

Quick-Start Protocol

QIAamp[®] DNA Host-Free Microbiome Kit

The components of the QIAamp DNA Host-Free Microbiome Kit are shipped in two boxes: Box 1 is shipped on wet ice at 4°C, and Box 2 is shipped at room temperature (15–25°C). Upon receipt, store the Host Depletion Solution and the QIAamp UCP Mini Spin Columns at 2–8°C, Benzonase at –20°C, and all other kit components dry at room temperature.

Further information

- QIAamp DNA Host-Free Microbiome Kit Handbook (www.qiagen.com/resources/kithandbook/qiaamp-dna-host-free-microbiome-kit-handbook)
- Safety Data Sheets: www.qiagen.com/safety
- Technical assistance: support.qiagen.com

Notes before starting

- Dissolve lyophilized Host Depletion Solution in 1.8 mL PBS. Mix gently by inverting. Do not spin down. For long-term storage, store single-use aliquots at –15°C to –25°C. Thawed aliquots can be stored at 2–8°C for 6 weeks. A precipitate may form during storage; however, this does not compromise product activity.
- If a precipitate has formed in Buffer ATL or Buffer APL2, dissolve by incubating at 56°C.
- Ensure that Buffers AW1 and AW2 have been prepared according to the instructions on the bottle.
- Add 100 µL Reagent DX to 14 mL Buffer ATL. If smaller amounts are needed, transfer 1.5 mL of Buffer ATL into a sterile 2 mL vial and add 10 µL Reagent DX. Mix well. After

preparation, the mixture is stable for 6 months at room temperature.

- Spin the PowerBead Pro Tube briefly to ensure that the beads have settled at the bottom.
- Tissue samples: mince or grind sample before starting the protocol (e.g., using TissueRuptor® II, cat. no. 990890). Spin down (10 min, 10,000 × g) and remove the supernatant.
- Swab samples: swirl or vortex the swab in 1 mL transport media or PBS for at least 20 s and dry off by pressing against the wall of the tube multiple times. Spin down (10 min, 10,000 × g) and remove the supernatant.
- Bodily fluids: spin down (10 min, 10,000 × g) and remove the supernatant.
- Viscous respiratory samples: refer to the *QIAamp Power Pro Quick-Start Protocol* (www.qiagen.com/HB-3575). Spin down (10 min, 10,000 × g) and remove the supernatant.

Equipment and reagents to be supplied by user

- Ethanol (96–100%)
- Phosphate-buffered saline (PBS)
- Microcentrifuge (with rotor for 2 mL tubes)
- Shaker–incubator
- Equipment for sample disruption and homogenization, one of the following:
 - Vortex-Genie® 2 and Vortex Adapter for 24 (1.5–2 mL) tubes (cat. no. 13000-V1-24)
 - TissueLyser III (cat. no. 9003240) with adapter sets for use with the PowerBead Pro Tubes (TissueLyser Adapter Set 2 × 24, cat. no. 69982), or 2 mL Tube Holder (cat. no. 11993), in conjunction with Plate Adapter Set (cat. no. 11990)

Procedure

1. Add, in the following order, 220 µL Buffer RDD, 3 µL Benzonase and 30 µL Host Depletion Solution to up to 100 mg of tissue or to the preprocessed swab sample in a 2 mL microcentrifuge tube. Mix well and incubate at 37°C for 45 min at 600 rpm in a heating

block or water bath.

2. Add 20 μ L Proteinase K and mix well. Incubate at 56°C for 20 min at 600 rpm in a heating block or water bath.

Note: If needed, the sample may be stored overnight at 2–8°C or at room temperature after this step. Do not store samples for longer periods. Before continuing with step 3, thaw the sample if required.

3. Add 200 μ L Buffer ATL (containing Reagent DX). Mix well to avoid loss of sample material, and transfer into a PowerBead Pro Tube.
4. Mechanically disrupt samples using one of the following methods:
 - a. Secure the PowerBead Pro Tube horizontally on a Vortex Adapter for 24 (1.5–2 mL) tubes (cat. no. 13000-V1-24). Orient the tube caps to point toward the center of the vortex adapter. Vortex at maximum (20,000 $\times$ g) for 10 min.
 - b. Use TissueLyser III. Place the PowerBead Pro Tube into the TissueLyser Adapter Set 2 $\times$ 24 (cat. no. 69982) or 2 mL Tube Holder (cat. no. 11993) and Plate Adapter Set (cat. no. 11990). Fasten the adapter into the instrument and shake for 5 min at 30 Hz. Reposition the adapter so that the side that was closest to the machine body is now furthest from it. Shake again for 5 min at 30 Hz.

5. Centrifuge the PowerBead Pro Tube at 10,000 $\times$ g for 1 min. Transfer the supernatant to a fresh microcentrifuge tube.

Note: If needed, the sample may be stored overnight at 2–8°C or at room temperature after this step. Do not store samples for longer periods. Before continuing with step 6, thaw the sample if required.

6. Add 200 μ L APL2. Mix by pulse vortexing for 30 s.
7. Incubate at 70°C for 10 min and briefly spin the tube to remove condensation.
8. Carefully apply up to 700 μ L of the mixture from step 7 to the QIAamp UCP Mini Column. Close the cap and centrifuge at 6000 $\times$ g for 1 min.
9. Discard the flow-through. Put the column back into the collection tube to repeat step 8 with any remaining mixture from step 7.

- 10. Transfer the QIAamp UCP Mini Column to a fresh collection tube. Carefully open the cap and add 500 µL Buffer AW1 without wetting the rim. Close the cap and centrifuge at 6000 × g for 1 min. Discard the flow-through.
- 11. Add 500 µL Buffer AW2 to the QIAamp UCP Mini Column without wetting the rim. Centrifuge at full speed (20,000 × g) for 3 min. Discard the flow-through.
- 12. Centrifuge at full speed (20,000 × g) for 1 min.
- 13. Place the QIAamp UCP Mini Column into a fresh 1.5 mL tube and apply 50 µL RNase-Free Water directly onto the center of the membrane. Close the lid and incubate at room temperature for 5 min.
- 14. Centrifuge at 6000 × g for 1 min to elute the DNA.

Document Revision History

Date	Changes
03/2025	Initial revision.
04/2026	Added sample preparation instructions for different sample type categories. Added information about stopping points.

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

Trademarks: QIAGEN®, Sample to Insight®, QIAamp®, TissueRuptor® (QIAGEN); Vortex-Genie® (Scientific Industries, Inc.). Registered names, trademarks, etc. used in this document, even when not specifically marked as such, are not to be considered unprotected by law.

1139783 04/2026 HB-3652-002 © 2026 QIAGEN, all rights reserved.