

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

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

KRT8 RNA or DNA/RNA in cells and tissues
AMPIVIEW® KRT8 (AS) Dig RNA Probes (Human) are digoxigenin-labeled 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 Notes	AMPIVIEW® 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 function 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 results that can be visualized with a light microscope.
Contents	2 x 1 mL AMPIVIEW® KRT8 (AS) Dig RNA Probes (Human) (2µg/mL) 1 x 2 mL AMPIVIEW® Hybridization Buffer (1X)
Technical Info / Product Notes	AMPIVIEW® 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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.