AMPIVIEW[®] Nestin (AS) Dig RNA Probes (Rat) Set

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

RNA probes targeting nestin for *in situ* hybridization in tissues and cells. The probes are formulated in a buffered formamide solution with hybridization enhancers.

Ordering Information

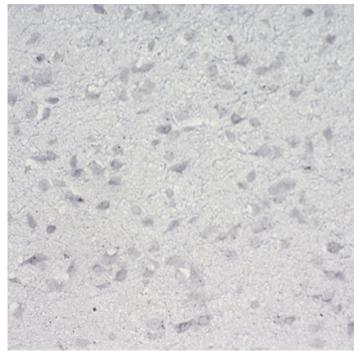
Order Online »

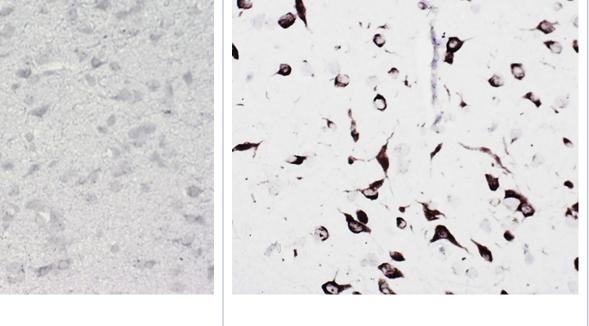
ENZ-GEN135-2000

2ml

Manuals, SDS & CofA

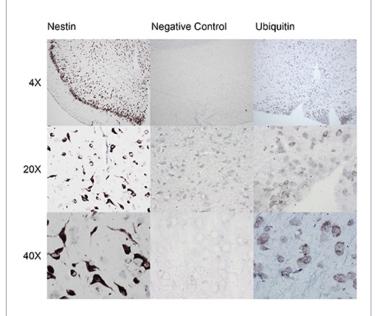
View Online »





Nestin (no signal) detection in rat brain tissue with no probes, amplified with DIGX® anti-digoxigenin linker, detected with POLYVIEW® PLUS HRP combined with HIGHDEF® DAB chromogen and counterstained with HIGHDEF® hematoxylin.

Nestin (brown) detection in rat brain tissue with AMPIVIEWâ,,¢ Nestin (AS) Dig RNA probes (rat) (ENZ-GEN135), amplified with DIGX® anti-digoxigenin linker, detected with POLYVIEW® PLUS HRP combined with HIGHDEF® DAB chromogen and counterstained with HIGHDEF® hematoxylin.



Nestin detection in rat brain tissue with AMPIVIEWâ,¢
Nestin (AS) Dig RNA probes (rat) (ENZ-GEN135)
(brown) (left column) and no probes (no signal) (middle column). Ubiquitin (brown) detection was used as a positive control with AMPIVIEWâ,¢ Ubiquitin (AS) Dig RNA probes (ENZ-GEN125) (right column). Images have been captured in 4X, 20X and 40X magnification.

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® Nestin (AS) Dig RNA probes have been designed to target rat neuroepithelial stem cell proteins (Nestin) in tissues and cells. Nestin is a type IV intermediate filament protein expressed in normal and diseased cells in different tissues and organs. Nestin is expressed by many types of cells during development. In adult cells, Nestin is reinduced during pathological situations, such as central nervous system injury and during regeneration of injured muscle tissue. Nestin is also used as a biomarker for neural stem cells and angiogenesis of proliferating endothelial cells in colorectal cancer.

AMPIVIEW[®] Nestin (AS) Dig RNA probes are conjugated with digoxigenin and have been optimized to produce clear results with Enzo's DIGX[®] anti-digoxigenin linker and nanopolymner-based detection systems, POLYVIEW[®] PLUS, combined with HIGHDEF[®] chromogens and counterstain to produce clear results that can be visualized with a light microscope.

Contents 2 x 1 mL AN

 $2 \times 1 \text{ mL AMPIVIEW}^{\textcircled{\$}} \text{ Nestin (AS) Dig RNA Probes (Rat)}$ (2 μ g/mL)

1 x 2 mL AMPIVIEW[®] Hybridization Buffer (1X)

Technical Info / Product Notes

Neuroepithelial stem cell protein (NESTIN) is a type IV intermediate filament (IF) protein expressed in normal and diseased cells in different tissues and organs. Among neural cells in the developing and adult CNS, nestin expression is thought to occur exclusively in uncommitted neural progenitor cells (NPCs). After NPCs differentiate, nestin expression typically is replaced by the expression of neuronal or glial specific markers.

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

Last modified: October 17, 2024



uk@enzolifesciences.com