AMPIVIEW® eGFP (AS) Dig RNA Probes Set

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

Probes targeting the nucleic acid (DNA/RNA or RNA) for eGFP 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-GEN352-2000

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[®] eGFP (AS) Dig RNA Probes have been

designed to target detect the nucleic acid that encodes the enhanced green fluorescent protein (eGFP). AMPIVIEW®

eGFP (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[®] eGFP (AS) Dig RNA Probes

 $(2\mu g/mL)$

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

Technical Info / Product NotesAMPIVIEW® HSP90 (AS) Dig RNA Probes (*L.giganteum*)

are optimized for detection of *L.giganteum* 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.

ENZO LIFE SCIENCE
INC.
Phone: 800.942.0430
info-

NCES, European Sales Office ENZO LIFE SCIENCES 430 (ELS) AG Phone: +41 61 926 8989 Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com