

Bioanalyzer System

Troubleshooting Section



Where and what to look for when analyzing a run file

- ➢ Gel like image overview
- Did the ladder run correctly?
- Are the markers present?
- Is this the profile I expected to see?





Data Analysis: Alignment

• For correct analysis: internal markers need to be identified correctly





Data Analysis: DNA Ladder

Check if all DNA ladder fragments have been identified correctly:



The standard curve is generated with a point to point fit using the migration time and sizes of the ladder peaks

The size of each sample peak is calculated from the standard curve



Data Analysis: Samples

- Specified DNA fragments are used as Lower and Upper Markers
- The 2100 Expert SW automatically detects the Markers
- Lower and Upper Marker are used for the alignment -> Sizing
- Upper Marker is used for the quantitation





DNA Chip Troubleshooting







Why is it not possible to toggle the x-axis between time and size?

Hints:

- Look for Markers present in every well
- Check sample concentrations loaded onto the chip

Solutions

- Manually set Marker(s)
- Get peak recognized as a peak, use Manual Integration
- If one sample is missing its Upper Marker, go to File > Save Selected Sample > choose ladder and all samples except the affected sample. This will salvage the remaining samples.





Upper Marker not automatically recognized

- Some samples show no sizing/quantitation information, red flags -

Issue:

Late migration - shifts the Upper Marker outside the expected time window

Solution:

Manually assign Upper Marker on Peak Table tab





Overloaded chip

- No sizing/quantitation information, red flags, error message: "optical signal too high" -

Issue:

Overloaded chip

Solution:

Adjust sample concentration by dilution





No detectable peaks, distorted baseline

Hints:

- Check for liquid spillage on chip surface
- Check chip priming station settings
- How much water was added to the cleaning chip
- When was maintanance for chip priming statior last performed
- When was maintenance for pin set last performed

Issues for this particular case:

- Chip priming station channel blocked
- Incorrect pressure was applied pressure not constant over time

Solution:

Prepare a new chip and re-run the samples





Low signal intensity (FU) with the HS DNA Ladder

HS DNA Ladder peaks are barely detectable

Solution:

Vortex ladder vial for 5 s before use





Broad dips in electropherogram (usually seen with HS DNA)

Broad dips in all/some samples and/or ladder

Explanation:

Residual RNaseZap or SDS on the electrodes

Solution:

Additional washes with cleaning chip with H₂O (2x)





RNA Chip Troubleshooting









Data Analysis: RNA Ladders

RNA Ladder is used for sizing and quantitation:

6 ladder peaks + Lower Marker

Ladder concentration:

RNA Nano: 150 ng/µl RNA Pico: 1000 pg/µl

Small RNA ladder: 1000pg/µl





Use of Lower Marker for Alignment





For Research Use Only. Not for use in diagnostic procedures.
Agilent Technologies
Rev. 1 [January 13, 2016]

Alignment

- A DNA Fragment (25 nt) is used as Lower Marker in RNA Nano/Pico Assay
- The 2100 Expert SW automatically detects the Lower Marker
- Lower Marker is used for the alignment





Why is RIN N/A?

Hints:

- Is the correct assay selected?
- Look for the red-flagged lanes on the gel-like image





Why is RIN N/A?

Solutions:

- Run new chip and select correct assay
- Manually assign peaks
- Clear critical error message(s) to get RIN displayed





2100 Expert Software

Setpoint Explorer

Troubleshooting example: RIN N/A – adjust threshold(s) in the Setpoint Explorer to clear critical error(s) \rightarrow RIN displayed



Relax 5S Region Anomaly Threshold to 1 (most relaxed) will clear the critical error



2100 Expert Software

Setpoint Explorer

Troubleshooting example: RIN N/A – adjust threshold(s) in the Setpoint Explorer to clear critical error(s) => RIN displayed





For Research Use Only. Not for use in diagnostic procedures. Rev. 1 [January 13, 2016]

Troubleshooting: additional ladder fragments

Solutions:

- Heat denature samples and ladder at 72°C for 2 min and immediately place on ice
- Aliquot the ladder after heat denaturation





Troubleshooting: additional RNA sample peak

Solution:

- DNAse treatment
- Heat denature samples at 72°C for 2 min to resolve secondary structures





Troubleshooting: low signal intensity

Solutions:

- Prepare a fresh geldye mix; protect dye and gel-dye mix from light
- Check kit (both chips and reagents) is within expiration
- Run full set of hardware diagnostic tests





Troubleshooting: degraded RNA Ladder

Improper handling, preparation, and/or storage

RNA Ladders are available separately



For Research Use Only. Not for use in diagnostic procedures.
Agilent Technologies
Rev. 1 [January 13, 2016]



Protein Chip Troubleshooting







Data Analysis: Protein Ladder

Check if all Protein Ladder fragments have been identified correctly:



The standard curve is generated with a point to point fit using the migration times and sizes of the ladder peaks

The size of each sample peak is calculated from the standard curve



Data Analysis: Protein Samples

- Lower and Upper Markers are included in the Sample Buffer
- The concentration of the Upper Marker is known [60 ng/ μ l]
- The ratio between sample and Sample Buffer as well as sample preparation are fixed in the protocol
- Comparing TIME-CORRECTED PEAK areas between Upper Marker and sample peak returns a relative sample peak concentration





Absolute Quantification Using the Calibration Curve

- Use internal standards (user defines number of wells to be used as standard)
- To select samples for generation of calibration curve, check "Use for Calibration" box next to these samples



For Research Use Only. Not for use in diagnostic procedures. Rev. 1 [January 13, 2016]

Troubleshooting: clogged priming station

Issue: Plastic adapter of priming station clogged

Symptom: Broad and fuzzy peaks

Solutions: Routine maintenance of priming station; avoid spillage of gel



Recommended after chip priming step:

- Quick view if residual gel is visible on the adapter/gasket;
- If yes: use Kimwipe to clean adapter, flush out adapter with syringe and water. If dried gel in hole, pick out with a needle.

29

Troubleshooting: peak not detected

Issue: A sample peak is not detected by the software automatically

Example:



Solutions:

- Use Setpoint Explorer to modify integration limits e.g. height threshold
- Activate Manual Integration and choose 'Add peak'



Want to know more?

- The 2100 Expert software has many features and a very extensive Help menu
- For the latest version of the Maintenance & Troubleshooting Guide: <u>http://www.chem.agilent.com/Library/usermanuals/Public/G2946-90003.pdf</u>





Need help?



Are you stuck? Do you need our help? Please see the following slides for details on the support process.



Need to get in contact with us?



- Email contact: <u>bioanalyzer@agilent.com</u>
- Phone: 1-800-227-9770 options 3x4x1



Need to get in contact with us?



Please provide the following information to help us with troubleshooting:

- Error description, in your words. Also, the expected result in your words.
- Kit used lot #s and expiration dates of chips and reagents used.
- Sample information:
 - DNA/RNA or Protein?
 - Which species, which tissue organ are the samples extracted from?
 - Extraction method, post-extraction processing (e.g. DNase digest), purification method
 - What kind of buffer are the samples in?
 - What is the expected concentration range (and sizing range, if applicable) of the samples?
- The data file (**.xad**) from the run that had problems. Also, any data file from a recent successful run.

Data file: e.g. 2100expert_EukaryoteTotalRNAPico_12345_2013-05-17_13-24-56.xad

Location: C:\Program Files (x86)\Agilent\2100 bioanalyzer\2100 expert\data\



Questions?





For Research Use Only. Not for use in diagnostic procedures.
Agilent Technologies
Rev. 1 [January 13, 2016]