

# AMPIVIEW® DLX6 (AS) DNP RNA Probes (Chicken) Set

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

DLX6 RNA or DNA/RNA in cells and tissues

AMPIVIEW® DLX6 (AS) DNP RNA Probes (Chicken) are DNP-labeled RNA probes targeting the nucleic acid (DNA/RNA or RNA) for DLX6 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-GEN350-2000	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	-20°C
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	ISH (in situ hybridization)
Application Notes	<p>AMPIVIEW® DLX6 (AS) DNP RNA probes (Chicken) have been designed to detect the nucleic acid for chicken DLX6 gene, part of the Distal-less homeobox (DLX) gene family, which plays a critical role in craniofacial development, limb formation, and chondrogenesis (cartilage formation).</p> <p>AMPIVIEW® DLX6 (AS) DNP RNA probes (Chicken) are conjugated with dinitrophenol (DNP) and have been optimized to produce clear results with AMPIVIEW® anti-DNP linker and nanopolymer-based detection systems, POLYVIEW® PLUS, combined with HIGHDEF® chromogens and counterstain to produce clear results that can be visualized with a light microscope.</p>
Contents	<p>2 x 1 mL AMPIVIEW® DLX6 (AS) DNP RNA Probes (Chicken) (2µg/mL)</p> <p>1 x 2 mL AMPIVIEW® Hybridization Buffer (1X)</p>
Technical Info / Product Notes	<p>AMPIVIEW® DLX6 (AS) DNP RNA Probes (Chicken) are optimized for detection of the nucleic acid (RNA/DNA or RNA) of chicken DLX6 in FFPE tissues or cells with AMPIVIEW® anti-DNP 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.</p>



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)