

AMPIVIEW[®] SARS-CoV-2 RNA Probes Set

A mixture of biotin-labeled RNA probes for the *in situ* hybridization detection SARS-CoV-2 virus.

The AMPIVIEW[®] SARS-CoV-2 RNA Probe is a mixture of biotin-labeled probes targeting the Spike gene (S) and the Nucleocapsid gene (N) of SARS-CoV-2 for *in situ* hybridization. The probes are formulated in a buffered formamide solution with hybridization enhancers.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-GEN159-2000	2ml
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored as recommended. Following dilution in appropriate buffer system, use product immediately. Recommended buffer ENZ-ACC152 AMPIVIEW® Hybridization Buffer (1X).
Handling	Allow contents to warm up to room temperature prior to use.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application ISH (in situ hybridization)

Application Notes AMPIVIEW® SARS-CoV-2 RNA probes have been designed to target the N and S protein sequences of SARS-CoV-2 for its detection in tissues and cells. SARS-CoV-2 RNA probes are conjugated with biotin and have been optimized to produce clear results with Enzo's biotin-based detection system, SAVIEW® Plus, Enzo's anti-biotin linker and nanopolymer-based detection systems, POLYVIEW® Plus, and high-quality chromogens, HIGHDEF® to produce clear results that can be visualized with a light microscope.

Contents AMPIVIEW® SARS-CoV-2 RNA Probes (ENZ-GEN157)
AMPIVIEW® Hybridization Buffer (ENZ-ACC152)
AMPIVIEW® NSpf RNA Probes (negative control)

Sensitivity AMPIVIEW® RNA probes uniquely combine the precision of targeted, sequence-specific RNA probes with the superior sensitivity of Enzo's novel loopRNA™ technology, thus making these probes one of the most sensitive means for the detection of nucleic acid targets in tissue samples and cells, while preserving the morphology of the sample.

Technical Info / Product Notes AMPIVIEW® SARS-CoV-2 RNA Probes 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, 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)