

AMPIVIEW[®] SARS-CoV-2 RNA Probes Kit

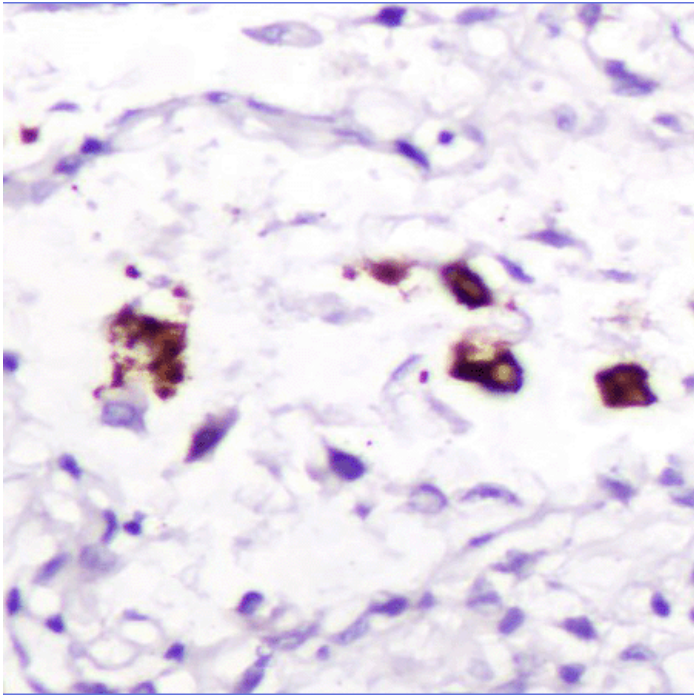
AMPIVIEW[®] SARS-CoV-2 RNA Probes kit comes, AMPIVIEW[®] SARS-CoV-2 RNA Probes, powered by Enzo's LoopRNA[™] ISH technology to deliver superior sensitivity with SAVIEW[®] PLUS detection system.

Citations: 1

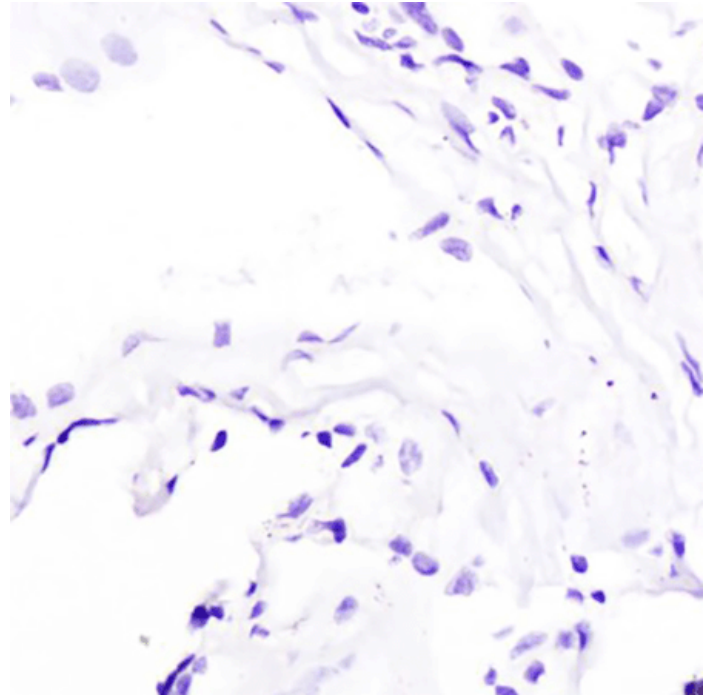
[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



SARS-CoV-2 detection with AMPIVIEW SARS-CoV-2 RNA probes (brown) in lung tissue of a COVID-19 pneumonia patient.



SARS-CoV-2 infection not detected with AMPIVIEW SARS-CoV-2 in normal lung tissue specimen

Handling & Storage

Use/Stability	Store the probes at -20°C to -80°C and the detection reagents at +4°C. Stable until expiration date when stored as recommended. Follow storage information provided in the package.
Handling	Allow contents to warm up to room temperature prior to use.
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ISH (in situ hybridization)
Contents	AMPIVIEW [®] SARS-CoV-2 RNA Probes (ENZ-GEN157) AMPIVIEW [®] Hybridization Buffer (ENZ-ACC152) Proteinase K (ENZ-33801) AMPIVIEW [®] Wash Buffer 1 AMPIVIEW [®] Wash Buffer 2 SAVIEW [®] PLUS AP (ENZ-ACC111) HIGHDEF [®] Red AP Chromogen/Substrate kit HIGHDEF [®] Hematoxylin (ENZ-ACC106)
Quantity	20 tests
Species Reactivity	Virus

Last modified: October 17, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com