

Protocol for Viral RNA Extraction Using the Aurum Total RNA Mini Kit or the Aurum Total RNA 96 Kit




Protocol

Nucleic Acid Sample Preparation

Bulletin 7331

The Aurum Total RNA Mini and Aurum Total RNA 96 Kits have been tested for recovery and purification of viral RNA using SARS-CoV-2 Standard (catalog #COV019). Bio-Rad's new SARS-CoV-2 Droplet Digital PCR (ddPCR) Kit was used to verify the presence of the purified viral RNA.

Product Information

Catalog #	Description	Manual
7326820 	Aurum Total RNA Mini Kit, 50 preps	4110133 
7326800 	Aurum Total RNA 96 Kit	4110131 

Kits can be ordered and manuals can be downloaded from [bio-rad.com](https://www.bio-rad.com).

Important Note about Aurum Vacuum Manifold

The Aurum Vacuum Manifold (#7326470) is no longer available and no replacement product is available from Bio-Rad at this time. Customers are encouraged to validate other commercially available vacuum manifolds. Please contact your Bio-Rad sales representative for sample availability.

Protocol

- Mix 100 μ l of a specimen (for example, a nasopharyngeal swab) in transport media with 350 μ l of lysis solution from the kit. Pipet up and down several times to mix thoroughly.
- Add 350 μ l of 70% ethanol (not supplied) to each tube or well, depending on the Aurum Kit used. Pipet up and down to mix thoroughly. Make sure that no bilayer is visible and that the viscosity is substantially reduced.
- Proceed with the extraction protocol as indicated below:
 - Aurum Total RNA Mini Kit**
 - Vacuum protocol (using a vacuum manifold): All Starting Sample Types ([page 13](#))
 - Spin protocol: All Starting Sample Types ([page 17](#))
 - Aurum Total RNA 96 Kit**
 - Vacuum protocol: All Starting Cell Types ([page 10](#))
 - Spin protocol (using a plate centrifuge): All Starting Cell Types ([page 13](#))

Related Products

Catalog #	Description
12010176	Reliance One-Step Multiplex Supermix, 200 x 20 μ l reactions
12010220	Reliance One-Step Multiplex Supermix, 1,000 x 20 μ l reactions
12010221	Reliance One-Step Multiplex Supermix, 2,000 x 20 μ l reactions
1725140	iTaq Universal Probes One-Step Kit, 100 x 20 μ l reactions
1725141	iTaq Universal Probes One-Step Kit, 500 x 20 μ l reactions

BIO-RAD, DDPCR, and DROPLET DIGITAL PCR are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

