



LASER  
PRECISION



For research use only. Not for use in diagnostic procedures.

**ProteinEXact™ Assay**  
For the LabChip® GXII Touch™ Protein Characterization System

  
**PerkinElmer®**  
*For the Better*

# LASER PRECISION: PROTEINEXACT™ ASSAY

Bring high sensitivity to your protein quantitation and sizing applications with the ProteinExact™ assay for the LabChip® GXII Touch™ protein characterization system. The ProteinExact™ assay is ideal for use in both upstream and downstream biotherapeutics workflows which require the utmost in precision. With the analytical power and speed of the LabChip® GXII Touch™ protein characterization system and IntelliChip™ assay optimization technology, you can accelerate your protein development workflows with high-precision reproducibility, sensitivity, and an expanded sizing range.

- Dependable reproducibility for quantitation (<10% CV) and sizing (<2% CV)
- Broad sizing and quantitation range: 6.5-250 kDa and 10-2000 ng/μL
- Highly sensitive: 0.2 ng/μL
- Flexible: Run in batches of 48 or 96 samples at a time
- Fast: Analysis in ~65 sec per sample

The LabChip® GXII Touch™ automated capillary electrophoresis platform delivers comparable results to traditional capillary electrophoresis with 46X faster run times without compromising sensitivity, resolution or reproducibility.

## PROTEINEXACT™ REPRODUCIBILITY, SENSITIVITY, & SIZING RANGE

### Reproducibility

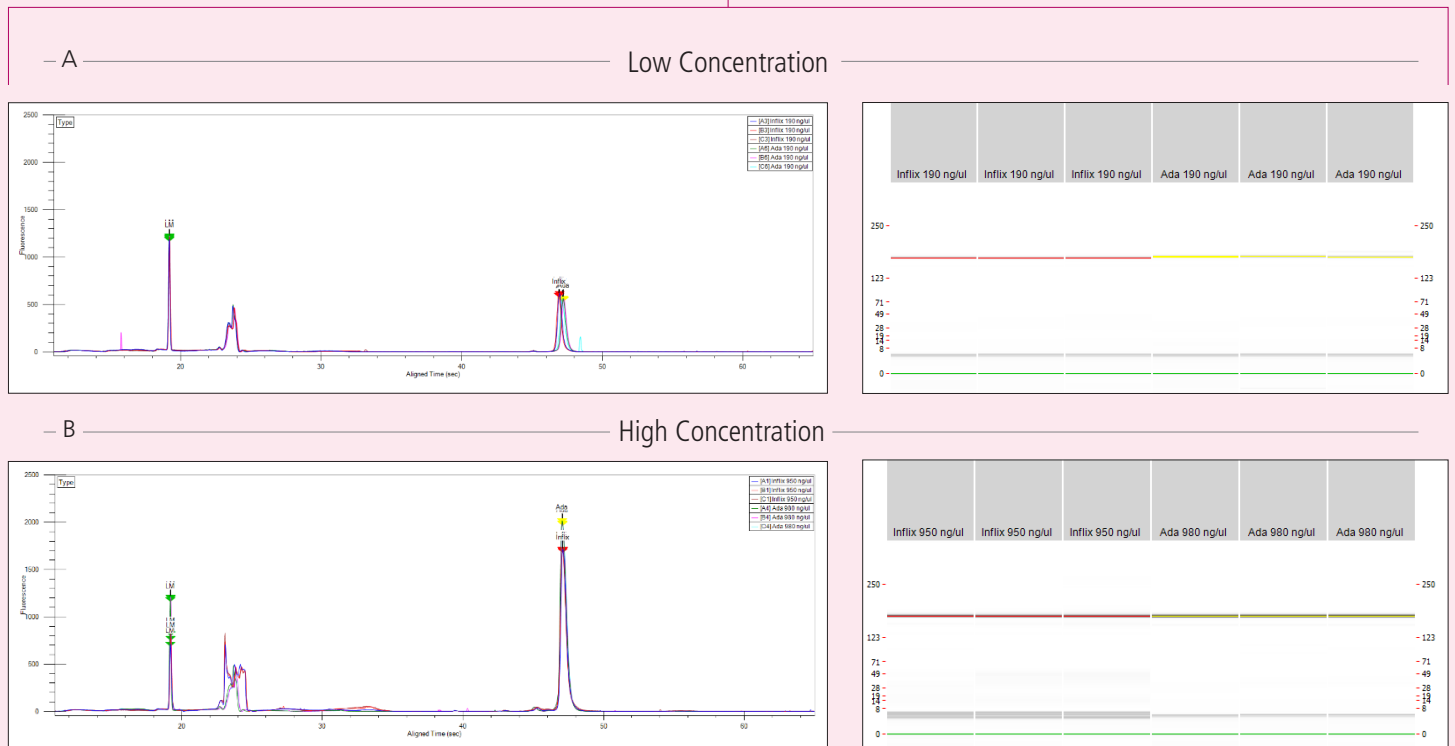
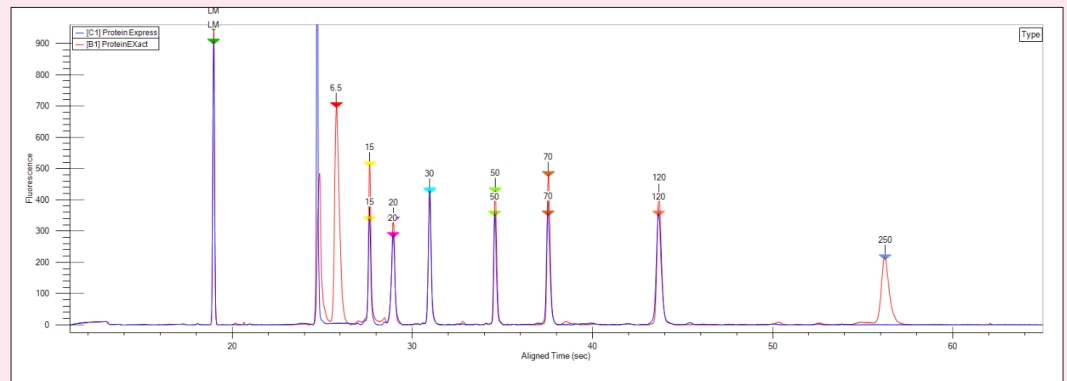
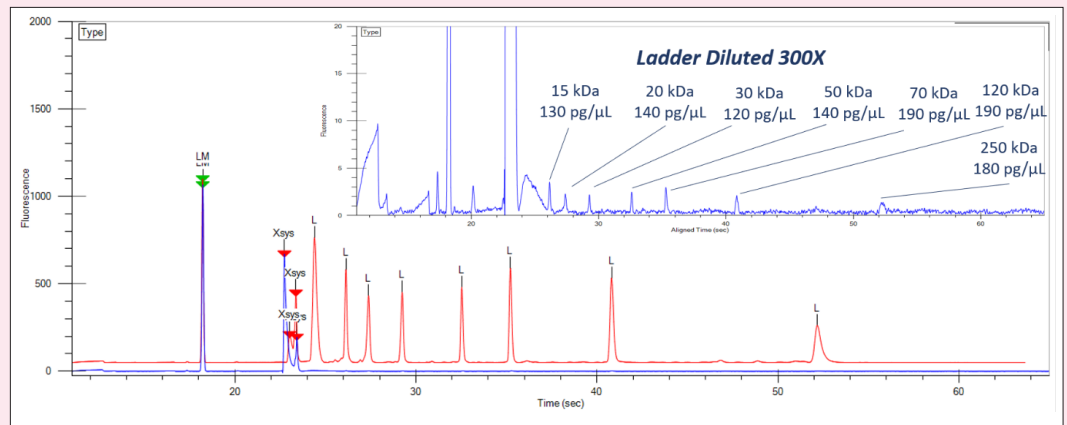
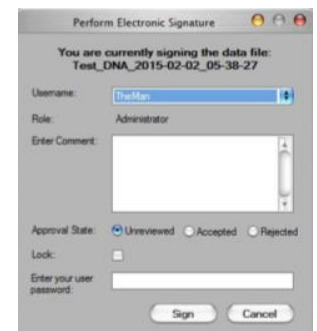


Figure 1. Electropherogram overlays and corresponding virtual gels for samples run 3 times at low (a) and high (b) concentrations.

## Expanded Sizing Range



The LabChip® GXII Touch™ System and Reviewer Software contains built-in technical controls and features specifically designed to be compatible with 21 CFR Part 11 requirements. 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. LabChip® Touch™ instruments generate data records in electronic form which are archived in a Central Data Repository. The LabChip® GX (GxP) Reviewer application can be run from any computer connected to the instrument network and allows modification and maintenance of the records with the ability to perform electronic signatures on the records generated from instruments in the lab system. The robust and comprehensive LabChip® GXP Software uses secured, computer-generated, time-stamped audit trails to independently record the date and time of operator entries and actions that create, modify, or delete electronic records (Figure 4). In addition, unique combinations of user ID and password for electronic signature are used. Once locked, a record cannot be further modified without a separate signed unlock action (Figure 5).



## ProteinEXact™ Assay Specifications

Sizing Range	6.5 kDa - 250 kDa
Sizing Precision RSD (CV)	<2%
Sizing Resolution	≤10%
Linear Dynamic Range	10-2000 ng/μL
Maximum Total Protein Concentration	2 mg/mL
Sensitivity Limit of Detection (LOD)	0.2 ng/μL
Quantitation Reproducibility	<10%
Maximum Salt Concentration	1M NaCl at pH 6.5 to 8.5
Chip Primes per Reagent Kit	10
Chip Sample Lifetime	400
Sample Analysis Time	65 sec

**System Requirements:** Running of the ProteinEXact™ assay requires a LabChip® GXII Touch™ protein characterization system with V1.7.819.0 software or newer and Data Reviewer V5.5.2312.0 or newer

## Ordering Information

Assay	LabChip® GXII Touch™ HT Chips	LabChip® GXII Touch™ 24 Chips	Reagent Kit
ProteinEXact™ Assay	CLS150337	CLS150338	CLS150466

Discover more at [perkinelmer-appliedgenomics.com/ProteinEXact](http://perkinelmer-appliedgenomics.com/ProteinEXact)

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



For more information, visit [www.applied-genomics.com](http://www.applied-genomics.com)

Copyright ©2018 PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.