

Cell-free DNA ScreenTape Assay for TapeStation Systems

Quick Guide

The Agilent 4150 (G2992AA) and 4200 (G2991AA and G2991BA) TapeStation systems are automated platforms for scalable, flexible, fast, and reliable electrophoresis of nucleic acids. This Quick Guide is intended for use with the Agilent 4150 and 4200 TapeStation systems only. The Cell-free DNA ScreenTape assay is designed for analyzing cell-free DNA samples and double-stranded DNA molecules from 50 to 800 basepairs.

Specifications

Analytical specifications	Cell-free DNA ScreenTape assay
Sizing range	50 – 800 bp
Sensitivity ¹	20 pg/μL
Sizing precision ²	10 % CV
Sizing accuracy ^{2,3}	±15 %
Quantitative precision ²	15 % CV
Quantitative accuracy ²	±20 %
Quantitative range	100 – 4000 pg/μL
%cfDNA functional range	100 – 5000 pg/μL
Maximum buffer concentration in sample	25 mM NaCl, 25 mM KCl, 3 mM EDTA, 0.1 % NaN ₃ , 5 mM phosphate buffer, 10 % ethanol, 10 % 2-propanol
Physical specifications	
Analysis time	16 samples: <25 min, 96 samples: <150 min
Samples per consumable	16
Sample volume required	2 μL
Kit stability	6 months
Kit size	112 samples

¹ Signal-to-noise >3 (single peak)

² Determined using the Cell-free DNA Ladder as sample





³ Sizing Accuracy for analysis with electronic ladder: ±20 %

Cell-free DNA ScreenTape Assay for TapeStation Systems

Storage Conditions

- Reagent vials and ScreenTape devices: 2 – 8 °C (36 – 46 °F).
- Store partially used ScreenTape devices upright at 2 – 8 °C (36 – 46 °F) for a maximum of 2 weeks.
- Never freeze ScreenTape devices. Discard any accidentally frozen ScreenTape devices.

Kit Components

Part Number	Name	Color	Amount
5067-5630	Cell-free DNA ScreenTape		7 ScreenTape devices
5067-5631	Cell-free DNA Reagents <ul style="list-style-type: none">• Cell-free DNA Ladder• Cell-free DNA Sample Buffer	 	2 vials 50 µL 300 µL
5067-5632	Cell-free DNA Ladder		1 vial, 50 µL
5067-5633	Cell-free DNA Sample Buffer		1 vial, 300 µL

Limited Use Label License

Some products within this assay contain SYBR™ Green I, which is licensed from Life Technologies Corporation for use in research and development only. SYBR™ is a registered trademark of Life Technologies Corporation.

For Research Use Only

Not for use in Diagnostic Procedures.

Additional Material Required for Analysis with the TapeStation Systems

- Loading tips (5067-5598, 1 pk or 5067-5599, 10 pk)
- Optical Tube 8x Strip (401428) and Optical Tube Cap 8x Strip (401425)
- Vortex mixer IKA MS3 with 96-well sample plate adapter
- 96-well sample plates (5042-8502) and 96-well Plate Foil Seal (5067-5154) (4200 TapeStation systems only)

Additional Equipment Required (Not Supplied)

- Volumetric micropipettes for handling volumes from 1 to 15 µL
- Centrifuges for tube strips and 96-well sample plates

WARNING

Toxic agents

- ✓ Refer to product material safety datasheets for further information.
- ✓ When working with the ScreenTape assay follow the appropriate safety procedures such as wearing safety goggles, laboratory gloves and protective clothing.

CAUTION

Damage to the TapeStation systems

- ✓ Only use the recommended consumables and reagents with the TapeStation systems.

Cell-free DNA ScreenTape Assay for TapeStation Systems

Essential Measurement Practices

Read about good measurement practices in the TapeStation Information Center and/or in the System Manual.

Environmental conditions	<ul style="list-style-type: none">• Operating temperature: 16 – 30 °C (61 – 86 °F)• Keep reagents during sample preparation at room temperature
Steps before sample preparation	<ul style="list-style-type: none">• Allow Sample Buffer to equilibrate at room temperature for 30 min prior to use• Vortex each vial and briefly spin down• Flick ScreenTape device to eliminate bubbles in the buffer chamber
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 Sample Buffer remains within or on the outside of the tip• Care must be taken due to viscosity of the Sample Buffer
Mixing and centrifugation recommendations	<ul style="list-style-type: none">• Apply foil seal to 96-well sample plate or cap the tube strips prior to mixing and centrifugation• Centrifuge to collect liquid at the base; then vortex using the IKA MS3 vortexer and adaptor at 2000 rpm for 1 min• 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• Run samples immediately after preparation

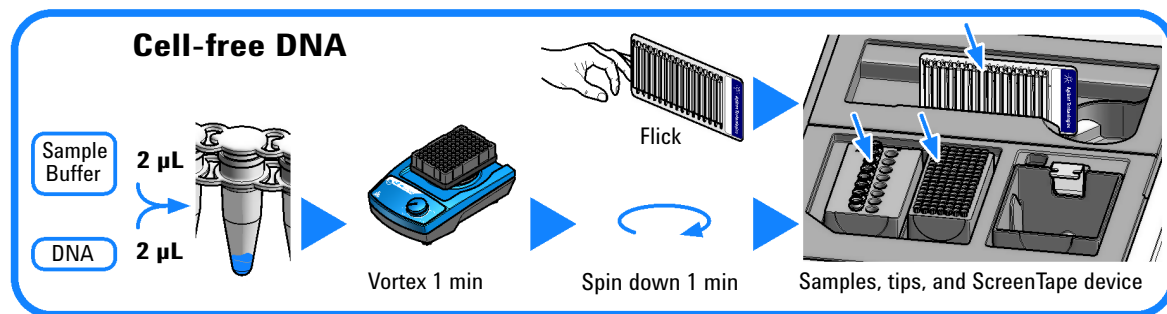
Assay Considerations

- All assay specifications are based on average values.
- Ladder is exclusively loaded from location A1 on the tube strip holder.
- Always use a complete tube strip when running ladder or samples from the tube strip holder.
- For best sizing and molarity quantification results, a ladder per ScreenTape device is recommended. Alternatively, an electronic ladder is available, which can be selected in the Agilent TapeStation Controller software.

Cell-free DNA ScreenTape Assay for TapeStation Systems

Cell-free DNA ScreenTape Assay Operating Procedure

- 1 Allow Cell-free DNA Reagents to equilibrate at room temperature for 30 minutes.
- 2 Launch the Agilent TapeStation Controller software.
- 3 Flick the Cell-free DNA ScreenTape device and insert it into the ScreenTape nest of the TapeStation instrument.
- 4 Select required sample positions in the TapeStation Controller software.
- 5 The required consumables (tips, further ScreenTape devices) are displayed in the TapeStation Controller software.
- 6 Vortex reagents and samples. Spin down before use.
- 7 Prepare ladder:
 - For 1 ScreenTape device: pipette 2 μ L Cell-free DNA Sample Buffer (●) and 2 μ L Cell-free DNA Ladder (●) at position A1 in a tube strip.
 - For 2 ScreenTape devices: pipette 4 μ L Cell-free DNA Sample Buffer (●) and 4 μ L Cell-free DNA Ladder (●) at position A1 in a tube strip.
 - For more than 2 ScreenTape devices¹: pipette 15 μ L Cell-free DNA Sample Buffer (●) and 15 μ L Cell-free DNA Ladder (●) at position A1 in a tube strip.
- 8 For each sample, pipette 2 μ L Cell-free DNA Sample Buffer (●) and 2 μ L DNA sample in a tube strip or 96-well sample plate¹.
- 9 Apply caps to tube strips and/or foil seals to 96-well sample plates.
- 10 Mix liquids using the IKA MS3 vortexer at 2000 rpm for 1 min.
- 11 Spin down samples and ladder for 1 min.



Sample Analysis

- 1 Load samples into the TapeStation instrument. Place ladder in position A1 on tube strip holder.
- 2 Carefully remove caps of tube strips. Visually confirm that liquid is positioned at the bottom.
- 3 Click **Start**.
- 4 The TapeStation Analysis software opens automatically after the run and displays results.

Technical Support and Further Information

For technical support, please visit www.agilent.com/chem/contactus. Visit Agilent Technologies' web site. It offers useful information, support and current developments about the products and technology: www.agilent.com/genomics/tapestation.

¹ Agilent 4200 TapeStation system only

www.agilent.com

© Agilent Technologies Inc. 2018-2025
Printed in Germany, Edition: 04/2025



Part No: G2991-90060 Rev. D
Document No: D0000319 Rev. D

