

AMPIVIEW® Exo 1,3-β Glucanase (AS) Bio RNA Probes (P. insidiosum) Set

AMPIVIEW® Exo 1,3-β Glucanase (AS) Bio RNA Probes (*P. insidiosum*)
A biotin-labeled RNA probe for the *in situ*
hybridization (ISH) detection of *P. insidiosum*
RNA or DNA/RNA in cells and tissues
are biotin-labeled RNA probes targeting *P. insidiosum* (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-GEN192-2000	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[®] Exo 1,3-β Glucanase (AS) Bio RNA probes (*P.insidiosum*) have been designed to target *Pythium insidiosum* Exo 1,3-β Glucanase nucleic acid to detect *P. insidiosum* infections in tissues and cells. AMPIVIEW[®] Exo 1,3-β Glucanase (AS) Bio RNA Probes (*P.insidiosum*) are conjugated with biotin and have been optimized to produce clear results with Enzo's anti-biotin 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[®] Exo 1,3-β Glucanase (AS) Bio RNA Probes (*P. insidiosum*) (2μg/mL)
1 x 2 mL AMPIVIEW[®] Hybridization Buffer (1X)

Technical Info / Product Notes AMPIVIEW[®] Exo 1,3-β Glucanase (AS) Bio RNA Probes (*P. insidiosum*) are optimized for detection with either SAVIEW[®] PLUS and POLYVIEW[®] PLUS detection reagents, HIGHDEF[®] chromogens and counterstains (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)