

HUMAN HEALTH

ENVIRONMENTAL HEALTH

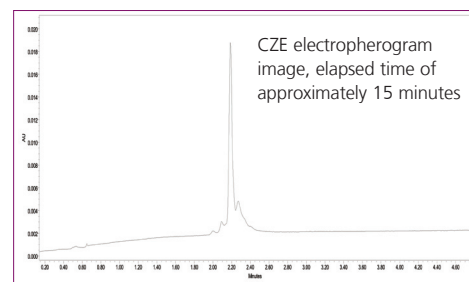
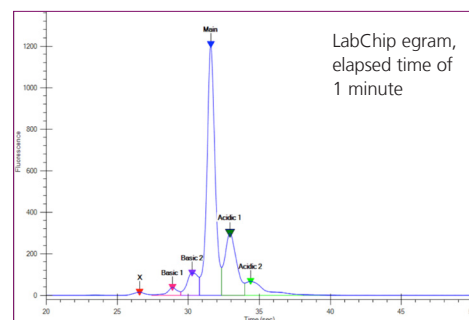
# MICROFLUIDICS CHARGE VARIANT ASSAY 68 SECONDS ANALYSIS TO RESULTS

## LabChip® GX II Charge Variant Assay: Fastest results, from clone selection to formulation

PerkinElmer LabChip®  
GX II technology  
delivers identification

of Basic and Acidic variants relative to the Main Peak of recombinant monoclonal antibodies (mAb ) with 15x faster results than conventional IEX and can calculate relative percentages of variants automatically. Greater throughput means you can optimize valuable research time to accelerate your drug to the clinic.

- Analysis run times selectable at either 68, 90, or 110 seconds
- Sample and chip preparation < 30min
- Sample pI range: ~7.0 - 9.5
- Resolution: comparable to IEX and conventional CZE
- Consistent, reproducible results: CV < 5% for fixed input sample concentration
- Easy assignment of Basic and Acidic variant peaks relative to the Main mAb peak

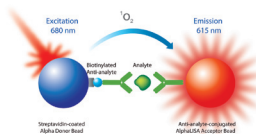


**Figure 1:** Comparison of LabChip HT Charge Variant Assay (top), and traditional CZE assay (bottom). Basic region migrates faster than Main Peak, Acidic region migrates more slowly. User names peaks using system callouts for Expected Peaks. Tabulate % relative amount of each variant.



Alpha Technology

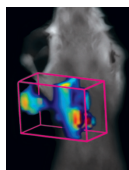
Operetta® High Content Imaging System



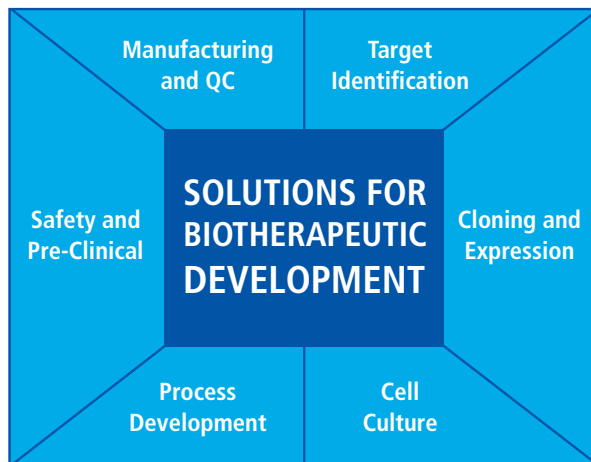
UltraView® VoX 3D Live Cell Imaging System



Fluorescent Agents, *in vivo*

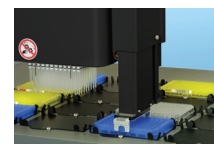


IVIS® and FMT™ *in vivo* imaging solutions

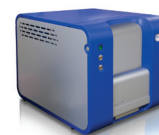


EnSpire® Multimode Plate Reader with label-free technology

Microplates



Sciclone® G3 Automated Workstation



LabChip GXII Protein Analysis and Quantitation System



JANUS Automated Workstations with EnVision® Multilabel Plate Reader

## SPECIFICATIONS

Sample Type	Monoclonal antibody (mAb)
pI Range	7.0 - 9.5
Amount of Sample Required	25 µl with concentration between 0.5 - 10 mg/mL (12.5 µg to 250 µg of mAb, total) Optimal concentration: 2 mg/mL
Resolution	Comparable to IEX and conventional CZE
Reproducibility	CV < 5% for varying concentration from 1 - 3 mg/mL CV < 3% at constant concentration
Carryover	< 0.5% of previous sample can be detected in following sample
Assay Run Time	1.8 - 3 hr for 96-well plate Three Assay durations, 68 s, 90 s, 110 s
Number of Samples per Kit	120 samples
Number of Samples per Chip Prep	96 samples
Chip Life Time	500 samples per chip

## ORDER INFORMATION

## PART NO.

LabChip 5K/RNA/CZE Chip	760435
HT Protein Charge Variant Reagent Kit	CLS760670

For more information please visit [www.perkinelmer.com/LabChip](http://www.perkinelmer.com/LabChip)

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)

  
**PerkinElmer®**  
For the Better

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

Copyright ©2012, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

010542\_01

Printed in USA