AMPIVIEW® NORAD (AS) Dig RNA Probes (Mouse) Set

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

AMDIVATIVE PROPERTY PROPERTY PROPERTY AND PROPERTY PROPERTY AND INCRNA for *in situ* hybridization in tissues and cells. The probes are formulated in a buffered formamide solution with hybridization enhancers.

Ordering Information

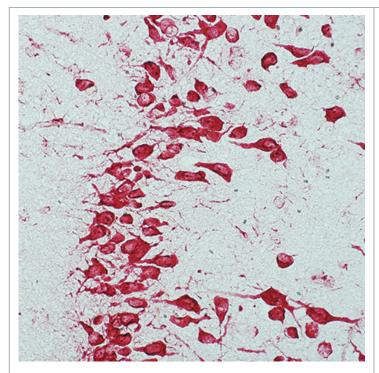
Order Online »

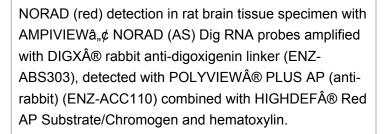
ENZ-GEN178-2000

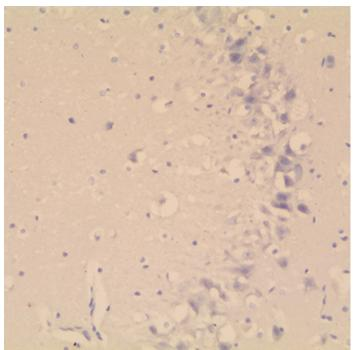
2ml

Manuals, SDS & CofA

View Online »







NORAD detection in rat brain tissue specimen AMPIVIEWâ,¢ NSP-Dig RNA probes (negative control), amplified with DIGX® rabbit anti-digoxigenin linker (ENZ-ABS303), detected with POLYVIEW® PLUS AP (anti-rabbit) (ENZ-ACC110) combined with HIGHDEF® Red AP Substrate/Chromogen and 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 NotesAMPIVIEW[®] NORAD (AS) Dig RNA probes have been

designed to target mouse or rat non-coding RNA activated by DNA damage (NORAD) in tissues and cells. NORAD is a highly abundant, highly conserved mammalian long noncoding RNA (IncRNA) that is induced after DNA damage.

AMPIVIEW® NORAD (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 AMPIVIEW[®] NORAD (AS) Dig RNA Probes

(2µg/mL)

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

Technical Info / Product Notes

NORAD is a long non-coding RNA that is induced after DNA damage. LncRNAs are identified as critical regulators in cerebral ischemia/reperfusion injury and are known to play an important role in the process of aging mammals.

AMPIVIEW[®] NORAD (AS) Dig RNA Probes are optimized for detection of NORAD 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

