

AMPIVIEW® NSP (AS) DNP RNA Probes Set

Non-specific negative control RNA probe

The AMPIVIEW® NSP (AS) DNP RNA Probes set is a ready-to-use in situ hybridization (ISH) probe, powered by Enzo's LoopRNA ISH™ technology, for chromogenic detection of a non-specific probe (NSP) sequence with no significant homology to human tissue RNA. NSP probes establish the baseline nonspecific background signal in ISH assays, enabling accurate interpretation of specific probe staining results and confirmation of assay specificity. in formalin-fixed paraffin-embedded (FFPE) tissue sections and cells. The probe is DNP-labeled and detected using an unconjugated anti-DNP linker followed by a polymer-based POLYVIEW® PLUS detection system combined with HIGHDEF® chromogens and counterstain, visualized under a standard light microscope. Supplied as 2 x 1.0 mL ready-to-use probe (ENZ-GEN199) with 1 x 2.0 mL AMPIVIEW® Hybridization Buffer (Fast Hyb Solution) (ENZ-ACC187).

AMPIVIEW® RNA probes are designed to detect the RNA or DNA/RNA of a non-specific probe (NSP) sequence with no significant homology to human tissue RNA. NSP probes establish the baseline nonspecific background signal in ISH assays, enabling accurate interpretation of specific probe staining results and confirmation of assay specificity. In situ hybridization (ISH) is a molecular technique that uses labeled nucleic acid probes to detect and localize specific RNA or DNA sequences directly within tissue sections or cells. Powered by Enzo's LoopRNA ISH™ technology, AMPIVIEW® RNA probes deliver superior sensitivity for chromogenic detection of RNA, DNA, or RNA/DNA targets on formalin-fixed paraffin-embedded (FFPE) tissue and cells, and are compatible with Enzo's complete immunohistochemistry detection workflow, from antigen retrieval to counterstain.

Ordering Information

[Order Online »](#)

ENZ-GEN199-2000	
-----------------	--

Manuals, SDS & CofA

[View Online »](#)

- Powered by Enzo's LoopRNA ISH™ technology for superior sensitivity
- Ready-to-use, DNP-labeled probe compatible with AMPIVIEW® Rabbit Anti-DNP Linker, POLYVIEW® PLUS detection and HIGHDEF® chromogens
- 1-hour hybridization with AMPIVIEW® Hybridization Buffer (Fast Hyb Solution) (ENZ-ACC187) for fast, same-day results
- Compatible with manual staining and automated platforms (LEICA BOND, Roche DISCOVERY/ULTRA) with prior optimization

Handling & Storage

Use/Stability	Stable under recommended storage conditions up to the expiration date indicated on the vial label. Avoid freeze-thaw cycles.
Handling	Protect from light and humidity. Avoid freeze-thaw cycles. Avoid RNases; wear gloves and work in a clean environment.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ISH
Contents	2 x 1.0 mL AMPIVIEW® NSP (AS) DNP RNA Probes set (ENZ-GEN199); 1 x 2.0 mL AMPIVIEW® Hybridization Buffer (ENZ-ACC187)
Quantity	2 x 1.0 mL probe + 1 x 2.0 mL hybridization buffer
Recommendation Dilutions/Conditions	Ready to use solution for manual use; Optimization may be required for automated staining systems.
Specificity	Non-specific sequence (antisense) — negative control probe
Technical Info / Product Notes	Ready-to-use probe for manual staining. Prior optimization required for automated staining systems. AMPIVIEW® RNA probes have been optimized with Enzo's AMPIVIEW® and POLYVIEW® Detection Systems combined with HIGHDEF® chromogens. Use of other linkers or polymer-based reagents will require prior optimization. Autostainers such as LEICA BOND or Roche DISCOVERY/ULTRA can be used but may require prior optimization



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)