

DV₂₀₀ RNA quality metric for FFPE analysis using LabChip GX Touch nucleic acid analyzer.

The process of fixing the tissue sample and embedding it in paraffin can cause severe degradation of the RNA. The fixation process and storage cause significant RNA degradation in some samples. The formalin fixation process causes cross linkage between nucleic acids and proteins, and the covalent modification of RNA by mono-methylol (-CH₂OH) addition to the bases¹. Thus, it is very important to assess RNA quality after FFPE (Formalin-Fixed, Paraffin-Embedded) extraction of RNA for downstream applications.

One measure of RNA quality is the RNA Quality Score, or 'RQS'. The RQS is a metric calculated by the LabChip™ RNA Assay and has a high degree of correlation to the Agilent RNA Integrity Number, or 'RIN'. While these are the most prominent available RNA metrics, they may not apply equally to all types of RNA applications.

The DV₂₀₀ metric has been promoted by Illumina®. As noted in the Illumina tech note² "Many researchers use the Agilent RNA Integrity Number (RIN) to determine RNA quality for gene expression analysis. However, we have found that RIN values from degraded FFPE samples are not a sensitive measure of RNA quality nor are they a reliable predictor of successful library preparation".

To improve the RNA integrity assessment, Illumina developed the DV₂₀₀ metric - the percentage of RNA fragments > 200 nucleotides to assess FFPE RNA quality. Following is the calculation of DV₂₀₀ on the LabChip™ GX Touch™ nucleic acid analyzer.

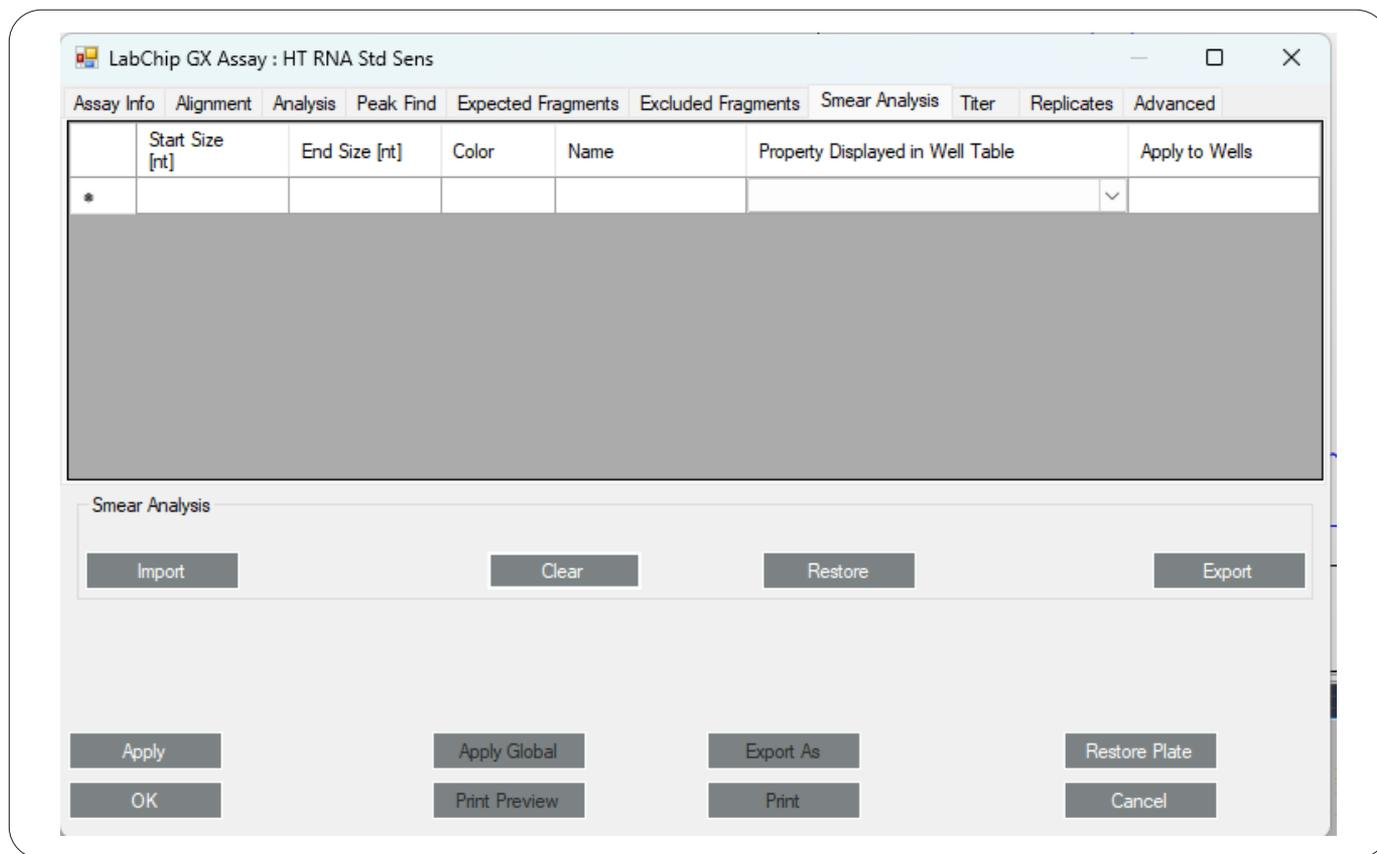


Evaluation of FFPE Samples for RNA Quality on LabChip GX Touch

Here we show how to obtain the DV₂₀₀ value by analyzing RNA FFPE samples on the LabChip GX Touch system. Prior to the DV₂₀₀ metric, the LabChip GX Touch software calculated a RQS (RNA Quality Score) value. It has been well demonstrated that there is a high correlation between the RQS value and Agilent RIN value for a large number of sample types, regardless of denaturation method used³.

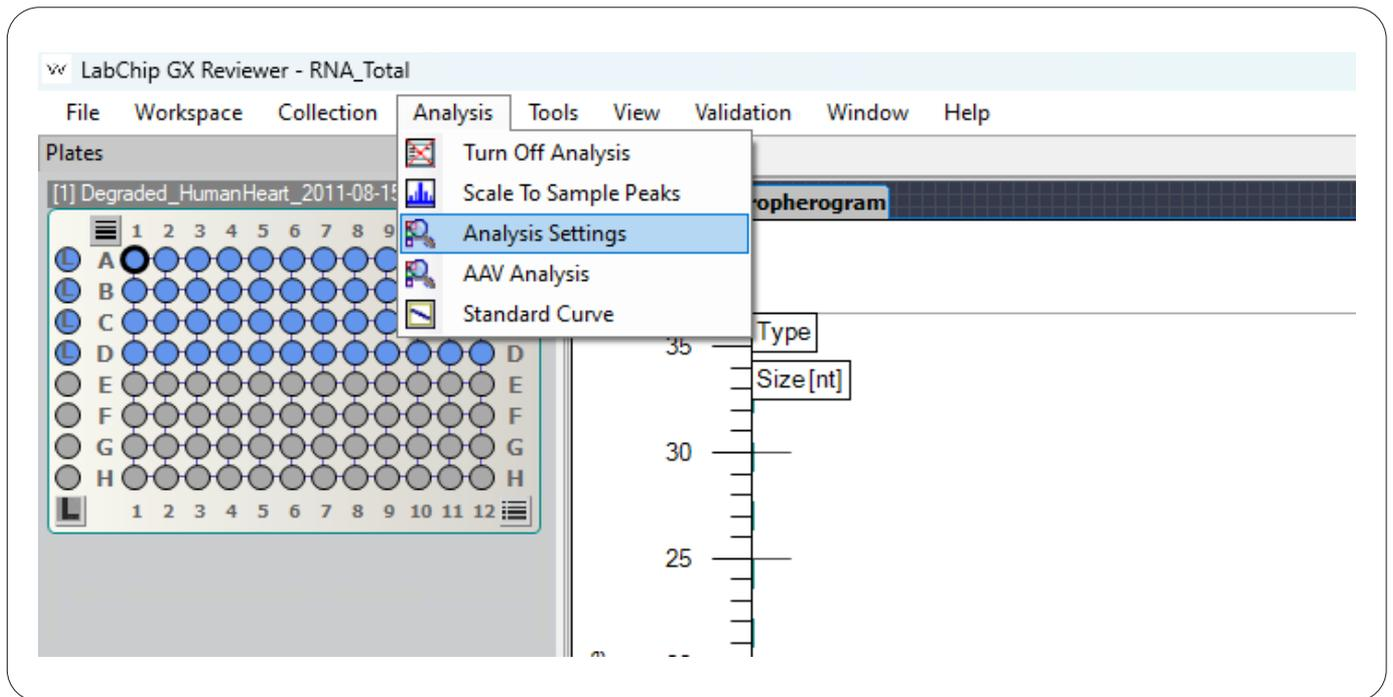
In addition to the RQS value, the LabChip Reviewer software can now calculate the DV₂₀₀ value. Both values can be exported and utilized to determine the RNA quality of FFPE samples.

1. Run samples on the LabChip GX Touch instrument. For further information on sample preparation and running the samples, refer to RNA Assay User Guide(s).
2. Open data files in the LabChip GX Reviewer Software.
3. Select 'Analysis Settings' under 'Analysis' tab.



4. Select 'Smear Analysis' tab.
5. Define region from 200 to 10000 nt. Enter the size were region starts under Start Size. Add the size were region ends under End Size
6. Under Name column, type in DV₂₀₀ or give any other name to the value being reported.
7. Select '% of Total Area' from the drop-down menu under 'Property Displayed' in Well Table.
8. Analysis can be applied to all or select samples under 'Apply to Wells' tab.
9. Click OK.

10. If multiple plates have been analyzed click 'Apply Global'.



11. To highlight the selected region in the electropherograms and gel image check 'Show Smears' in the properties.

