

Quick-Start Protocol

QIAmini Pathogen Kit

For use with the QIAmini instrument

The QIAmini Pathogen Kit (48) (cat. no. 590205) is shipped at ambient temperature. All kit components should be stored dry at room temperature (15–25°C).

Further information

- *QIAmini Pathogen Kit Handbook*: www.qiagen.com/HB-3809
- *QIAmini User Manual*: www.qiagen.com/HB-3799
- Safety Data Sheets: www.qiagen.com/safety
- Technical assistance: support.qiagen.com

Equipment and reagents to be supplied by the user

- Microcentrifuge tubes
- Thermal shaker

Notes before starting

- Buffer ACL and the QIAmini Pathogen cartridge contain guanidine salt and are therefore not compatible with disinfecting reagents containing bleach.
- Dissolve the lyophilized carrier RNA in RNase-free water to a final concentration of 1 µg/µL. See the *QIAmini Pathogen Kit Handbook* (www.qiagen.com/HB-3809) for carrier RNA storage details.
- For a 50 µL elution, add 75 µL RNase-free water into well 5 of the cartridge. For 100 µL elution, add 125 µL into well 5 of the cartridge. Select the appropriate protocol.
- Equilibrate buffers and cartridges to room temperature.
- Perform all steps at room temperature.
- Up to eight cartridges can be processed in one run.
- Download the appropriate protocol file from www.qiagen.com/QIAmini or from QIAmini Suite (www.qiagen.com/QIAminiSuite).
- If necessary, pretreat samples according to the appropriate procedure, as described in the *QIAmini Pathogen Kit Handbook* (www.qiagen.com/HB-3809).

Sample preparation

1. Pipet 20 µL QIAGEN Proteinase K into the bottom of a microcentrifuge tube.
Important: It is possible to add QIAGEN Proteinase K after dispensing samples, but it is best added beforehand.
Note: Do not add QIAGEN Proteinase K directly to the Buffer ACL.
2. Add 200 µL of the sample.
3. Add 240 µL Buffer ACL and mix thoroughly either by pulse vortexing for 20 s or by pipetting up and down at least 10 times. To ensure efficient lysis, it is essential that the sample and Buffer ACL are mixed thoroughly to yield a homogeneous solution.
4. Add 1 µL reconstituted Carrier RNA.
5. Incubate at room temperature for 15 min with shaking at 800 rpm (e.g., thermal shaker).

Cartridge preparation

6. Invert pre-filled cartridge(s) to resuspend magnetic beads. Flick down to remove reagents and bead suspension from under the foil.
7. Place cartridge(s) into the Cartridge Frame. Make sure to keep the pull tab accessible. Ensure cartridge(s) are fully inserted into the Cartridge Frame.
8. **Important:** Carefully remove the sealing foil. Avoid splashing of reagents.
9. Pipet RNase-free water for the required elution volume (75 μL for 50 μL eluate/125 μL for 100 μL eluate) in well 5 of the cartridge.
10. Transfer the lysate from step 5 into well 1 of the cartridge.

Procedure on the QIAmini instrument

11. Turn on the QIAmini instrument.
12. Select the desired protocol.
Note: Find a list of all imported protocols in the list view of the run menu.
Note: For details on the protocol selection, refer to the *QIAmini Pathogen Kit Handbook* (www.qiagen.com/HB-3809).
13. Open the instrument door. Load the Cartridge Frame holding the cartridge(s) into the instrument.
Note: Ensure that the Cartridge Frame is placed in the correct orientation. Press the cartridges down to fully engage with the heater block.
14. Slide the Rod Cover(s) into the rod cover holder of the instrument.
Note: Press the clip down to allow full insertion of the rod cover holder.
15. Close the instrument door, press  and confirm that consumables have been installed correctly to initiate the protocol run.
16. The display shows “End of run” when the run is completed. Select **Confirm** and press . Open the instrument door. Remove and discard the Rod Cover(s). Then remove the Cartridge Frame holding the cartridge(s) and transfer the purified nucleic acids from well 5 of each cartridge into a new 1.5 mL tube. Discard the cartridge(s).
Recommended: Clean the instrument and perform a UV decontamination of the instrument chamber.

Document Revision History

Date	Changes
08/2026	Initial release.



Scan QR code for handbook.
For up-to-date licensing information and product-specific disclaimers, see the respective QIAGEN kit handbook or user manual.

Trademarks: QIAGEN®, Sample to Insight®, (QIAGEN Group). Registered names, trademarks, etc. used in this document, even when not specifically marked as such, are not to be considered unprotected by law.
1139666 08/2026 HB-3810-001 © 2026 QIAGEN, all rights reserved.