



Agilent DNF-910 dsDNA Reagent Kit (35-1500 bp) Quick Guide for the Fragment Analyzer Systems

The Agilent Fragment Analyzer systems are automated capillary electrophoresis platforms for scalable, flexible, fast, and reliable electrophoresis of nucleic acids.

This Quick Guide is intended for use with the Agilent 5200, 5300, and 5400 Fragment Analyzer systems only. The dsDNA 910 Reagent Kit from Agilent is for the analysis of dsDNA fragments between 35 bp and 1,500 bp. Sizing and relative quantification between samples can be obtained using this kit. Example applications include general PCR fragment sizing and QC, and genotyping.

Specifications

| Analytical specifications | dsDNA 910 assay |
|---|--|
| DNA Sizing Range | 35 bp – 1,500 bp |
| DNA Sizing Accuracy ¹ | + 5% or better |
| DNA Fragment Concentration Range ¹ | 0.5 ng/μL – 50 ng/μL input DNA (adjustable by dilution sample) |
| Separation Resolution | 35 bp – 100 bp ≤ 10%; 100 bp – 700 bp ≤ 5%; 700 -1,500 bp ≤ 5% (22-47 Array) 35 bp – 100 bp ≤ 10%; 100 bp – 1,500 bp ≤ 5% (33-55 Array) 35 bp – 100 bp ≤ 10%; 100 bp – 1,500 bp ≤ 5% (55-80 Array) |
| DNA Sizing Precision ¹ | 2% CV |

Physical Specifications

| | |
|--------------------------------|--|
| Total electrophoresis run time | 22cm ² : 22 minutes, 33cm: 45 minutes, 55cm: 70 minutes |
| Samples per run | 12, 48 or 96; depending on the instrument type |
| Sample volume required | 2 μL (adjustable depending on sample concentration) |
| Kit stability | 4 months |

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

² The 22 cm effective, 47 cm total length capillary is only available for 12-capillary Fragment Analyzer instruments

Kit Components – 500 Sample Kit

| Kit Component Number | Part Number (Re-order Number) | Description | Quantity Per Kit |
|----------------------|-------------------------------|--|------------------|
| 5191-6595* | | dsDNA 910/915, 500, 4°C | |
| | DNF-810-0240 | dsDNA 810 Gel, 240 mL | 1 |
| | DNF-355-0125 | 5x 930 dsDNA Inlet Buffer, 125 mL <ul style="list-style-type: none"> Dilute with sub-micron filtered water prior to use | 1 |
| | DNF-495-0060 | Dilution Buffer 1X TE, 60mL | 1 |
| DNF-910-FR* | | dsDNA 910 (35-1500), FR | |
| | DNF-600-U030 | Intercalating Dye, 30 µL | 1 |
| | FS-SLR910-0001 | 100bp DNA Ladder, 1mL | 1 |
| | FS-SMK910-0003 | Markers, 35bp & 1500bp, 3.2 mL | 1 |
| 5191-6614* | | Qualitative DNA, 500, RT | |
| | FS-SM015 | Mineral Oil Dropper Bottle, 15mL | 1 |
| | DNF-475-0050 | 5x Capillary Conditioning Soln, 50 mL | 1 |

*not orderable

WARNING

- Refer to product safety data sheets for further information
- When working with the Fragment Analyzer kit components follow the appropriate safety procedures such as wearing goggles, safety gloves and protective clothing.

Kit Components – 1000 Sample Kit

| Kit Component Number | Part Number (Re-order Number) | Description | Quantity Per Kit |
|----------------------|-------------------------------|--|------------------|
| 5191-6596* | | dsDNA 910/915, 1000, 4°C | |
| | DNF-810-0500 | dsDNA 910 Gel, 500 mL | 1 |
| | DNF-355-0300 | 5x 930 dsDNA Inlet Buffer, 300 mL <ul style="list-style-type: none"> Dilute with sub-micron filtered water prior to use | 1 |
| | DNF-495-0125 | Dilution Buffer 1X TE, 125mL | 1 |
| DNF-910-FR* | | dsDNA 910 (35bp-1500bp), FR | |
| | DNF-600-U030 | Intercalating Dye, 30 µL | 2 |
| | FS-SLR910-0001 | 100bp DNA Ladder, 1mL | 2 |
| | FS-SMK910-0003 | Markers, 35bp & 1500bp, 3.2 mL | 2 |
| 5191-6615* | | Qualitative DNA, 1000-5000, RT | |
| | FS-SM015 | Mineral Oil Dropper Bottle, 15mL | 1 |
| | DNF-475-0100 | 5x Capillary Conditioning Soln, 100 mL | 1 |

*not orderable

WARNING

- Refer to product safety data sheets for further information
- When working with the Fragment Analyzer kit components follow the appropriate safety procedures such as wearing goggles, safety gloves and protective clothing.

Additional Material Required for Analysis with the Fragment Analyzer Systems

- Fragment Analyzer systems with LED fluorescence detection:
 - 5200 Fragment Analyzer system (p/n M5310AA)
 - FA 12-Capillary Array Ultrashort, 22 cm (p/n A2300-1250-2247) OR
 - FA 12-Capillary Array Short, 33 cm (p/n A2300-1250-3355) OR
 - FA 12-Capillary Array Long, 55 cm (p/n A2300-1250-5580)
 - 5300 Fragment Analyzer system (p/n M5311AA)
 - FA 48-Capillary Array Short, 33 cm (p/n A2300-4850-3355) OR
 - FA/ZAG 96-Capillary Array Short, 33 cm (p/n A2300-9650-3355) OR
 - FA/ZAG 96-Capillary Array Long, 55 cm (p/n A2300-9650-5580)
 - 5400 Fragment Analyzer system (p/n M5312AA)
 - FA 48-Capillary Array Short, 33 cm (p/n A2300-4850-3355) OR
 - FA/ZAG 96-Capillary Array Short, 33 cm (p/n A2300-9650-3355) OR
 - FA/ZAG 96-Capillary Array Long, 55 cm (p/n A2300-9650-5580):
 - Agilent Fragment Analyzer controller software (Version 1.1.0.11 or higher)
 - Agilent ProSize data analysis software (Version 2.0.0.61 or higher)

Additional equipment/reagents required (not supplied)

- 96-well PCR sample plates. Please refer to Appendix – Fragment Analyzer Compatible Plates and Tubes in the Fragment Analyzer System User Manual for a complete approved sample plate list
- Multichannel pipettor(s) and/or liquid handling device capable of dispensing 1 – 100 µL volumes (sample plates) and 1,000 µL 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 Solution)
- 96-deepwell 1mL plate: Fisher Scientific #12-566-120 (inlet buffer and/or waste plate)
- Reagent reservoir, 50 mL (VWR #89094-680 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
 - 50 mL (for 5200 Fragment Analyzer system or 50 mL volumes): BD Falcon #352070, available from Fisher Scientific #14-432-22 or VWR #21008-940
 - 250 mL (for 5300 and 5400 Fragment Analyzer systems or larger volumes): Corning #430776, available from Fisher Scientific #05-538-53 or VWR #21008-771
- Vortexer (for mixing of samples, ladders, and/or markers in tubes and/or plates)
- Capillary Storage Solution (p/n GP-440-0100)

Essential Measurement Practices

| | |
|---|--|
| Environmental conditions | <ul style="list-style-type: none"> • Ambient operating temperature: 19 – 25 °C (66 – 77 °F) • Keep reagents during sample preparation at room temperature |
| Steps before sample preparation | <ul style="list-style-type: none"> • Allow reagents to equilibrate at room temperature for 30 min prior to use |
| Pipetting practice | <ul style="list-style-type: none"> • Pipette reagents carefully against the side of the 96-well sample plate or sample tube • Ensure that no sample or Diluent Marker remains within or on the outside of the tip |
| Mixing and centrifugation recommendations | <ul style="list-style-type: none"> • Apply a new seal to 96-well sample plate prior to mixing and centrifugation • When mixing, 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. After mixing, briefly centrifuge and visually confirm that all liquid is collected at the bottom of the 96-well sample plate or tube strips and any air bubble is removed <ul style="list-style-type: none"> • Place a plate seal on the sample plate and vortex the sample plate at 3,000 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. • Use a separate pipette tip set to a larger 20 µL volume, and pipette each well up/down to further mix. • Run samples immediately after preparation, or within a day with oil overlay. If not using right away, cover and keep at 4°C, warm to RT and centrifuge before running plate |

35 bp/1,500 bp Marker Preparation

1. Store the 35 bp and 1,500 bp Marker solution at -20°C upon arrival.
2. Bring the 35 bp and 1,500 bp 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 a 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), Rows A-D (48-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 the plate is loaded with well A1 toward the back left

100 bp DNA Ladder

1. Store the 100 bp DNA Ladder solution at -20°C upon arrival.
2. Bring the 100 bp 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 100 bp DNA Ladder solution is supplied as a ready-to-use solution, containing approximately 2.0 ng/μL total DNA concentration in a 1x TE buffer solution. It is used for calibrating the size of analyzed DNA fragments, and is typically added to a well of the sample plate and run in parallel with the samples:

12-Capillary System: Well 12 of each row to be analyzed

48-Capillary System:

- a) Well D12 if samples are in Row A to Row D, or
- b) Well H12 if samples are in Row E to Row H

96-Capillary System: Well H12

4. Alternatively, once the 100 bp 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, 48 or 96 wells of the sample plate. Refer to the ProSize software User Manual for information on exporting and importing calibration ladders.

Gel preparation

Prepare gel/dye mixture for 5200, 5300, and 5400 Fragment Analyzer Systems. To ensure the gel/dye mixture is mixed homogeneously without generating bubbles, gently invert the centrifuge tube 5 to 10 times, depending on the volume of the mixture. **NOTE:** Centrifuge dye prior to opening the vial to reduce risk of leaking.

5200 Fragment Analyzer system volume specifications

| # of Samples to be Analyzed ¹ | Volume of Intercalating Dye | Volume of Separation Gel ² | Volume of 1x Conditioning Solution ² |
|--|-----------------------------|---------------------------------------|---|
| 12 | 1.0 µL | 10 mL | 10 mL |
| 24 | 1.5 µL | 15 mL | 15 mL |
| 36 | 2.0 µL | 20 mL | 20 mL |
| 48 | 2.5 µL | 25 mL | 25 mL |
| 96 | 4.5 µL | 45 mL | 45 mL |

¹ One sample well per separation is dedicated to the ladder.

² A 5 mL minimum volume in the tube is included.

5300 Fragment Analyzer system volume specifications with 48-capillary array

| # of Samples to be Analyzed ¹ | Volume of Intercalating Dye | Volume of Separation Gel ² | Volume of 1x Conditioning Solution ² |
|--|-----------------------------|---------------------------------------|---|
| 48 | 2.5 µL | 25 mL | 25 mL |
| 96 | 4.0 µL | 40 mL | 40 mL |
| 144 | 5.5 µL | 55 mL | 55 mL |
| 192 | 7.0 µL | 70 mL | 70 mL |
| 240 | 8.5 µL | 85 mL | 85 mL |
| 288 | 10.0 µL | 100 mL | 100 mL |

¹ One sample well per separation is dedicated to the ladder.

² A 5 mL minimum volume in the tube is included.

5300 and 5400 Fragment Analyzer systems volume specifications with 96-capillary arrays

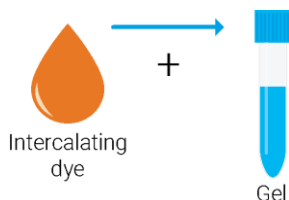
| # of Samples to be Analyzed ¹ | Volume of Intercalating Dye | Volume of Separation Gel ² | Volume of 1x Conditioning Solution ² |
|--|-----------------------------|---------------------------------------|---|
| 96 | 4.0 µL | 40 mL | 40 mL |
| 192 | 8.0 µL | 80 mL | 80 mL |
| 288 | 12.0 µL | 120 mL | 120 mL |
| 384 | 16.0 µL | 160 mL | 160 mL |
| 480 | 20.0 µL | 200 mL | 200 mL |

¹ One sample well per separation is dedicated to the ladder.

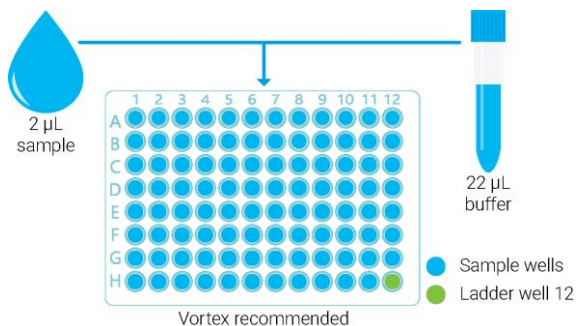
² A 5 mL minimum volume in the tube is included.

Agilent dsDNA 910 (35-1500bp) DNF-910 assay operating procedure

- Mix fresh gel and dye according to the volumes in the Gel preparation tables. Refill 1x Capillary Conditioning Solution as needed.



- Place a fresh 1x 930 dsDNA Inlet Buffer in drawer 'B' on the system, 1.0 mL/well. Replace daily.
 - 5200 system; Fill row A of buffer plate
 - 5300 system - 48 capillary; Fill rows A-D of buffer plate
 - 5300/5400 system - 96 capillary; Fill all rows of buffer plate
- Prepare Capillary Storage Solution plate. Replace every 2-4 weeks for optimal results.
 - 5200 system; Fill row H of buffer plate with 1.0mL/well, place in drawer "B"
 - 5300 system - 48 capillary; Fill rows A-D of a sample plate with 100 μ L/well, place in drawer '3'
 - 5300/5400 system - 96 capillary; Fill all rows of a sample plate with 100 μ L/well, place in drawer '3'
 - 5400 system; place in drawer "S"
- Prepare Marker plate and place in drawer 'M' on the system, 30 μ L/well. Add 1 drop or \sim 30 μ L of mineral oil to each well. The marker plate should last for 30+ injections or \sim 1 month.
 - 5200 system; Fill row A of sample plate
 - 5300 system - 48 capillary; Fill rows A-D of sample plate
 - 5300/5400 system - 96 capillary; Fill all rows of sample plate
- Mix samples with Diluent Buffer 1x TE in sample plate. Add ready to use ladder in corresponding well, dependent on the capillary size.



5200 system; Ladder – well 12, depending on which row is chosen

5300 system - 48 capillary; Ladder – well D12 or H12, depending on which group is chosen

5300/5400 system - 96 capillary; Ladder – well H12


WARNING

Working with Chemicals
The handling of reagents and chemicals might hold health risks.

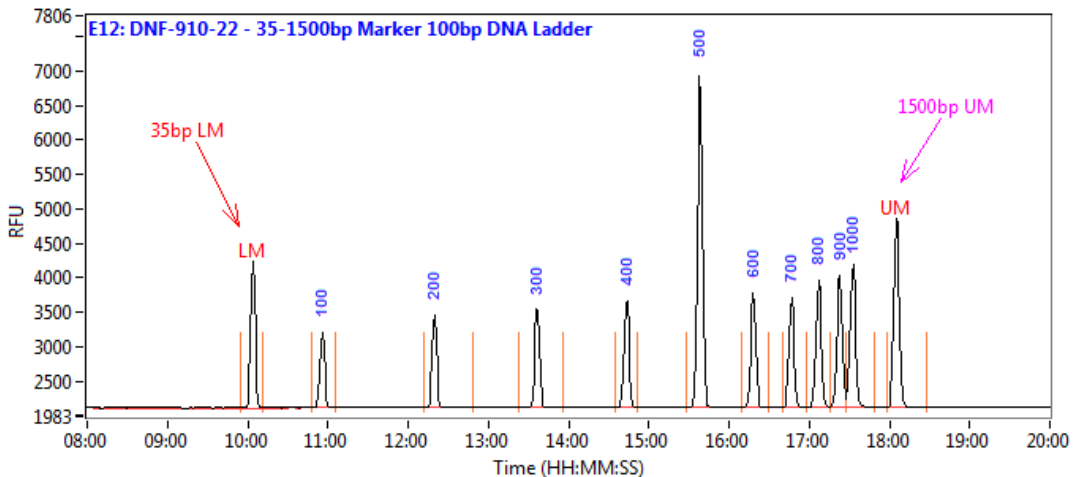
- Refer to product material safety datasheets for further chemical and biological safety information.

- Follow the appropriate safety procedures such as wearing goggles, safety gloves and protective clothing.

Agilent Fragment Analyzer software operating procedure

1. Select Row, Group or Tray to run.
2. Enter **sample ID** and **Tray ID**(optional).
3. Select **Add to Queue**, from the dropdown menus select the corresponding method based on your capillary length;
 - 3.1 DNF-910-22 – DNA 35-1500 bp
 - 3.2 DNF-910-33 – DNA 35-1500 bp **OR** DNF-910-33 – DNA 35-1500 bp-3kv
 - 3.3 DNF-910-55 – DNA 35-1500 bp
4. Enter **Tray Name**, **Folder Prefix**, and **Notes** (optional).
5. Select **OK** to add method to the queue.
6. Select  to start the separation.

100 bp DNA Ladder result



Representative 100 bp DNA Ladder result double injected with 35 bp lower marker and 1,500 bp upper marker, using the Fragment Analyzer system with the dsDNA 910 Reagent Kit. Method: DNF-910-22 (“ultra-short” array).

Troubleshooting

The following table lists several potential assay specific issues which may be encountered when using the dsDNA 910 Reagent Kit and suggested remedies. Contact Agilent technical support if you have any additional troubleshooting or maintenance questions.

| 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. | 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. |
| Sample peak(s) migrate before or co-migrate with 35 bp Lower Marker. | 1 Excess primer-dimer species in sample. | 1 Further dilute input DNA sample concentration with 1x TE buffer to minimize primer-dimer interference and repeat experiment. |
| Sample peak(s) migrate after or co-migrate with 1,500 bp Upper Marker. | 1 DNA sample size out of range of assay. | 1 Analyze samples with dsDNA 915 Reagent Kit, 35 bp – 5,000 bp (DNF-915), dsDNA 920 Reagent Kit, 75 bp – 15,000 bp (DNF-920), or dsDNA 930 Reagent Kit, 75 bp – 20,000 bp (DNF-930). |
| No peak observed for DNA sample when expected. Lower/Upper Marker peaks observed. | 1 Sample concentration too low and out of range. 2 Sample was not added to 1x TE diluent or not mixed well. | 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. |
| No sample peak or marker peak observed for individual sample. | 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. | 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 the Appendix – Capillary Array Cleaning of the Fragment Analyzer User Manual for |

unclogging a capillary array.

For Research Use Only

Not for use in Diagnostic Procedures.

Technical Support and Further Information

For technical support, please visit www.agilent.com. It offers useful information and support about the products and technology.

www.agilent.com

© Agilent Technologies, Inc. 2021

Edition 08/22

SD-AT000121

