

# AMPIVIEW<sup>®</sup> Myosin-like Gene (AS) Dig RNA Probes (P. karlingii) Set

AMPIVIEW<sup>®</sup> Myosin-like Gene (AS) Dig RNA Probes (P. karlingii) are digoxigenin-labeled RNA probes targeting the *P. karlingii* (DNA/RNA or RNA) for *in situ* hybridization in tissues and cells. The probes are formulated in a buffered formaldehyde solution with hybridization enhancers.

tissues

## Ordering Information

[Order Online »](#)

ENZ-GEN194-2000	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<sup>®</sup> Myosin-like Gene (AS) Dig RNA Probes (*P. karlingii*) have been designed to target *Paralagenidium karlingii* myosin-like gene nucleic acid to detect *P. karlingii* infections in tissues and cells. AMPIVIEW<sup>®</sup> Myosin-like Gene (AS) Dig RNA Probes (*P. karlingii*) are conjugated with digoxigenin and have been optimized to produce clear results with Enzo's DIGX<sup>®</sup> anti-digoxigenin linker and nanopolymer-based detection systems, POLYVIEW<sup>®</sup> PLUS, combined with HIGHDEF<sup>®</sup> chromogens and counterstain to produce clear results that can be visualized with a light microscope.

**Contents** 2 x 1 mL AMPIVIEW<sup>®</sup> Myosin-like Gene (AS) Dig RNA Probes (*P. karlingii*) (2µg/mL)  
1 x 2 mL AMPIVIEW<sup>®</sup> Hybridization Buffer (1X)

**Technical Info / Product Notes** AMPIVIEW<sup>®</sup> Myosin-like Gene (AS) Dig RNA Probes (*P. karlingii*) are optimized for detection of *P. karlingii* DNA or RNA/DNA in FFPE tissues or cells with DIGX<sup>®</sup> anti-digoxigenin linker, POLYVIEW<sup>®</sup> PLUS detection reagents combined with HIGHDEF<sup>®</sup> chromogens and counterstains (linker and detection solutions not included). Dilutions and concentration optimization can be done with AMPIVIEW<sup>®</sup> Hybridization Buffer (1X), included in this set.



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-  
fr@enzolifesciences.com](mailto:info-fr@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)