AMPIVIEW® PLV (AS) Dig RNA Probes (Kingsnake) Set

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

FINAL ASIDIS RNAP TO BEST (Kingsnake) are digoxigenin-labeled RNA probes targeting the picorna-like virus nucleic acid (DNA/RNA tissues of kingsnakes or RNA) 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-GEN188-2000

2ml

Manuals, SDS & CofA

View Online »

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® PLV (AS) Dig RNA probes (Kingsnake) have

been designed to target picorna-like virus infections in tissues and cells from kingsnakes. AMPIVIEW[®] PLV (AS)

Dig RNA probes (Kingsnake) 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[®] PLV (AS) Dig RNA Probes

(Kingsnake) (2µg/mL)

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

Technical Info / Product NotesAMPIVIEW® PLV (AS) Dig RNA Probes (Kingsnake) are

optimized for detection of picorna-like virus DNA 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.



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com