AMPIVIEW® KRT8 (AS) Dig RNA Probes (Human) Set

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

RNA probes targeting the nucleic acid (DNA/RNA or RNA) for KRT8 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-GEN182-2000

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® KRT8 (AS) Dig RNA probes (Human) have

been designed to target the nucelic acid that encodes for Keratin 8 (KRT8), a type II intermediate filament protein that plays a vital role in the structure and functioin of simple epithelial cells. Overexpression of KRT8 has been linked to tumor progression and metastasis, especially in cancer like clear cells renal carcinoma (ccRCC), breast cancer, colorectal and liver. AMPIVIEW[®] KRT8 (AS) Dig RNA probes (Human) 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

Contents 2 x 1 mL AMPIVIEW[®] KRT8 (AS) Dig RNA Probes

(Human) (2µg/mL)

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

results that can be visualized with a light microscope.

Technical Info / Product NotesAMPIVIEW[®] KRT8 (AS) Dig RNA Probes (Human) are

optimized for detection of the nucleic acid (RNA/DNA or RNA) encoding KRT8 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: September 30, 2025



eu@enzolifesciences.com