# AMPIVIEW<sup>®</sup> SARS-CoV-2 RNA Probes Set

## A mitxture of biotin-labeled RNA probes for the *in situ* hybridization detection SARS-

The MP VIEW<sup>®</sup> 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**

Stable for at least one year after receipt when stored as recommended. Following **Use/Stability** 

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)

AMPIVIEW® SARS-CoV-2 RNA probes have been **Application Notes** 

> 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 biotinbased detection system, SAVIEW® Plus, Enzo's anti-biotin

linker and nanopolymner-based detection systems, POLYVIEW® Plus, and high-quality chromogens,

HIGHDEF<sup>®</sup> to produce clear results that can be visualized

with a light microscope.

AMPIVIEW<sup>®</sup> SARS-CoV-2 RNA Probes (ENZ-GEN157) Contents

> AMPIVIEW® Hybridization Buffer (ENZ-ACC152) AMPIVIEW® NSpf RNA Probes (negative control)

AMPIVIEW® RNA probes uniquely combine the precision Sensitivity

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.

AMPIVIEW® SARS-CoV-2 RNA Probes are optimized for **Technical Info / Product Notes** 

detection with either SAVIEW® PLUS and POLYVIEW® PLUS detection reagents, HIGHDEF<sup>®</sup> chromogens and counterstains (not inlcuded). Dilutions and concentration optimization can be done with AMPIVIEW® hybridization

buffer, included in this set.

Last modified: October 17, 2024



info-

eu@enzolifesciences.com