



Agilent dsDNA 930 Reagent Kit (75–20000 bp)

Kit Guide

For Research Use Only.

Not for use in diagnostic procedures.

DNF-930 dsDNA Reagent kit, 75 bp – 20,000 bp, 500 Samples (Part # DNF-930-K0500)

DNF-930 dsDNA Reagent kit, 75 bp – 20,000 bp, 1,000 Samples (Part # DNF-930-K1000)

Notices

Manual Part Number

M5310-91930

Edition 12/2018

Copyright

© Agilent Technologies, Inc. 2018

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.

Agilent Technologies, Inc.
5301 Stevens Creek Blvd.
Santa Clara, CA 95051
USA

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	Agilent dsDNA 930 Reagent Kit	4
2	Additional Material and Equipment Required	7
	Material and Equipment Required for Analysis with the Fragment Analyzer	7
	Additional Equipment/Reagents Required (Not Supplied)	8
3	Agilent dsDNA 930 Reagent Kit Protocol	9
	Gel Preparation	9
	Inlet Buffer Preparation	12
	Capillary Conditioning Solution Preparation	12
	Instrument Preparation	13
	Marker/Ladder/Sample Preparation	14
	General Information	14
	75 bp/20,000 bp Marker Preparation	14
	1 kb Plus DNA Ladder Preparation	15
	Sample Plate Preparation	16
	Performing Experiments	18
	Running an Experiment	18
	Viewing and Editing Experimental Methods	22
	Processing Experimental Data	25
	Fragment Analyzer Shut Down/Storage	26
	Instrument Shut Down/Storage	26
4	Checking Your Separation Results	27
	1 kb Plus DNA Ladder	27
5	Troubleshooting	29

1

Agilent dsDNA 930 Reagent Kit

The DNF-930 Reagent kit from Agilent is for the analysis of dsDNA fragments between 75 bp and 20 kb. Sizing and relative quantification between samples can be obtained using this kit. Example applications include PCR fragment sizing, and restriction digest analysis.

Table 1 Physical Specifications

Type	Specifications
Sample Volume Required	2 μ L (adjustable depending upon sample concentration)
Number of Samples per Run	12-Capillary: 11 (+ 1 well DNA Ladder) or 12 (Imported DNA Ladder) ² 48-Capillary: 47 (+ 1 well DNA Ladder) or 48 (Imported DNA Ladder) ² 96-Capillary: 95 (+ 1 well DNA Ladder) or 96 (Imported DNA Ladder) ²
Total Electrophoresis Run Time	18 min (22-47 Array) ³ 30 min (33-55 Array) 75 min (55-80 Array)

Table 2 Analytical Specifications

Type	Specifications
DNA Sizing Range	75 bp – 20,000 bp
DNA Sizing Precision ¹	5% CV
DNA Sizing Accuracy ¹	\pm 10% or better
Separation Resolution	75 bp – 1,500 bp \leq 10%; 1,500 bp – 20,000 bp \leq 15%
DNA Fragment Concentration Range ¹	0.5 ng/ μ L – 50 ng/ μ L input DNA (adjustable by dilution of sample)

¹ Results using DNA Ladder or DNA Fragment standards initially prepared in 1x TE buffer.

² Refer to Marker/Ladder/Sample Preparation section for information on using an imported DNA Ladder.

³The 22 cm effective, 47 cm total length capillary is only available for 12-capillary Fragment Analyzer instruments.

Agilent dsDNA 930 Reagent Kit

Table 3 Storage Conditions

Store at –20°C:	Store at 4°C (DO NOT FREEZE):	Store at Room Temperature (DO NOT FREEZE):
Intercalating Dye	dsDNA 930 Gel	5x Capillary Conditioning Solution
Markers, 75 bp & 20 kb	5x 930 dsDNA Inlet Buffer	Mineral Oil Dropper Bottle
1 kb Plus DNA Ladder	Dilution Buffer 1x TE	

Ensure all reagents are completely warmed to room temperature prior to use.

dsDNA 930 Reagent kit, 75 bp – 20,000 bp, 500 Samples (Part # DNF-930-K0500)

Table 4 kit Components

Part Number	Name	Amount
DNF-930-0240	dsDNA 930 Gel	240 mL
DNF-600-U030	Intercalating Dye	30 µL
DNF-355-0125	5x 930 dsDNA Inlet Buffer	125 mL (dilute with sub-micron filtered water prior to use)
DNF-475-0050	5x Capillary Conditioning Solution	50 mL (dilute with sub-micron filtered water prior to use)
FS-SMK930-0003	Markers, 75 bp and 20 kb <ul style="list-style-type: none"> • 0.5 ng/µL concentration each in 1x TE buffer 	3.2 mL
FS-SLR930-U100	1 kb Plus DNA Ladder <ul style="list-style-type: none"> • 75 bp – 20,000 bp; 50 ng/µL total DNA concentration in 1x TE buffer 	100 µL
FS-SM015	Mineral Oil Dropper Bottle	15 mL
DNF-495-0060	Dilution Buffer 1x TE	60 mL

Agilent dsDNA 930 Reagent Kit

dsDNA 930 Reagent kit, 75 bp – 20,000 bp, 1,000 Samples (Part # DNF-930-K1000)

Table 5 kit Components

Part Number	Name	Amount
DNF-930-0500	dsDNA 930 Gel	500 mL
DNF-600-U030	Intercalating Dye	30 µL x 2
DNF-355-0300	5x 930 dsDNA Inlet Buffer	300 mL (dilute with sub-micron filtered water prior to use)
DNF-475-0100	5x Capillary Conditioning Solution	100 mL (dilute with sub-micron filtered water prior to use)
FS-SMK930-0003	Markers, 75 bp and 20 kb <ul style="list-style-type: none">0.5 ng/µL concentration each in 1x TE buffer	3.2 mL
FS-SLR930-U100	1 kb Plus DNA Ladder Lower Marker <ul style="list-style-type: none">75 bp – 20,000 bp; 50 ng/µL total DNA concentration in 1x TE buffer	100 µL x 2
FS-SM015	Mineral Oil Dropper Bottle	15 mL
DNF-495-0125	Dilution Buffer 1x TE	125 mL

WARNING

Working with Chemicals

The handling of reagents and chemicals might hold health risks.

- Refer to product material safety datasheets for further chemical and biological safety information.
- Follow the appropriate safety procedures such as wearing goggles, safety gloves and protective clothing.

2 Additional Material and Equipment Required

Material and Equipment Required for Analysis with the Fragment Analyzer

Hardware:

- Fragment Analyzer with LED fluorescence detection:
 - 5200 Fragment Analyzer (Part # M5310AA)
 - 5300 Fragment Analyzer (Part # M5311AA)
 - 5400 Fragment Analyzer (Part # M5312AA)
- FA 12-Capillary Array Ultrashort, 22cm (Part # A2300-1250-2247), OR
- FA 12-Capillary Array Short, 33cm (Part # A2300-1250-3355), OR
- FA 12-Capillary Array Long, 55cm (Part # A2300-1250-5580), OR
- FA/ZAG 96-Capillary Array Short, 33cm (Part # A2300-9650-3355), OR
- FA/ZAG 96-Capillary Array Long, 55cm (Part # A2300-9650-5580)

Software:

- Fragment Analyzer instrument control software (Version 1.1.0.11 or higher)
- ProSize data analysis software (Version 2.0.0.51 or higher)

Reagents:

- Capillary Storage Solution, 100 mL (Part #GP-440-0100)

Additional Equipment/Reagents Required (Not Supplied)

- 96-well PCR sample plates. Please refer to Appendix 3 – Fragment Analyzer Compatible Plates and Tubes in the Fragment Analyzer User Manual for a complete approved sample plate list.
- Multichannel pipettor(s) and/or liquid handling device capable of dispensing 1 – 100 μ L volumes (sample plates) and 1,000 μ L volumes (Inlet Buffer plate)
- Pipette tips
- 96-well plate centrifuge (for spinning down bubbles from sample plates)
- Sub-micron filtered DI water system (for diluting the 5x 930 dsDNA Inlet Buffer and 5x Capillary Conditioning Solutions)
- Fisherbrand 96 DeepWell 1 mL Plate, Natural Polypropylene, part # 12-566-120 (Inlet Buffer and waste plate)
- Reagent Reservoir, 50 mL (VWR 82026-355 or similar) (for use in pipetting Inlet Buffer plates/sample trays)
- Conical centrifuge tubes for prepared Separation Gel/Dye mixture and/or 1x Capillary Conditioning Solution
 - 250 mL (for 96-Capillary instruments or larger volumes): Corning #430776, available from Fisher #05-538-53 or VWR #21008-771
 - 50 mL (for 12-Capillary instruments or 50 mL volumes): BD Falcon #352070, available from Fisher #14-432-22 or VWR #21008-940
- Clean graduated cylinder (for measurement of dsDNA Gel volume and dilution of 5x 930 dsDNA Inlet Buffer and 5x Capillary Conditioning Solution)

3 Agilent dsDNA 930 Reagent Kit Protocol

Gel Preparation

- 1 Store the dsDNA 930 Gel at 4°C upon arrival.
- 2 The Intercalating Dye is supplied as a 20,000x concentrate in DMSO and should be stored at -20°C.

NOTE

For this assay, it is recommended to use the Intercalating Dye at 2x normal concentration (1:10,000 dilution).

- 3 Bring the separation gel and Intercalating Dye to room temperature prior to mixing.
- 4 Mix appropriate volumes of Intercalating Dye and Separation Gel necessary for less than two weeks of operation. Use the supplied 50 mL conical centrifuge tube to allow a small minimum working volume. For larger volumes, use a 250 mL conical centrifuge tube and remove the collar of the tube holder in the instrument reagent compartment. For maximum accuracy, it is recommended to dispense Separation Gel into a clean glass graduated cylinder for volume measurement and transfer to the working tube prior to adding Intercalating Dye.

NOTE

Some loss of detection sensitivity will be observed over a two-week period after the gel/dye mixture has been prepared. For best results, it is recommended to prepare gel/dye mixture daily. It is not recommended to use gel/dye mixture that is more than two weeks old.

- 5 The volume of Separation Gel required per run varies between 12-capillary, and 96-capillary Fragment Analyzer systems. The volumes required are summarized below.

Agilent dsDNA 930 Reagent Kit Protocol

Table 6 Volume Specifications for 12-Capillary Fragment Analyzer Systems

# of Samples to be Analyzed ¹	Volume of Intercalating Dye	Volume of Separation Gel
12	1.0 µL	10 mL ¹
24	1.5 µL	15 mL
36	2.0 µL	20 mL
48	2.5 µL	25 mL
96	4.5 µL	45 mL

¹ A 5 mL minimum volume should be initially added to the tube. One sample well per separation is dedicated to the ladder.

Table 7 Volume Specifications for 96-Capillary Fragment Analyzer Systems

# of Samples to be Analyzed	Volume of Intercalating Dye	Volume of Separation Gel
96	4.0 µL	40 mL
192	8.0 µL	80 mL
288	12.0 µL	120 mL
384	16.0 µL	160 mL
480	20.0 µL	200 mL

- 6 Place the prepared Separation Gel/Intercalating Dye mixture onto the instrument and insert into the desired gel fluid line (Gel 1 or Gel 2 pump position). Ensure the fluid line is positioned at the bottom of the conical tube to avoid introducing air bubbles, which can cause pressurization errors.

Agilent dsDNA 930 Reagent Kit Protocol

- When adding Separation Gel to the instrument, update the solution levels in the Fragment Analyzer instrument control software. From the main menu, select **Utilities > Solution Levels**. A menu will be displayed to enter in the updated fluid levels (Figure 1).

Check the fluid volumes before proceeding. Ensure that the waste is empty and that the gel and conditioning solutions are full.

Record the solution volumes here:

	Volume (mL)	Solutions
Gel 1	50.0	DNF-930
Gel 2	43.3	NaOH
Conditioning Solution	41.6	
Waste	99.8	

OK Cancel

Figure 1 Solution Levels menu

- When switching applications (e.g., between kits), prime the appropriate gel fluid line after loading fresh Gel solution. From the main menu of the Fragment Analyzer instrument control software, select **Utilities > Prime...** Select the desired fluid line(s) (Conditioning, Gel 1, or Gel 2) and press **OK** to purge the fluid line with fresh gel (Figure 2).

Prime...

Fluid Selected

- Conditioning
- Gel1
- Gel2

Cycles: 1

Fill Rate: 300 $\mu\text{L/s}$

Empty Rate: 300 $\mu\text{L/s}$

OK Cancel

Figure 2. Prime menu

Inlet Buffer Preparation

- 1 Store the 5x 930 dsDNA Inlet Buffer at 4°C upon arrival. Do not freeze.
- 2 Bring the 5x 930 dsDNA Inlet Buffer to room temperature prior to mixing and use.
- 3 In a clean container, add 20 mL of the 5x 930 dsDNA Inlet Buffer per 80 mL of deionized sub-micron filtered water. Agitate to mix. The entire bottle can be mixed to 1x concentration and stored at 4°C if desired.

Capillary Conditioning Solution Preparation

- 1 Store the 5x Capillary Conditioning Solution at room temperature upon arrival. Do not freeze.
- 2 In a clean container (e.g. 50 mL or 250 mL conical centrifuge tube), add 20 mL of the 5x Capillary Conditioning Solution per 80 mL of deionized sub-micron filtered water. Agitate to mix. The entire bottle can be mixed to 1x concentration and stored at room temperature if desired.
- 3 Once mixed, place the 1x Capillary Conditioning Solution onto the instrument and insert the conditioning fluid line (Conditioning Solution pump position). Ensure the fluid line is positioned at the bottom of the conical tube to avoid introducing air bubbles, which can cause pressurization errors.
- 4 The 1x Capillary Conditioning Solution should be added to the system as use demands. A typical 12-capillary experiment cycle consumes less than 4 mL; a typical 96-capillary experiment consumes less than 35 mL.
- 5 When adding fresh 1x Capillary Conditioning Solution to the instrument, update the solution levels in the Fragment Analyzer instrument control software. From the main menu, select **Utilities > Solution Levels**. A menu will be displayed to enter in the updated fluid levels (Figure 1).

Instrument Preparation

- 1 Check the fluid level of the waste bottle and waste tray daily and empty as needed.
- 2 Prepare a fresh 96 DeepWell 1mL Plate filled with 1.0 mL/well of 1x 930 dsDNA Inlet Buffer daily.
 - 12-capillary system: row A only
 - 96-capillary system: all rowsDo not overfill the wells of the inlet buffer plate.

12-Capillary Systems:

- a In Row H of the same prepared buffer plate, place 1.1 mL/well of Capillary Storage Solution (Part # GP-440-0100). Row H of the buffer plate is used for the Store location, and the array moves to this position at the end of the experimental sequence.

96-Capillary Systems:

- a In the Sample 3 drawer, place a sample plate filled with 100 μ L/well of Capillary Storage Solution (Part# GP-440-0100). Sample 3 is used for the Store location, and the array moves to this position at the end of the experimental sequence.

NOTE

Ensure Row H of the buffer tray (12-capillary systems) or Sample 3 (96-capillary systems) is always filled with Capillary Storage Solution, and the capillary array is placed against Storage Solution when not in use, to prevent the capillary tips from drying out and potentially plugging.

- 3 Place the prepared inlet buffer plate into Drawer "B" (top drawer) of the Fragment Analyzer. Ensure that the plate is loaded with well A1 toward the back left on the tray.
- 4 Place an empty 96 DeepWell 1mL Plate into Drawer "W" (second from top) of the Fragment Analyzer. This plate serves as the capillary waste tray, and should be emptied daily. Alternatively, the supplied open reservoir waste plate may be used.

Marker/Ladder/Sample Preparation

General Information

- 1 The recommended 96-well sample plate for use with the Fragment Analyzer system is a semi-skirted PCR plate from Eppendorf (Part #951020303). Please refer to Appendix 3 – Fragment Analyzer Compatible Plates and Tubes in the Fragment Analyzer User Manual for a complete approved sample plate list. The system has been designed to operate using these dimensions/styles of PCR plates.

NOTE

The use of PCR plates with different dimensions to the above recommended plate could lead to decreased injection quality and consistency. Damage to the capillary array cartridge tips is also possible.

75 bp/20,000 bp Marker Preparation

- 1 Store the 75 bp and 20 kb Marker solution at -20°C upon arrival.
- 2 Bring the 75 bp and 20 kb Marker solution to room temperature prior to use; agitate solution to ensure it is properly mixed and centrifuge vial prior to dispensing.
- 3 The Marker solution is supplied as a ready-to-use solution, containing 0.5 ng/μL of each fragment in 1x TE buffer solution. It is intended for use as an external standard marker plate.
- 4 Prepare the Marker solution plate by dispensing 30 μL/well into Row A only (12-Capillary) or every well (96-Capillary) of a separate sample plate. Cover the wells with 20 μL/well of the supplied mineral oil to allow reuse for at least 30+ injections.
- 5 The prepared Marker solution plate should be placed into Drawer “M” (third from top) of the Fragment Analyzer. Ensure that the plate is loaded with well A1 toward the back left on the tray.

1 kb Plus DNA Ladder Preparation

- 1 Store the 1 kb Plus DNA Ladder solution at -20°C upon arrival.
- 2 Bring the 1 kb Plus DNA Ladder solution to room temperature prior to use; agitate solution to ensure it is properly mixed and centrifuge vial prior to dispensing.
- 3 The 1 kb Plus DNA Ladder solution is supplied as a **concentrate**. This enables the solution to be diluted with either 1x TE or 0.1x TE depending upon the available sample concentration and matrix (see section Sample Plate Preparation). The solution contains 50 ng/μL total DNA concentration in a 1x TE buffer solution. It is intended for use as a sizing standard for calibration of DNA size, and is typically loaded into 1 well of the sample plate:

12-Capillary System:

- a Pipette 1 kb Plus DNA Ladder solution into well 12 of each row of the sample plate.

96-Capillary System:

- a Pipette 1 kb Plus DNA Ladder solution into well H12 of the sample plate.
- 4 Prepare the working 1 kb Plus DNA Ladder solution by diluting with either 1x TE buffer or 0.1x TE buffer. Suggested dilutions are:
 - **When working with higher sample concentrations** (total input concentration > 10 ng/ μL):
Dilute 1 kb Plus DNA Ladder solution 12x with 1x TE buffer in sample well (2 μL 1 kb Plus DNA Ladder + 22 μL 1x TE buffer).
 - **When working with lower sample concentrations** (total input concentration < 10 ng/ μL):
Dilute 1 kb Plus DNA Ladder solution 50x with 0.1x TE buffer in sample well (1 μL 1 kb Plus DNA Ladder + 49 μL 0.1x TE buffer).
- 5 The highest level of sizing accuracy is obtained when the 1 kb Plus DNA Ladder is diluted to a similar concentration range and with a similar diluent (1x or 0.1x TE) to the samples being analyzed.
- 6 Alternatively, once the 1 kb Plus DNA Ladder has been run under the experimental method and additional samples are to be run under the same experimental conditions, the ladder can be imported in the ProSize software, enabling use of all 12 wells per row or all 96 wells of the sample plate

Sample Plate Preparation

Some suggested sample preparation guidelines are presented below. It may be necessary to adjust the sample dilution and diluent concentration (1x TE or 0.1x TE) depending upon initial sample concentration and sample matrix. For best results, the 1 kb Plus DNA Ladder should be prepared with a similar concentration and diluent to the samples.

- 1** If total initial sample concentration is > 10 ng/ μ L (e.g., PCR products):
 - a** Using a clean 96-well sample plate, pipette 22 μ L of supplied 1x TE buffer solution to each well including any ladder well(s).
 - b** Pipette 2 μ L of each DNA sample into the respective wells of the sample plate; mix the contents of the well using the pipette by aspiration/expulsion in the pipette tip.
 - c** 1 kb Plus DNA Ladder: Pipette 2 μ L of 1 kb Plus DNA Ladder solution into the respective well(s) of the sample plate to contain ladder; mix the contents of the well using the pipette by aspiration/expulsion in the pipette tip.
- 2** If total initial sample concentration is < 10 ng/ μ L (e.g., restriction digests):
 - a** Prepare a 0.1x TE solution by diluting the supplied 1x TE buffer 10x with deionized water.
 - b** Using a clean 96-well sample plate, pipette 20 μ L of the 0.1x TE buffer solution to each well to contain sample. Pipette 49 μ L of the 0.1x TE buffer solution to any well(s) to contain 1 kb Plus DNA Ladder.
 - c** Pipette 4 μ L of each DNA sample into the respective wells of the sample plate; mix the contents of the well using the pipette by aspiration/expulsion in the pipette tip.
 - d** 1 kb Plus DNA Ladder: Pipette 1 μ L of 1 kb Plus DNA Ladder solution into the respective well(s) of the sample plate to contain ladder; mix the contents of the well using the pipette by aspiration/expulsion in the pipette tip.

NOTE

Important Sample Mixing Information

When mixing sample with diluent solution, it is important to mix the contents of the well thoroughly to achieve the most accurate quantification. It is highly suggested to perform one of the following methods to ensure complete mixing:

- When adding 2 μL of sample to the 22 μL of diluent, **swirl the pipette tip** while pipetting up/down to further mix.
- After adding 2 μL of sample to the 22 μL of diluent, place a plate seal on the sample plate and **vortex the sample plate at 3000 rpm for 2 min**. Any suitable benchtop plate vortexer can be used. Ensure that there is no well-to-well transfer of samples when vortexing. The plate should be spun via a centrifuge after vortexing to ensure there are no trapped air bubbles in the wells.
- After adding 2 μL of sample to the 22 μL of diluent, **use a separate pipette tip set to a larger 20 μL volume**, and pipette each well up/down to further mix.
- Use an **electronic pipettor capable of mixing a 10 μL volume in the tip** after dispensing the 2 μL sample volume. Some models enable using the pipette tip for both adding and mixing.

- 3 After mixing the samples with 1x TE dilution buffer in each well, check the wells of the sample plate to ensure there are no air bubbles trapped in the bottom of the wells. If necessary, centrifuge the plate to remove any air bubbles. The presence of trapped air bubbles can lead to injection failures.
- 4 Run the sample plate immediately once prepared, or cover the sample plate with a cover film, store at 4°C, and use as soon as possible. Alternatively, to prevent evaporation, place a mineral oil overlay on each sample (20 μL /well).
- 5 To run the samples, place the plate in one of the three sample plate trays (Drawers 4-6 from the top) of the Fragment Analyzer instrument. Load or create the experimental method as described in the following sections.
- 6 The CCD detection system of the Fragment Analyzer system provides a high dynamic range for detection. An ideal injection range would yield peak heights from 100 – 20,000 RFUs. Overloading of sample can decrease separation resolution and saturate the detector, leading to mismatched lower/upper marker peak heights and poor results. It is important to optimize sample dilution and concentration, and use experimental parameters to work with within the specified RFU range. The highest level of sizing accuracy is obtained when the sample and DNA Ladder peak heights are of similar RFU peak heights.

TIP

If the above methods yield peak heights consistently above 20,000 RFUs, decrease the marker/sample injection time or reduce the sample volume to 1 μ L sample + 23 μ L 1x TE.

If low signals are encountered, increase the marker/sample injection time, or alternatively add 4 μ L of sample + 20 μ L of DI water in each well. When making adjustments to the sample dilution, the total volume should be maintained to at least 24 μ L..

NOTE

Whenever making adjustments to the sample dilution, ensure the **Dilution Factor** of the ProSize software is adjusted accordingly when processing the data.

Performing Experiments

Running an Experiment

- 1 To set up an experiment, from the main menu of the Fragment Analyzer instrument control software, select the **Operation** tab (Figure 3). Select the sample tray location to be analyzed (1, 2, or 3) by left clicking the **Sample Tray #** dropdown or by clicking the appropriate sample plate tab (alternate plate view) and choosing the appropriate location.

NOTE

For 96-Capillary Systems: Sample 3 is typically assigned to the Capillary Storage Solution.

- 2 Left click a well of the desired sample plate row with the mouse. The selected row will be highlighted in the plate map (e.g., Row A in Figure 3). Enter the sample name if desired into the respective **Sample ID** cell by left clicking the cell and typing in the name. Alternatively, sample information can be imported from .txt or .csv file by selecting the **Load from File...** option.

Agilent dsDNA 930 Reagent Kit Protocol

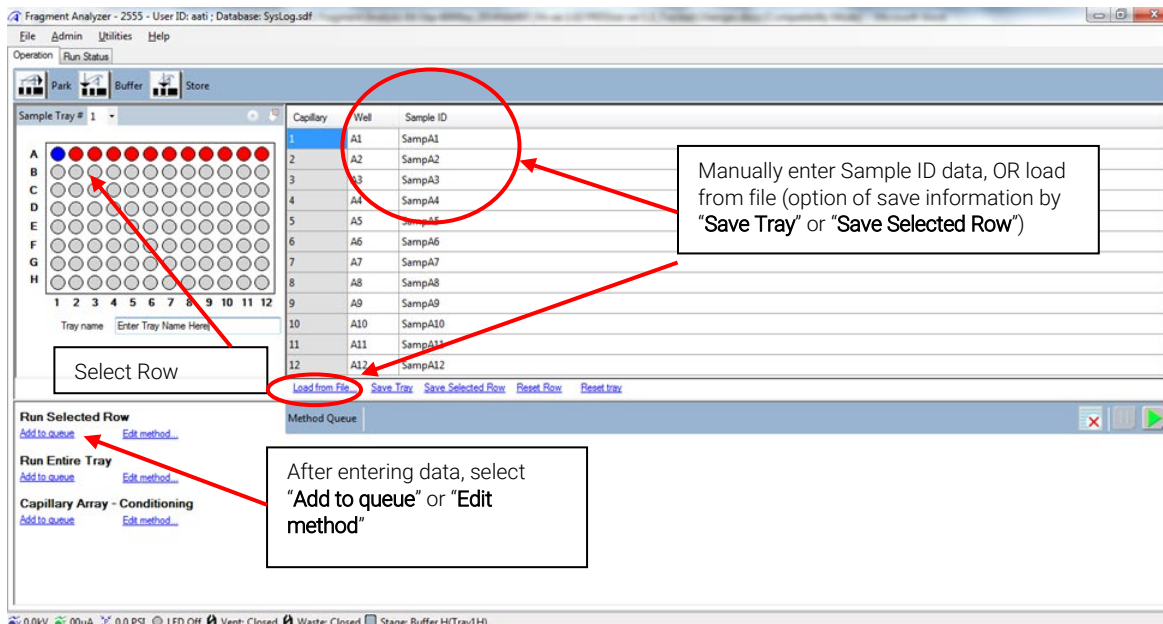


Figure 3. Main Screen showing selection of sample row and entering sample information

- 3 After sample information for the row or plate has been entered, under the **Run Selected Group** field press **Add to queue**. The **Separation Setup** form will be displayed enabling the user to select the experimental method and enter additional information (Figure 4).

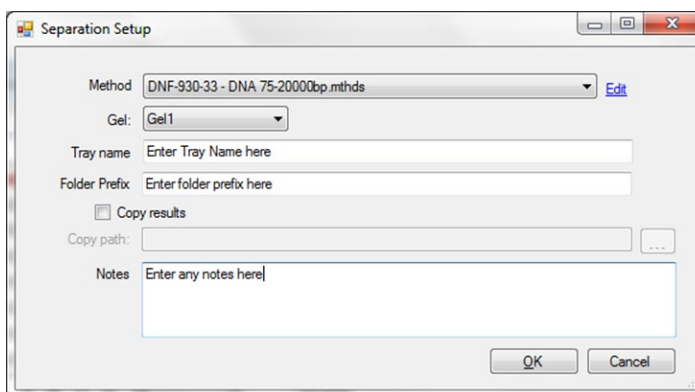





Figure 4. Separation Setup form to select experimental Method and enter tray/folder information

Agilent dsDNA 930 Reagent Kit Protocol

- 4 In the **Separation Setup** pop-up form, left click the dropdown and select the appropriate preloaded experimental **Method** file. The available methods are sorted by kit number and are linked to the directory containing methods for the currently installed capillary array length (e.g., 33cm or 55cm). Select the following method:
 - Select **DNF-930-33 – DNA 75-20000bp.mthds** when the 33 cm effective, 55 cm total “short” capillary array is installed.
 - Select **DNF-930-55 – DNA 75-20000bp.mthds** when the 55 cm effective, 80 cm total “long” capillary array is installed.
- 5 Select the appropriate **Gel** line being used for the experiment (Gel 1 or Gel 2) using the dropdown.
- 6 The **Tray Name** can be entered to identify the sample plate. The **Folder Prefix** if entered will amend the folder name (normally a time stamp of HH-MM-SS from the start of the CE run).
- 7 To copy the experimental results to another directory location in addition to the default save directory, check the **Copy results** box and select the desired **Copy path:** directory by clicking the ... button and navigating to the desired save directory.
- 8 Any **Notes** can be entered regarding the experiment; they will be saved and displayed in the final PDF report generated by the ProSize software.
- 9 Once all information has been entered, press **OK** to add the method to the instrument queue (press **Cancel** to abort adding the method).
- 10 Repeat Steps 3-9 for any remaining sample rows to be analyzed.
- 11 On 96-capillary systems, or in 12-capillary systems if the entire 96-well sample tray is to be run using the same experimental method, under the **Run Entire Tray** field press **Add to queue**. A form similar to Figure 4 will be displayed for entering information and adding the run to the instrument queue for the entire 96-well sample tray.
- 12 After a row or tray has been added to the queue, the method(s) will be listed on the main screen under the **Method Queue** field (Figure 5).
- 13 Prior to starting the experiment, verify all trays (buffer/storage, waste, marker, sample, etc.) have been loaded into their respective drawer locations.
- 14 Press the **Play** icon () to start the sequence loaded into the queue. To **Pause** the queue after the currently running experiment is completed, press the  button. To **Clear** the run queue of all loaded experiments, press the  button.

Agilent dsDNA 930 Reagent Kit Protocol

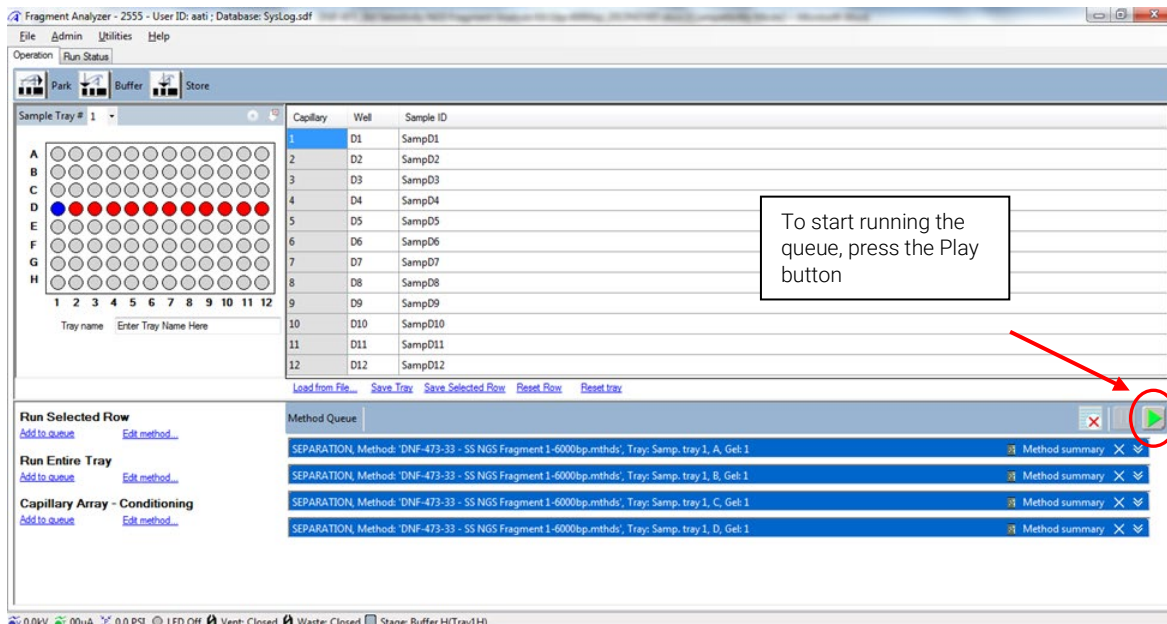


Figure 5. Main Screen after selection of samples to the run queue

- 15 Once an experiment has been loaded onto the queue, the user can view or edit the method (Administrator level only can edit a method) by pressing the **Method Summary** field. To remove the method from the queue, press the **x** button; to view the stepwise details of the method press the double down arrow icon.
- 16 The user may add a Pause or Prime step into the queue by right clicking the mouse while over the queue and selecting **Insert Pause** or **Insert Prime**.
- 17 The order of the experimental queue can be rearranged by dragging down individual entries. Further information regarding the **Method Queue** operation is provided in the Fragment Analyzer User Manual.
- 18 Once started, the instrument will perform all the programmed experiments in the **Method Queue** uninterrupted unless a pause step is present. Note that additional experiments can be programmed and added to the **Method Queue** at any time while the instrument is running if desired. After completion of the last queued experiment, the instrument stage will automatically move to the **Store** location (12-Capillary Systems: Row H of the inlet buffer tray containing the Capillary Storage Solution; 96-Capillary Systems: Sample 3 location)

Viewing and Editing Experimental Methods

- 1 A user level operator can **View** the steps of the experimental method by pressing the **View** link on the **Separation Setup** screen, or by pressing the **Method Summary** option once a method has been loaded onto the experimental queue. User level operators cannot edit any steps of a queued separation method.
- 2 Administrator level operators can **Edit** certain steps of the experimental method. To open the method editor screen, press the **Edit** link from the **Separation Setup** screen (Figure 4). The method editor screen is displayed, showing the steps of the method (Figure 6).
- 3 The preloaded, optimized steps for the **DNF-930-33** (Figure 6), **DNF-930-55** (Figure 7) methods are shown below. The general steps of the method are as follows:
 - a **Full Conditioning** flushing method (Automatically enabled).
Gel Selection: Gel 1.
 - b Perform Prerun (ENABLED) (8 kV, 30 sec)
 - c Rinse (DISABLED)
 - d Marker Injection (ENABLED) Voltage Injection (5 kV, 5 sec). This step injects the 75 bp/20,000 bp marker plate.
 - e Rinse (DISABLED)
 - f Sample Injection (ENABLED) Voltage Injection (5 kV, 5 sec). This step injects the prepared sample plate.
 - g Separation (ENABLED) Voltage (8 kV, 30 min). This step performs the CE Separation.

Agilent dsDNA 930 Reagent Kit Protocol

Separation Method:

Full Condition Gel prime to buffer Gel Selection: **Gel 1** ▼
 Gel Prime

Perform Prerun Voltage: **8.0** kV Time: **30** Sec

Rinse Tray: **Buffer** ▼ Row: **A** ▼ # Dips: **1**

Marker Injection Row: **A** ▼

Voltage Injection Voltage: **5.00** kV Time: **5** Sec
 Vacuum Injection Pressure: **-2.0** PSI

Rinse Tray: **Buffer** ▼ Row: **A** ▼ # Dips: **1**

Sample Injection

Voltage Injection Voltage: **5.00** kV Time: **5** Sec
 Vacuum Injection Pressure: **-2.0** PSI

Separation Voltage: **8.0** kV Time: **30.0** Min

OK **Cancel**

Figure 6. DNF-930-33 dsDNA Reagent kit (75 bp – 20,000 bp) method

Agilent dsDNA 930 Reagent Kit Protocol

- Figure 7 shows the preloaded method for the 55 cm effective, 80 cm total length “long” array. The Prerun and Separation voltage is set to 8 kV, the Injection voltage to 7.5 kV, and the Separation time to 75 min.

Separation Method:

Full Condition Gel prime to buffer Gel Selection: Gel 1

Gel Prime

Perform Prerun Voltage: 8.0 kV Time: 30 Sec

Rinse Tray: Buffer Row: A # Dips: 1

Marker Injection Row: A

Voltage Injection Voltage: 7.50 kV Time: 5 Sec

Vacuum Injection Pressure: -2.0 PSI

Rinse Tray: Buffer Row: A # Dips: 1

Sample Injection

Voltage Injection Voltage: 7.50 kV Time: 5 Sec

Vacuum Injection Pressure: -2.0 PSI

Separation Voltage: 8.0 kV Time: 75.0 Min

OK Cancel

Figure 7. DNF-930-55 dsDNA Reagent kit (75 p – 20,000 bp) method

- An Administrator level user has the option to adjust the **Gel Selection**; **Prerun** settings; **Rinse** settings including **Tray**, **Row** and **# Dips**; **Marker Injection** settings including **Row**; **Sample Injection** settings; and the **Separation** settings. For example, if the marker solution is loaded into a row other than Row A on a 12-capillary instrument, this can be adjusted prior to or while the method is loaded on the experimental queue
- To apply any adjustments to the method being placed on the experimental queue, press the **OK** button. To exit the editor screen without applying any changes press the **Cancel** button.

NOTE

Any edits made to the experimental method from the **Separation Setup** or **Method Summary** screen will only apply to the currently loaded experiment in the queue. No changes are made to the original separation method file.

Processing Experimental Data

- 1 When processing data, the ProSize software (Version 1.3 and higher) will automatically recognize the separation method performed and apply the appropriate matching configuration file from the *C:\ProSize 2.0\Configurations* directory:
 - The **DNF-930-33** separation method will be processed using the **DNF-930-33 – DNA 75-20000bp** configuration file.
 - The **DNF-930-55** separation method will be processed using the **DNF-930-55 – DNA 75-20000bp** configuration file.

NOTE

If the preloaded ProSize software configuration files “**DNF-930-33 – DNA 75-20000bp**” and “**DNF-930-55 – DNA 75-20000bp**” are not located in the *C:\ProSize 2.0\Configurations* directory, contact Agilent Technical Support to obtain these files.

- 2 The data is normalized to the lower marker (set to 75 bp) and upper marker (set to 20,000 bp), and calibrated to the 1 kb Plus DNA Ladder run in parallel to the samples. Figures 8-9 show examples of the 75 bp and 15,000 bp markers injected with the 1 kb Plus DNA Ladder. A total of 15 peaks should be observed.
- 3 When processing data, the ProSize software is set to the **DNA** mode in the **Advanced Settings**. The **Quantification** settings are set to **Use Lower Marker** for quantification with a **Final Conc. (ng/μL)** of 0.5 and a **Dilution Factor** of 12 (2 μL sample + 22 μL Diluent Marker). Note that if a pre-dilution was performed prior to the experiment, the **Dilution Factor** setting should be changed to reflect the estimated final sample concentration.

NOTE

UPDATE (June 25th 2013): The **Quantification** settings should now be set to **Use Lower Marker** for quantification from the previous **Use Upper Marker** setting.

- 4 For full information on processing data, refer to the ProSize User Manual.

Fragment Analyzer Shut Down/Storage

Instrument Shut Down/Storage

After each experiment, the instrument automatically places the capillary array in the **Store** position against Capillary Storage Solution:

- 12-capillary systems: row H of the buffer tray
- 96-capillary systems: sample 3

No further action is required.

If the instrument is to be idle for more than one day, it is recommended to turn off the power to the system.

4 Checking Your Separation Results

1 kb Plus DNA Ladder

Figure 8 shows the typical expected results for the 1 kb Plus DNA Ladder diluted 12x with 1x TE diluent, co-injected with the 75 bp lower marker and 20,000 bp upper marker using a 33 cm effective, 55 cm total length capillary array. A total of 15 peaks should be observed, with the sizes annotated as in Figure 8. All fragments in the ladder should be resolved.

Figure 9 shows the same separation performed on a 55 cm effective, 80 cm total length array.

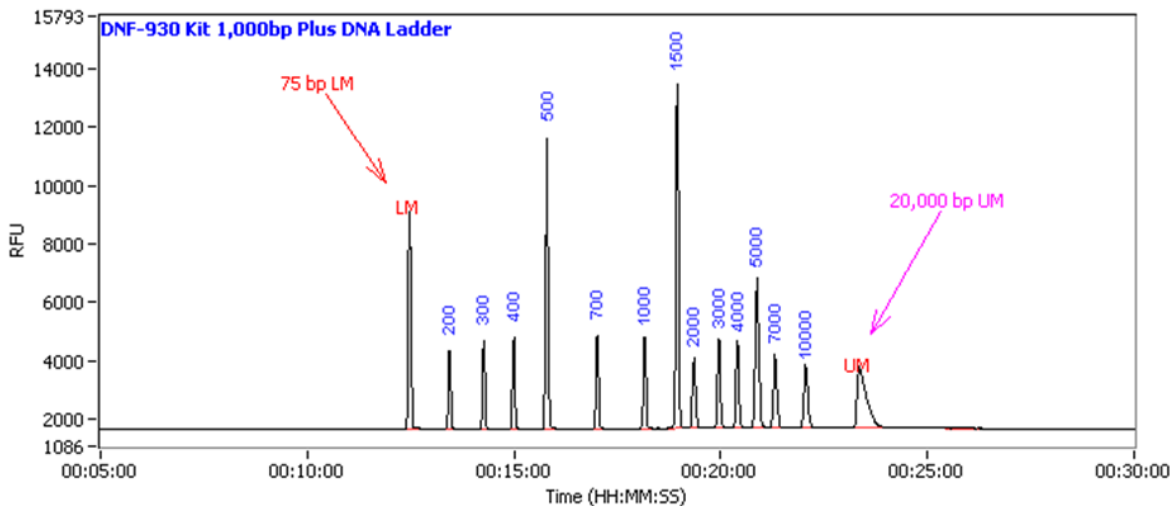


Figure 8. Representative 1 kb Plus DNA Ladder result injected with 75 bp lower marker and 20,000 bp upper marker, using the Fragment Analyzer system with the DNF-930 reagent kit. Array: 33 cm effective, 55 cm total length "short" array. Separation: 8 kV, 30 min.

Checking Your Separation Results

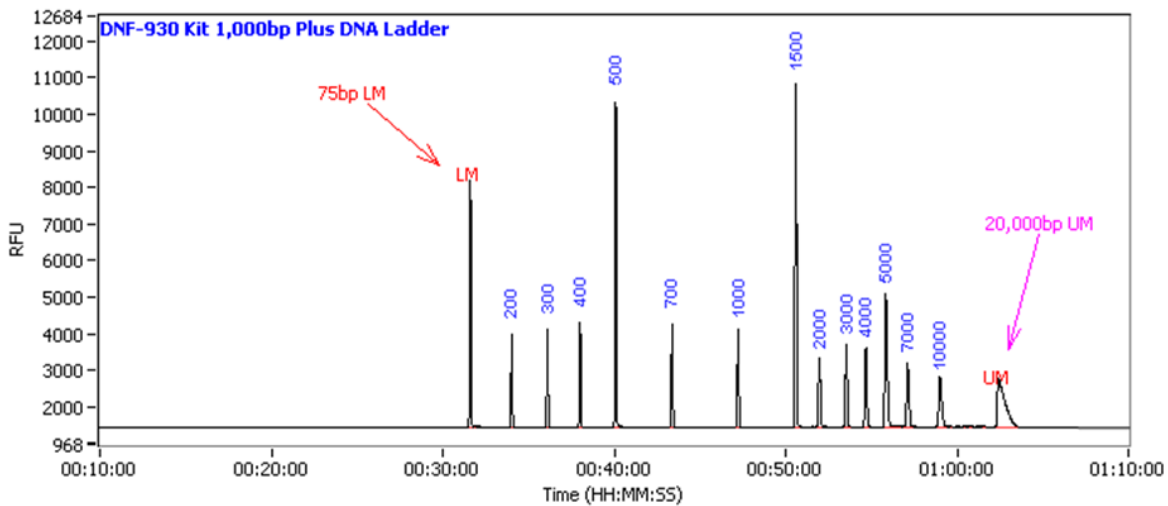


Figure 9. Representative 1 kb Plus DNA Ladder result injected with 75 bp lower marker and 20,000 bp upper marker, using the Fragment Analyzer system with the DNF-930 reagent kit. Array: 55 cm effective, 80 cm total length “short” array. Separation: 8 kV, 75 min.

5 Troubleshooting

The following table lists several potential assay specific issues which may be encountered when using the DNF-930 Reagent kit and suggested remedies. For a full list of instrument specific troubleshooting information, refer to the Troubleshooting and Maintenance Guide for the Fragment Analyzer system.

Table 8 Troubleshooting actions for assay specific issues

Issue	Cause	Corrective Action
The peak signal is >> 20,000 RFU; upper marker peak is low or not detected relative to lower marker.	1 Input DNA sample concentration is too high.	<ol style="list-style-type: none"> 1 Further dilute input DNA sample concentration with 1x TE buffer and repeat experiment. 2 Reduce injection time and/or injection voltage, and repeat experiment. Use the same injection voltage/time settings for the Marker Plate and Sample Plate to maximize quantification accuracy.
No peak observed for DNA sample when expected. Lower/Upper Marker peaks observed.	<ol style="list-style-type: none"> 1 Sample concentration too low and out of range. 2 Sample was not added to 1x TE diluent or not mixed well. 	<ol style="list-style-type: none"> 1 Prepare more concentrated sample and repeat experiment (e.g. 4 µL sample + 20 µL DI water); OR Repeat experiment using increased injection time and/or injection voltage for Marker Plate and Sample Plate. 2 Verify sample was correctly added and mixed to sample well.
Sample peak(s) migrate before or co-migrate with 75 bp Lower Marker.	1 Excess primer-dimer species in sample.	<ol style="list-style-type: none"> 1 Further dilute input DNA sample concentration with 1x TE buffer to minimize primer-dimer interference, and repeat experiment. 2 If fragment size is below 1,500 bp or 5,000 bp, analyze using DNF-910 Reagent kit (DNF-910-K0500; 35 bp – 1,500 bp range) or DNF-915 Reagent kit (DNF-915-K0500; 35 bp – 5,000 bp) to better resolve primer-dimer species.
Sample peak(s) migrate after or co-migrate with 20,000 bp Upper Marker.	1 DNA sample size out of range of assay.	1 Analyze samples with Genomic DNA Analysis kit (no upper marker), or the Large Fragment DNA Analysis kits, which have a large upper size limit.).
No sample peak or Lower Marker peak observed for individual sample.	<ol style="list-style-type: none"> 1 Air trapped at the bottom of sample plate well, or bubbles present in sample well. 2 Insufficient sample volume. A minimum of 20 µL is required. 3 Capillary is plugged. 	<ol style="list-style-type: none"> 1 Check sample plate wells for trapped air bubbles. Centrifuge plate. 2 Verify proper volume of solution was added to sample well. 3 Check waste plate for liquid in the capillary well. If no liquid is observed, follow the steps outlined in <i>Appendix G – Capillary Array Cleaning</i> of the Fragment Analyzer User Manual for unclogging a capillary array.

In This Book

This Kit Guide describes the following:

- Agilent dsDNA 930 Reagent Kit
- Additional Material and Equipment Required
- Agilent dsDNA 930 Reagent Kit Protocol
- Checking Your Separation Results
- Troubleshooting

www.agilent.com

© Agilent Technologies, Inc. 2018

Edition 12/2018



M5310-91930

