

RNA 6000 Nano Kit for 2100 Bioanalyzer Systems

Kit Guide

For Research Use Only - Not for use in diagnostic procedures



Notices

Document Information

Document No: SD-UF0000075
Rev. E.00
Edition: 09/2023

Copyright

© Agilent Technologies, Inc. 2000-2023

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Warranty

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

Safety Notices

CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

Contents

1	RNA 6000 Nano Kit	4
2	Required Equipment for RNA 6000 Nano Assay	6
3	Setting up the Assay Equipment and Bioanalyzer	7
	Setting up the Chip Priming Station	8
	Setting up the Bioanalyzer	9
	Vortex Mixer	9
	Starting the 2100 Expert Software	10
4	Essential Measurement Practices	11
5	Preparing the RNA Ladder after Arrival	12
6	RNA 6000 Nano Assay Protocol	13
	Decontaminating the Electrodes	14
	Preparing the Gel	15
	Preparing the Gel-Dye Mix	16
	Loading the Gel-Dye Mix	17
	Loading the RNA 6000 Nano Marker	18
	Loading the Ladder and Samples	19
	Inserting a Chip in the 2100 Bioanalyzer Instrument	20
	Starting the Chip Run	21
	Cleaning Electrodes after a RNA 6000 Nano Chip Run	23
7	Checking Your RNA 6000 Nano Assay Results	24
	RNA 6000 Nano Ladder Well Results	24
	RNA 6000 Nano Sample Well Results	25



1 RNA 6000 Nano Kit

Agilent RNA 6000 Nano Kit (5067-1511)	
Agilent RNA 6000 Nano Chips	Agilent RNA 6000 Nano Reagents (5067-1512) & Supplies
25 RNA Nano Chips	● (yellow) RNA 6000 NanoLadder (1 vial, 5067-1529)
2 Electrode Cleaners	● (blue) RNA 6000 Nano Dye Concentrate (1 vial)
	● (green) RNA 6000 Nano Marker (2 vials)
Syringe Kit	● (red) RNA 6000 Nano Gel Matrix (2 vials)
1 Syringe	4 Spin Filters (5185-5990)
Tubes for Gel-Dye Mix	
30 Safe-Lock Eppendorf Tubes PCR clean (DNase/RNase free) for gel-dye mix	

Table 1 Physical Specifications

Type	Specification
Analysis run time	30 min
Number of samples	12 samples/chip
Sample volume	1 µL
Kit stability	4 months
Kit size	25chips 12 samples/chip = 300 samples/kit

Table 2 Analytical Specifications

Specification	Total RNA Assay	mRNA Assay
Quantitative range	25 – 500 ng/µL	25 – 250 ng/µL
Qualitative range	5 – 500 ng/µL	5 – 250 ng/µL
Sensitivity (S/N>3)	5 ng/µL in water	25 ng/µL in water
Quantitative precision (within a chip)	10 % CV	10 % CV
Quantitative accuracy ¹	20 %	20 %
Maximum buffer concentration in sample	100 mM Tris 0.1 mM EDTA or 125 mM NaCl 15 mM MgCl ₂	100 mM Tris 0.1 mM EDTA or 125 mM NaCl 15 mM MgCl ₂

¹ Determined analyzing the RNA ladder as sample

2

Required Equipment for RNA 6000 Nano Assay

Equipment Supplied with the Agilent 2100 Bioanalyzer System

- Chip priming station (5065-4401)
- IKA vortex mixer
- 16-pin bayonet electrode cartridge (5065-4413)

Additional Material Required (Not Supplied)

- RNaseZAP® recommended for electrode decontamination
- RNase-free water
- Pipettes (10 µL and 1000 µL) with compatible tips (RNase-free, no filter tips, no autoclaved tips)
- 0.5 mL microcentrifuge tubes (RNase-free). Eppendorf Safe-lock PCR clean or Eppendorf DNA LoBind microcentrifuge tubes are highly recommended.
- Microcentrifuge (≥ 1300 g)
- Heating block or water bath for ladder/sample denaturation

Further Information

Visit the Agilent website. It offers useful information, support, and current developments about the products and technology:

www.agilent.com/en/product/automated-electrophoresis/bioanalyzer-systems.

3

Setting up the Assay Equipment and Bioanalyzer

Before beginning the chip preparation protocol, ensure that the chip priming station and the Bioanalyzer are set up and ready to use.

You have to

- replace the syringe at the chip priming station with each new kit
- adjust the base plate of the chip priming station
- adjust the syringe clip at the chip priming station
- set up the vortex mixer
- finally, make sure that you start the software before you load the chip.

NOTE

Please read this guide carefully and follow all instructions to guarantee satisfactory results.

Setting up the Chip Priming Station

NOTE

Replace the syringe with each new reagent kit.

- 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.



Setting up the Bioanalyzer

- 1 Open the lid of the Bioanalyzer and make sure that the electrode cartridge is inserted in the instrument. If not, open the latch and insert the electrode cartridge.

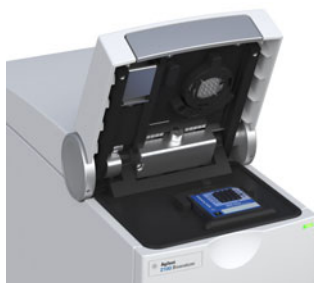


Figure 1 Electrode cartridge inserted in the instrument (graphic shows an example).

- 2 Remove any remaining chip.

Vortex Mixer

IKA - Model MS3

- 1 To set up the vortex mixer, adjust the speed knob to 2400 rpm.



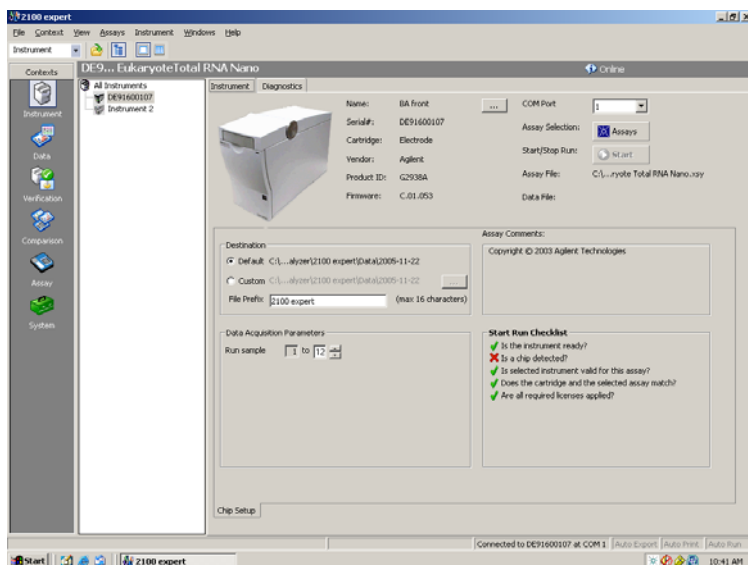
Starting the 2100 Expert Software

To start the software:

- 1 Go to your desktop and double-click the following icon.



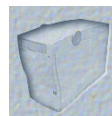
The screen of the software appears in the **Instrument** context. The icon in the upper part of the screen represents the current instrument/PC communication status:



Lid closed, no chip or chip empty



Lid open



Dimmed icon: no communication



Lid closed, chip inserted, RNA or demo assay selected

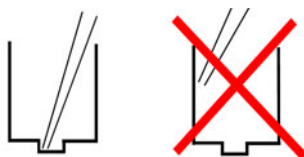
- 2 If more than one instrument is connected to your PC, select the instrument you want to use in the tree view.



4

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 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.
- Perform the RNase decontamination procedure for the electrodes daily before running any assays. Refer to the kit guide for details on electrode cleaning and decontamination.

5

Preparing the RNA Ladder after Arrival

For proper handling of the ladder, following steps are necessary:

- 1 After reagent kit arrival, spin ladder down. The ladder can be ordered separately (5067-1529).
- 2 Heat denature the ladder for 2 min at 70 °C.
- 3 Immediately cool the vial on ice.
- 4 Prepare aliquots in RNase- free vials with the required amount for a typical daily use.
- 5 Store aliquots at -28 – -15 °C (-18 – 5 °F).
- 6 Before use, thaw ladder aliquots and keep them on ice (avoid extensive warming upon thawing process).

6

RNA 6000 Nano Assay Protocol

After completing the initial steps in “Setting up the Assay Equipment and Bioanalyzer” on page 7, you can prepare the assay, load the chip, and run the assay, as described in the following procedures.

NOTE

If you use the RNA 6000 Nano kit for the first time, you must read these detailed instructions. If you have some experience, you might want to use the *RNA 6000 Nano Kit for 2100 Bioanalyzer Systems Quick Guide*.

Decontaminating the Electrodes

To avoid decomposition of your RNA sample, follow this electrode decontamination procedure on a daily basis before running any RNA Nano assays.

NOTE

Perform the following RNase decontamination procedure on a daily basis before running any assays.

- 1 Slowly fill one of the wells of an electrode cleaner with 350 μ L RNaseZAP.
- 2 Open the lid and place the electrode cleaner in the 2100 Bioanalyzer instrument.
- 3 Close the lid and leave it closed for about 1 minute.
- 4 Open the lid and remove the electrode cleaner. Label the electrode cleaner and keep it for future use. You can reuse the electrode cleaner for all 25 chips in the kit.
- 5 Slowly fill one of the wells of another electrode cleaner with 350 μ L RNase-free water.
- 6 Place electrode cleaner in the 2100 Bioanalyzer instrument.
- 7 Close the lid and leave it closed for about 10 seconds.
- 8 Open the lid and remove the electrode cleaner. Label the electrode cleaner and keep it for future use.
- 9 Wait another 10 seconds to allow the water on the electrodes to evaporate before closing the lid.

Preparing the Gel

- 1 Allow all reagents to equilibrate to room temperature for 30 minutes before use.
- 2 Place 550 μL of RNA 6000 Nano gel matrix (red ●) into the top receptacle of a spin filter.
- 3 Place the spin filter in a microcentrifuge and spin for 10 minutes at $1500\text{ g} \pm 20\%$ (for Eppendorf microcentrifuge, this corresponds to 4000 rpm).
- 4 Aliquot 65 μL filtered gel into 0.5 mL RNase-free microcentrifuge tubes that are included in the kit. Store the aliquots at $2 - 8\text{ }^{\circ}\text{C}$ ($36 - 46\text{ }^{\circ}\text{F}$) and use them within one month of preparation.

Preparing the Gel-Dye Mix

WARNING**Handling Reagents**

The dye can cause eye irritation. Because the dye binds to nucleic acids, it should be treated as a potential mutagen.

Kit components contain DMSO. DMSO is skin-permeable and can elevate the permeability of other substances through the skin.

- ✓ Follow the appropriate safety procedures and wear personal protective equipment including protective gloves and clothes as well as eye protection.
- ✓ Follow good laboratory practices when preparing and handling reagents and samples.
- ✓ Always use reagents with appropriate care.
- ✓ For more information, refer to the material safety data sheet (MSDS) on www.agilent.com.

- 1 Allow all reagents to equilibrate to room temperature for 30 minutes before use. Protect the dye concentrate from light while bringing it to room temperature.
- 2 Vortex RNA 6000 Nano dye concentrate (blue ●) for 10 seconds and spin down.
- 3 Add 1 μ L of blue-capped RNA 6000 Nano dye concentrate (blue ●) to a 65 μ L aliquot of filtered gel (prepared as described in "Preparing the Gel" on page 15).
- 4 Cap the tube, vortex thoroughly and visually inspect proper mixing of gel and dye. Store the dye concentrate at 2 – 8 °C (36 – 46 °F) in the dark again.
- 5 Spin tube for 10 minutes at room temperature at 13000 g (for Eppendorf microcentrifuge, this corresponds to 14000 rpm). Use prepared gel-dye mix within one day.


**NOTE**

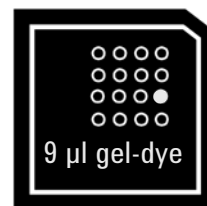
A larger volume of gel-dye mix can be prepared in multiples of the 65+1 ratio, if more than one chip will be used within one day. Always re-spin the gel-dye mix at 13000 g for 10 minutes before each use.

Loading the Gel-Dye Mix

NOTE

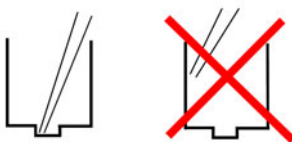
Before loading the gel-dye mix, make sure that the base plate of the chip priming station is in position (C) and the adjustable clip is set to the top position. Refer to [“Setting up the Chip Priming Station”](#) on page 8 for details.

- 1 Allow the gel-dye mix to equilibrate to room temperature for 30 minutes before use. Protect the gel-dye mix from light during this time.
- 2 Take a new RNA chip out of its sealed bag.
- 3 Place the chip on the chip priming station.
- 4 Pipette 9.0 μ L of the gel-dye mix at the bottom of the well marked  and dispense the gel-dye mix.

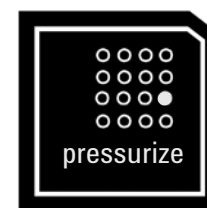


NOTE

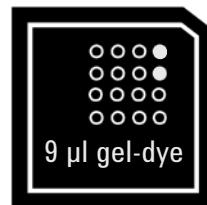
When pipetting the gel-dye mix, make sure not to draw up particles that may sit at the bottom of the gel-dye mix vial. Insert the tip of the pipette to the bottom of the chip well when dispensing. This prevents a large air bubble forming under the gel-dye mix. Placing the pipette at the edge of the well may lead to poor results.



- 5 Set the timer to 30 seconds, make sure that the plunger is positioned at 1 mL and then close the chip priming station. The lock of the latch will click when the chip priming station is closed correctly.
- 6 Press the plunger of the syringe down until it is held by the clip.
- 7 Wait for exactly 30 seconds and then release the plunger with the clip release mechanism.
- 8 Visually inspect that the plunger moves back at least to the 0.3 mL mark.
- 9 Wait for 5 seconds, then slowly pull back the plunger to the 1 mL position.




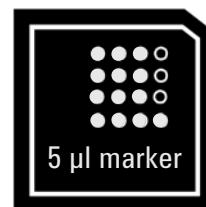
- 10 Open the chip priming station.
- 11 Pipette 9.0 μ L of the gel-dye mix in each of the wells marked **G**

**NOTE**

Please discard the remaining vial with gel-dye mix.

Loading the RNA 6000 Nano Marker

- 1 Pipette 5 μ L of the green-capped RNA 6000 Nano marker (green ●) into the well marked with the ladder symbol  and each of the 12 sample wells.


**NOTE**

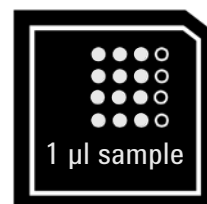
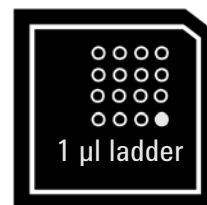
Do not leave any wells empty or the chip will not run properly. Unused wells must be filled with 5 μ L of the green-capped RNA 6000 Nano marker (green ●) plus 1 μ L of the buffer in which the samples are diluted.

Loading the Ladder and Samples

NOTE

Always use RNase-free microcentrifuge tubes, pipette tips and water.

- 1 Before use, thaw ladder aliquots and keep them on ice (avoid extensive warming upon thawing process).
- 2 To minimize secondary structure, heat denature (70 °C, 2 minutes) the samples before loading on the chip.
- 3 Pipette 1 μ L of the RNA ladder into the well marked with the ladder symbol .
- 4 Pipette 1 μ L of each sample into each of the 12 sample wells.
- 5 Set the timer to 60 seconds.
- 6 Place the chip horizontally in the adapter of the IKA vortex mixer and make sure not to damage the buldge that fixes the chip during vortexing.
If there is liquid spill at the top of the chip, carefully remove it with a tissue.



CAUTION

Wrong vortexing speed

If the vortexing speed is too high, liquid spill that disturbs the analysis may occur for samples generated with detergent containing buffers.

- ✓ **Reduce vortexing speed to 2000 rpm!**

- 7 Vortex for 60 seconds at 2400 rpm.
- 8 Refer to the next topic on how to insert the chip in the 2100 Bioanalyzer instrument. Make sure that the run is started within 5 minutes.

NOTE

Depending on the RNA isolation protocol, varying results can be expected. Known dependencies include: salt content, cell fixation method and tissue stain. Best results are achieved for RNA samples which are dissolved in deionized and RNase-free water. Avoid genomic DNA contamination by including DNase treatment in the preparation protocol.

Inserting a Chip in the 2100 Bioanalyzer Instrument

- 1 Open the lid of the 2100 Bioanalyzer instrument.
- 2 Check that the electrode cartridge is inserted properly. Refer to “Setting up the Bioanalyzer” on page 9 for details.
- 3 Place the chip carefully into the receptacle. The chip fits only one way.

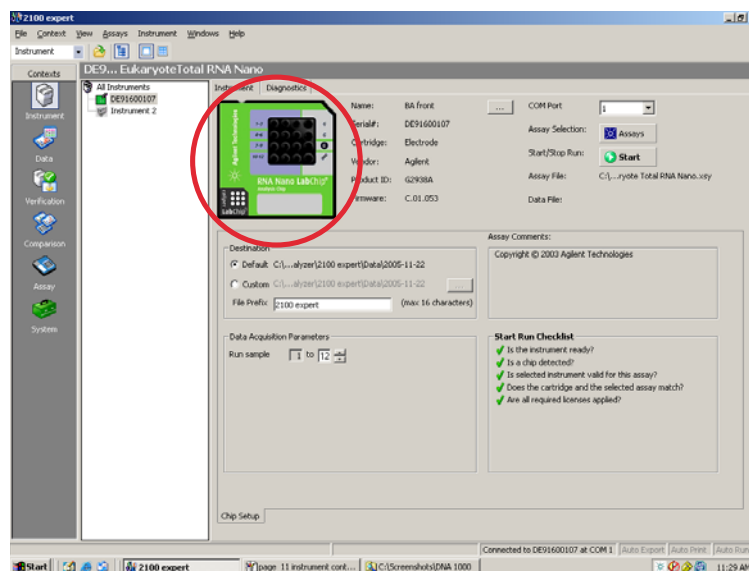
CAUTION

Sensitive electrodes and liquid spills

Forced closing of the lid may damage the electrodes and dropping the lid may cause liquid spills resulting in bad results.

- ✓ Do not use force to close the lid and do not drop the lid onto the inserted chip.

- 4 Carefully close the lid. The electrodes in the cartridge fit into the wells of the chip.
- 5 The 2100 Expert software screen shows that you have inserted a chip and closed the lid by displaying the chip icon at the top left of the **Instrument** context.

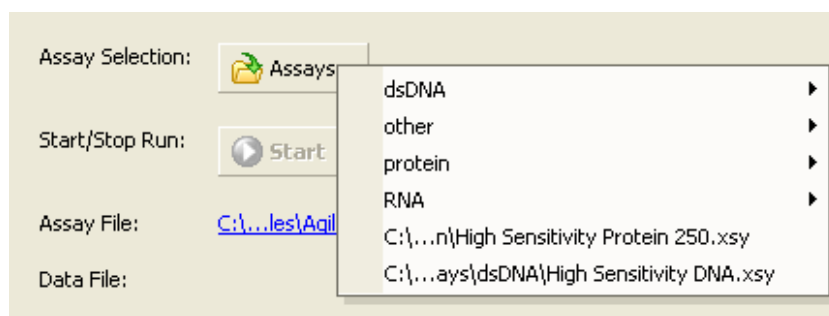


Starting the Chip Run

NOTE

Please note that the order of executing the chip run may change if the Agilent Security Pack software (only applicable for 2100 Expert software revision B.02.02 and higher) is installed. For more details please read the 'User's Guide' which is part of the Online Help of your 2100 Expert software.

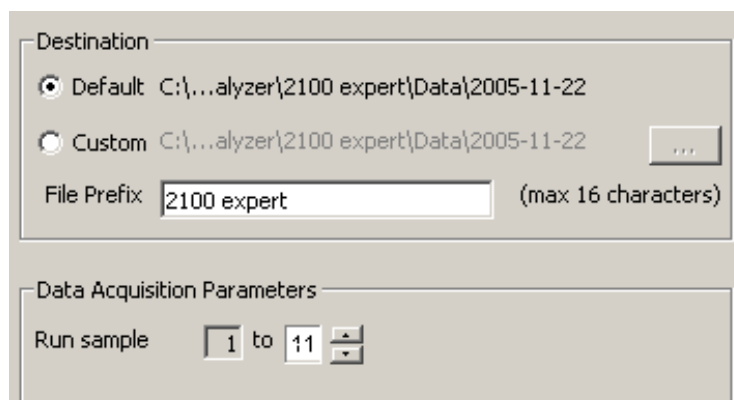
- 1 In the **Instrument** context, select the appropriate assay from the Assay menu.



- 2 Accept the current **File Prefix** or modify it.

NOTE

Run sample numbers can vary between assays.



Data will be saved automatically to a file with a name using the prefix you have just entered. At this time you can also customize the file storage location and the number of samples that will be analyzed.

- To enter sample information like sample names and comments, complete the sample name table.

	Sample Name	Sample Comment	Status	Observation	Result Label	Result Color
1	Sample 1					
2	Sample 2					
3	Sample 3					
4	Sample 4					
5	Sample 5					
6	Sample 6					
7	Sample 7					
8	Sample 8					
9	Sample 9					
10	Sample 10					
11	Sample 11					
12	Sample 12					

Chip Comments :

Sample Information

Study Information

- Click the **Start** button in the upper right of the window to start the chip run. The incoming raw signals are displayed in the **Instrument** context.



CAUTION

Contamination of electrodes

Leaving the chip for a period longer than 1 hour (e.g. over night) in the Bioanalyzer may cause contamination of the electrodes.

- ✓ Immediately remove the chip after a run.

- After the chip run is finished, remove the chip from the receptacle of the Bioanalyzer and dispose of it according to good laboratory practices.

Cleaning Electrodes after a RNA 6000 Nano Chip Run

When the assay is complete, *immediately* remove the used chip from the 2100 Bioanalyzer instrument and dispose of it according to good laboratory practice. After a chip run, perform the following procedure to ensure that the electrodes are clean (no residues are left over from the previous assay).

NOTE

Use a new electrode cleaner with each new kit.

CAUTION

Leak currents between electrodes

Liquid spill may cause leak currents between the electrodes.

✓ **Never fill more than 350 µL of water in the electrode cleaner.**

- 1 Slowly fill one of the wells of the electrode cleaner with 350 µL of fresh RNase-free water.
- 2 Open the lid and place the electrode cleaner in the 2100 Bioanalyzer instrument.
- 3 Close the lid and leave it closed for about 10 seconds.
- 4 Open the lid and remove the electrode cleaner.
- 5 Wait another 10 seconds to allow the water on the electrodes to evaporate before closing the lid.

NOTE

After 5 chip runs, empty and refill the electrode cleaner.

After 25 chip runs, replace the used electrode cleaner by a new one.

NOTE

Remove the RNase-free water out of the electrode cleaner at the end of the day. For a more thorough cleaning of the electrodes, refer to the *2100 Expert Maintenance and Troubleshooting Guide* which is part of the Online Help of the 2100 Bioanalyzer software.

7

Checking Your RNA 6000 Nano Assay Results

RNA 6000 Nano Ladder Well Results

To check the results of your run, select the Gel or Electropherogram tab in the **Data** context. The electropherogram of the ladder well window should resemble those shown below.

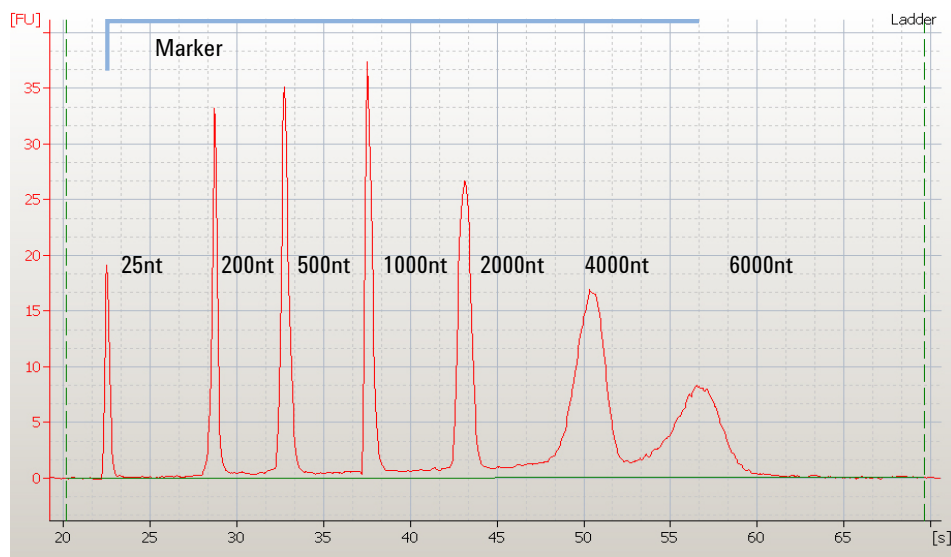


Figure 2 RNA 6000 Nano ladder

Major features of a successful ladder run are:

- 1 marker peak
- 6 RNA peaks (2100 Expert software calls for 5 first ladder peaks only)
- All 7 peaks are well resolved
- Correct peak size assignment in the electropherogram

If the electropherogram of the ladder well window does not resemble the one shown above, refer to the *2100 Expert Maintenance and Troubleshooting Guide* for assistance.

RNA 6000 Nano Sample Well Results

To review the results of a specific sample, select the sample name in the tree view and highlight the *Results* sub-tab. The electropherogram of the sample well window for total RNA (eukaryotic) should resemble the one shown here.

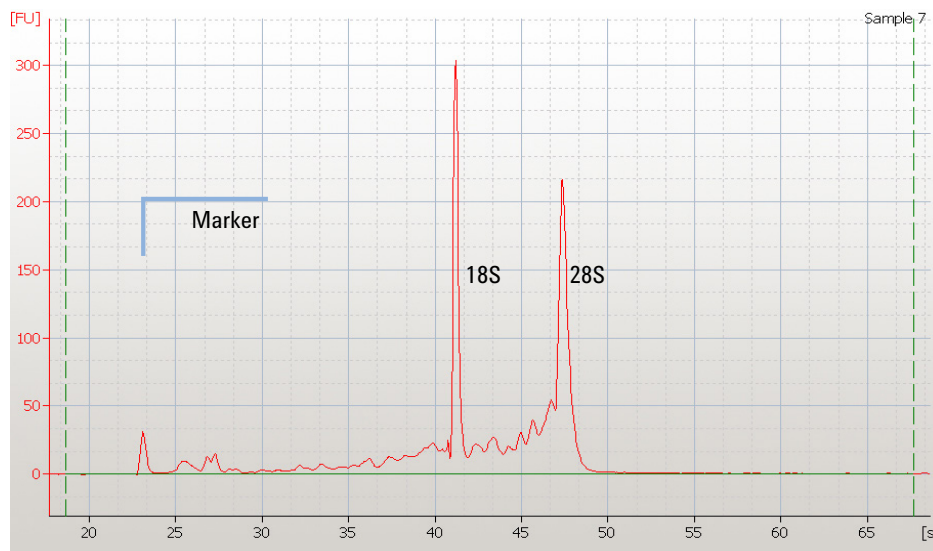


Figure 3 RNA peaks of a successful sample run

Major features of a successful total RNA run are:

- 1 marker peak
- 2 ribosomal peaks (with successful sample preparation)

By selecting the *Results* sub-tab, values for the calculated RNA concentration, the ribosomal ratio and the RNA Integrity Number (RIN), implemented with 2100 Expert software version B.02.02 or higher, are displayed.

NOTE

For information on the RNA data analysis and the RNA integrity number (RIN) please refer to the *2100 Expert User's Guide or Online Help*.

The electropherogram of the sample well window for mRNA should resemble the one shown here.

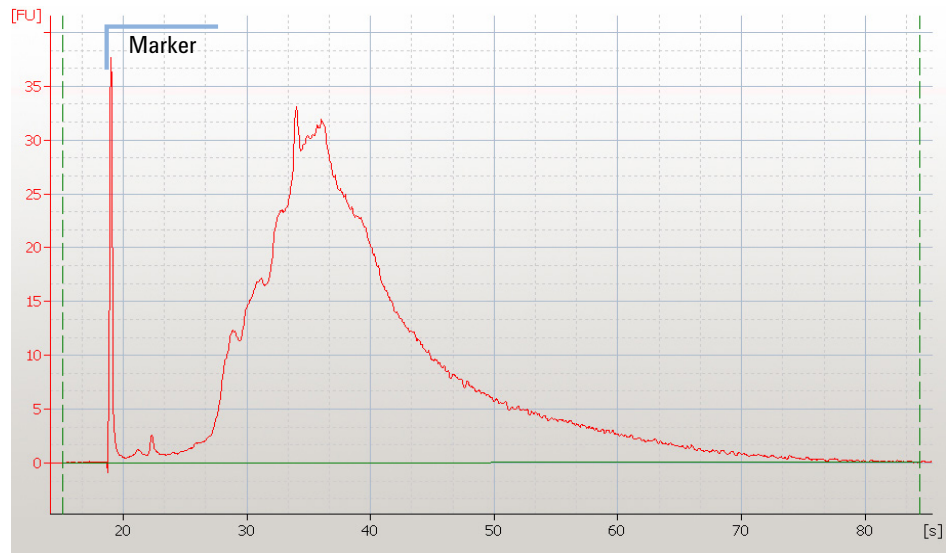


Figure 4 mRNA peaks of a successful sample run

Major features of a successful mRNA run are:

- 1 marker peak
- Broad hump (with successful sample preparation)
- Contamination with ribosomal RNA shown as 2 overlaid peaks (if present)

In This Book

you find the procedure to analyze RNA samples with the Agilent RNA 6000 Nano reagent kit and the Agilent 2100 Bioanalyzer instrument.

www.agilent.com

© Agilent Technologies Inc. 2000-2023
Edition: 09/2023

Document No: SD-UF0000075 Rev. E.00

