

## Automated Electrophoresis for Protein Characterization



Figure 1. LabChip GXII Touch

The LabChip® GXII Touch System offers researchers an automated alternative to traditional methods by streamlining the multiple, manual steps of slab gel electrophoresis, while also providing the throughput and data quality essential in the biopharmaceuticals workflow (Figure 1).

The platform supports multiple assays for characterizing proteins in reduced and non-reduced samples including:

- Purity
- Titre
- Glycan screening
- mAb charge heterogeneity
- Fragmentation

Analysis can be performed in as few as forty seconds per sample that deliver comparable data to traditional capillary electrophoresis with as much as a 70X increase in throughput. Choose 96-well or up to 384-well platforms depending upon throughput needs. With an easy to use touch screen interface, even occasional users get up and running samples quickly. TIBCO SpotFire® data visualization further enhances data output.

PerkinElmer offers solutions to ensure the consistent, reproducible results your biopharmaceuticals research and patient safety considerations demand. The LabChip GxP Security Software provides users with electronic signature, electronic record retention, and access management to streamline meeting requirements for compliance.

## LabChip GXII Touch Screen Simplifies Sample Analysis

### Touch – User friendly operation

- Load sample plate and chip
- Select samples (up to 384 in a run)
- Select assay type
- Touch 'Run' to start
- You can even have the system automatically export data directly to your network or LIMS system

### Run – Observe runs in real time

- Sample analysis in as few as 40 seconds
- View electropherogram in real time during data collection
- Overlay collected data to compare sample profiles within software
- Select from various run time analytical feature annotations

### Review –See data in real time or export for later analysis

- Choose display in E-gram, virtual gel or data table format (Figure 2)
- Enhance results with TIBCO SpotFire® data visualization
- Pull multiple archived plates into data review or analytical comparisons
- Apply data mining filter functions on key attributes
- Highlight expected peaks
- Track relevant user access and data history parameters with 21 CFR Part 11 compliance software

The LabChip GXII Touch operator controls are designed to allow users to easily setup and execute a run with as few as three easy steps. Run templates can include such things as well selections, sample names, expected peak tables, and more, which facilitate operator ease of use. Data can be automatically exported to network or LIMS directories for subsequent analysis. Every instrument also comes with a full software package for data review, allowing analysis from current or archived data sets.

LabChip GX Touch and Reviewer Software contain built-in technical controls and features specifically designed to support the 21 CFR Part 11 compliance. These features include a shared user account database, access controls, device check, enforced sequencing of run steps, audit trails, record copying, record retention, system documentation, and electronic signature controls (Figure 3).

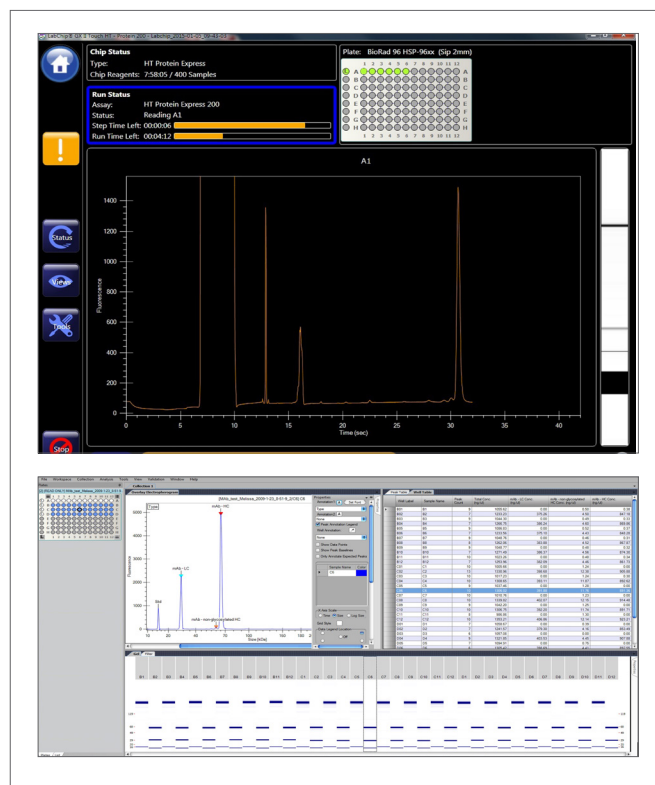


Figure 2. The Touch interface offers simple instructions for project initiation with real-time run sample evaluation (top) and complete sample analysis (bottom).

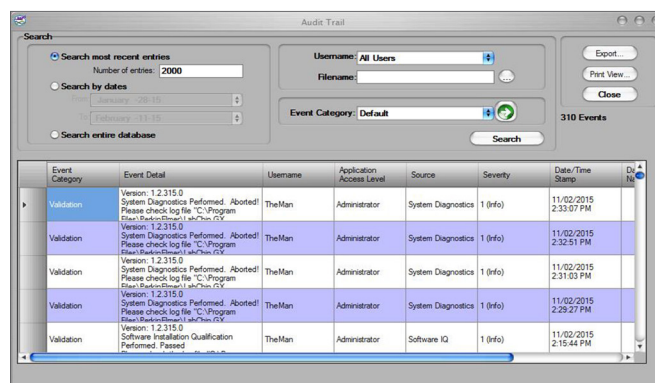


Figure 3. 21 CFR Part 11 Software, part of a suite of digital governance, helps ensure process compliance.

## Rapid Analysis Throughout the Protein Workflow



The LabChip GXII Touch offers rapid quantification and quality control throughout the biotherapeutics workflow. For example, automating the characterization process allows multiple, critical quality attributes to be obtained significantly faster. Researchers can now screen for optimal protein characteristics earlier in the process, and integrate Quality by Design initiatives into their biotherapeutics development workflow (Figure 4).

### Small-Scale High Throughput Process Development

#### Process Parameters

#### Purification & Sample Prep

JANUS G3 BioTx  
Pro Plus Workstation



#### Determination of Critical Quality Attributes

LabChip GXII  
Touch



- ✓ Purity & Fragmentation
- ✓ N-Glycan profiling
- ✓ Change Variant

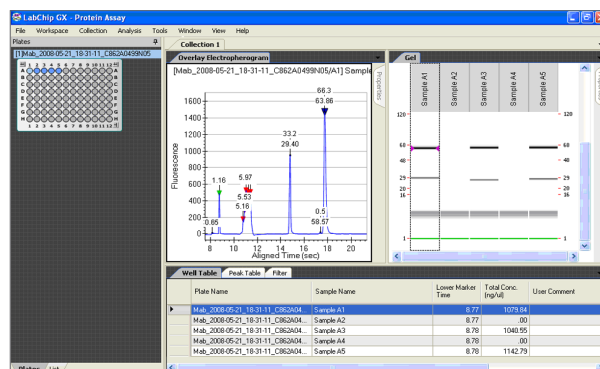


LabChip DS

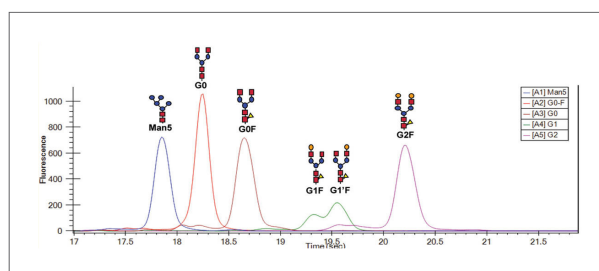
- ✓ Recovery & Yield

#### Informatics

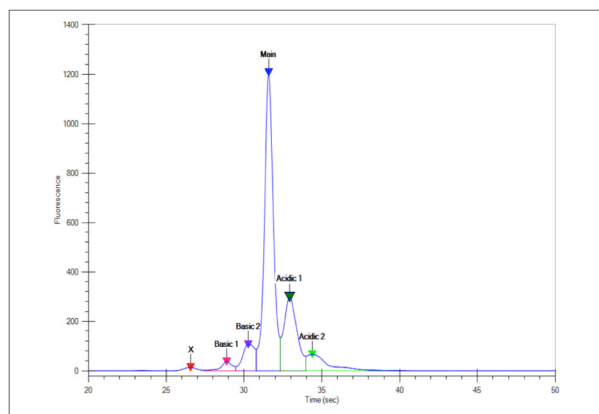
#### Process Outcomes



Standard and Pico Sensitivity Protein Assay



Glycan Profiling



Charge Heterogeneity Assay

Figure 4. With LabChip microfluidics technology and the JANUS® G3 BioTx Workstation, you can purify and analyze in one day what typically takes weeks with alternative methods. Explore a broader range of experimental conditions and save in both the development time and cost of your biotherapeutics research (above). LabChip Touch reagents are available for multiple protein attribute analysis including standard and pico protein, glycan profiling and charge heterogeneity (right).

## LabChip Electrophoresis

### How Does it Work?

LabChip electrophoresis is performed on a small, microfluidic chip. Prior to analysis, reagents are loaded into the individual wells of the chip. These wells are connected to tiny microchannels about the size of a human hair etched within the quartz microfluidic chip. (Figure 5).

When the chip is loaded into the LabChip GXII Touch system, the chip's wells interface with platinum electrodes that provide voltage and current control. The system robot moves the microtiter plate wells directly under the chip's capillary 'sipper', and approximately 150 nL of sample is aspirated onto the chip. Sample staining and destaining are performed automatically on the instrument platform.

Individual sample analytes are separated electrophoretically and the bands are detected via laser induced fluorescence. Sizing and concentration for each band are determined using ladder and internal markers. Because the sipper is rinsed between samples, cross-contamination or carryover is eliminated.

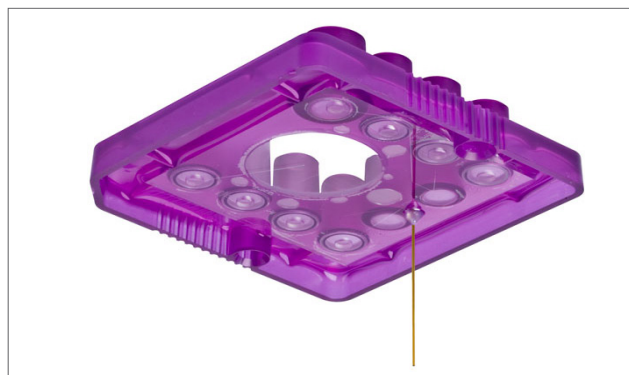


Figure 5. LabChip for protein research

### Ordering Information

|   | LabChip GXII | LabChip GXII Touch HT | LabChip GXII Touch 24 |
|---|--------------|-----------------------|-----------------------|
| <b>Protein Express LabChip</b>                  | 760499       | 760499                | CLS138950             |
| <b>Protein Express LabChip (Economy 4 pack)</b> | 760528       | 760528                | N/A                   |
| Protein Express Reagent Kit                     | CLS960008    | CLS960008             | CLS960008             |
| Bulk Protein Express Sample Buffer              | 760518       | 760518                | N/A                   |
| Pico Protein Reagent Kit                        | 760498       | 760498                | 760498                |
| Bulk Pico Protein Sample Buffer                 | 760414       | 760414                | N/A                   |
| <b>High Resolution Protein LabChip</b>          | 760524       | 760524                | CLS138951             |
| Low Molecular Weight Reagent Kit                | 760573       | 760573                | 760573                |
| Glycan Screening Reagent Kit                    | 760525       | 760525                | 760525                |
| Glycan Release & Labeling Kit                   | 760523       | 760523                | 760523                |
| <b>DNA 5K/RNA/CZE LabChip</b>                   | 760435       | 760435                | CLS138949             |
| Charge Variant Reagent Kit                      | CLS760670    | CLS760670             | CLS760670             |

| LabChip GXII Touch Specifications |                  |                            |                 |
|-----------------------------------|------------------|----------------------------|-----------------|
| <b>Height</b>                     | 25.75 in         | <b>Power Requirements</b>  | 100-240 Vac     |
| <b>Width</b>                      | 19.25 in         | <b>Power Consumption</b>   | N/A             |
| <b>Depth</b>                      | 18.25 in         | <b>Plate Formats</b>       | 96- or 384-well |
| <b>Weight</b>                     | 54 lbs (24.5 kg) | <b>Excitation/Emission</b> | 635 and 700 nm  |
| <b>Temperature Range</b>          | 18-26 deg C      | <b>Humidity Range</b>      | 20% - 80% RH    |

PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2014-2016, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.