

QIAmini Protocol Overview

Available QIAmini protocol files for dedicated QIAmini kits.

Important note before starting

QIAmini instruments are delivered without pre-installed protocols. Before starting a run, download the protocol file that matches the kit, starting material, and analyte required for your application, then transfer the file to the instrument.

How to use this overview

- Use this document together with the relevant QIAmini kit instructions for use to identify the appropriate protocol file for each kit and sample type.
- Confirm that the selected protocol file matches the kit, starting material, analyte, and elution volume stated in the table.
- After importing the protocol file, use the instrument touchscreen display to select and start the required protocol.

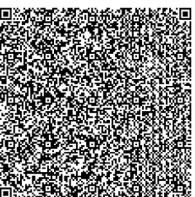
Download and transfer of protocol files

- **Download:** Please download the appropriate protocol file from www.qiagen.com/QIAmini or from QIAmini Suite at www.qiagen.com/QIAminiSuite to make sure you use the current file for your kit and application.
- **Transfer:** Transfer the protocol file to QIAmini using a USB drive, Bluetooth or QIAmini barcode scanner.
- **Select and run:** A list of all imported protocols is available in the Run menu list view. Select the appropriate protocol on the touchscreen display and start the run.

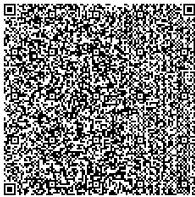
Available QIAmini protocol files

These tables list protocol file names with the associated kit, starting material, analyte, and protocol- specific notes. Always use the latest available protocol file from the download locations listed above.

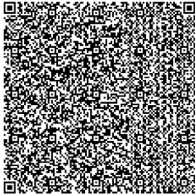
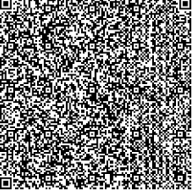
QIAmini DNA Tissue/Cells Kit

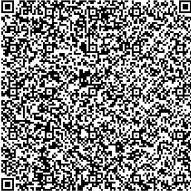
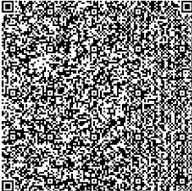
Protocol file	Starting material	Analyte(s)	Protocol notes
DNA_TC_200_Q1 	cells and tissue	gDNA	Extraction of gDNA from tissues and cultured cells using 200 µL lysate and 200 µL elution volume.
DNA_TC_50_Q1 	Swabs and low biomass samples	gDNA	Extraction of gDNA from buccal swabs and low biomass samples using approx. 250 µL lysate and 50 µL elution volume.

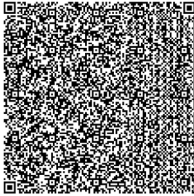
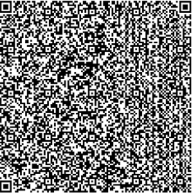
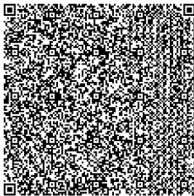
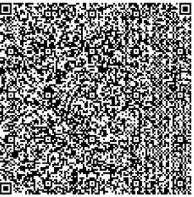
QIAmini DNA Blood Kit

Protocol file	Starting material	Analyte(s)	Protocol notes
DNA_BI_80_Q2 	whole blood	gDNA	Extraction of gDNA from whole blood using 415 µL lysate. The protocol yields approximately 80 µL of eluate when using an elution input volume of 100 µL.

QIAmini PowerFecal Pro DNA/RNA Kit

Protocol file	Starting material	Analyte(s)	Protocol notes
NA_PF_100_Q1 	feces	total nucleic acids, DNA	For DNA-only purification, mix 420 µL sample supernatant with 4 µL RNase A for RNA digestion before loading into the cartridge. Prefilled elution volume: 100 µL RNase-free water.
RNA_PF_100_Q2 - part 1 of 2  RNA_PF_100_Q2 - part 2 of 2 	feces	RNA	For RNA-only purification, pipet 75 µL Buffer RDD, 150 µL isopropanol, and 10 µL resuspended DNase I to the bead suspension in well 6 of the cartridge. Prefilled elution volume: 100 µL RNase-free water.

Protocol file	Starting material	Analyte(s)	Protocol notes
NA_TC_80_Q1 	cells and tissue	total nucleic acids	Extraction of total nucleic acids, including miRNA, from cell and tissue samples. The protocol yields approximately 80 µL of eluate when using an elution input volume of 100 µL.
RNA_TC_80_Q2 - part 1 of 2  RNA_TC_80_Q2 - part 2 of 2 	cells and tissue	RNA	Extraction of total RNA, including miRNA, from cell and tissue samples. The protocol yields approximately 80 µL of eluate when using an elution input volume of 100 µL.

Protocol file	Starting material	Analyte(s)	Protocol notes
Path_100_Q1 	Serum; Plasma; Whole blood; Urine; Transport media; Swabs; Respiratory samples	viral total nucleic acids, bacterial DNA	Extraction of viral nucleic acids and bacterial DNA from biofluids and transport media using 200 µL sample input and 100 µL elution volume.
Path_50_Q1 	Serum; Plasma; Whole blood; Urine; Transport media; Swabs; Respiratory samples	viral total nucleic acids, bacterial DNA	Extraction of viral nucleic acids and bacterial DNA from biofluids and transport media using 200 µL sample input and 50 µL elution volume.
PathXL_100_Q1 	Serum; Plasma	viral total nucleic acids, bacterial DNA	Extraction of viral nucleic acids and bacterial DNA from serum or plasma using 300 µL sample input and 100 µL elution volume.
PathXL_50_Q1 	Serum; Plasma	viral total nucleic acids, bacterial DNA	Extraction of viral nucleic acids and bacterial DNA from serum or plasma using 300 µL sample input and 50 µL elution volume.

Notes and disclaimers

This overview supports protocol selection and download. It does not replace the kit instructions for use, user manual, or any applicable safety regulatory and instrument instructions. Refer to the relevant kit instructions for use, for complete protocol details and sample preparation requirements.



Scan QR code for user manual.

For up-to-date licensing information and product-specific disclaimers, see the respective QIAGEN instructions for use or user manual. QIAGEN kit instructions for use and user manuals are available at www.qiagen.com or can be requested from QIAGEN Technical Service or your local distributor.

Document Revision History

Date	Changes
08/2026	Initial release
09/2026	Corrections and adjustments

Trademarks: QIAGEN®, Sample to Insight® Registered names, trademarks, etc. used in this document, even when not specifically marked as such, are not to be considered unprotected by law.

09/2026 HB-3870-002 © 2026 QIAGEN, all rights reserved.