

AMPIVIEW[®] HSP90 (AS) Dig RNA Probes (*L. giganteum*) Set

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

giganteum form *caninum* RNA or DNA/RNA in cells and tissues. AMPIVIEW[®] HSP90 (AS) Dig RNA Probes (*L. giganteum*) are digoxigenin-labeled RNA probes targeting the *L. giganteum* nucleic acid (DNA/RNA or RNA) 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-GEN196-2000	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 [®] HSP90 (AS) Dig RNA Probes (<i>L. giganteum</i>) have been designed to target <i>Lagenidium giganteum forma caninum</i> HSP90 nucleic acid to detect <i>L. giganteum</i> infections in tissues and cells. AMPIVIEW [®] HSP90 (AS) Dig RNA Probes (<i>L. giganteum</i>) are conjugated with digoxigenin and have been optimized to produce clear results with Enzo's DIGX [®] anti-digoxigenin 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.
Contents	2 x 1 mL AMPIVIEW [®] HSP90 (AS) Dig RNA Probes (<i>L. giganteum</i>) (2µg/mL) 1 x 2 mL AMPIVIEW [®] Hybridization Buffer (1X)
Technical Info / Product Notes	AMPIVIEW [®] HSP90 (AS) Dig RNA Probes (<i>L. giganteum</i>) are optimized for detection of <i>L. giganteum</i> 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 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)