# AMPIVIEW® Wnt5a (AS) Dig RNA Probes (Rat) Set

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

RNA probes targeting Wnt5a (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**

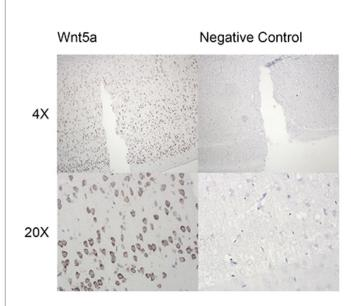
Order Online »

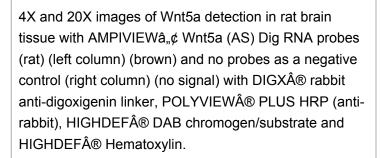
ENZ-GEN137-2000

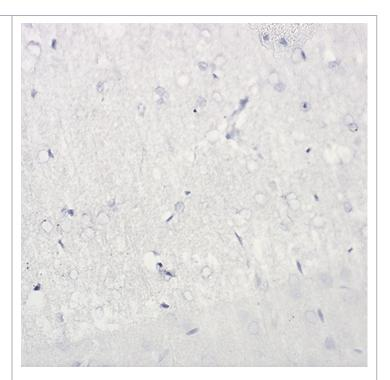
2ml

Manuals, SDS & CofA

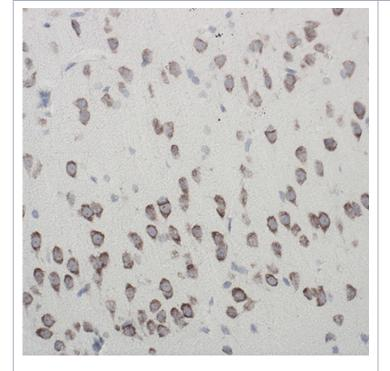
**View Online** »







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



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

#### **Handling & Storage**

**Use/Stability** Aliquot and store at -20°C or -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<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 nanopolymner-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)

 $(2\mu g/mL)$ 

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.

Last modified: October 17, 2024