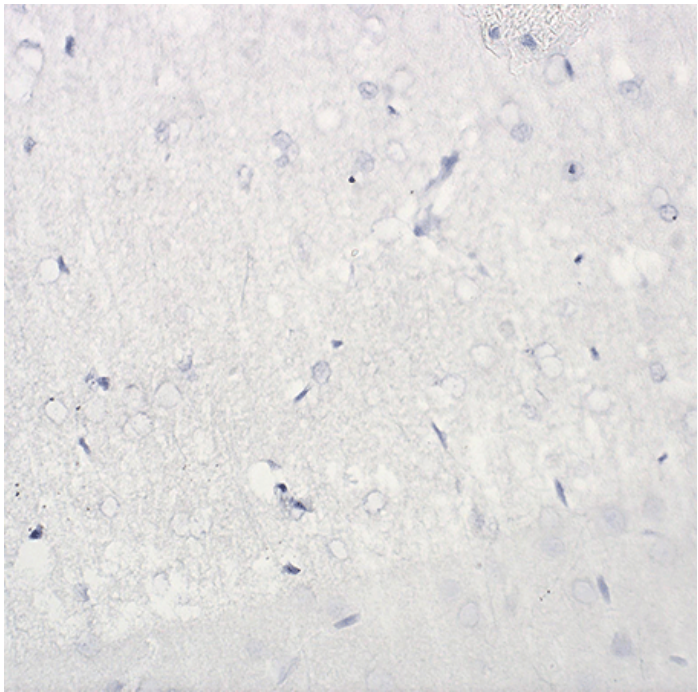
ENZ-GEN137-2000	2ml
-----------------	-----

## Manuals, SDS & CofA

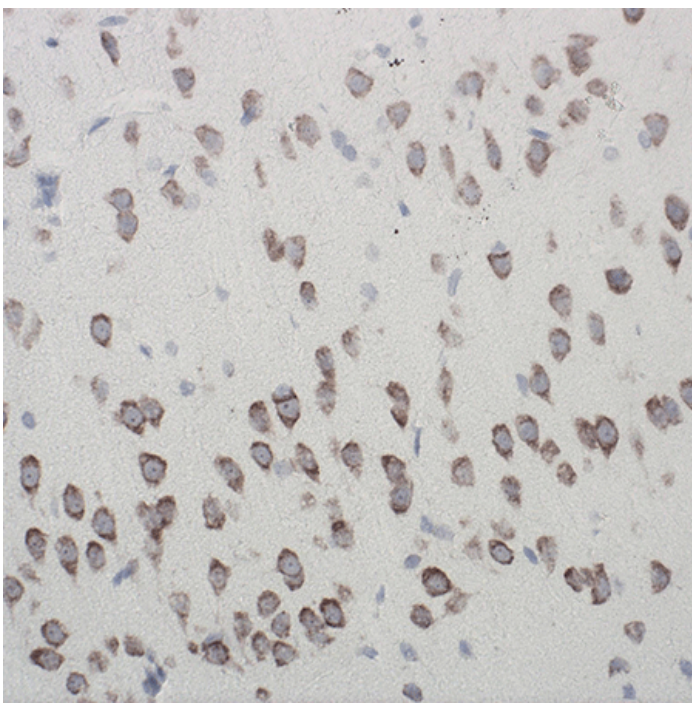
[View Online »](#)



4X and 20X images of Wnt5a detection in rat brain tissue with AMPIVIEW<sup>®</sup> Wnt5a (AS) Dig RNA probes (rat) (left column) (brown) and no probes as a negative control (right column) (no signal) with DIGX<sup>®</sup> rabbit anti-digoxigenin linker, POLYVIEW<sup>®</sup> PLUS HRP (anti-rabbit), HIGHDEF<sup>®</sup> DAB chromogen/substrate and HIGHDEF<sup>®</sup> Hematoxylin.



20X image of Wnt5a detection (no signal) in rat brain tissue with no probes as a negative control with DIGX<sup>®</sup> rabbit anti-digoxigenin linker, POLYVIEW<sup>®</sup> PLUS HRP (anti-rabbit), HIGHDEF<sup>®</sup> DAB chromogen/substrate and HIGHDEF<sup>®</sup> Hematoxylin.



20 X image of Wnt5a detection (brown) in rat brain tissue with AMPIVIEW<sup>®</sup> Wnt5a (AS) Dig RNA probes (rat) with DIGX<sup>®</sup> rabbit anti-digoxigenin linker, POLYVIEW<sup>®</sup> PLUS HRP (anti-rabbit), HIGHDEF<sup>®</sup> DAB chromogen/substrate and HIGHDEF<sup>®</sup> Hematoxylin.

## Handling & Storage

<b>Use/Stability</b>	Aliquot and store at -20°C or -80°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>	-80°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) Dig RNA probes have been designed to target rat Wnt5a, a signaling molecule expressed embryonically during development. In addition to various aspects of development, Wnt5a has been implicated in many different cancers, cardiovascular and neurodegenerative diseases.

AMPIVIEW<sup>®</sup> Wnt5a (AS) Dig RNA probes are conjugated with digoxigenin and have been optimized to produce clear results with Enzo's DIGX<sup>®</sup> anti-digoxigenin linker and nanopolymer-based detection systems, 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** 2 x 1 mL AMPIVIEW<sup>®</sup> Wnt5a (AS) Dig RNA Probes (Rat) (2µg/mL)  
1 x 2 mL AMPIVIEW<sup>®</sup> Hybridization Buffer (1X)

**Technical Info / Product Notes** One of the key signaling pathways that are known to play a crucial role in regulating various aspects of neurogenesis is Wnt signaling. Wnt5a is a signaling molecule expressed embryonically during gastrulation in various developing body regions. The role of Wnt5a has been extensively explored in various aspects of development. Loss of Wnt5a results in cerebellar hypoplasia and depletion of GABAergic and glutamatergic neurons. In addition to development, Wnt5a is implicated in many different types of cancer and cardiovascular disease, such as atherosclerosis.

AMPIVIEW<sup>®</sup> Wnt5a (AS) Dig RNA Probes (Rat) are optimized for detection of nestin RNA or RNA/DNA in FFPE tissues or cells with DIGX<sup>®</sup> anti-digoxigenin linker, POLYVIEW<sup>®</sup> PLUS detection reagents combined with HIGHDEF<sup>®</sup> chromogens and counterstains (linker and detection solutions not included). Dilutions and concentration optimization can be done with AMPIVIEW<sup>®</sup> Hybridization Buffer (1X), included in this set.



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

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 90 26 8888  
info-  
usa@enzolifesciences.com

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
info-  
be@enzolifesciences.com

France  
Phone: +33 1 72 44 633  
info-  
fr@enzolifesciences.com

Germany  
Phone: +49 7621 5500 622  
info-  
de@enzolifesciences.com

UK & Ireland  
Phone: +44 1552 224949  
0845 601 1488  
Phone: +44 1552 224949  
info-  
uk@enzolifesciences.com