

Quick-Start Protocol

QIAmini DNA Blood Kit

For use with the QIAmini instrument

The QIAmini DNA Blood Kit (48) (cat. no. 590105) is shipped at ambient temperature. All reagents and kit components should be stored at room temperature (15–25°C).

Further information

- *QIAmini DNA Blood Kit Handbook*: www.qiagen.com/HB-3803
- *QIAmini User Manual*: www.qiagen.com/HB-3799
- Safety Data Sheets: www.qiagen.com/safety
- Technical assistance: support.qiagen.com

Equipment and reagents supplied by the user

- 2 mL tubes
- Elution tubes

Notes before starting

- Buffer AL and buffers in cartridge contain guanidine salt and are therefore not compatible with disinfecting reagents containing bleach.
- If necessary, redissolve any precipitate in Buffer AL by warming.
- Equilibrate cartridges to room temperature.
- Thaw and equilibrate whole blood samples at room temperature (15–25°C).
- Perform all steps at room temperature.
- Up to eight cartridges can be processed in one run.

Sample preparation

1. Pipet 15 µL QIAGEN Proteinase K into the bottom of a 2 mL tube (not provided). Add 200 µL blood to the Proteinase K.
2. Add 200 µL Buffer AL and mix thoroughly either by vortexing for 20 s or by vigorously pipetting up and down at least 10 times. To ensure efficient lysis, it is essential that the sample and Buffer AL are mixed thoroughly to yield a homogenous solution.

Note: Do not add QIAGEN Proteinase K directly to Buffer AL.

Cartridge preparation

3. Invert pre-filled cartridge(s) to resuspend magnetic beads. Flick down to remove reagents and bead suspension from under the foil.
4. 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.
5. **Important:** Carefully remove the sealing foil. Avoid splashing of reagents.
6. Transfer RNase-free water into well 5 of the cartridges for elution with the respective volume according to the selected protocol.
7. Add 415 µL lysate to the well 1 of the cartridge.

Procedure on the QIAmini instrument

8. Turn on the QIAmini instrument.

9. 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 DNA Blood Kit Handbook* (www.qiagen.com/HB-3803).

10. 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.

11. 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.

12. Close the instrument door, press , and confirm that consumables have been installed correctly to initiate the protocol run.

13. 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.

1139663 08/2026 HB-3804-001 © 2026 QIAGEN, all rights reserved.