

Agilent RNA 6000 Pico Kit Quick Start Guide

The complete RNA 6000 Pico Guide can be found in the online help of the 2100 Expert software.

| Agilent RNA 6000 Pico Kit (reorder-no 5067-1513) | |
|---|---|
| <i>Agilent RNA 6000 Pico Chips</i> | <i>Agilent RNA 6000 Pico Reagents (reorder-no 5067-1514) & Supplies</i> |
| 25 RNA Pico Chips | ● (blue) RNA 6000 Pico Dye Concentrate ¹ |
| 3 Electrode Cleaners | ● (green) RNA 6000 Pico Marker (4 vials) ○ (white) RNA 6000 Pico Conditioning Solution |
| <i>Syringe Kit</i> | ● (red) RNA 6000 Pico Gel Matrix (2 vials) |
| 1 Syringe | ● (yellow) RNA 6000 Pico Ladder (reorder-no 5067-1535) (1 vial, 10x concentrate) 4 Spin Filters (reorder-no 5185-5990) |
| <i>Tubes for Gel-Dye Mix</i> | |
| 30 Safe-Lock Eppendorf Tubes PCR clean (DNase/RNase free) for gel-dye mix | |

¹ “This product is provided under a license by Life Technologies Corporation to Agilent Technologies. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product only as described in accompanying product literature. The sale of this product is expressly conditioned on the buyer not using the product or its components (1) in manufacturing; (2) to provide a service, information, or data to an unaffiliated third party for payment; (3) for therapeutic, diagnostic or prophylactic purposes; (4) to resell, sell or otherwise transfer this product or its components to any third party, or use for any use other than use in the subfields of research and development, quality control, forensics, environmental analysis, biodefense or food safety testing. For information on purchasing a license to this product for purposes other than described above contact Life Technologies Corporation, Cell Analysis Business Unit, Business Development, 29851 Willow Creek Road, Eugene, OR 97402, Tel: (541) 465-8300. Fax: (541) 335-0354.”

Assay Principles

Agilent RNA kits contain chips and reagents designed for analysis of RNA fragments. Each RNA chip contains an interconnected set of microchannels that is used for separation of nucleic acid fragments based on their size as they are driven through it electrophoretically. Agilent RNA kits are designed for use with the Agilent 2100 Bioanalyzer instrument only.

Applications and Kits

Agilent RNA kits are designed for the analysis of total RNA (eukaryotic, prokaryotic, and plant) and mRNA samples.

Agilent RNA kits: RNA 6000 Nano Kit (reorder-no 5067-1511), RNA 6000 Pico Kit (reorder-no 5067-1513), and Small RNA Kit (reorder-no 5067-1548)

Storage Conditions

- Freeze unopened RNA ladder at -20 °C. Prepared ladder aliquots need to be stored at -70 °C. Keep all other reagents and reagent mixes refrigerated at 4 °C when not in use to avoid poor results caused by reagent decomposition.
- Protect dye and dye mixtures from light. Remove light covers only when pipetting. Dye decomposes when exposed to light.



Agilent Technologies

Agilent RNA 6000 Pico Kit Quick Start Guide

Equipment Supplied with the Agilent 2100 Bioanalyzer System

- Chip priming station (reorder-no 5065-4401)
- IKA vortex mixer

Additional Material Required (Not Supplied)

- RNaseZAP® recommended for electrode decontamination *only* (Ambion, Inc. cat. no. 9780)
- RNase-free water recommended for routine electrode cleaning
- Pipettes (10 µL and 1000 µL) with compatible tips (RNase-free, no filter tips, no autoclaved tips)
- 0.5 mL and 1.5 mL microcentrifuge tubes (RNase-free Eppendorf Safe-lock PCR clean or Eppendorf DNA LoBind microcentrifuge tubes are highly recommended)
- Microcentrifuge (> 13000g)
- Heating block or water bath for ladder/sample preparation
- Mandatory: Dedicated bayonet electrode cartridge (reorder-no 5065-4413) for RNA 6000 Pico and Small RNA assays.

Sample Preparation

- Prepare RNA samples in deionized water. For estimation of RNA concentration, the total RNA concentration in the sample must be between 200 – 5000pg/µL. The mRNA concentration must be between 500 – 5000pg/µL. If concentration of a sample is above this range, dilute with RNase-free water.

Specifications

| Physical Specifications | | Analytical Specifications | | |
|-------------------------|--|--|---|---|
| Type | Specification | Specification | Total RNA Assay | mRNA Assay |
| Analysis time | 30 min | Qualitative range | 50 – 5000 pg/µL in water | 250 – 5000 pg/µL in water |
| Samples per chip | 11 | Sensitivity (S/N>3) | 50 pg/µL in water or 200 pg/µL in TE | 250 pg/µL in water or 500 pg/µL in TE |
| Sample volume | 1 µL | Quantitation reproducibility (within a chip) | 20 % CV | 20 % CV |
| Kit stability | = 4 months at 4 °C | Quantitation accuracy ¹ | 30 % | - |
| Kit size | 25 chips 11 samples/chip = 275 samples/kit | Maximum sample buffer strength | 50 mM Tris 0.1 mM EDTA or 50 mM NaCl 15 mM MgCl ₂ | 50 mM Tris 0.1 mM EDTA or 50 mM NaCl 15 mM MgCl ₂ |

¹ Determined analyzing the RNA ladder as sample

Setting up the Chip Priming Station

- 1 Replace the syringe:
 - a Unscrew the old syringe from the lid of the chip priming station.
 - b Release the old syringe from the clip. Discard the old syringe.
 - c Remove the plastic cap of the new syringe and insert it into the clip.
 - d Slide it into the hole of the luer lock adapter and screw it tightly to the chip priming station.
- 2 Adjust the base plate:
 - a Open the chip priming station by pulling the latch.
 - b Using a screwdriver, open the screw at the underside of the base plate.
 - c Lift the base plate and insert it again in position C. Retighten the screw.
- 3 Adjust the syringe clip:
 - a Release the lever of the clip and slide it up to the top position.



Essential Measurement Practices

- Handle and store all reagents according to the instructions on the label of the individual box.
- Avoid sources of dust or other contaminants. Foreign matter in reagents and samples or in the wells of the chip will interfere with assay results.
- Allow all reagents to equilibrate to room temperature for 30 min before use. Thaw samples on ice.
- Protect dye and dye mixtures from light. Remove light covers only when pipetting. The dye decomposes when exposed to light and this reduces the signal intensity.
- Always insert the pipette tip to the bottom of the well when dispensing the liquid. Placing the pipette at the edge of the well may lead to poor results.
- Always wear gloves when handling RNA and use RNase-free tips, microcentrifuge tubes and water.
- It is recommended to heat denature all RNA samples and RNA ladder before use for 2 min and 70 °C (once) and keep them on ice.
- Do not touch the Agilent 2100 Bioanalyzer instrument during analysis and never place it on vibrating surface.
- Always vortex the dye concentrate for 10 s before preparing the gel-dye mix and spin down afterwards.
- Use a new syringe and electrode cleaners with each new kit.
- Use loaded chips within 5 min after preparation. Reagents might evaporate, leading to poor results.
- To prevent contamination (e.g. RNase), it is strongly recommended to use a dedicated electrode cartridge for RNA assays. For running the RNA 6000 Pico assay, the 16 pin bayonet cartridge is mandatory.



Agilent RNA 6000 Pico Assay Protocol

WARNING

Handling DMSO

Kit components contain DMSO. Because the dye binds to nucleic acids, it should be treated as a potential mutagen and used with appropriate care.

- Wear hand and eye protection and follow good laboratory practices when preparing and handling reagents and samples.
- Handle solutions with particular caution as DMSO is known to facilitate the entry of organic molecules into tissues.

Preparing the RNA Ladder

- 1 Spin the ladder down and pipette in an RNase-free vial.
- 2 Heat denature the ladder for 2 min at 70 °C.
- 3 Immediately cool the vial on ice.
- 4 Add 90 µL of RNase-free water and mix thoroughly.
- 5 Prepare aliquots in recommended 0.5 mL RNase-free vials with the required amount for typical daily use.
- 6 Store aliquots at -70 °C. After initial heat denaturation, the frozen aliquots should not require repeated heat denaturation.
- 7 Before use, thaw ladder aliquots on ice (avoid extensive warming).

Preparing the Gel

- 1 Pipette 550 µL of RNA gel matrix (red ●) into a spin filter.
- 2 Centrifuge at 1500g ± 20 % for 10 min at room temperature.
- 3 Aliquot 65 µL filtered gel into 0.5 mL RNase-free microcentrifuge tubes. Use filtered gel within 4 weeks. Store at 4 °C.



Agilent RNA 6000 Pico Kit Quick Start Guide

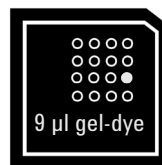
Preparing the Gel-Dye Mix

- 1 Allow the RNA dye concentrate (blue ●) to equilibrate to room temperature for 30 min.
- 2 Vortex RNA dye concentrate (blue ●) for 10 s, spin down and add 1 μL of dye into a 65 μL aliquot of filtered gel.
- 3 Vortex solution well. Spin tube at 13000g for 10 min at room temperature. Use prepared gel-dye mix within one day.




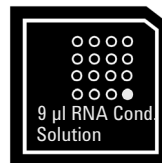
Loading the Gel-Dye Mix

- 1 Put a new RNA chip on the chip priming station.
- 2 Pipette 9 μL of gel-dye mix in the well marked .
- 3 Make sure that the plunger is positioned at 1 mL and then close the chip priming station.
- 4 Press plunger until it is held by the clip.
- 5 Wait for exactly 30 s then release clip.
- 6 Wait for 5 s. Slowly pull back plunger to 1 mL position.
- 7 Open the chip priming station and pipette 9 μL of gel-dye mix in the wells marked .
- 8 Discard the remaining gel-dye mix.




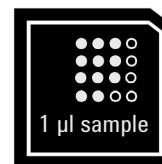
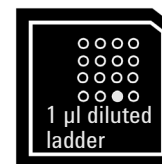
Loading the Conditioning Solution and Marker

- 1 Pipette 9 μL of the RNA conditioning solution (white ○) into the well marked CS.
- 2 Pipette 5 μL of RNA marker (green ●) in all 11 sample wells and in the well marked .



Loading the Diluted Ladder and Samples

- 1 Pipette 1 μL of the heat denatured and aliquoted ladder in the well marked .
- 2 Pipette 1 μL of sample in each of the 11 sample wells. Pipette 1 μL of RNA Marker (green ●) in each unused sample well.
- 3 Put the chip horizontally in the IKA vortexer and vortex for 1 min at 2400rpm.
- 4 Run the chip in the Agilent 2100 Bioanalyzer instrument within 5 min.



Technical Support

Please visit our support web page <http://www.agilent.com/genomics/contactus> to find information on your local Contact Center.

Further Information

Visit the 2100 Bioanalyzer site at <http://www.agilent.com/genomics/bioanalyzer>. You can find useful information, support and current developments about the products and the technology.



G2938-90049 Rev. D

Edition 01/2017

For Research Use Only
Not for use in diagnostic procedures

Printed in Germany

© Agilent Technologies, Inc. 2002-2016, 2017

Agilent Technologies
Hewlett-Packard-Straße 8
76337 Waldbronn, Germany