

# AcroMetrix™ Coronavirus 2019 (COVID-19) RNA Control

**RUO** For Research Use Only

**REF** 954519 AcroMetrix Coronavirus 2019 (COVID-19) RNA Control

## Summary and Explanation

AcroMetrix™ Coronavirus 2019 (COVID-19) RNA Control is prepared by formulating genomic RNA from Severe Acute Respiratory Syndrome-Related Coronavirus 2 (SARS-CoV-2) into a proprietary buffer. The RNA is ready for reverse transcription, PCR amplification and detection, as appropriate to the test. The kit contains two vials of SARS-CoV-2 specific RNA at the concentrations that will result low positive and ultra-low positive levels in most commonly used PCR based Coronavirus 2019 (COVID-19) nucleic acid testing methods.

AcroMetrix Coronavirus 2019 (COVID-19) RNA Control is quantified using Bio-Rad Droplet Digital™ PCR (ddPCR™).

AcroMetrix Coronavirus 2019 (COVID-19) RNA Control sequence design is compatible with, but not limited to, the following assays:

- Thermo Fisher Scientific™ TaqPath™ COVID-19 Combo kit (A47814)
- USA CDC assay
- China CDC assay
- Japan National Institute of Infectious Diseases assay
- Thailand National Institute of Health assay
- WHO Assay from Charité Institute of Virology

## Content

Catalog Number	Control Name	Copies/μL	Quantity
954519	AcroMetrix COVID-19 Low Positive Control	500	1 x 20 μL
	AcroMetrix COVID-19 Ultra-Low Positive Control	100	1 x 20 μL

## Warnings and Precautions

AcroMetrix Coronavirus 2019 (COVID-19) RNA Control contains genomic RNA. Although the contents are not known to be infectious, it is recommended that the product is handled as potentially hazardous to health. It should be used and discarded according to your own laboratory's safety procedures. Such safety procedures should include the wearing of protective gloves and avoiding the generation of aerosols.

AcroMetrix Coronavirus 2019 (COVID-19) RNA Control is packaged in polypropylene vials. Care should be taken during handling to avoid breakage and injuries inflicted upon by broken plastics.

## Instruction For Use

1. Thaw the control on ice, vortex gently, pulse-spin in a microcentrifuge to collect the material at the bottom of vial.
2. RNA reverse transcription should be performed as though the control is extracted RNA from patient samples.
3. Up to three freeze-thaw cycles is allowed for each vial.
4. The material may be stored at 2-8°C for 7 days with up to 5 uses.
5. The input amount of control for one-step or two-step RT-qPCR should be determined by sensitivity and limit of detection of the assay.

## Storage Condition

All vials should be stored immediately at -20°C or lower upon receipt.

See Safety Data Sheet for safety information.

## Glossary:

<http://www.thermofisher.com/symbols-glossary>



Microgenics Corporation  
46500 Kato Road  
Fremont, CA 94538 USA  
Technical Support:  
1-800-232-3342



For insert updates go to:  
[www.thermofisher.com/Acrometrix](http://www.thermofisher.com/Acrometrix)

© 2020 Thermo Fisher Scientific Inc. All rights reserved.