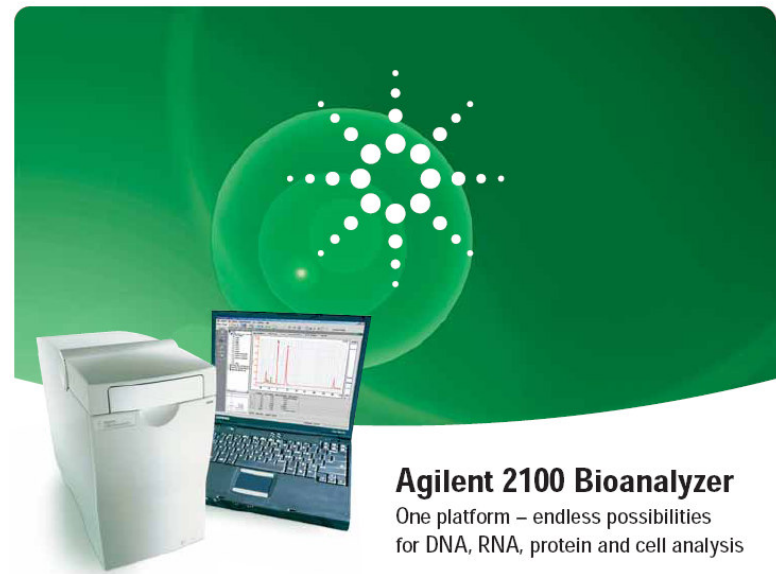


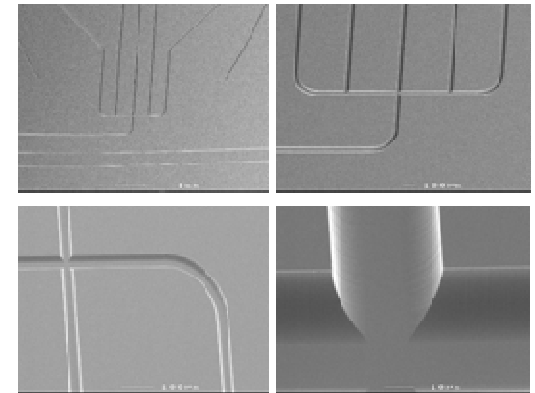
2100 Bioanalyzer

Iman Kishawi, PhD
Senior Application Scientist

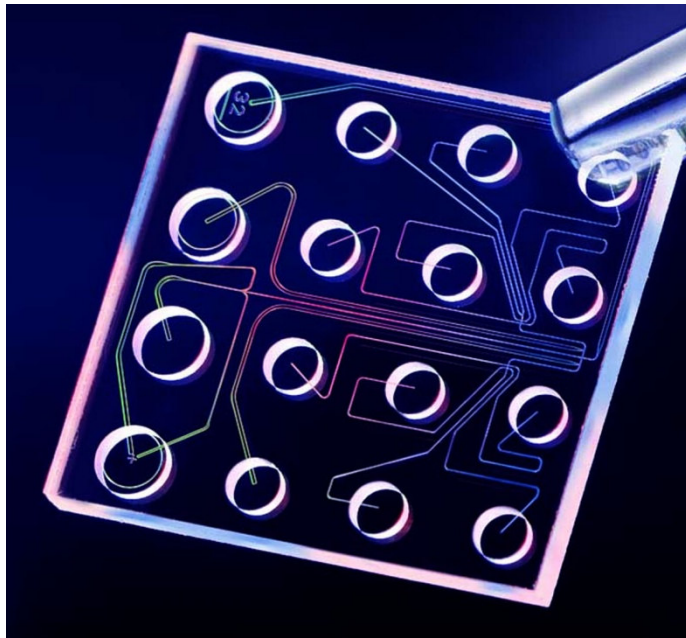


Agilent 2100 Bioanalyzer
One platform – endless possibilities
for DNA, RNA, protein and cell analysis

The Lab-on-a-Chip Approach



Increasing quality and speed of gel electrophoresis



Sample volumes 1 - 5 μ l

10 -12 samples depending on Assay

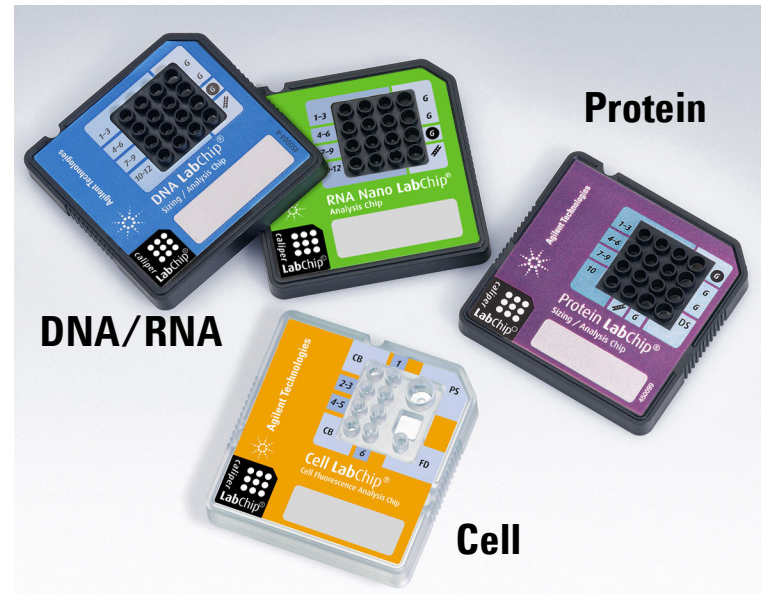
Separation, staining, detection of samples

Results in 5-30 minutes available

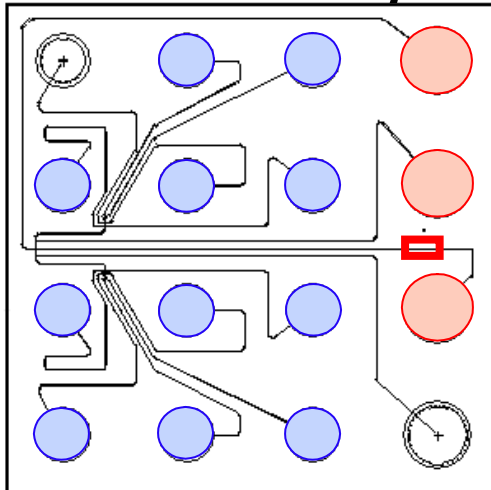
No extra waste removal needed

Disposable Chip, no crosscontamination

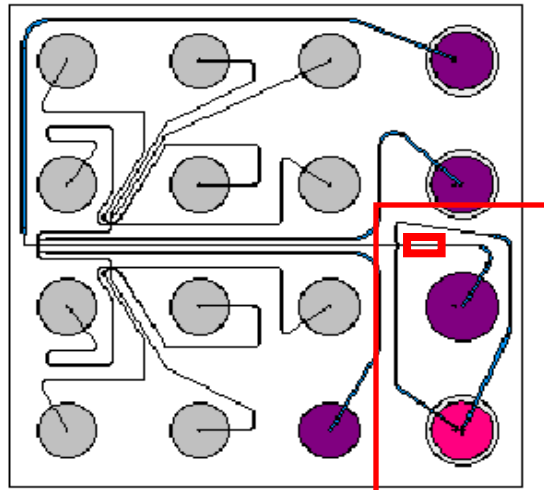
Three Chip Types



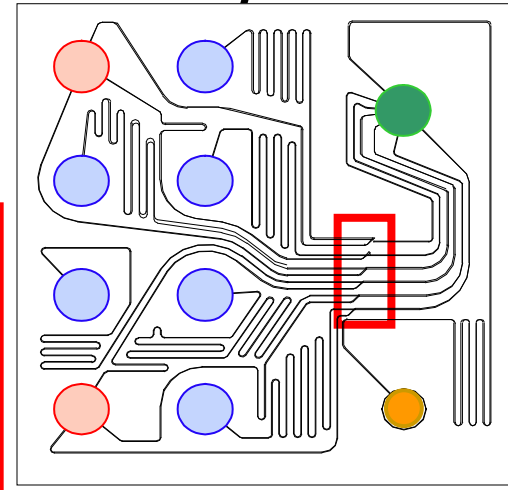
DNA/RNA analysis



Protein analysis

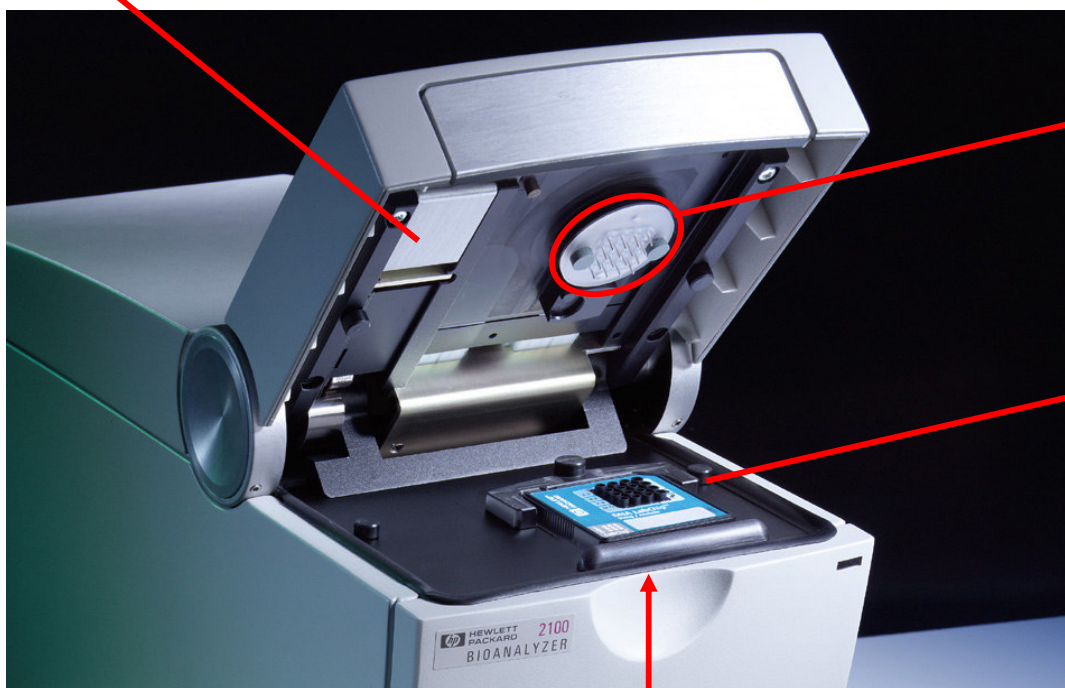
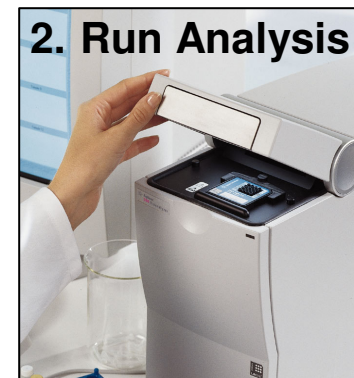


Cell analysis



Agilent 2100 Bioanalyzer

Exchangeable cartridge for different assays



**16 pin electrodes
connected to
HV-sources**

**Chip holder with
heater plate**

Optics for detection



Agilent Technologies

Current 2100 Analysis Kits

Electrophoretic Separations

DNA Assays:

1000, 7500, 12000

- Sizing
- Quantitation
- PCR products, digests, larger DNA fragments
- 12 samples in 30 min.

RNA Assays:

nano, pico, Small RNA

- Quantitation (Sizing in Small RNA)
- total RNA, mRNA
- purity & integrity determination
- 10 samples in 30 min.

Protein Assays:

P80, P230, HSP-250

- Sizing
- Quantitation
- cell lysates, column fractions, purified proteins, antibodies etc.
- 10 samples in 40 min.

Flow Cytometry

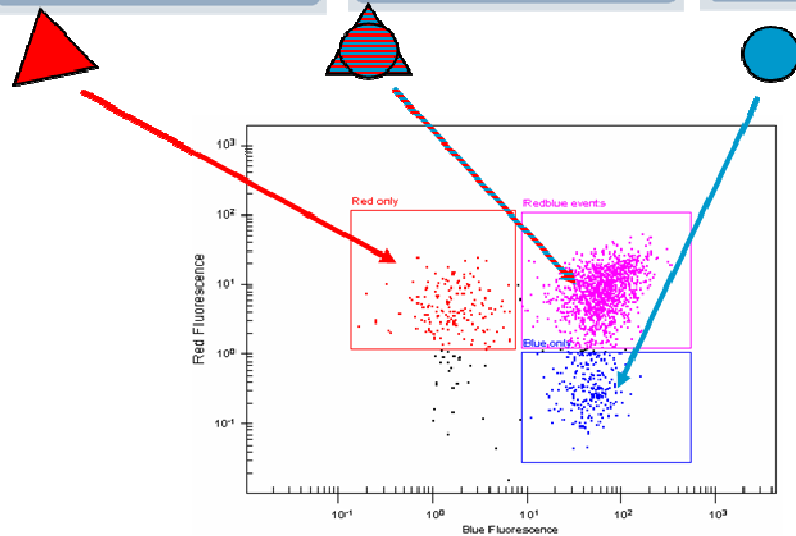
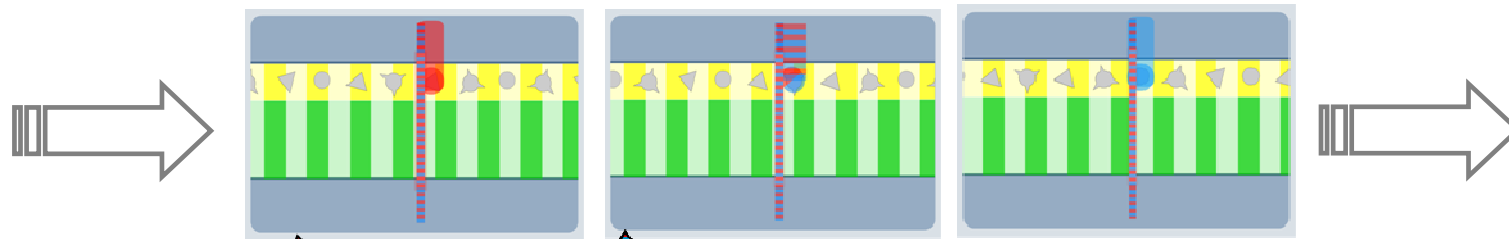
Cell Assays:

Flexible use

- Analysis of 6 samples
- Two color detection
- Analysis of protein expression in cells



Cell Applications



Dot plot view for easy data evaluation

[Video](#)

Cell Assay



Apoptosis

Transfection Efficiency Monitoring

- Detection of GFP-transfected cells
- Antibody staining: Detection of transfected cells expressing the encoded protein

Protein Expression Monitoring

- Extracellular and Intracellular Antibody staining for detection of protein expressed on the cell surface, in the cytoplasm, or in the nucleus

Gene silencing

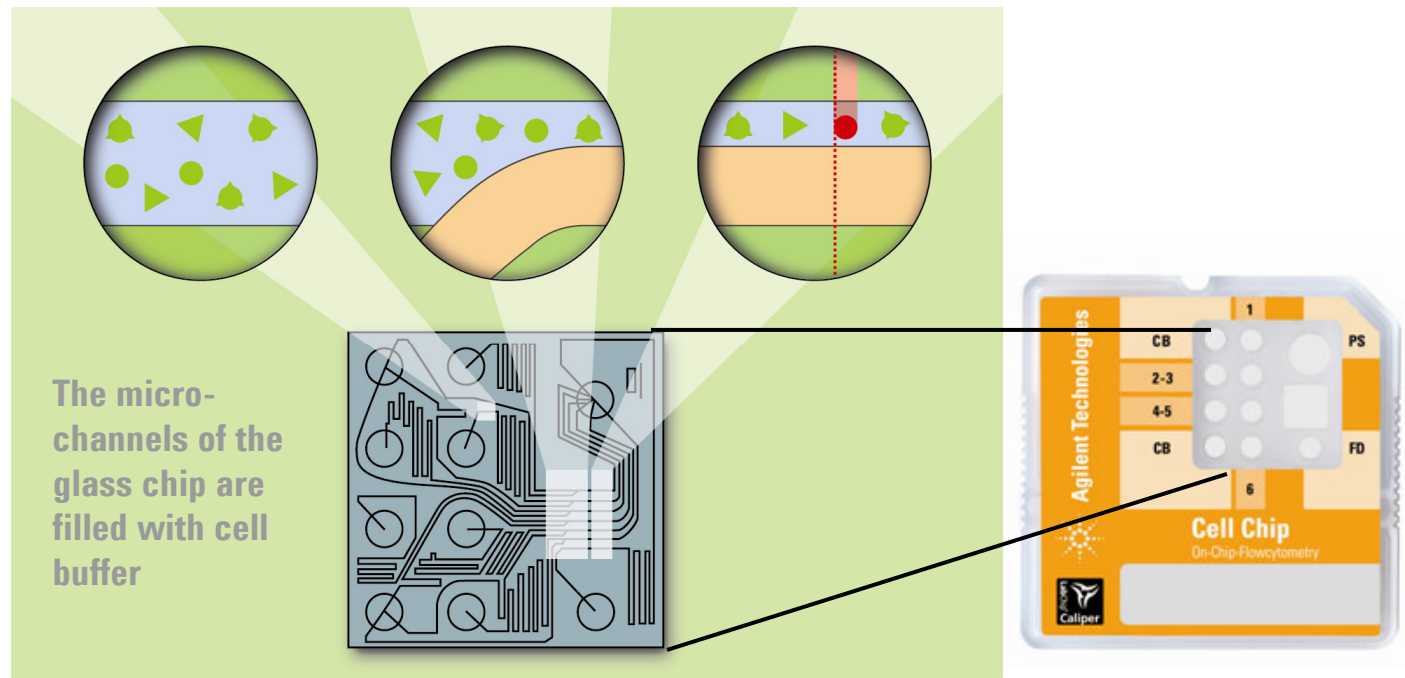
Principle of Pressure-Driven Flow

For cell assays (analysis of cell fluorescence parameters)
On-chip simple flow cytometric studies

Pressure driven flow is used to move cells in a controlled manner through the micro-channels

Cells are hydrodynamically focused to a portion of the channel by a side stream of buffer

Cells pass the fluorescence detector in single file and each event is monitored in a histogram or dot plot



The Bioanalyzer Lab-on-a-Chip Approach

Separation on disposable, μ -fabricated glass chips

- made of two glass layers:
 - one with micro-channels ($\times 10\mu\text{m}$, etched),
 - one with through-holes
- glued into a plastic caddy which accommodates wells for gel, sample, standard (ladder), buffer and other reagents
- for handling nl-amounts of liquids
- one separation channel for ladder and sample
- microfluidic sample movement with fluorescence detection

Setup

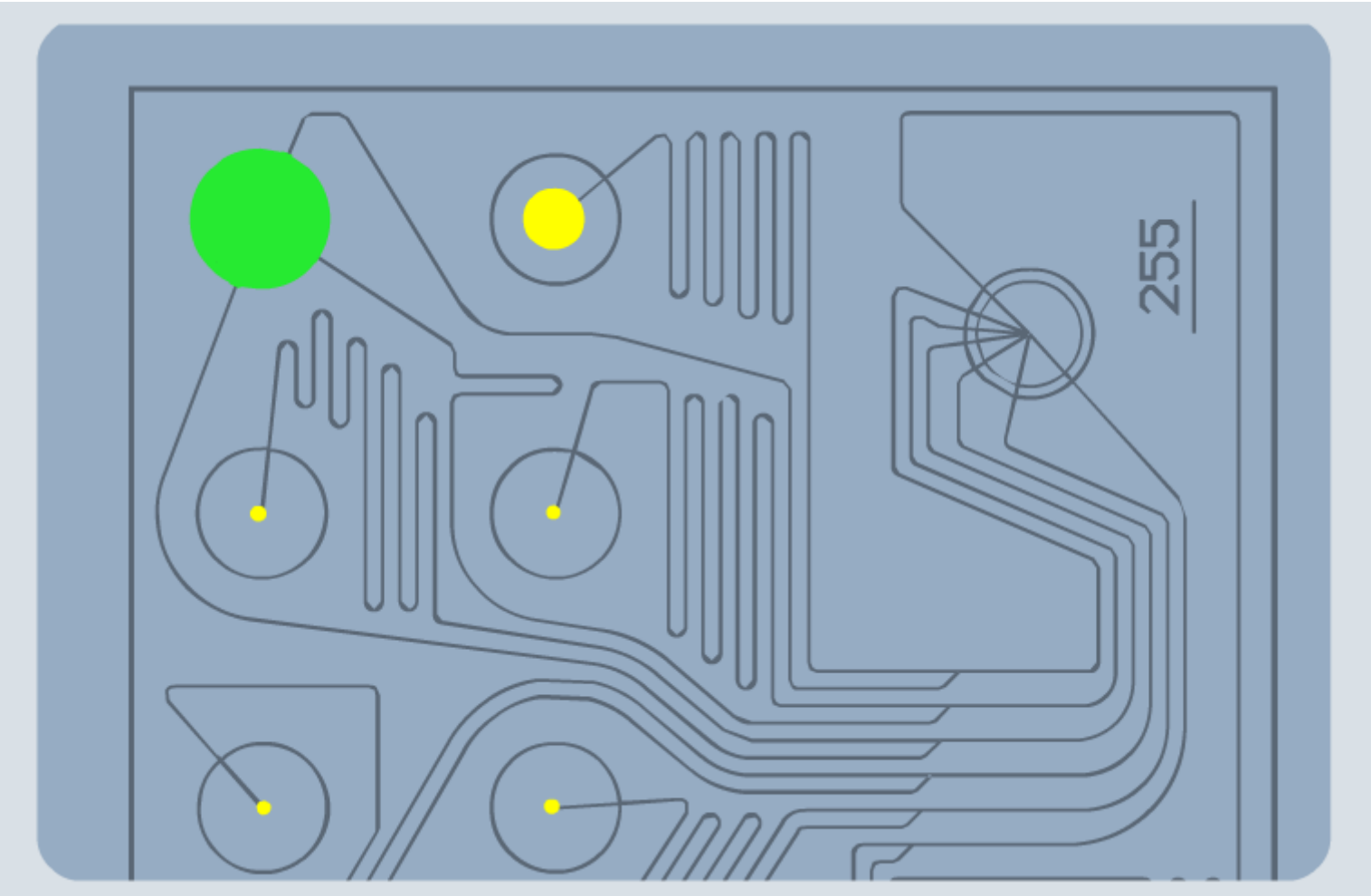
- micro-channels are filled with gel or buffer
- sample, ladder and reagents are filled into the respective wells
- chip preparation in less than 5 minutes

Benefits

- convenient handling
- minimized risk of cross-contamination
- versatile design for multiple experiments on one platform



Flow cytometry on the chip - Detection Animation



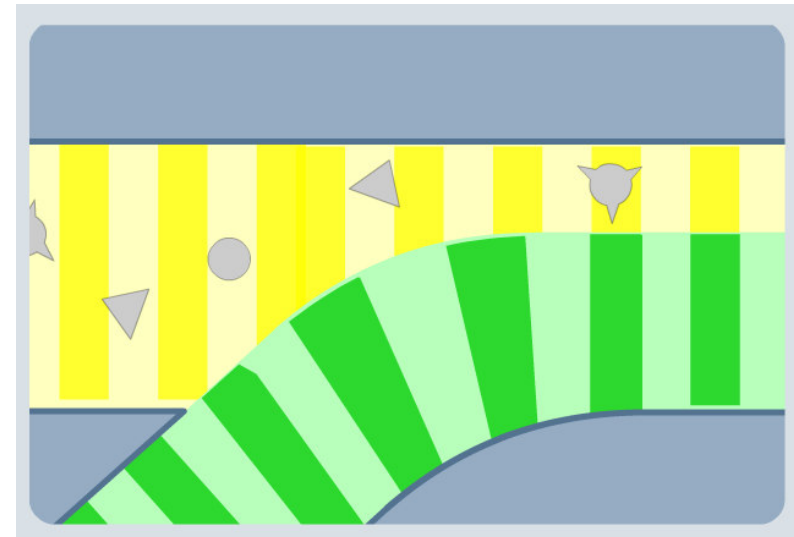
Flow Cytometry on a Chip - Hydrodynamic Focusing

All six cell samples are hydrodynamicly focused to one side of the micro channel

At each of the six pinch points the cell stream is joined by a buffer stream from one of the two buffer wells

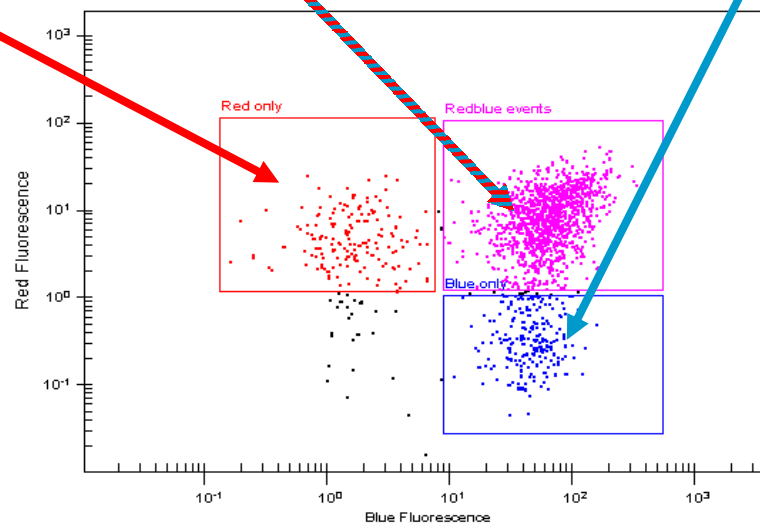
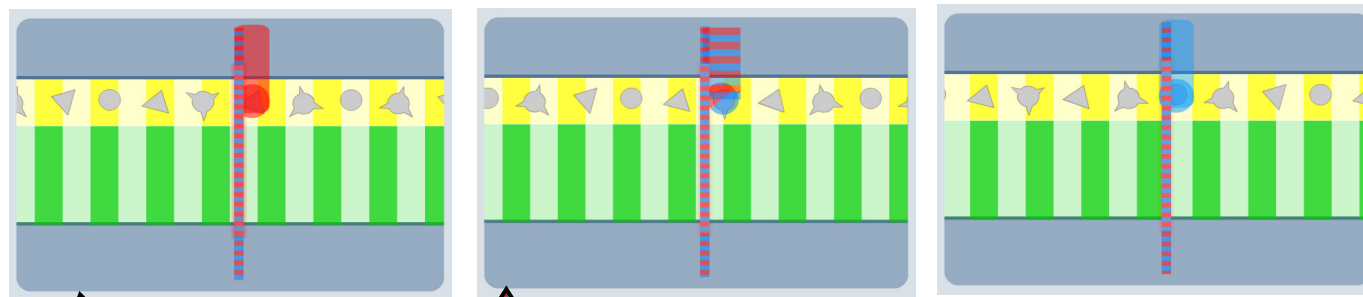
The two liquids do not mix immediately

The cells then move towards the detector in single file



Flow Cytometry on a Chip

- Two Color Detection- Three Types of Events



Dot plot view for easy data evaluation

Some Target applications



Apoptosis:

- Annexin V** **Detection of phosphatidylserine on the cell surface**
- Caspase-3** **Detection of activated caspase-3 in the cytoplasm**

Transfection Efficiency Monitoring:

- GFP:** **Detection of GFP-transfected cells**
- Antibody staining:** **Detection of transfected cells expressing the encoded protein**

Protein Expression Monitoring:

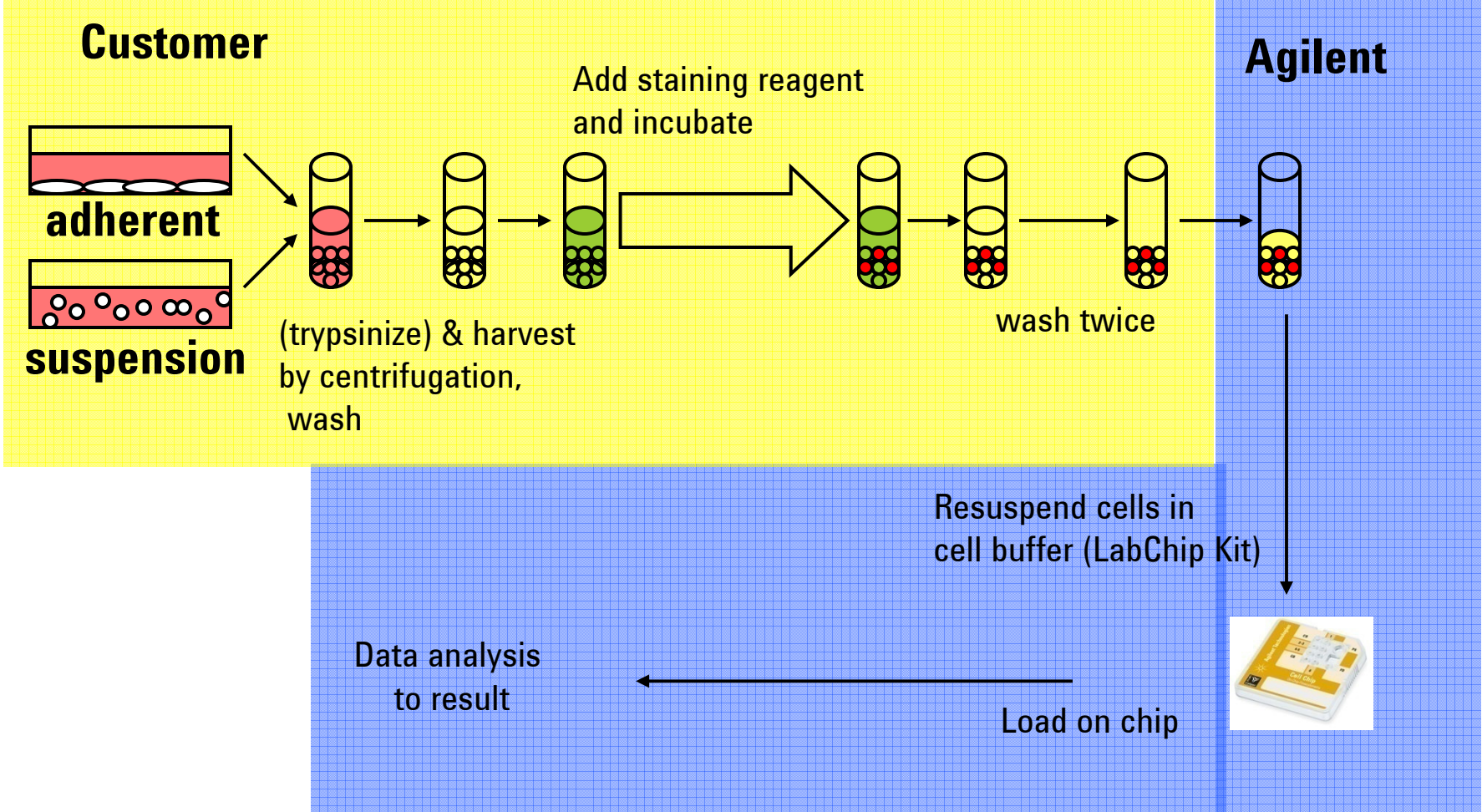
Extracellular and Intracellular Antibody staining for detection of protein expressed on the cell surface, in the cytoplasm, or in the nucleus

Gene silencing:

- Optimization of siRNA transfection procedure**
- Verify silencing by cellular protein expression measurement**
- Correlation of siRNA uptake and gene knockdown**

Cell assays: sample preparation

Typical workflow:

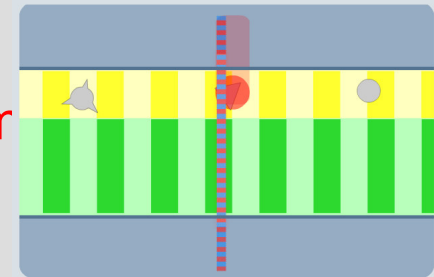


Flow Cytometry on a Chip - Optics & Detection

2100 Bioanalyzer

Red detection channel:

- 620-645 nm excitation with Laser (Maximum 630 nm)
- 674-696 nm detection range (Maximum 680 nm)

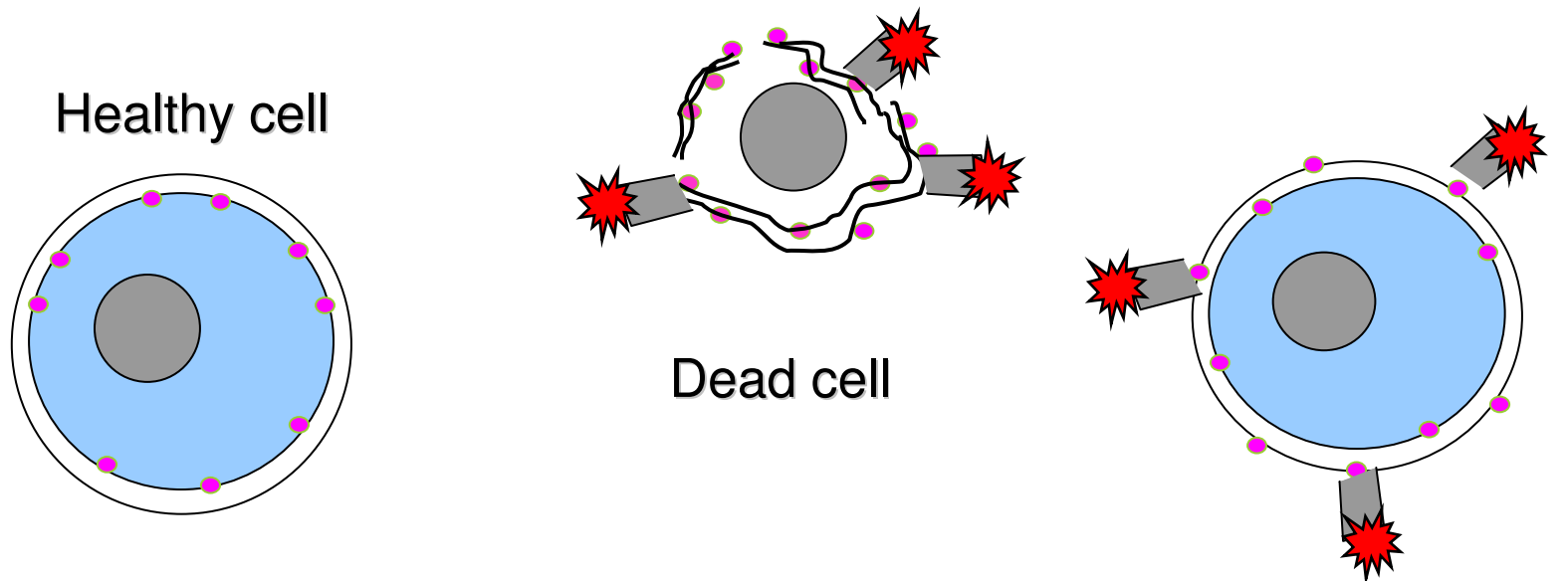


Blue detection channel:

- 458-482 nm excitation with LED (Maximum 470 nm)
- 510-540 nm detection range (Maximum 525 nm)

Cell Assays - Applications: Apoptosis

Annexin Binding

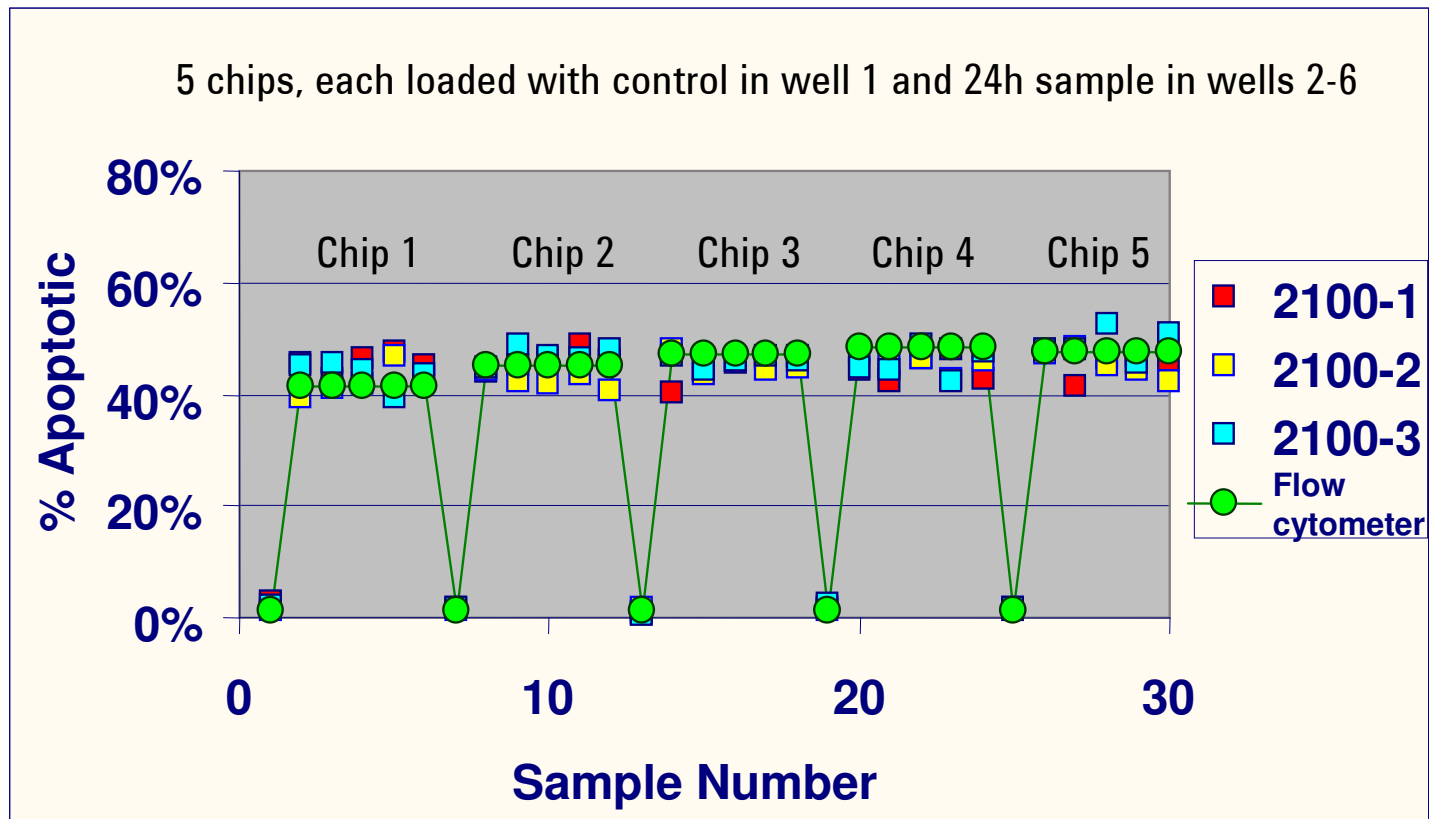


Live dye: Calcein
biotin-Annexin+ Cy5-streptavidin

Phosphatidyl-serine from inner leaflet flips to outer membrane during apoptosis and can be labeled by Annexin V

Annexin V Assay (24h Induction)

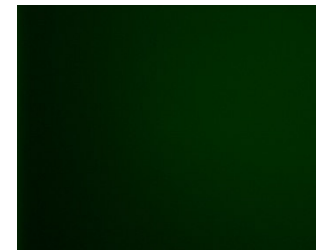
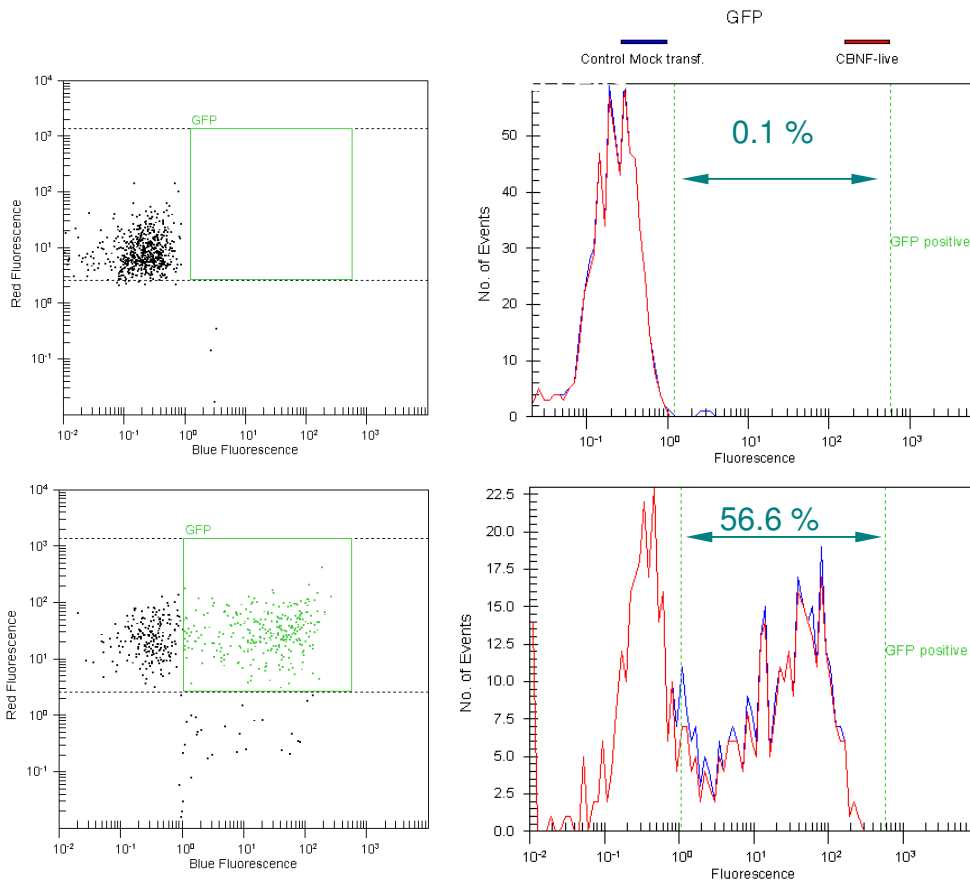
Three Bioanalyzer instruments vs a flow cytometer reference instrument



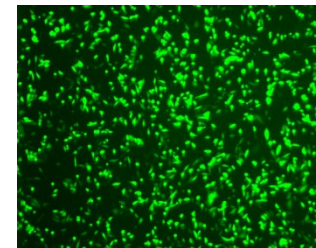
Applications: Protein Expression Analysis

GFP Transfection Efficiency Control

CHO-K1 cells were transfected with EGFP DNA and Lipofectamine.

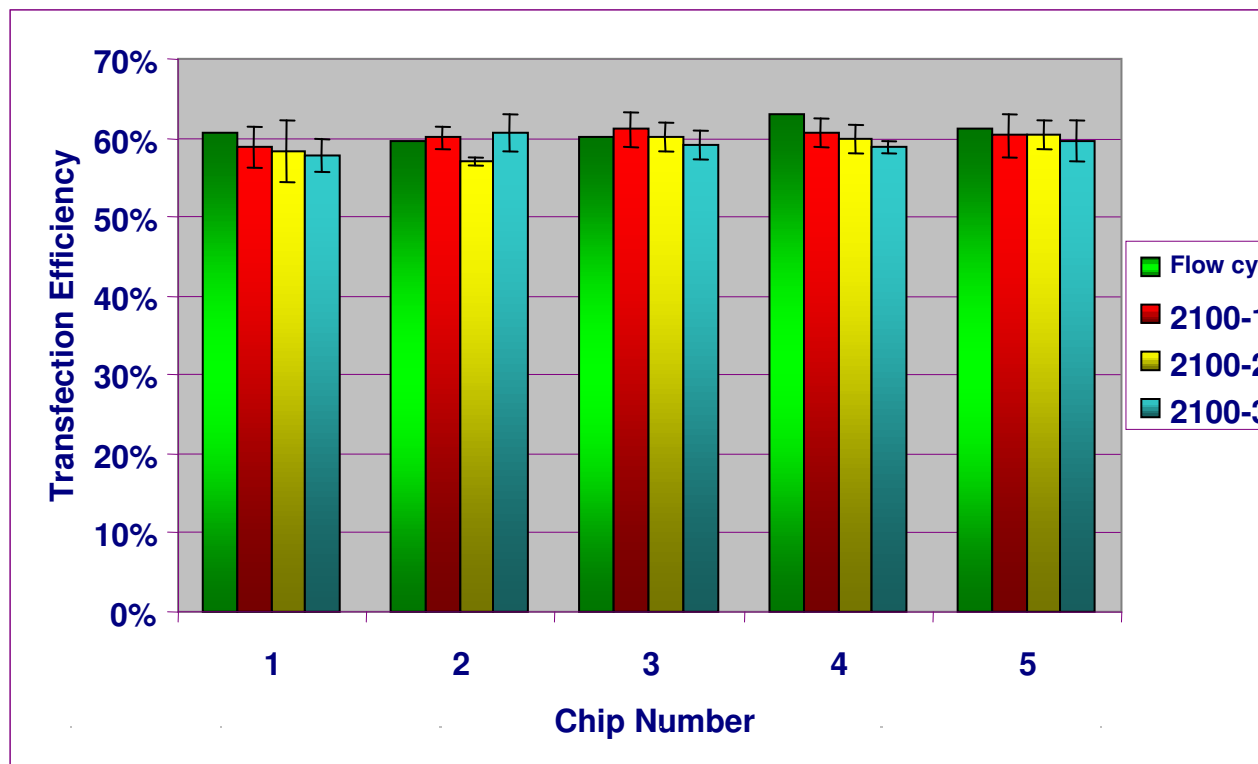


Control



EGFP transfected

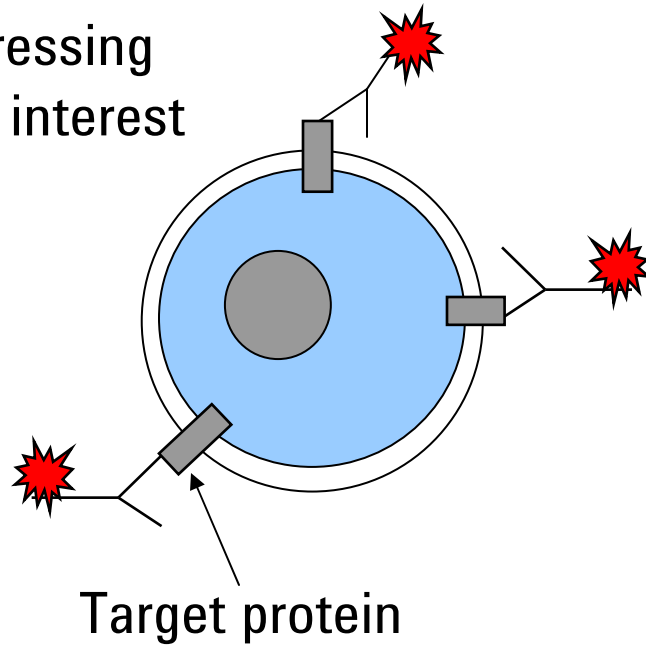
GFP Transfection Efficiency



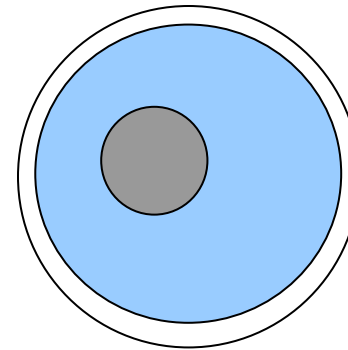
		2100-1	2100-2	2100-3	All	Flow cyt.
ctrl	mean	0.46	0.31	0.47	0.40	0.16
	SD	0.08	0.29	0.43	0.29	0.12
GFP	mean	60.19	59.15	59.26	59.53	60.90
	SD	2.13	2.48	2.10	2.26	1.22
	%CV	3.54	4.19	3.54	3.80	2.01

Flow Cytometry Assays Applications - Cell surface Antibody staining

Cell expressing protein of interest

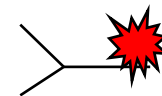


Cell not expressing protein of interest



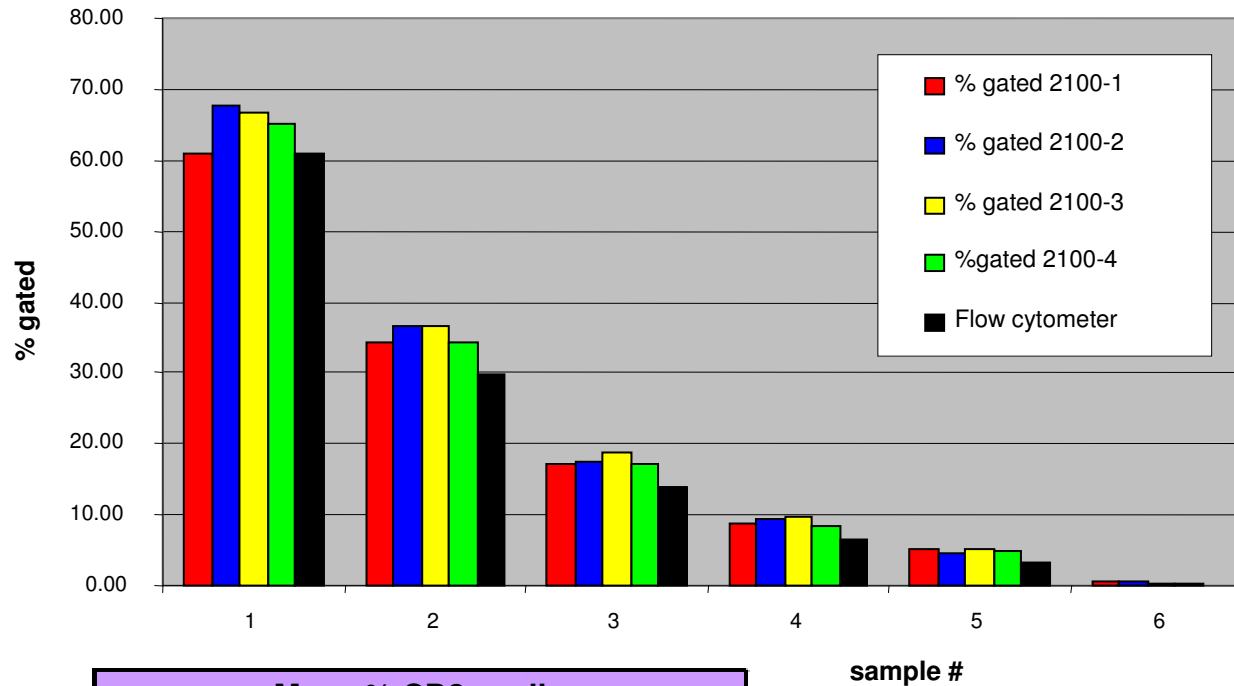
Live dye: Calcein

Cy5 or APC-labeled Antibody



Extracellular Antibody Staining

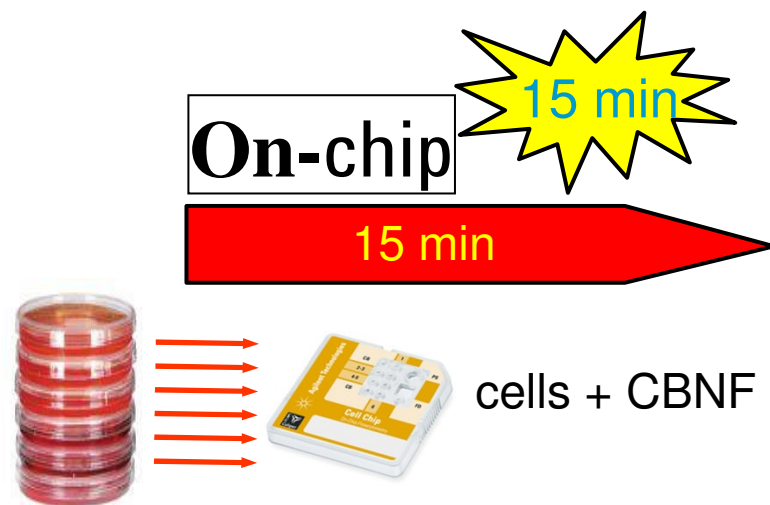
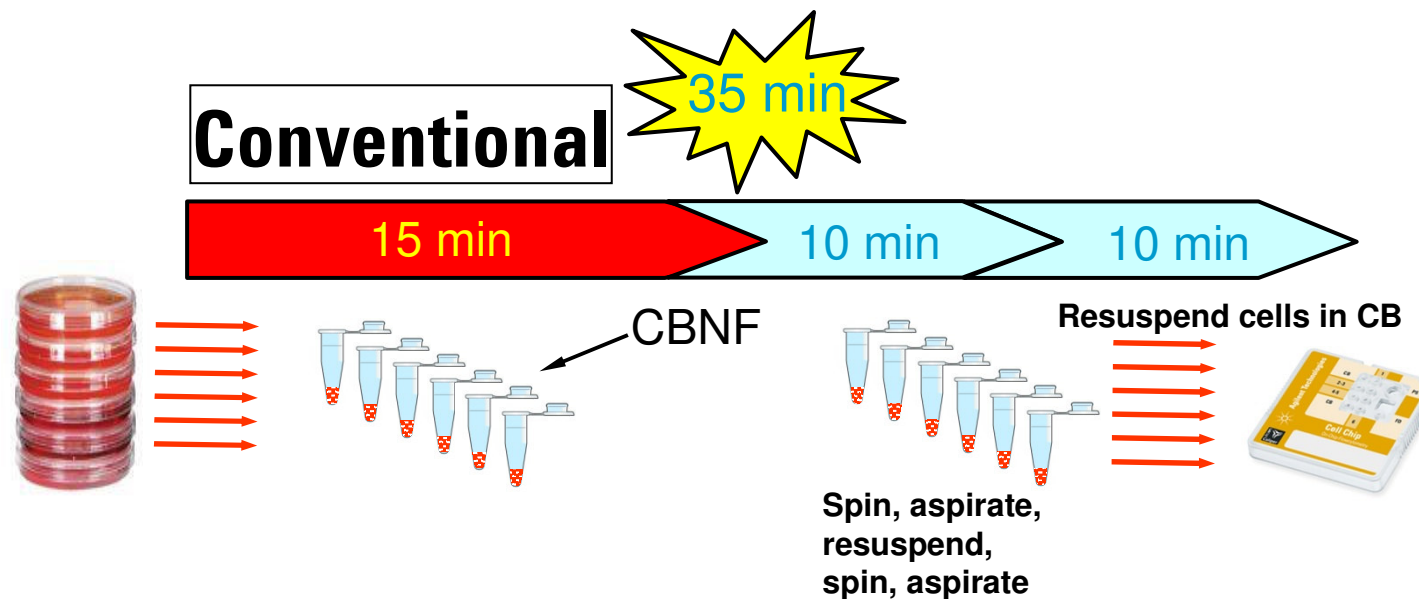
Averaged data per instrument



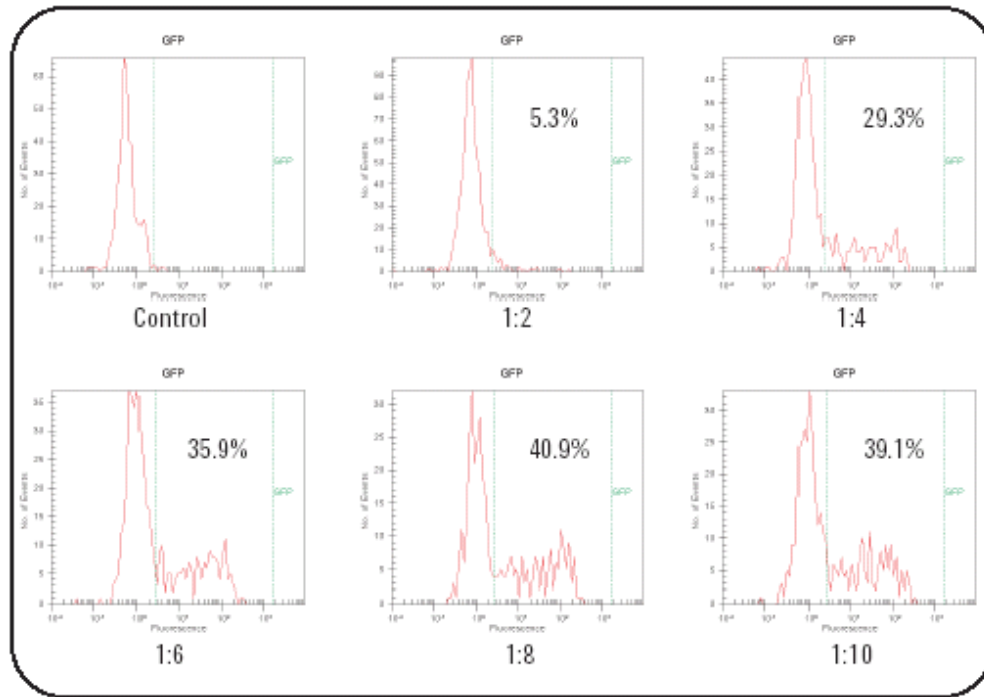
Mean % CD3+ cells				
2100-1	2100-2	2100-3	2100-4	Flow cyt.
60.9	67.8	66.6	65.0	60.9
34.4	36.7	36.7	34.3	29.8
17.3	17.6	18.7	17.2	13.8
8.9	9.4	9.9	8.3	6.5
5.1	4.4	5.3	4.9	3.2
0.8	0.6	0.3	0.3	0.0

Jurkat cells were stained with calcein alone or with calcein and APC-labeled anti-CD3 antibody. Mixtures of both populations were prepared at various ratios. Samples were analyzed with four 2100 instruments on 5 chips and compared to a flow cytometer reference instrument

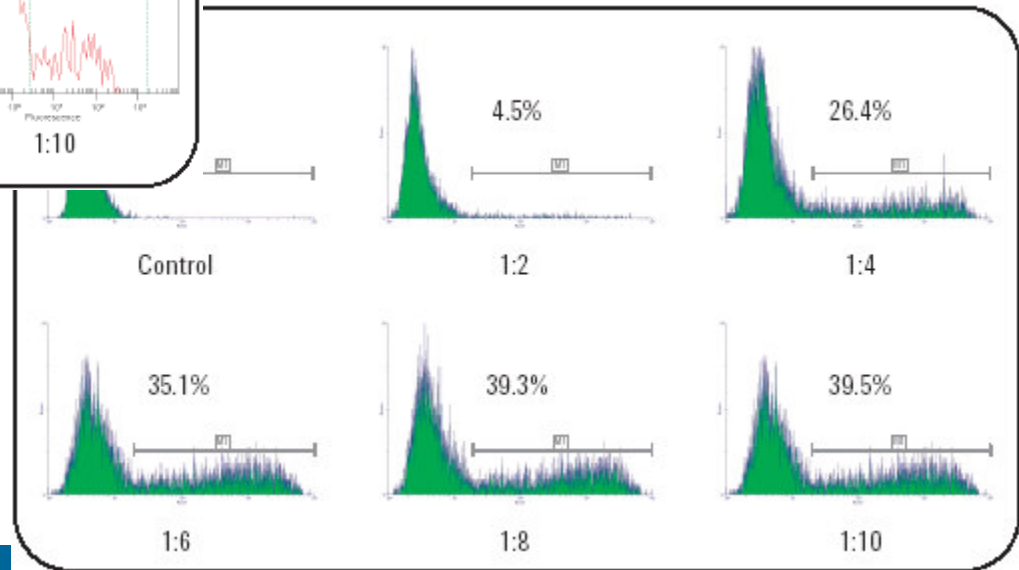
GFP On-Chip Staining - Workflow



GFP On-Chip Staining - Histogram Quality



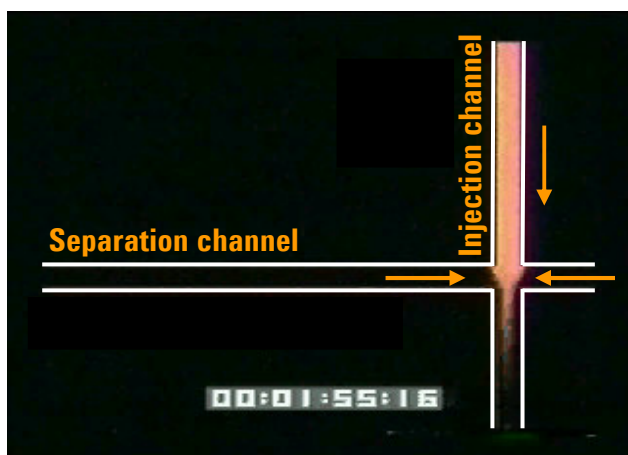
2100 bioanalyzer



Flow
Cytometer

Lab-on-a-Chip - Principle of Injection & Separation

1



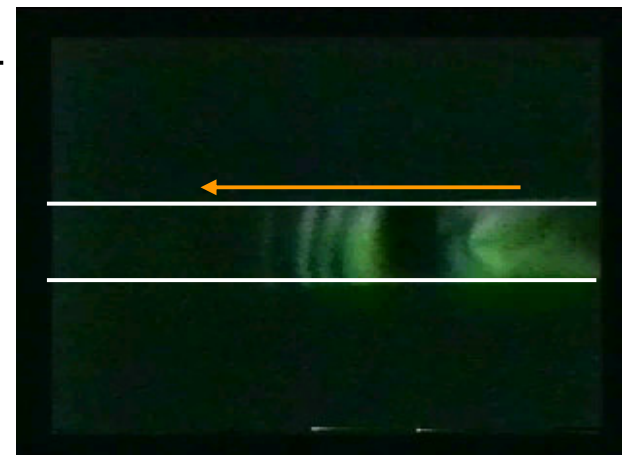
2



3



4



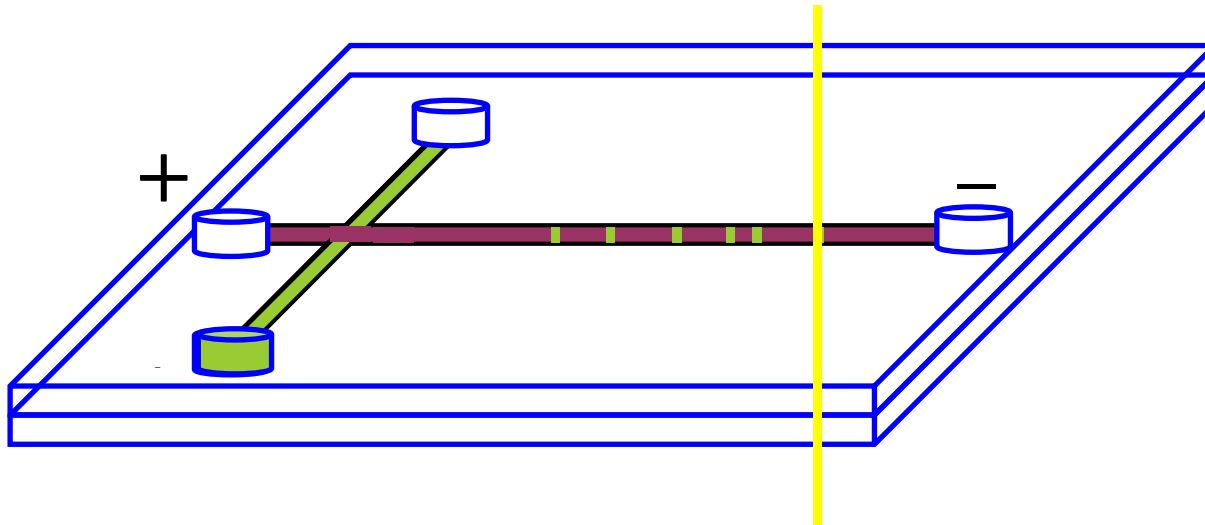
← Direction of electrodriven movement of liquids and molecules within liquids



Agilent Technologies

The LabChip® Approach - Simplified Model

(see chip animation.ppt)



DNA Applications



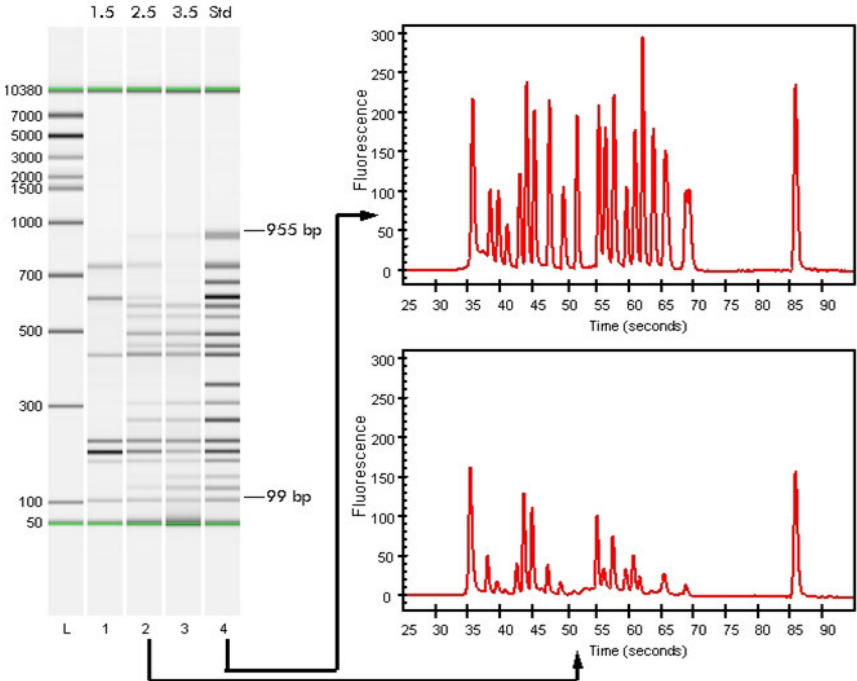
mPCR validation,
impurity check

Gene
Expression

Restriction
Digest Analysis

Food
Analysis

Forensic
Testing



Application Areas for the DNA Assays

PCR product purity

Multiplex PCR Applications

Gene expression analysis via RT-PCR (target validation)

GMO testing

Pathogen detection (homeland defense, hospitals, environmental)

Genotyping applications

- **Duplications/ deletions**
- **Allele frequency**
- **Bacterial sub-typing**
- **Forensics**

Cancer diagnostics

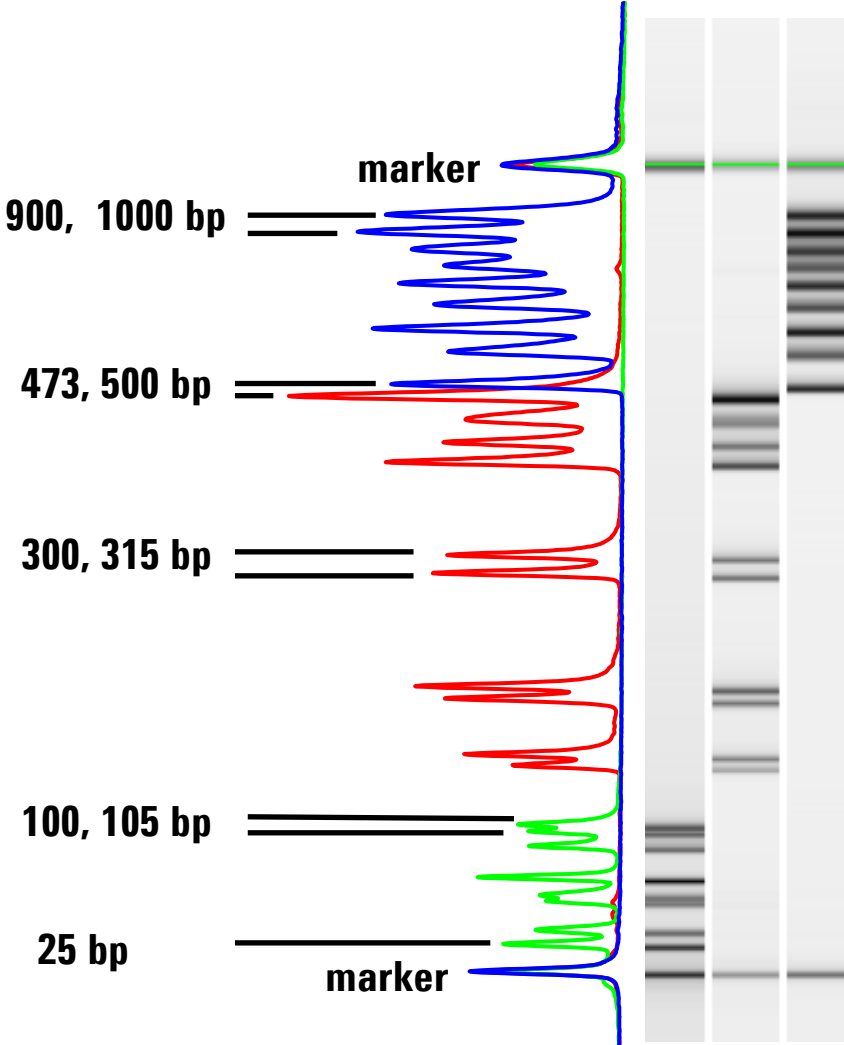


DNA Kit Specification

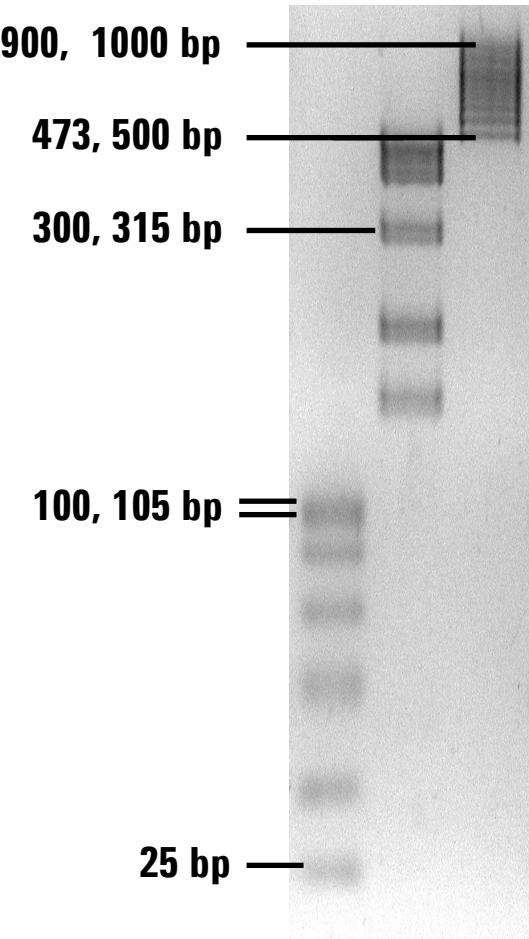
	DNA 1000 Assay	DNA 7500 Assay	DNA 12000 Assay
Sizing range	25–1000 bp	100–7500 bp	100–12000 bp
Sizing resolution	25–100 bp: 5 bp 100–500 bp: 5 % CV 500–1000 bp: 10 % CV	100–1000 bp: 5 % 1000–7500 bp: 15 %	100–1000 bp: 5 % 1000–12000 bp: 10 %
Sizing accuracy	± 10 %	± 10 % CV *	±15 %
Sizing reproducibility	5% CV *	5% CV *	5% CV *
Quantitation accuracy	20% CV *	20% CV *	25% CV *
Quantitation reproducibility	25–500 bp: 15 % CV * 500–1000 bp: 5 % CV *	100–1000 bp: 10 % CV * 1000–7500 bp: 5 % CV *	100–1000 bp: 15 % CV * 1000–12000 bp: 10 % CV *
Quantitative range	0.1 - 50 ng/μL *	0.1 - 50 ng/μL *	0.1 - 50 ng/μL *
Maximum salt concentration in sample	250 mM for KCl, 15 mM for MgCl ₂ 250 mM NaCl	250 mM for KCl 15 mM for MgCl ₂ 250 mM NaCl	250 mM for KCl 15 mM for MgCl ₂ 250 mM NaCl
25 chips per kit	DNA 12/chip = 300 samples/kit		

* Respective DNA ladder as sample

Data Format - Gel-Like Image c/w Agarose Gel

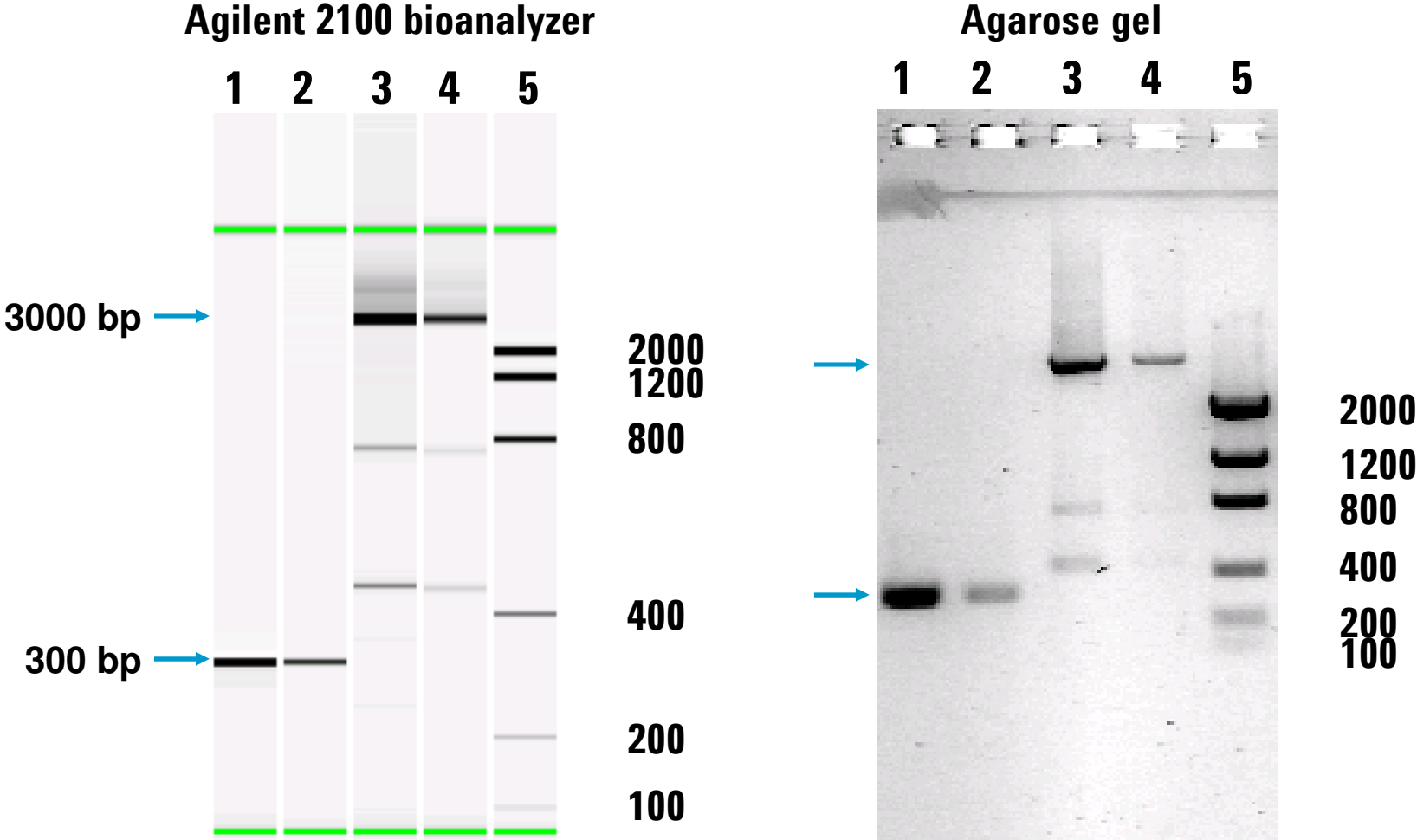


2100 bioanalyzer data
Gel-like image

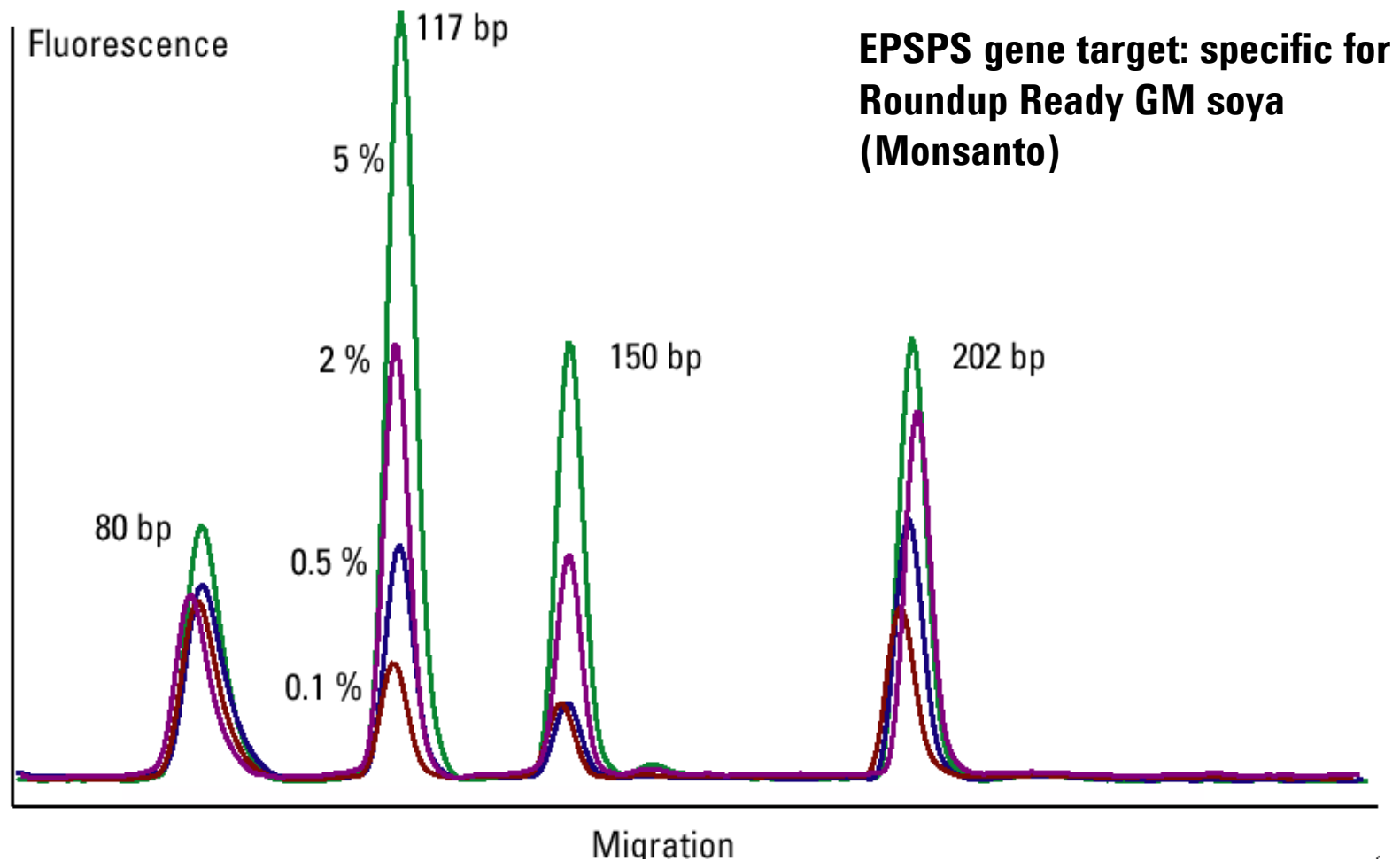


2 % agarose gel stained
with Ethidiumbromide

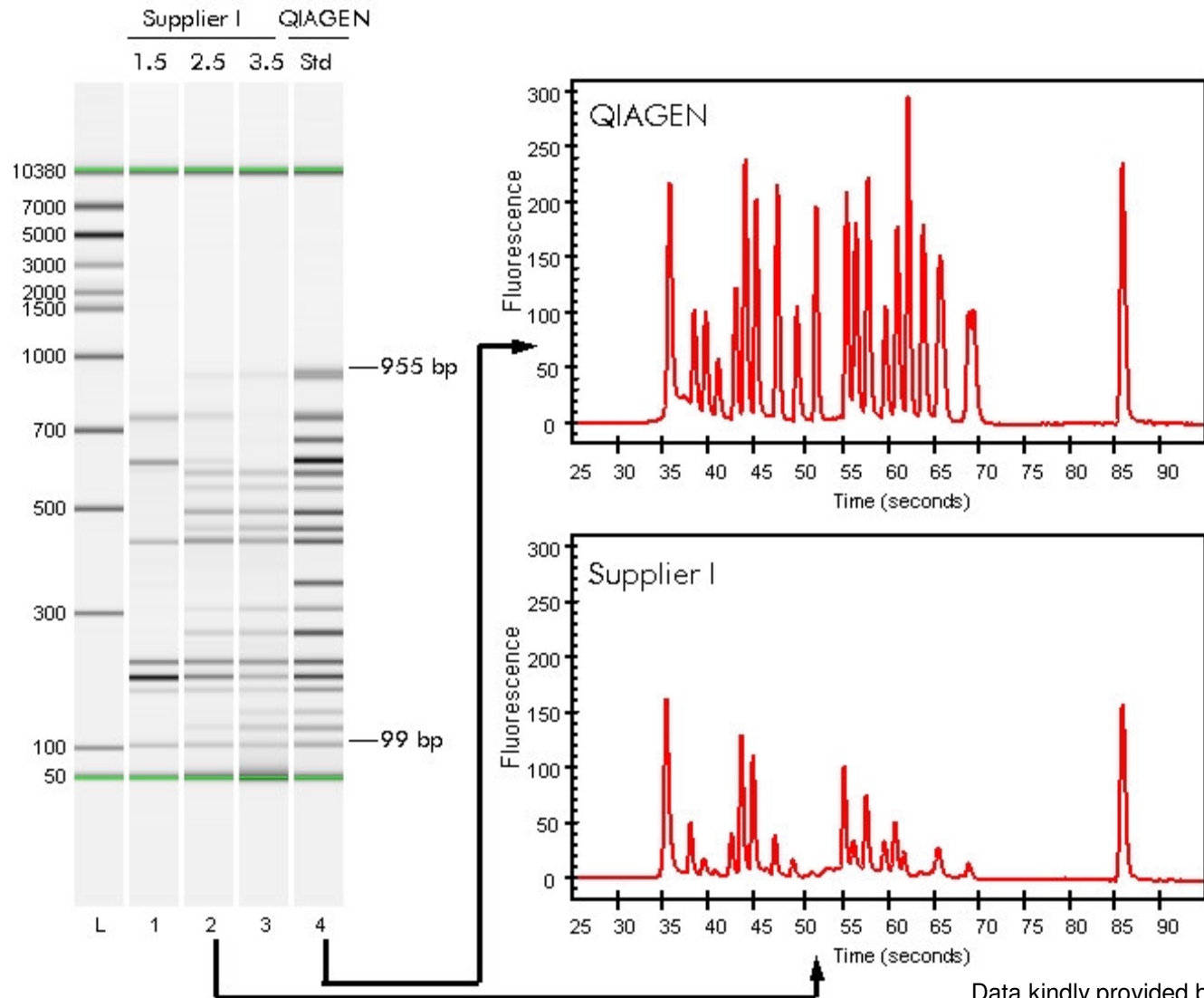
Determination of PCR Product Impurity



GMO Detection: Determination of GM Soya Percentage

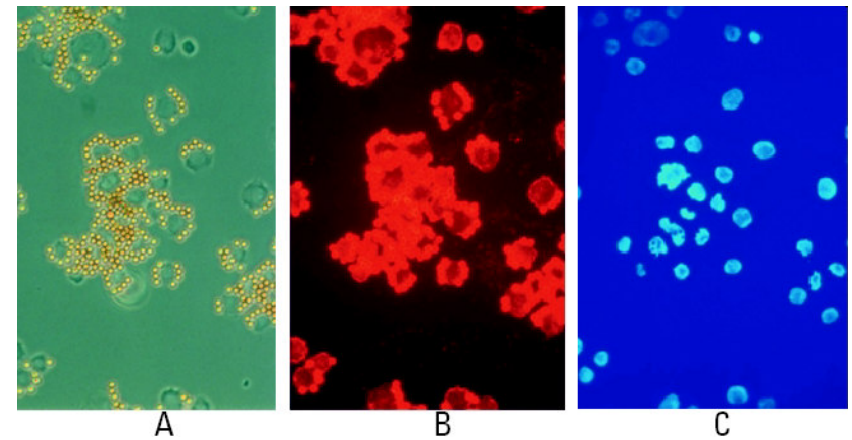
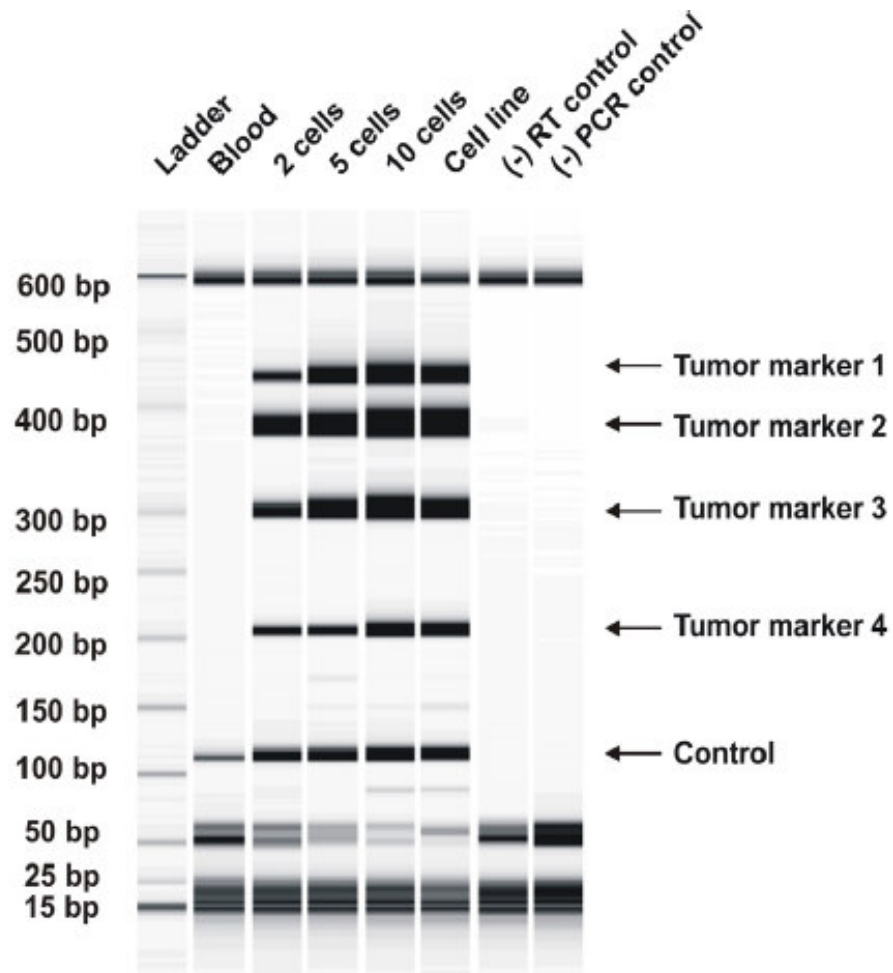


Optimization of Multiplex PCR on a 19-plex PCR



Data kindly provided by QIAGEN GmbH, Germany

Tumor Diagnostics



1. Spiking experiment with given amount of cancer cells
2. Enrichment with AdneGen Cancer Select kit (antibody based immunomagnetic enrichment.)
3. Multiplex Amplification with AdnaGen CancerDetect kit
4. Detection with Agilent 2100 Bioanalyzer and DNA 500 LabChip kit

Data kindly provided by Adnagen

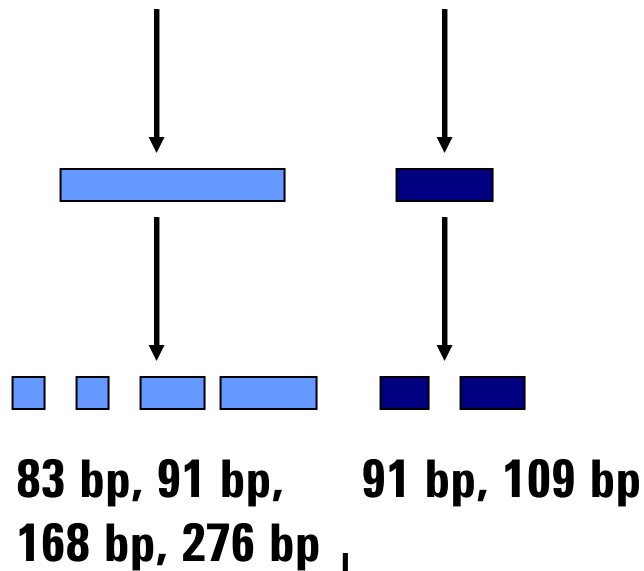
Detection of Single Base Mutations (1)

in Exons 7 and 8 of the Human p53 Gene by RFLP Mapping using the DNA 7500 kit



exon 7

exon 8



**83 bp, 91 bp,
168 bp, 276 bp**

91 bp, 109 bp



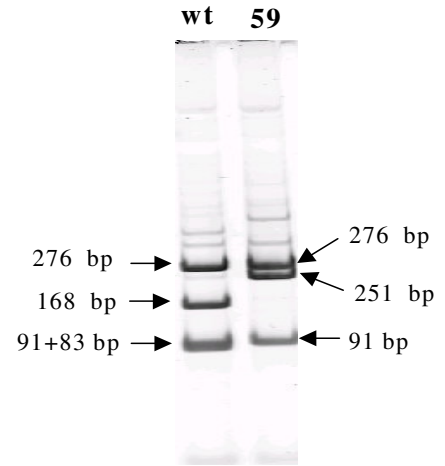
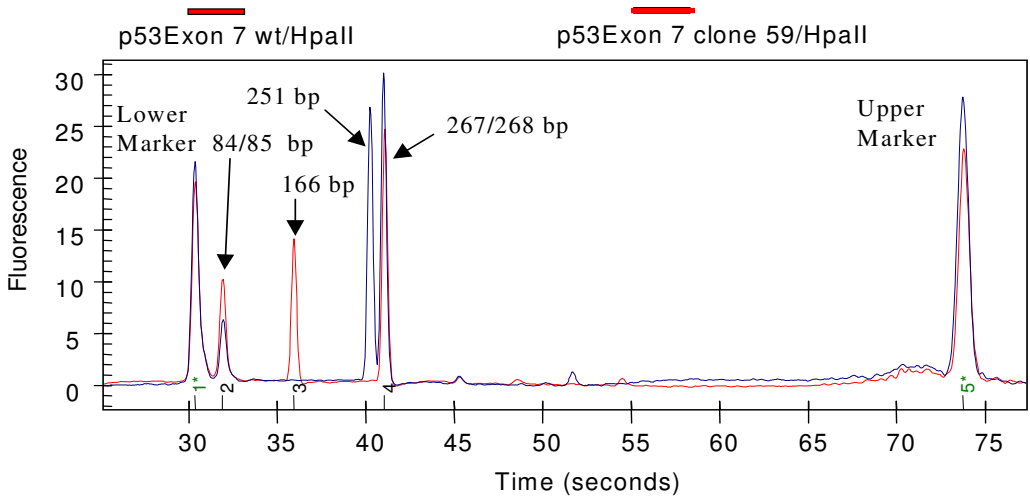
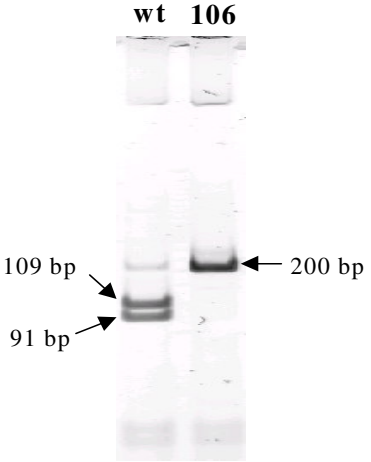
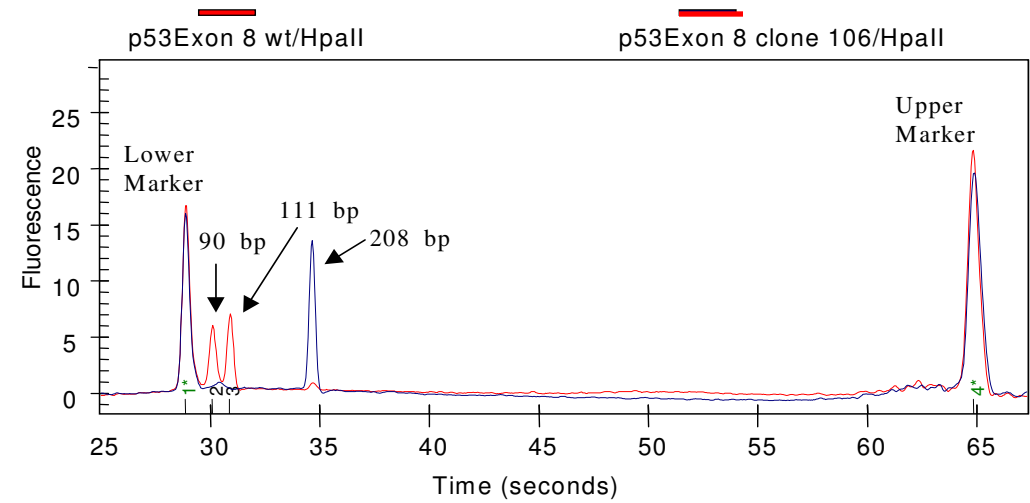
Amplify exons 7 and 8 (resulting products:
618 bp fragment and 200 bp fragment)

Digest with Hpa II

In each example one of the restriction sites can be deleted by a point mutation

Analyze using Agilent 2100 bioanalyzer and 4-20 % acrylamide gel

Detection of Single Base Mutations (2)



Label free Analysis of Microsatellite Instabilities

Clinical Diagnostics and Molecular Diagnostics of Cancer

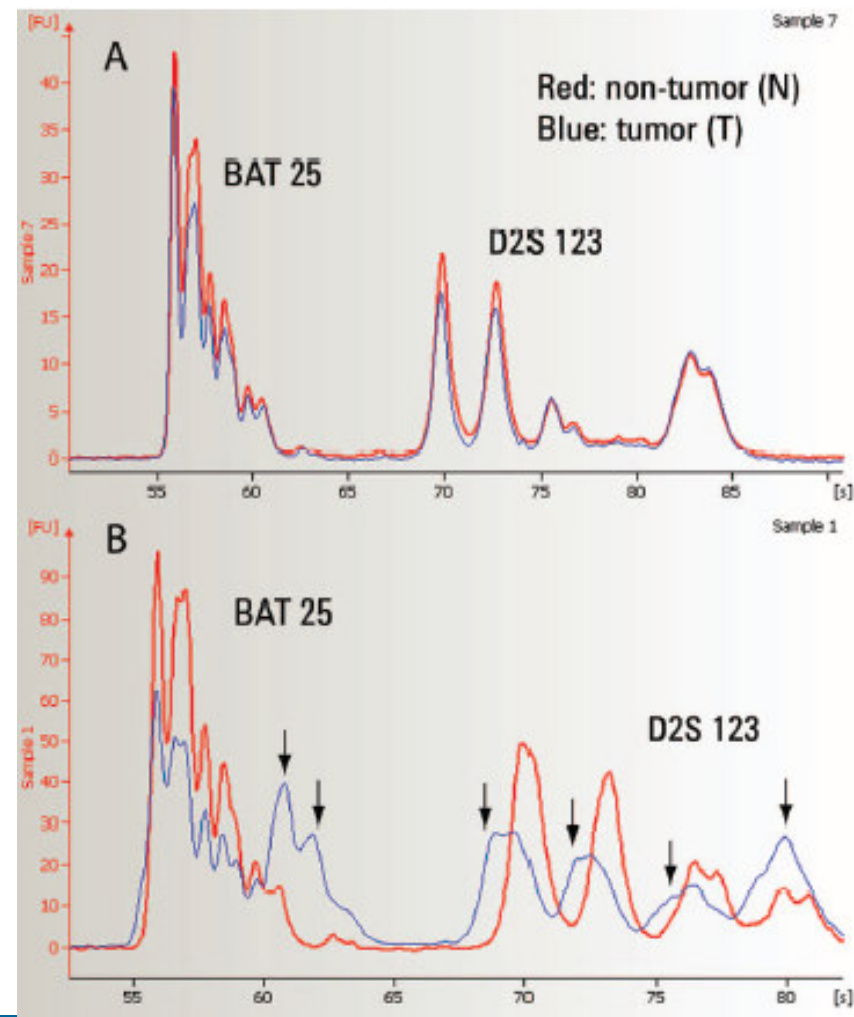
Microsatellite instabilities present in 10-15% of colon and gastric carcinomas

Study: 40 cases of colon carcinoma

5 microsatellite loci investigated

Results compared with traditional PAGE:

95% concordance rate



RNA Applications



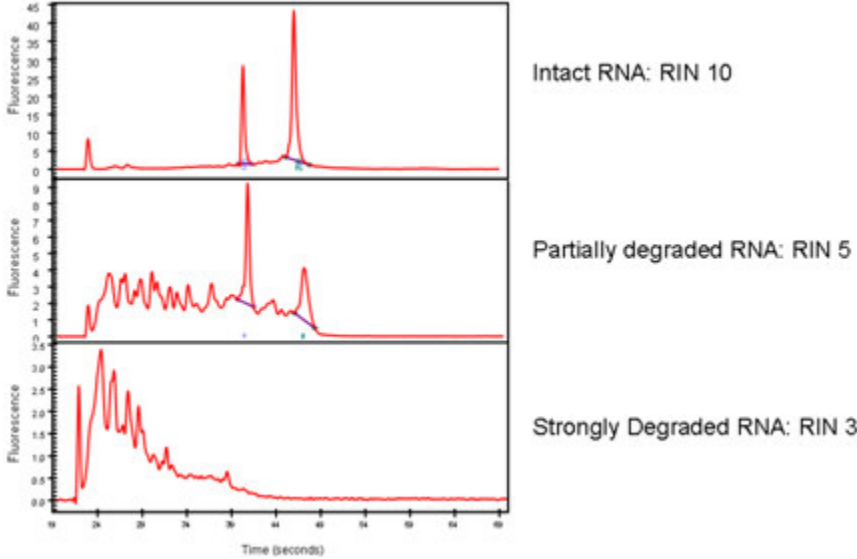
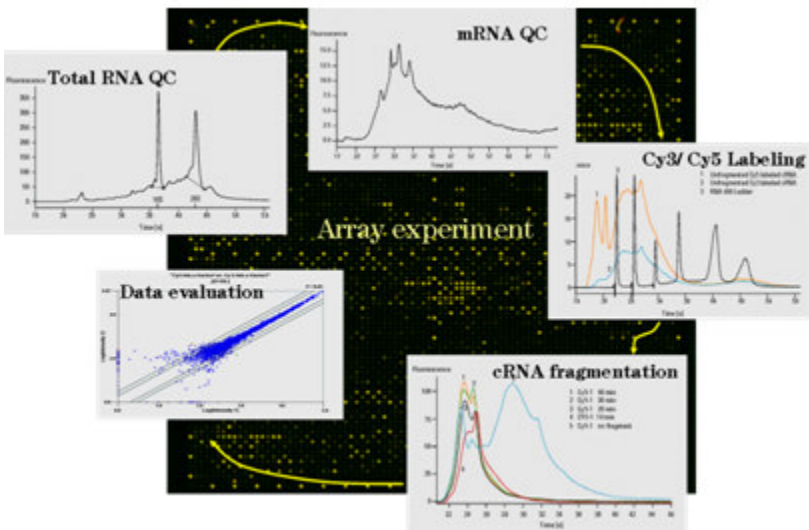
RNA QA/QC for
Microarrays

Gene
Expression

RNA QA/QC for
qPCR

RNA QA/QC for
mPCR

smallRNA
QA/QC



Agilent 2100 bioanalyzer: the industry standard in RNA QC

Electrophoretic sizing, quantitation and QC of XNA and Proteins on a small glass Chip as done traditionally on slab gels (Agarose or SDS-PAGE)

First commercially available Lab-on-a-Chip product (since October 1999)

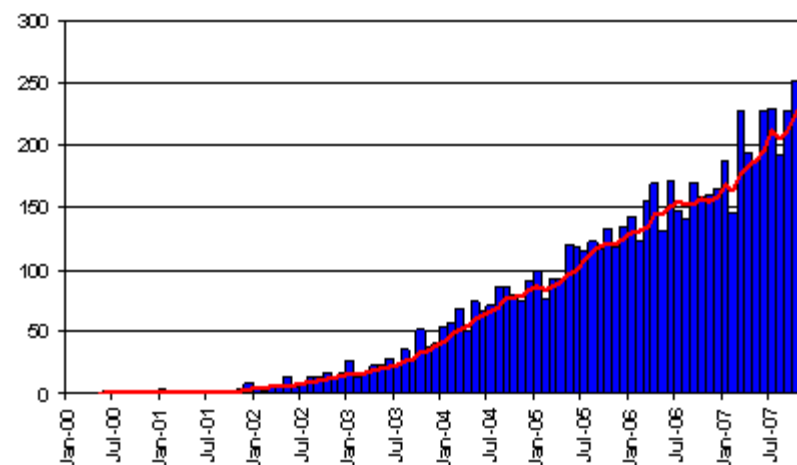
Analysis of totalRNA, mRNA and small RNA samples in ng and pg concentration range

Standardized RNA integrity assessment with **RIN*** algorithm

Multi-analysis capabilities: DNA, RNA, Proteins and Flow Cytometry

RIN = RNA Integrity Number, an Agilent patented algorithm to Determine RNA quality in a normalized way

Number of BioA publications/month



January 2008: ~ 7200 citations

RNA Kit Specifications

Kit	RNA 6000 Nano total RNA	RNA 6000 Nano mRNA	RNA 6000 Pico total RNA	RNA 6000 Pico mRNA	Small RNA Assay
Analytical Specifications					
Quantitative range	25–500 ng/μL	25–250 ng/μL			50-2000 pg/μL of purified miRNA in water
Qualitative range	5–500 ng/μL	25–250 ng/μL	50–5000 pg/μL in water	250–5000 pg/μL in water	50-2000 pg/μL of purified miRNA in water
Sensitivity (S/N>3)	5 ng/μL in water	25 ng/μL in water	50 pg/μL in water 200 pg/μL in TE	250 pg/μL in water 500 pg/μL in TE	50 pg/μL in water**
Quantitation reproducibility	10% CV (within a chip)	10% CV (within a chip)	20 % CV (within a chip)	20 % CV (within a chip)	25 % CV (within a chip)
Quantitation accuracy	20 % CV*	20 % CV*	30 % CV*	30 % CV*	
Maximum sample buffer strength	100 mM Tris, 0.1 mM EDTA or 125 mM NaCl or 15 mM MgCl ₂	100 mM Tris, 0.1 mM EDTA or 125 mM NaCl or 15 mM MgCl ₂	50 mM Tris, 0.1 mM EDTA or 50 mM NaCl or 15 mM MgCl ₂	50 mM Tris, 0.1 mM EDTA or 50 mM NaCl or 15 mM MgCl ₂	10 mM Tris, 0.1 mM EDTA
Physical Specifications					
Analysis time	30 minutes	30 minutes	30 minutes	30 minutes	30 minutes
Samples per chip	12	12	11	11	11
Sample volume	1 μL	1 μL	1 μL	1 μL	1 μL
Kit stability	≥4 months at 4 °C	≥4 months at 4 °C	≥4 months at 4 °C	≥4 months at 4 °C	≥4 months at 4 °C
25 chips per kit	RNA Nano 12/chip = 300 samples/kit		RNA Pico 11/chip = 275 samples/kit		

* Determined analyzing the RNA ladder as sample

** Measured for the 40 nt fragment of the Small RNA ladder

Features of the RNA 6000 Assays

total RNA

determine integrity and quality of total RNA

determination of RNA concentration

identify ribosomal peaks

calculate the ratio of ribosomal peaks (18S/28S or 16S/23S)

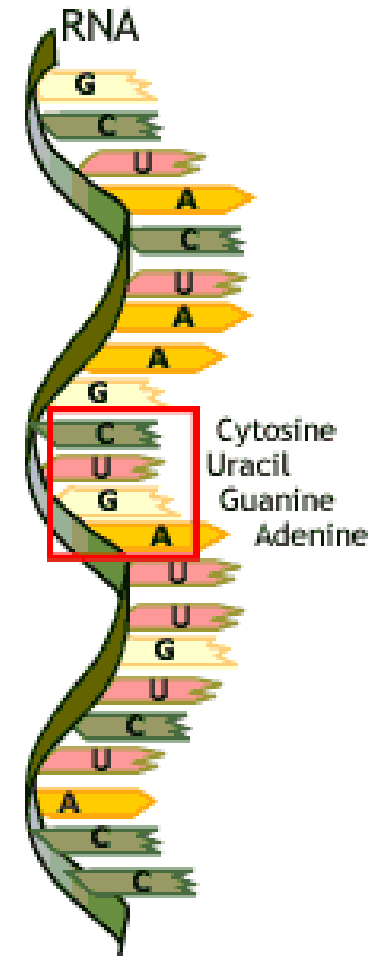
RNA integrity number (RIN)

mRNA

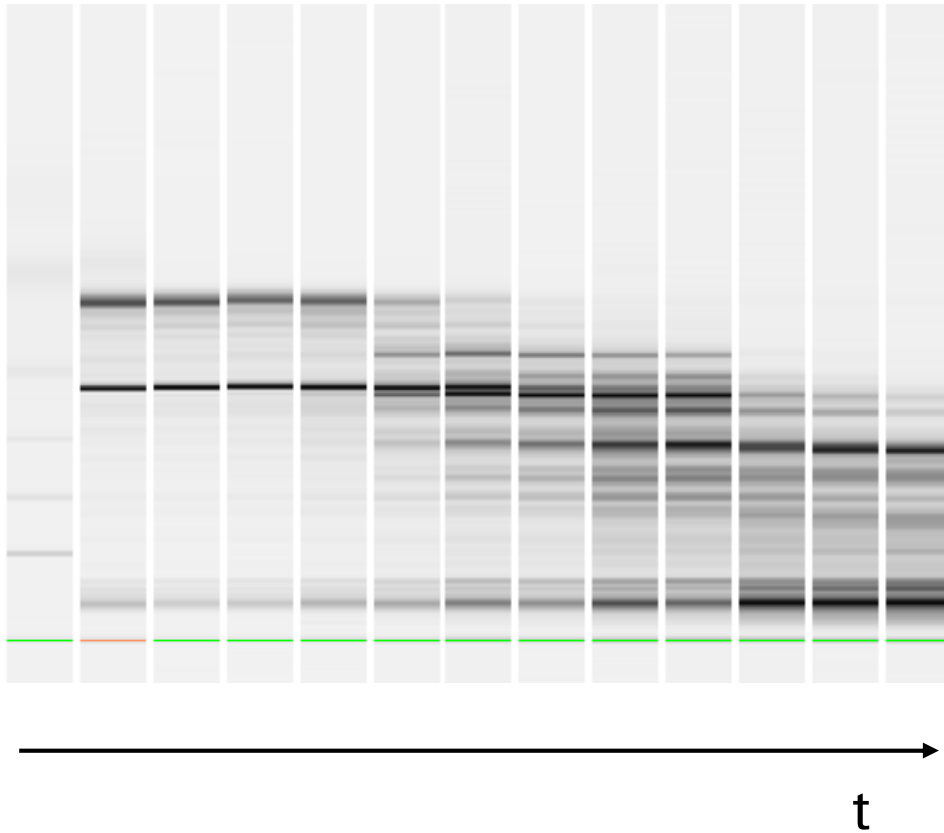
determine integrity and quality of mRNA samples

Determination of mRNA concentration

calculate % ribosomal RNA in mRNA samples



Problem Description



The ratio of ribosomal bands is not sufficient to describe RNA integrity!

RNA degradation is a gradual process.

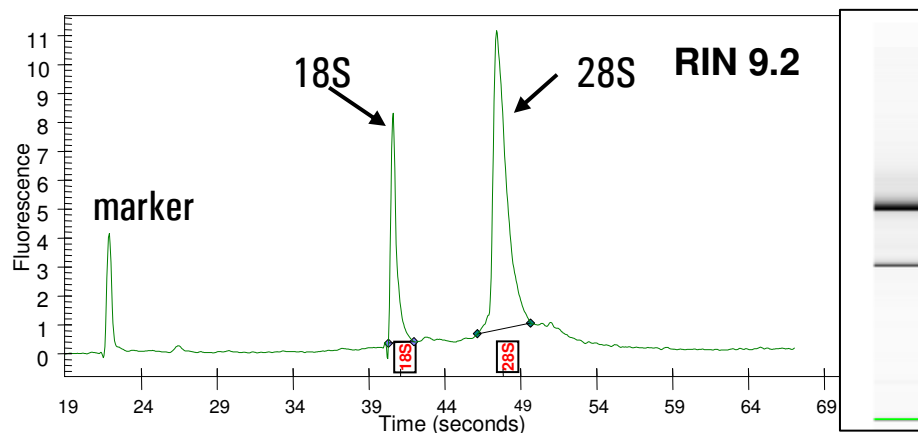
Results have to be interpreted by visual inspection.

Overlay of electropherograms only works well for samples with the same concentration.

Instrument dependency in signal height

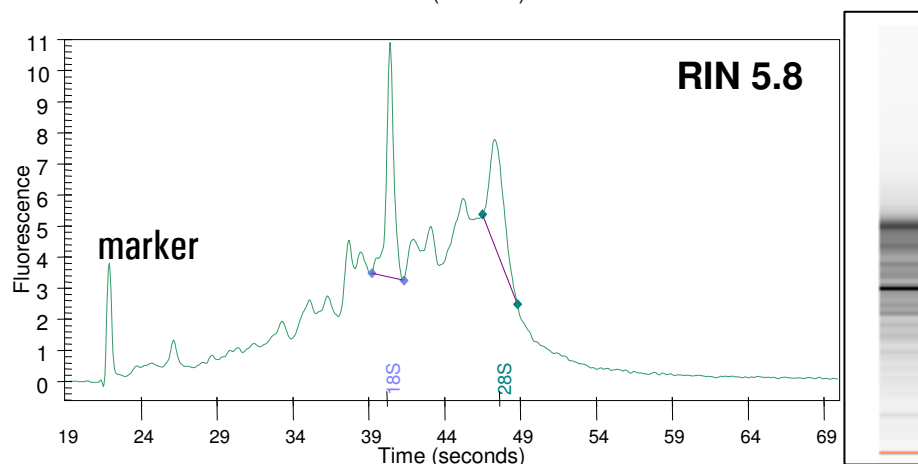


RNA Quality Control: Assessing Total RNA Integrity



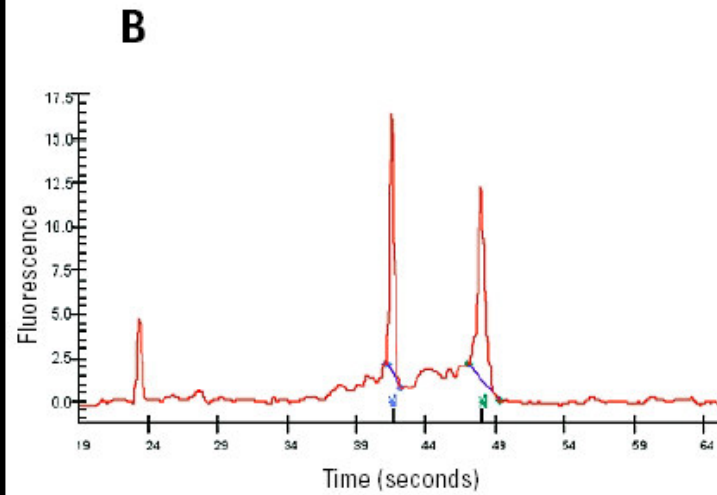
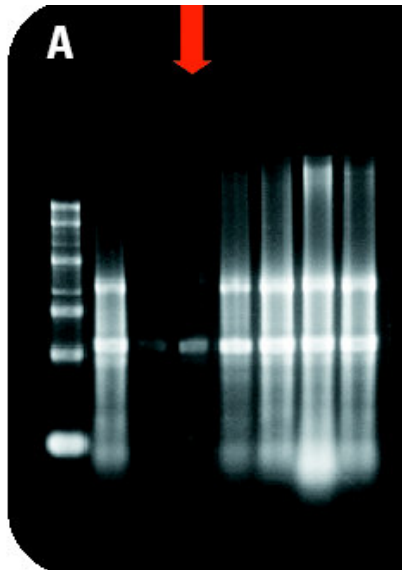
Typical first QC step after RNA sample prep prior to microarrays or real-time PCR

High quality total RNA

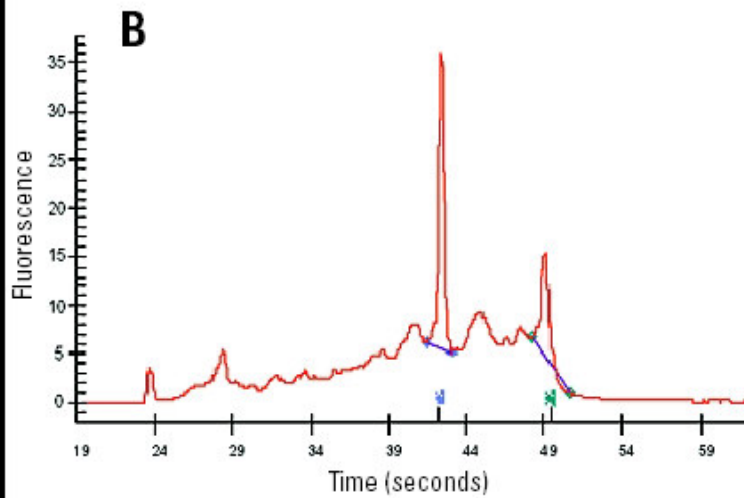
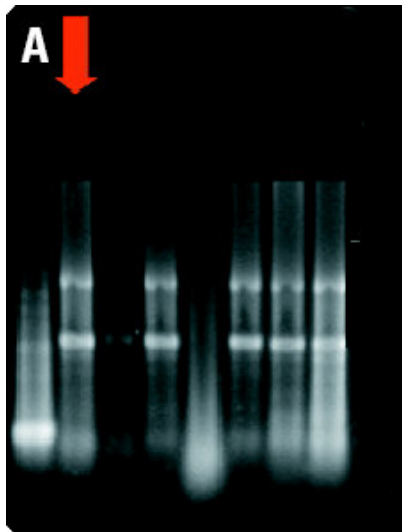


Partially degraded total RNA

Gel Chip Comparison

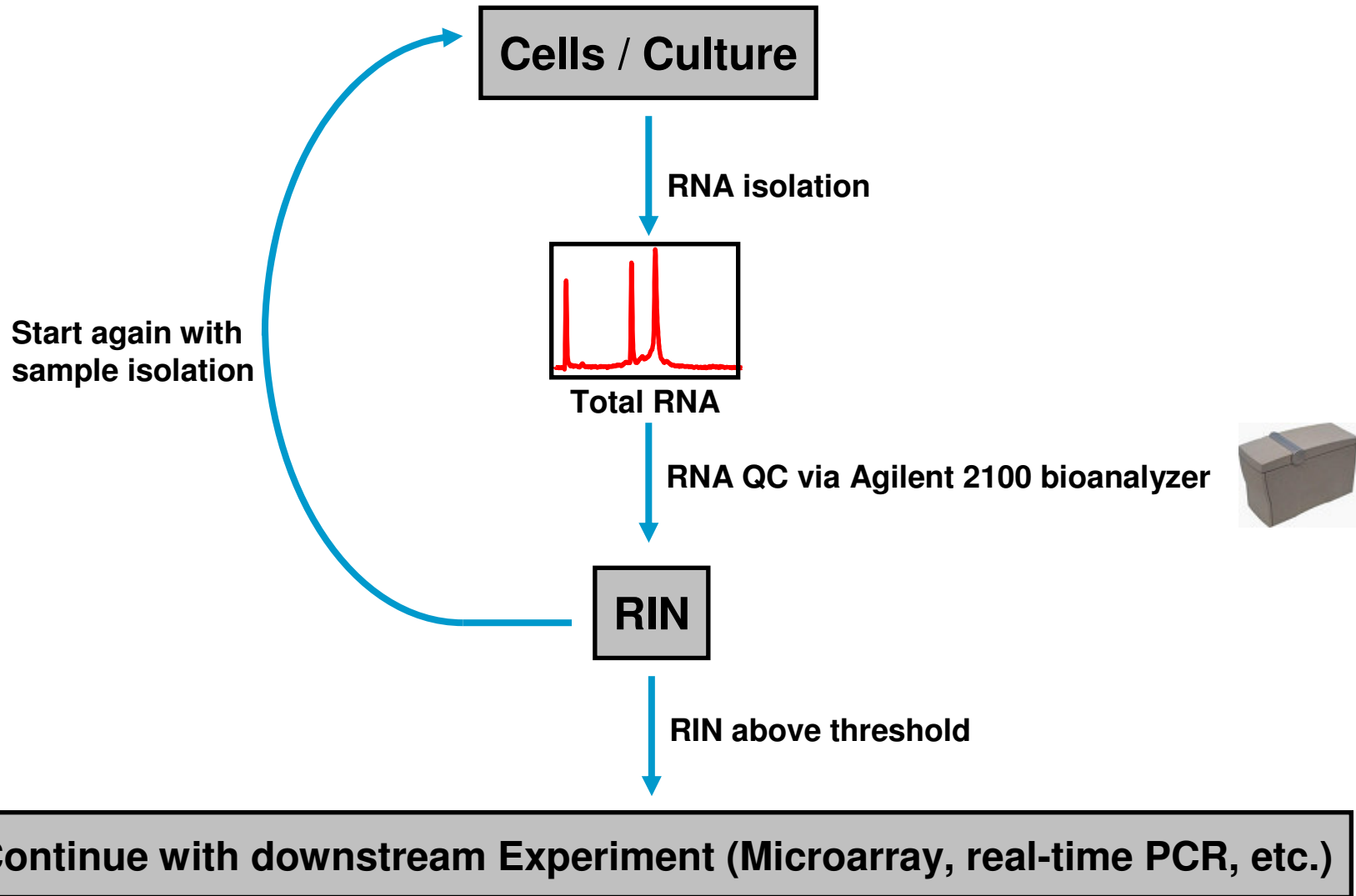


False Negative

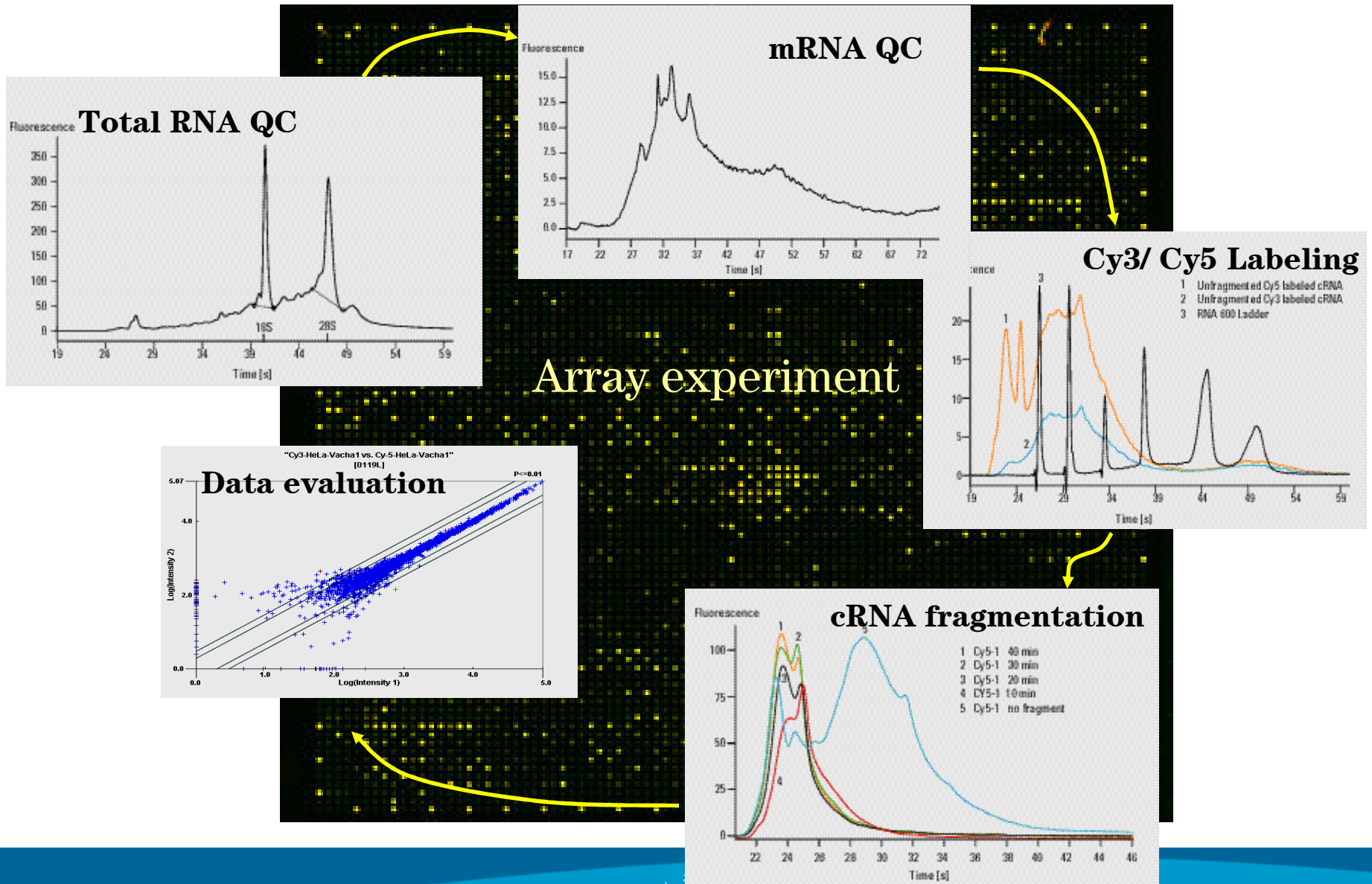


False Positive

RNA QC in Routine Gene Expression Workflow

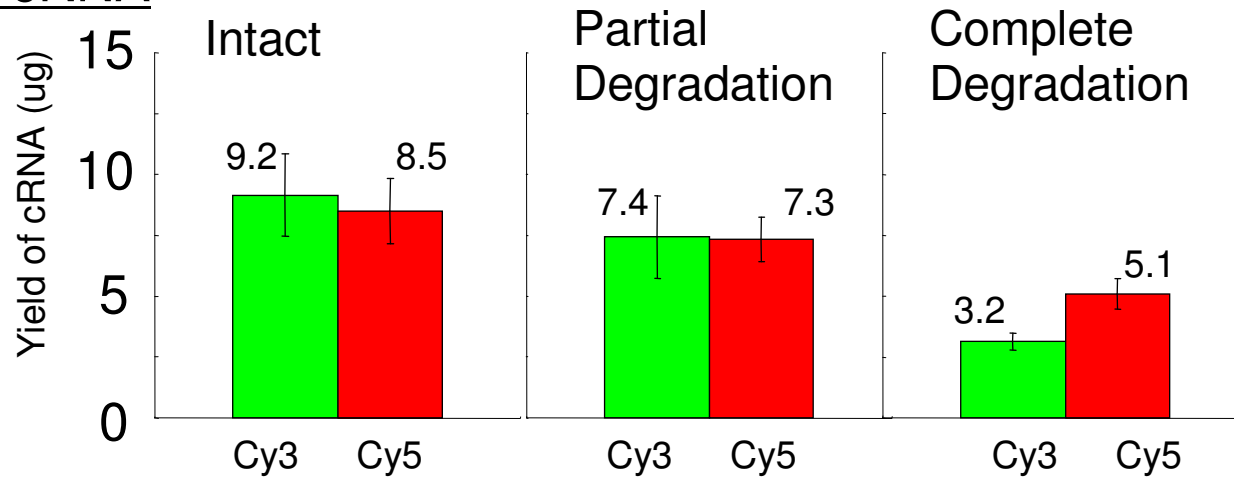


cRNA Hybridization - Workflow

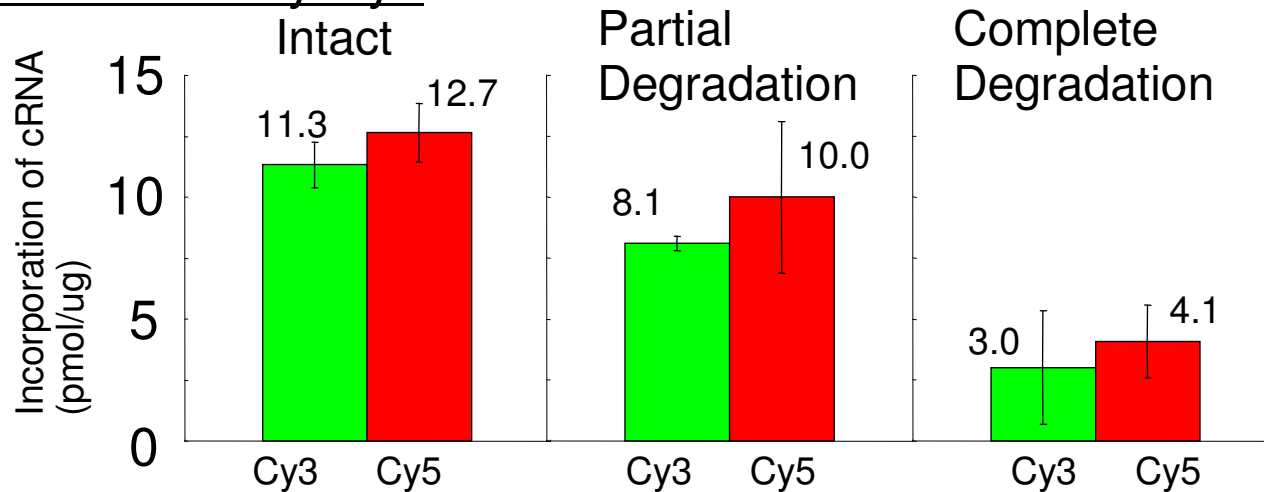


Labeled cRNA Quality Check by NanoDrop

Yield of cRNA

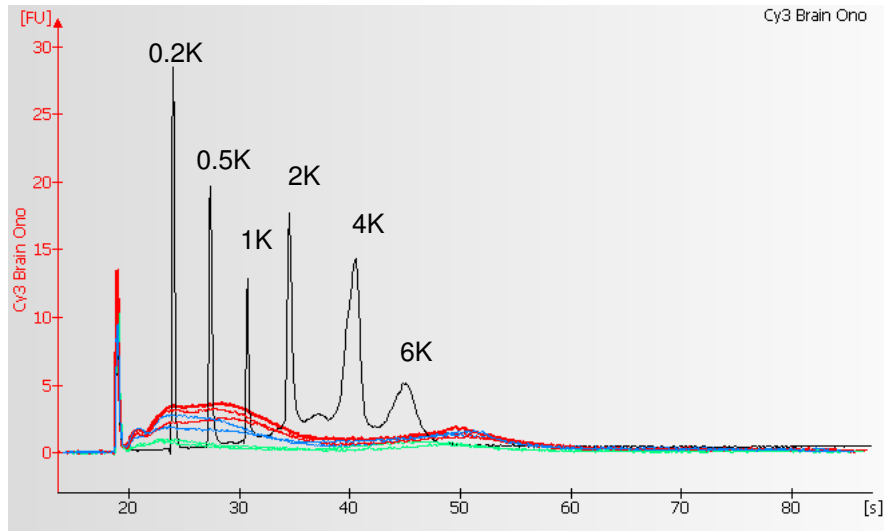


Incorporation of Cy Dye

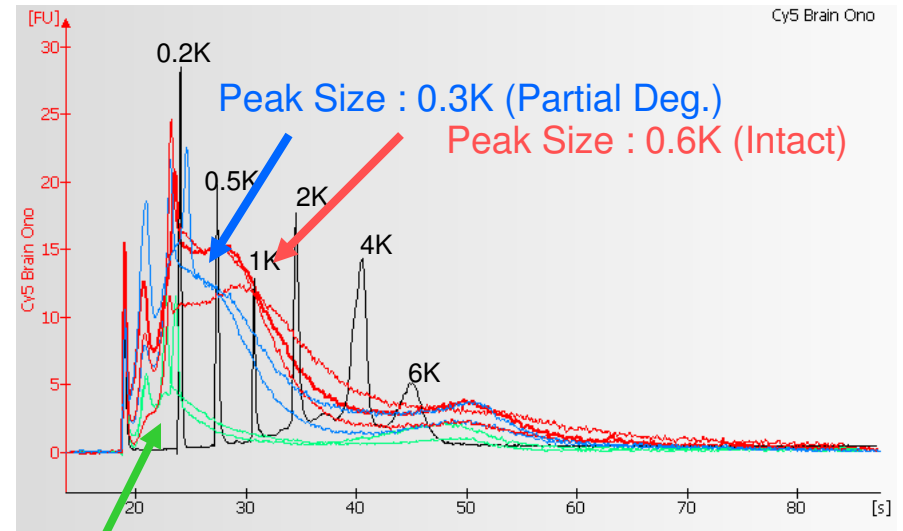


Labeled cRNA Quality Check by BioAnalyzer

Cy3 labeled cRNA



Cy5 labeled cRNA

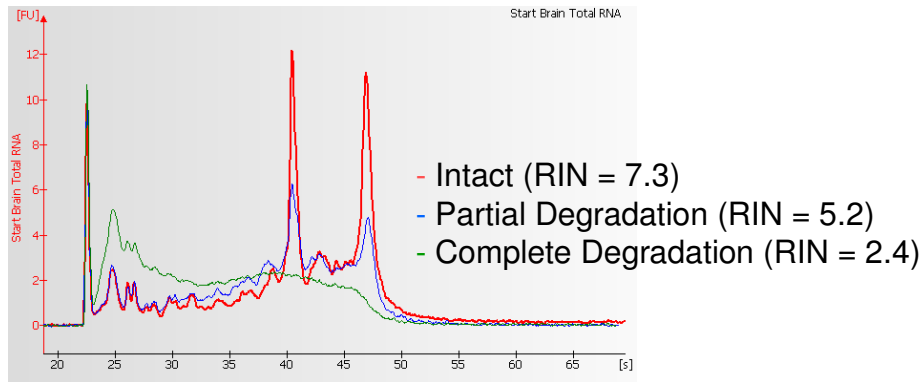


Peak Size : 0.1K (Complete Deg.)

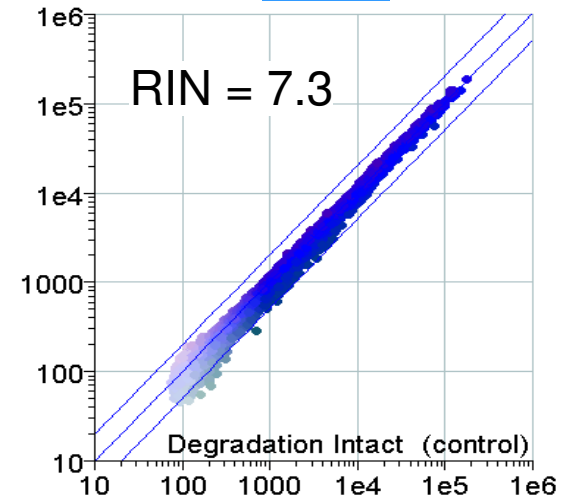
- Ladder
- Intact (RIN = 7.3)
- Partial Degradation (RIN = 5.2)
- Complete Degradation (RIN = 2.4)



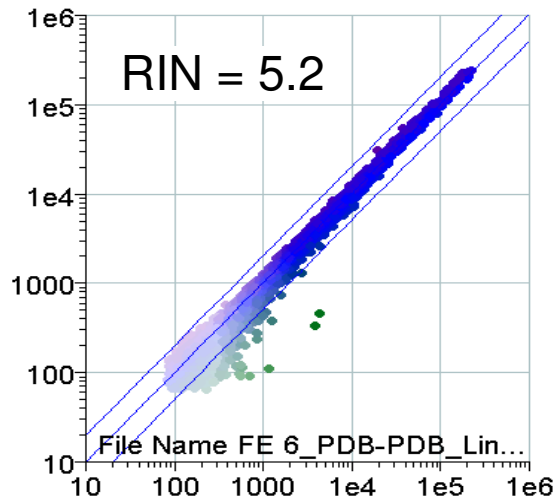
Scatter Plots : Self vs. Self



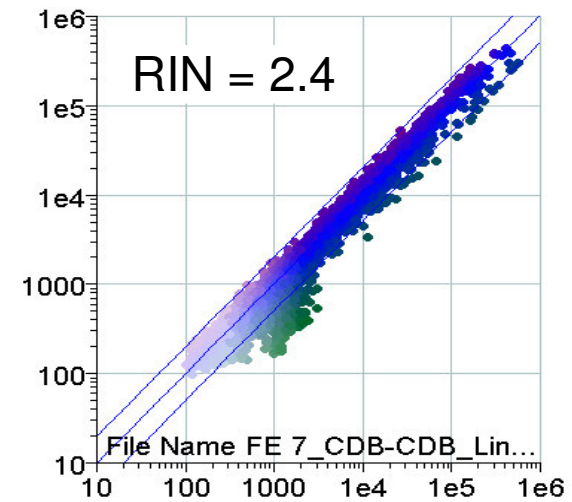
Intact



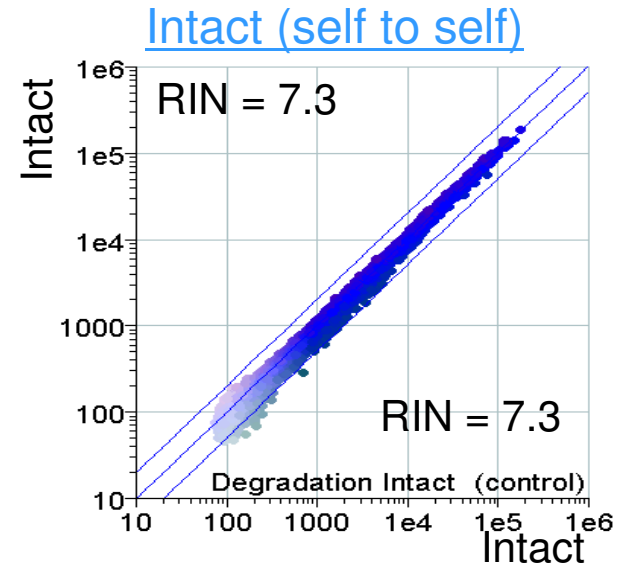
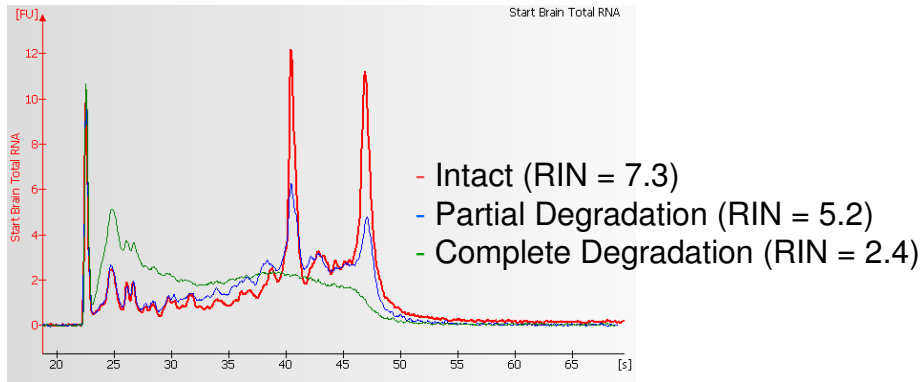
Partial Degradation



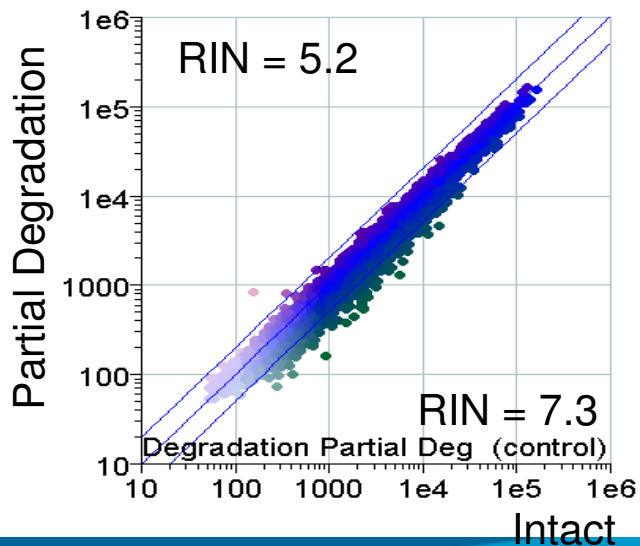
Complete Degradation



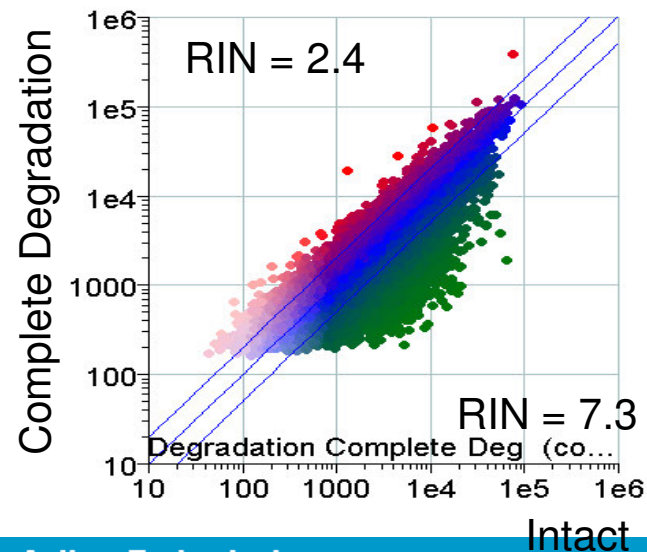
Scatter Plots : Intact vs. Degradation



Intact vs. Partial Degradation



Intact vs. Complete Degradation

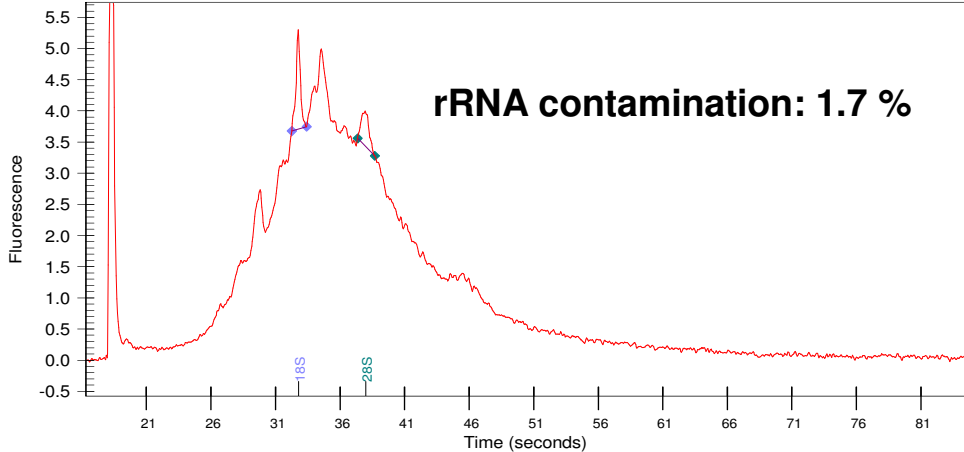
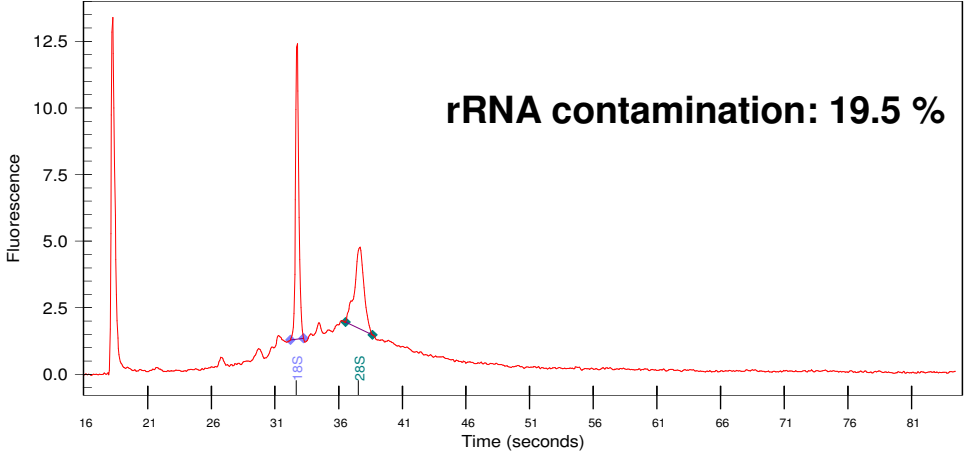
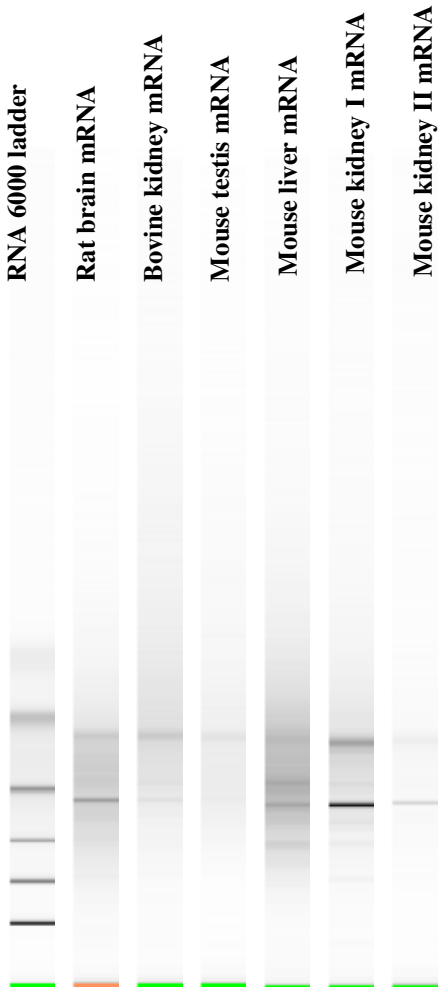


Conclusions

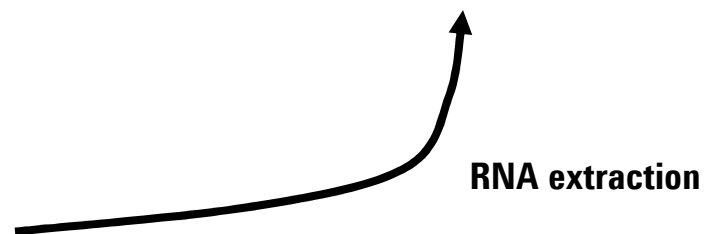
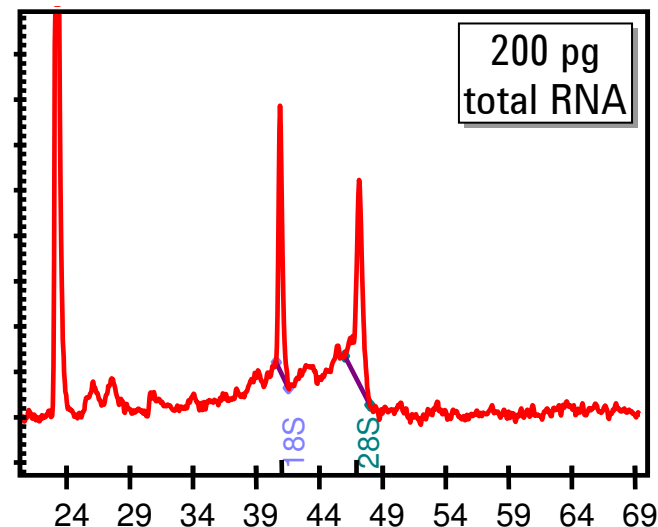
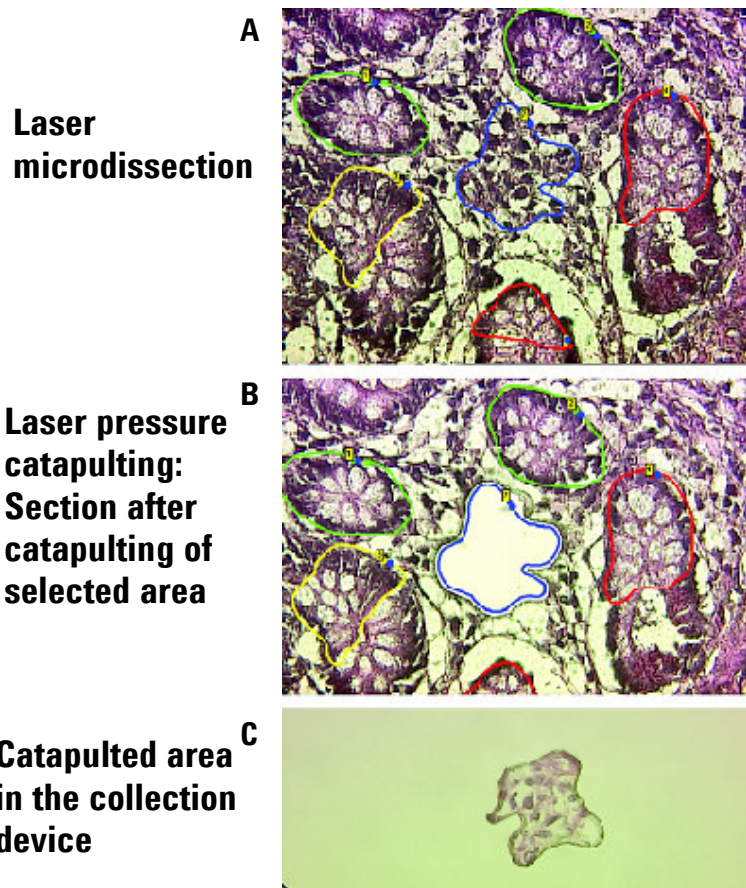
- **Labeled cRNA from different levels of RNA degradation results in low Cy Dye incorporation and low yield of cRNA**
- **RNA integrity levels of starting material had serious impact on downstream gene expression microarray results**
- **The RIN is an effective tool that can be used to evaluate RNA integrity objectively**



Ribosomal RNA contamination in mRNA samples



Laser Microdissection – PALM MicroBeam System and RNA Pico kit

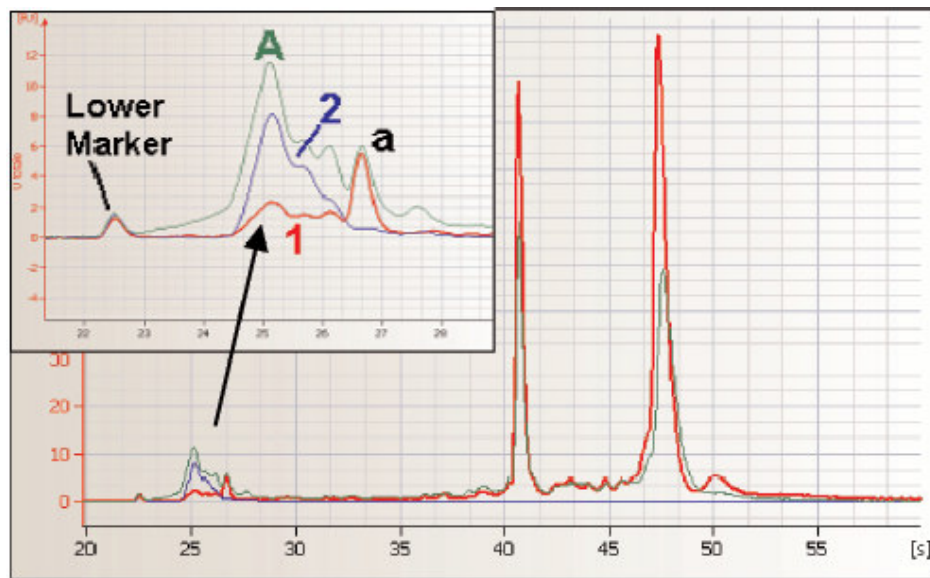
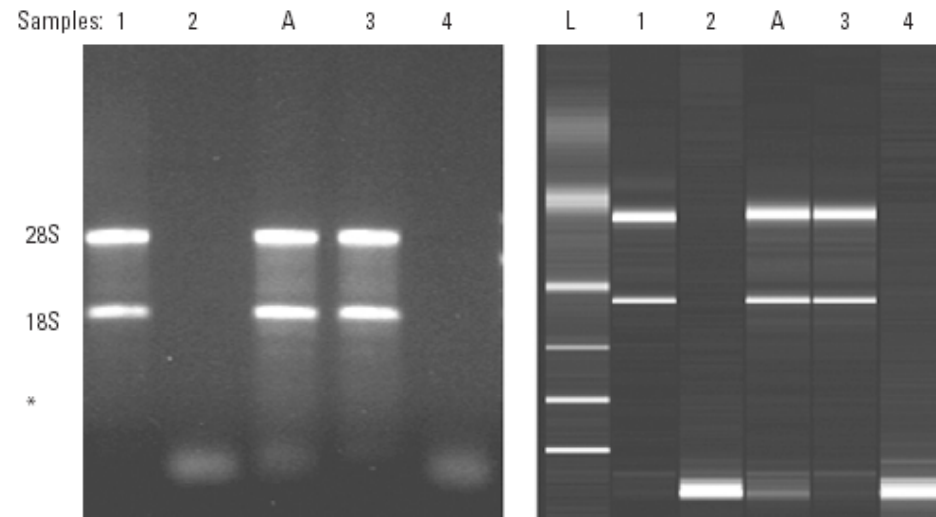


Laser Microdissection and Pressure Catapulting (LMPC)

RNA sample QC using the Agilent 2100 bioanalyzer and the RNA 6000 Pico LabChip kit

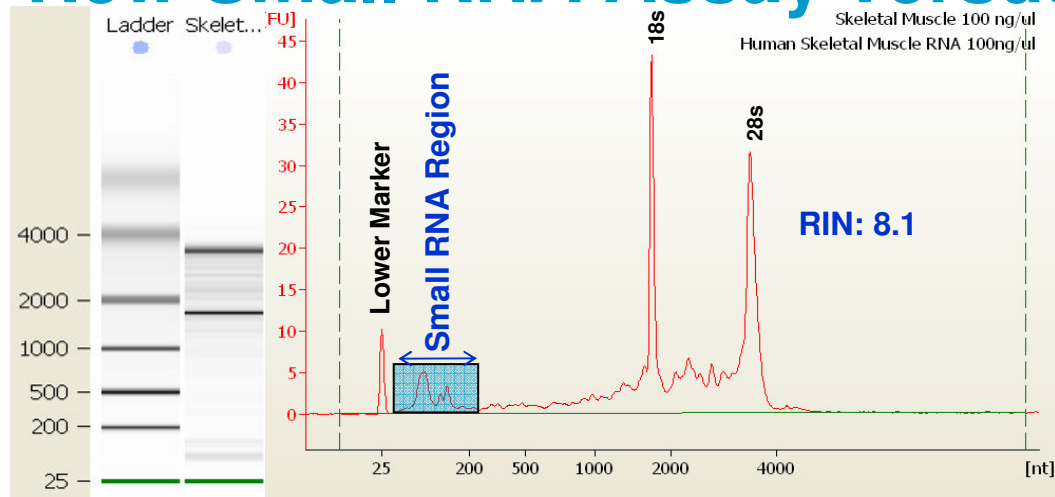
Analysis of Small RNA (using RNA 6000 Assay)

Small RNA
fraction: < 200 nts
e.g. miRNA,
siRNA, snRNA,
tRNA, 5S RNA



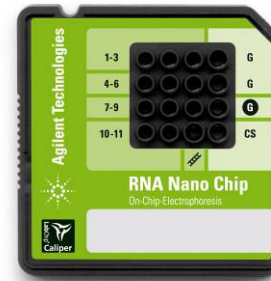
Bioanalyzer
allows
discrimination of
different profiles

New Small RNA Assay versus existing RNA Assay



RNA 6000Nano kit

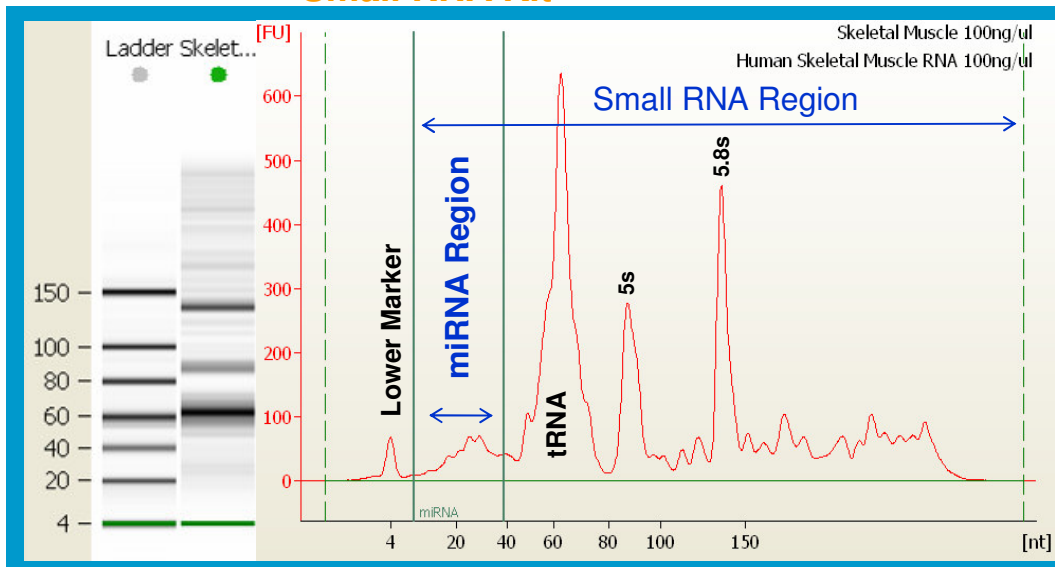
Small RNA Kit



RNA 6000Nano

Size range: 25-6000nt

Results: Integrity, Total RNA amount, gDNA contamination



NEW! Small RNA

Size range: 6-150nt

Results: miRNA amount, Ratio and amount of other Small RNA

Applications

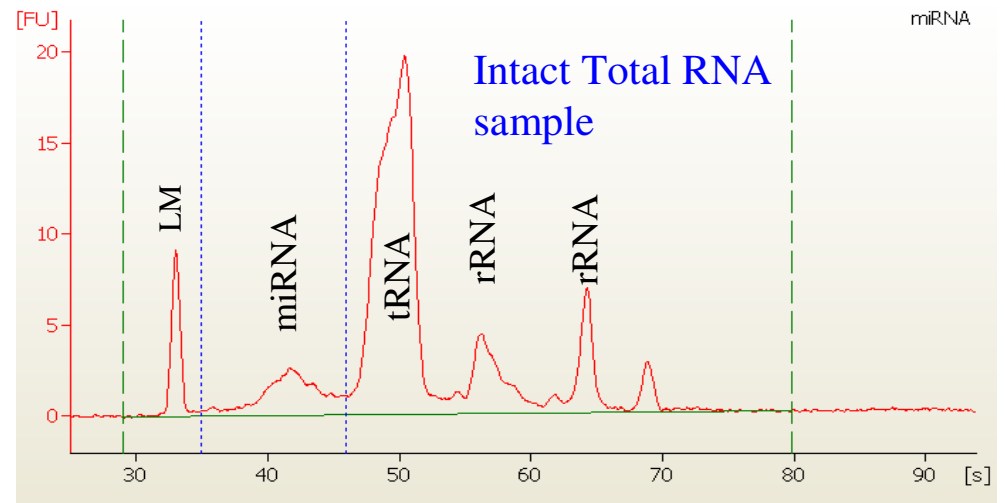
The new small RNA Assay as a tool for:

Verification, comparison and optimization in the small RNA region:

- High sensitivity to detect low abundant fragments
- High resolution for ss oligos, miRNA, pre-, t-, 5S-RNA's
- compatible with Total RNA samples or purified small RNAs.
- Semi-quantitative for single stranded RNA.
- semi- Denaturing
- Analysis up to 150nt

Plus: Qualitative assessment of dsDNA, siRNA or other hairpin RNA up to 150bp

(size separation and relative amount estimation)



Small RNA Assay specifications

Analytical Range 6 -150 nt (to avoid overlap)

Sensitivity 50 pg/ μ l

(diluted Ladder - 40 nt fragment; S/N > 3:1)

Quantitative range 50 pg/ μ l – 2000 pg/ μ l

(purified miRNA in water after extraction ~<200nt)

Quantitation Reproducibility 25 % CV

(defined on Ladder)

Max amount total RNA 100 ng/ μ l total RNA

Carryover Below detection limit

Protein Applications



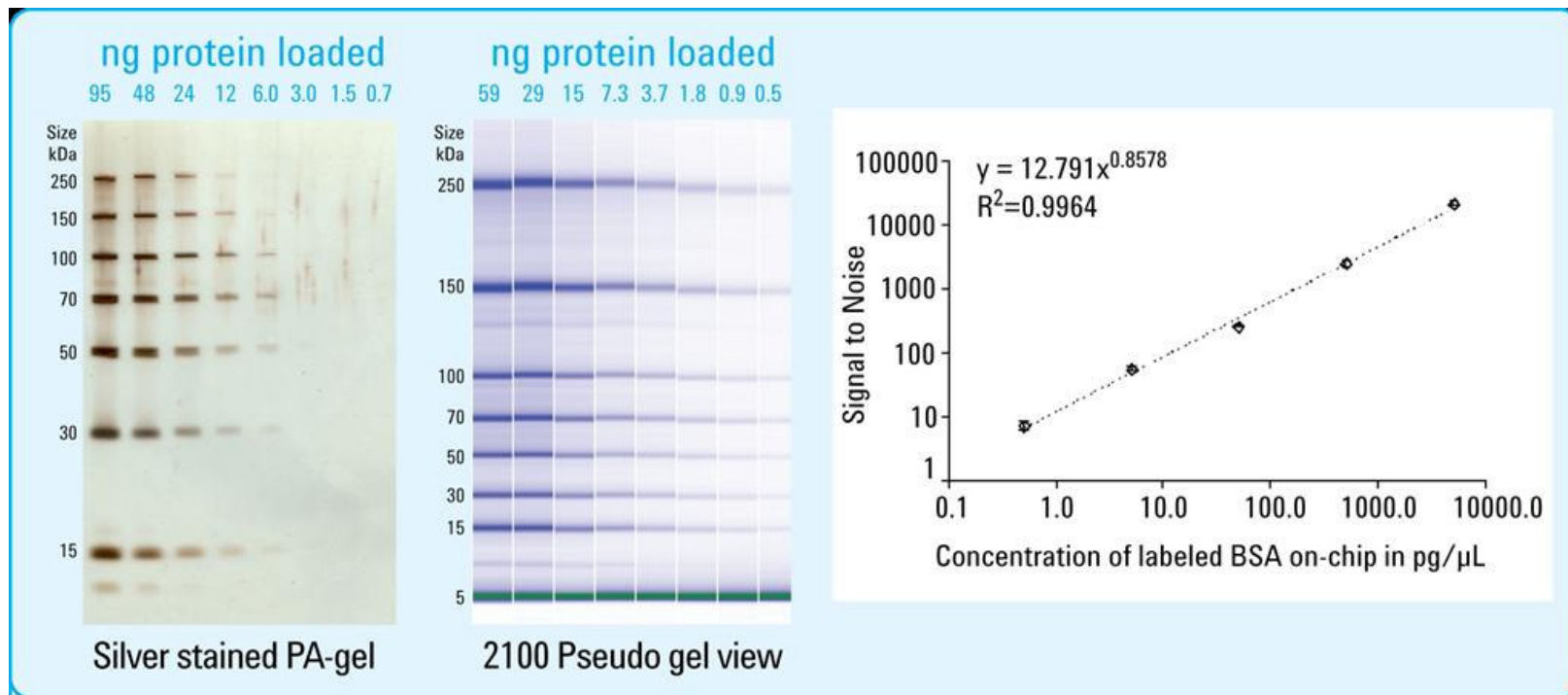
Protein Purification

Protein Expression

Protein Production

Food Analysis

Purity and QA/QC



Bioanalyzer Protein Kit portfolio

Agilent Protein 80 kit	Prod Number	5067-1515
Agilent Protein 230 kit	Prod Number	5067-1517
Agilent High Sensitivity Protein 250 kit	Prod Number	5067-1575



P 80

Range: 5 - 80 kDa
 Sensitivity: Coomassie
 Samples: 10

Samples
 -Antibodies (reduced)
 -Small Proteins

Coomassie Range (5 ng/μL BSA)



P 230

Range: 14 - 230 kDa
 Sensitivity: Coomassie
 Samples: 10

Samples
 -Antibodies (all types)
 -Standard Proteins



HSP 250

Range: 10 - 250 kDa
 Sensitivity: 1 pg/μl BSA on Chip
 Samples #: 10 per Chip
 Chips #: 10 per Kit
 Labeling Conc: 1 ng – 1 μg /μl

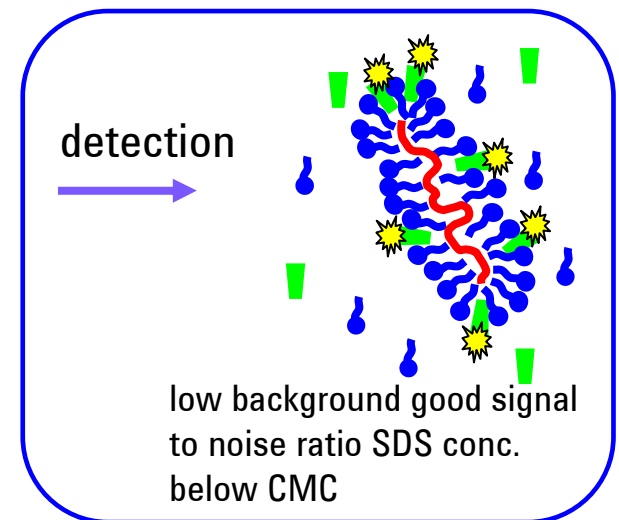
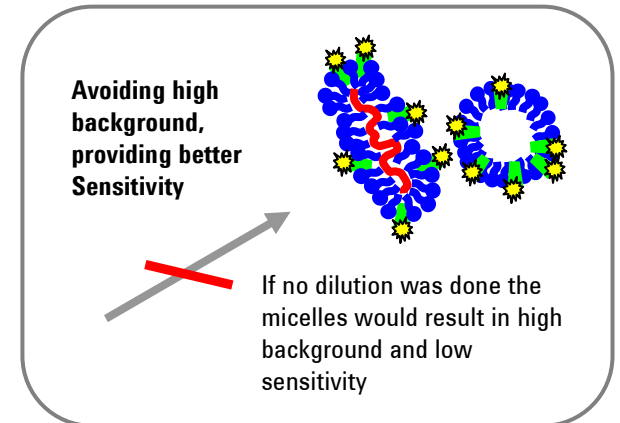
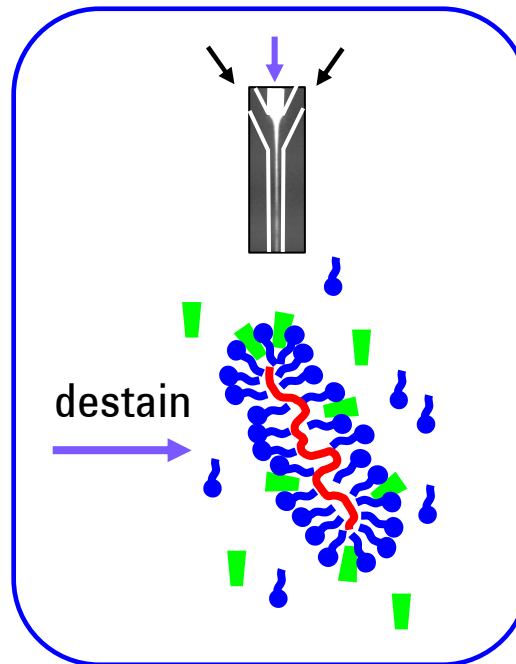
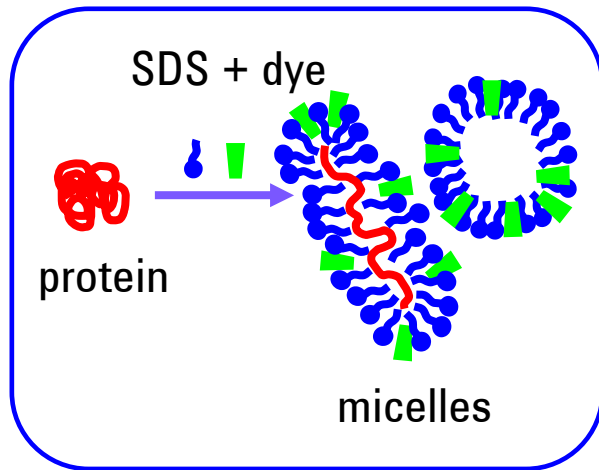
Silver stain Range (200 pg/μL BSA)

Protein Kit Specifications

	Product No. 5067-1515	Product No. 5067-1517	Product No. 5067-1575
	Protein 80 Assay	Protein 230 Assay	High Sensitivity Protein 250 Assay
Analytical specifications			
Sizing range	5-80 kDa	14-230 kDa	10-250 kDa
Typical sizing resolution	10%	10%	10%
Typical sizing accuracy	10% CV (CAII, BLG)	10% CV (BSA, CAII)	10% CV (BSA)
Sizing reproducibility	3% CV (CAII, BLG)	3% CV (BSA, CAII)	3% CV (BSA)
Sensitivity (Signal/Noise > 3)	6 ng/μL CAII (15 ng/μL BSA) in PBS, 10 ng/μL (CAII) in 0.5 M NaCl (30 ng/μL BSA in 0.5 M NaCl)	6 ng/μL CAII (15 ng/μL BSA) in PBS 30 ng/μL (BSA) in 0.5 M NaCl	1 pg/μL (labeled BSA) in water on chip 5 pg/μL (labeled BSA) in PBS on chip Labeling reaction at 1 ng/μL of total protein *
Quantitative range	60-2000 ng/μL CAII in PBS	15-2000 ng/μL CAII, 30-2000 ng/μL BSA in PBS	up to 4 orders of magnitude (0.3 to 3000 ng/μL BSA)
Qualitative range	6-4000 ng/μL CAII and BLG	6-5000 ng/μL CAII, 15-5000 ng/μL BSA in PBS	-
Quantitation reproducibility	20% CV (CAII, BLG)	20% CV (BSA, CAII)	20% CV (BSA)
Physical specifications			
Analysis run time	30 minutes	25 minutes	30 minutes
Number of samples	10 samples/chip	10 samples/chip	10 samples/chip
Sample volume	4 μL	4 μL	5 μL
Kit stability	4 months (for storage temperature see individual box)	4 months (for storage temperature see individual box)	6 months at -20° C
Compatible buffers	List of compatible buffers	List of compatible buffers	List of compatible buffers
CAII = Carbonic Anhydrase, BSA = Bovine Serum Albumin, BLG = beta-Lactoglobulin			

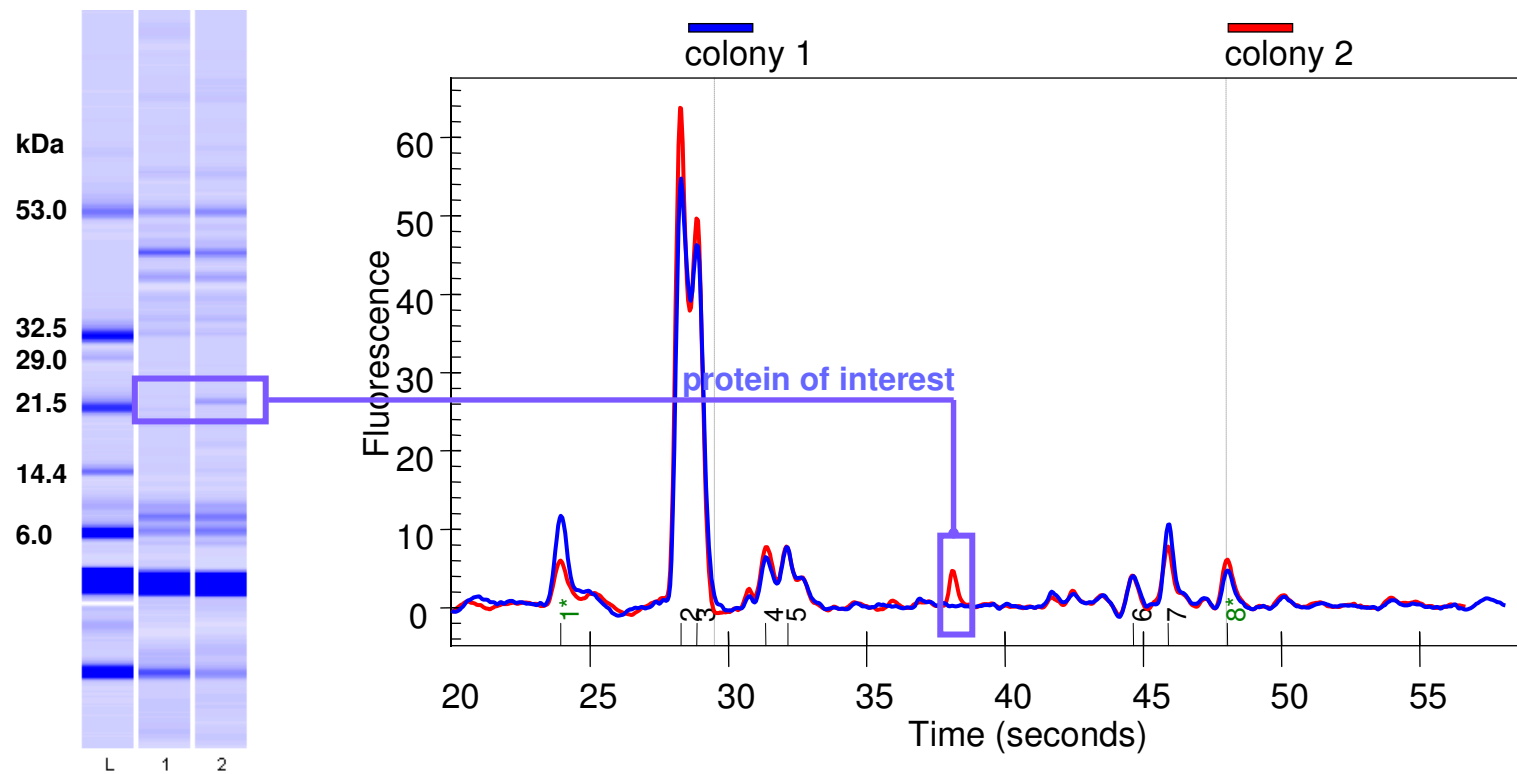
* Prior to measurement on Chip we recommend within the High Sensitivity Protein 250 labeling protocol to dilute the labeled sample by a factor of 200.

Staining, Destaining and Detection (P-80 and P-230)



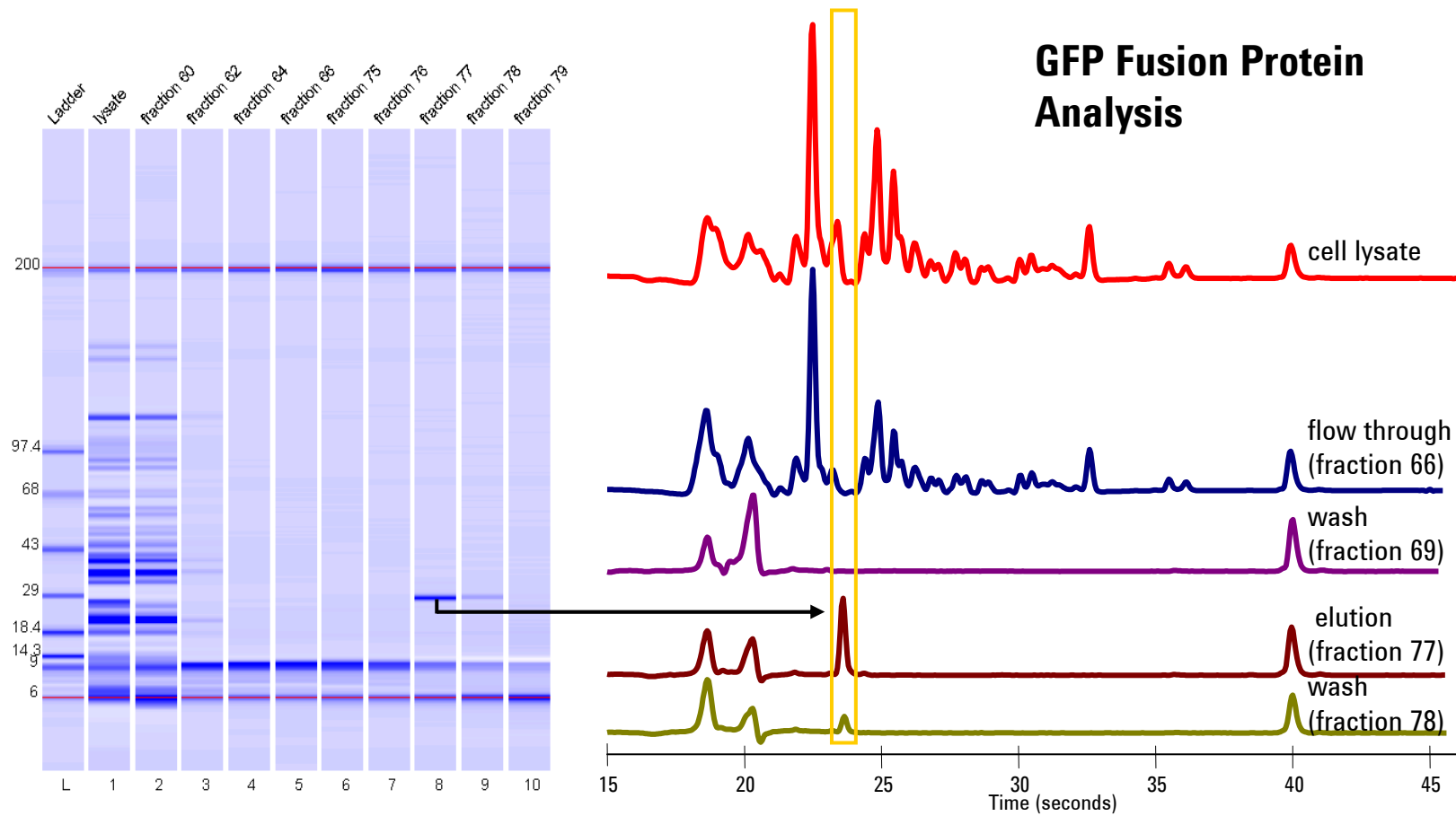
Clone Selection based on Protein Expression

Example measured with Protein 50 kit



Monitoring of Protein Purification Process

Example measured with Protein 200 plus kit



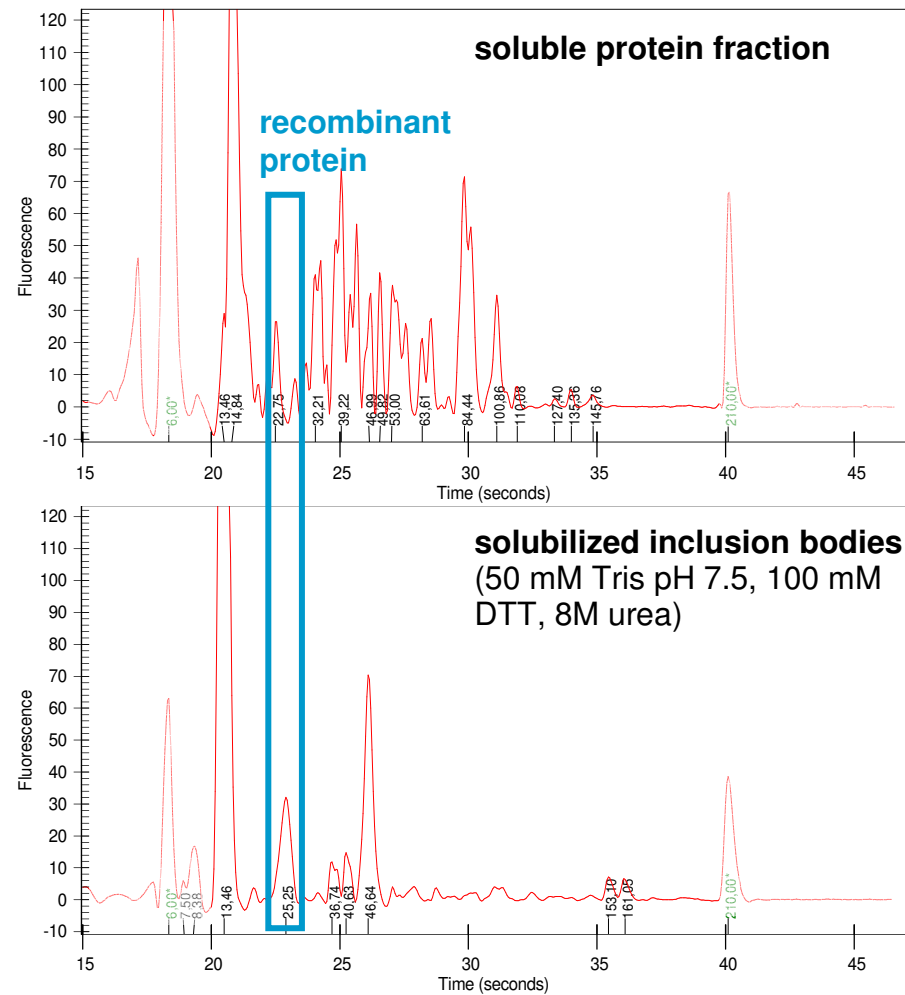
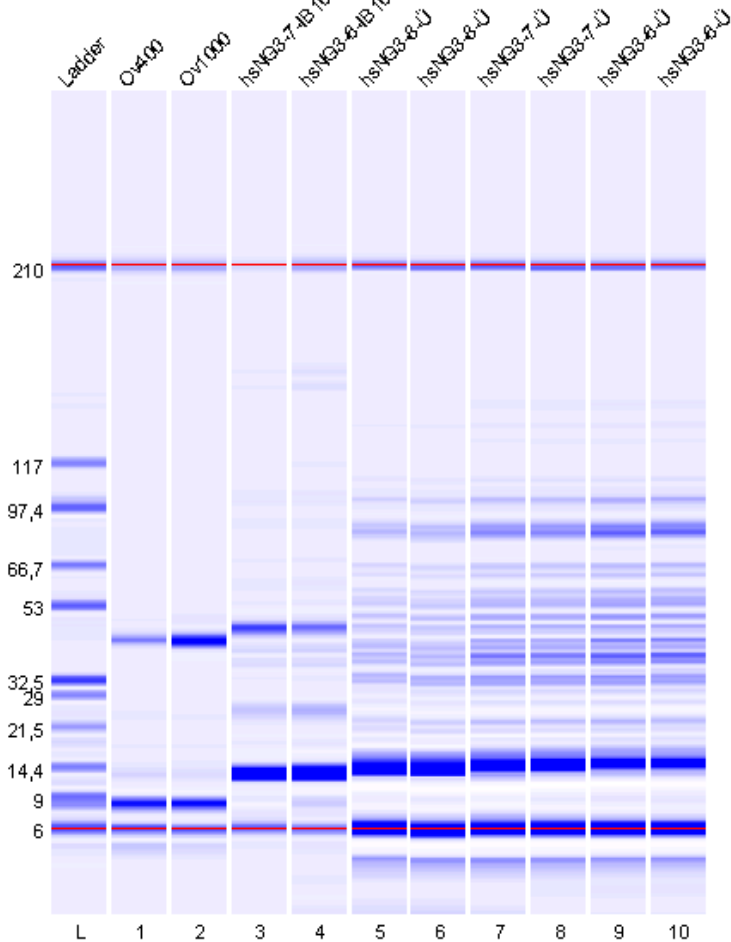
↑ 2100 bioanalyzer: gel-like image

↑ 2100 bioanalyzer: electropherogram

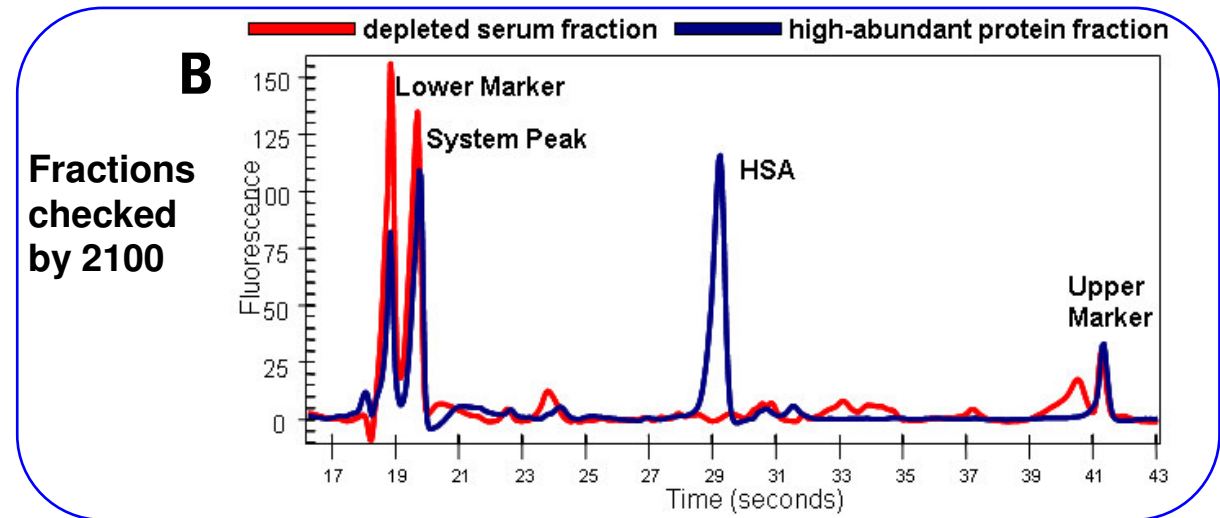
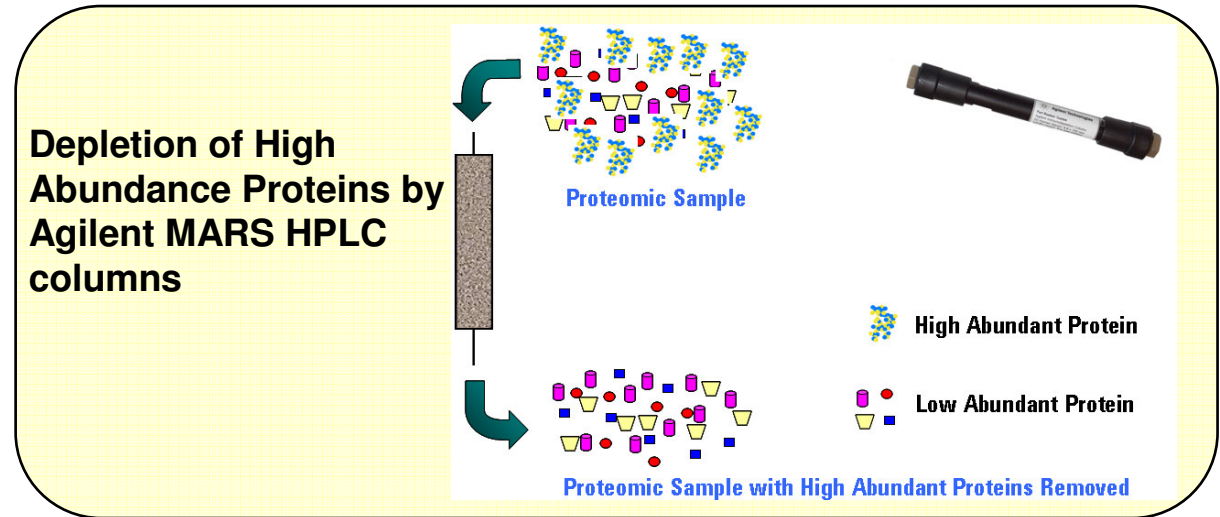
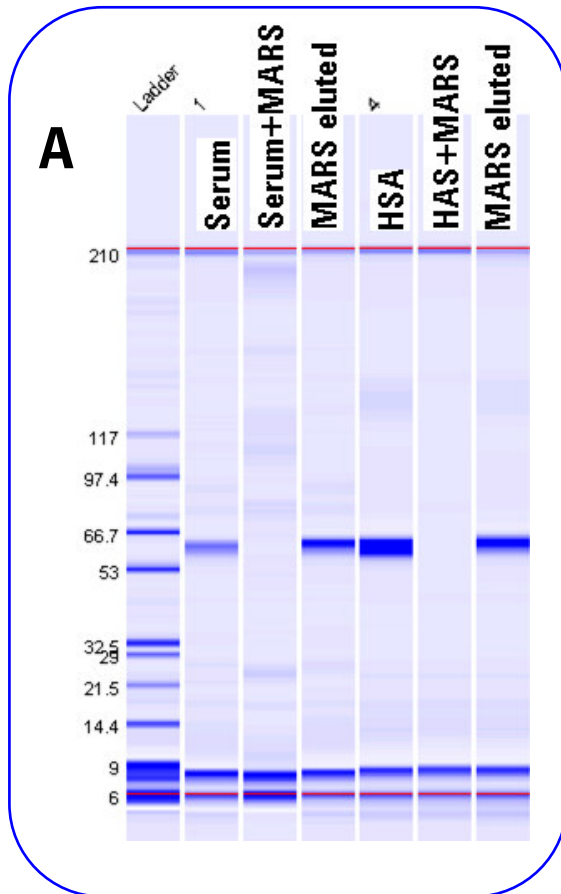
Courtesy of P. Sebastian and S.R. Schmidt
GPC-Biotech AG, Martinsried, Germany

Expression of a Recombinant Protein in *E.coli*

- Optimization of Fermentation and Induction Conditions

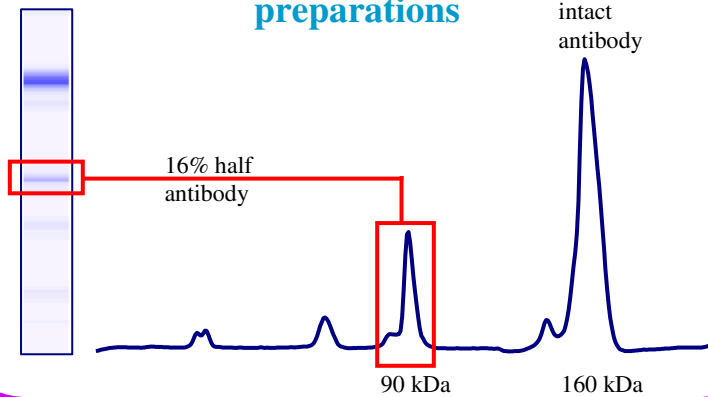


Quality control of the Depletion of High Abundance Proteins in Human Serum

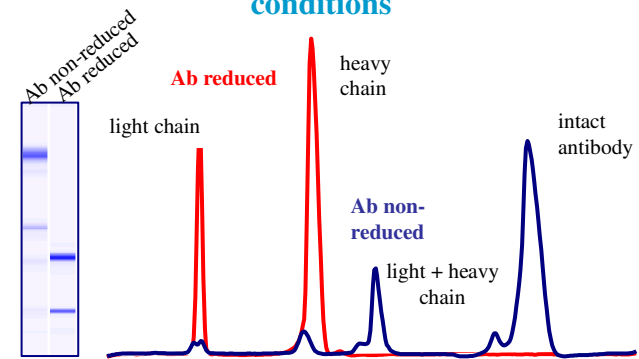


Quality Control of Antibodies

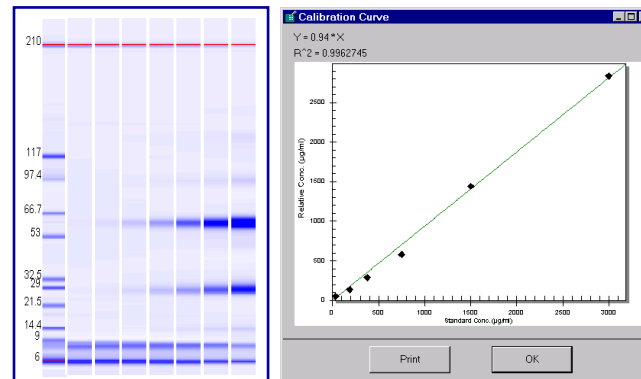
Determine the half antibody content in IgG preparations



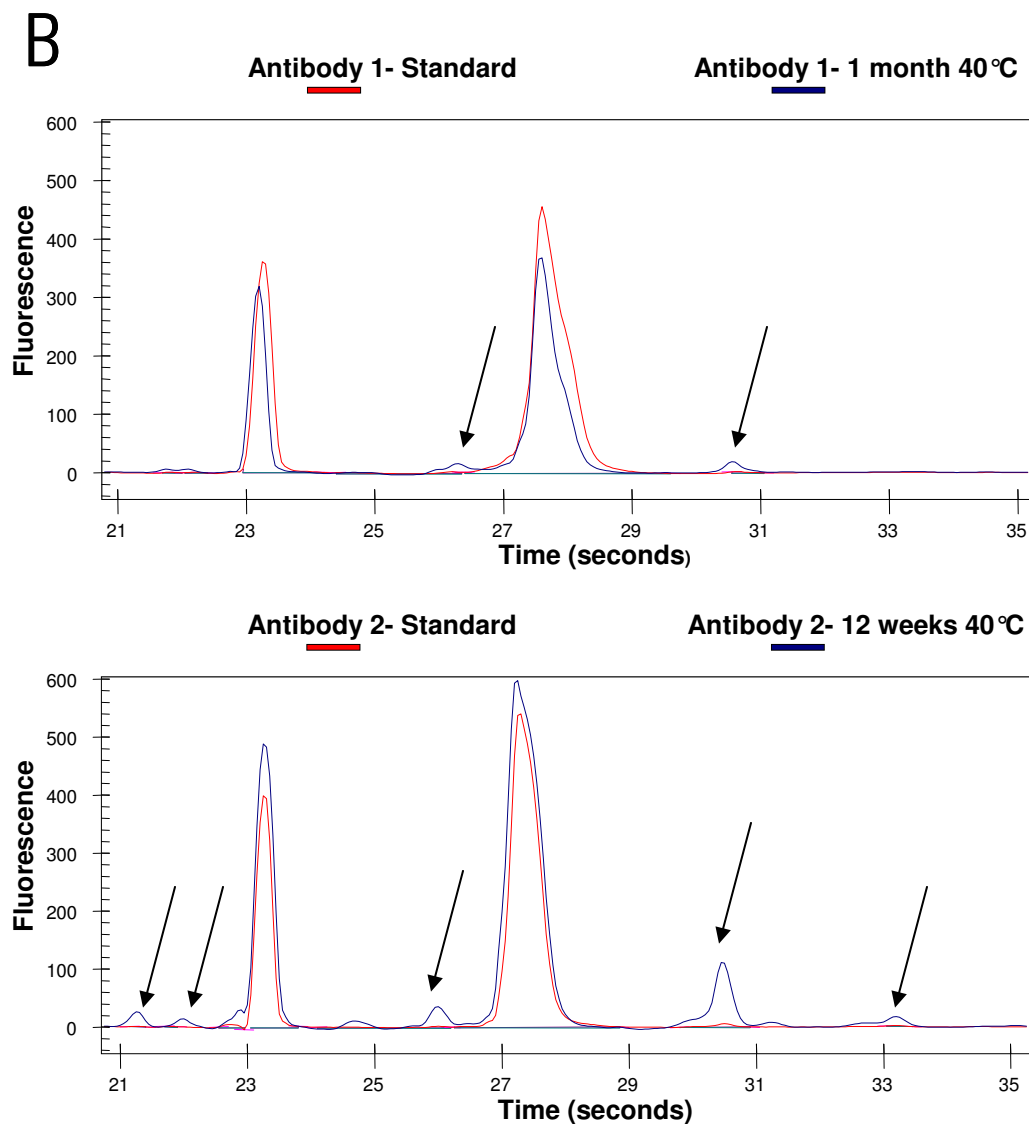
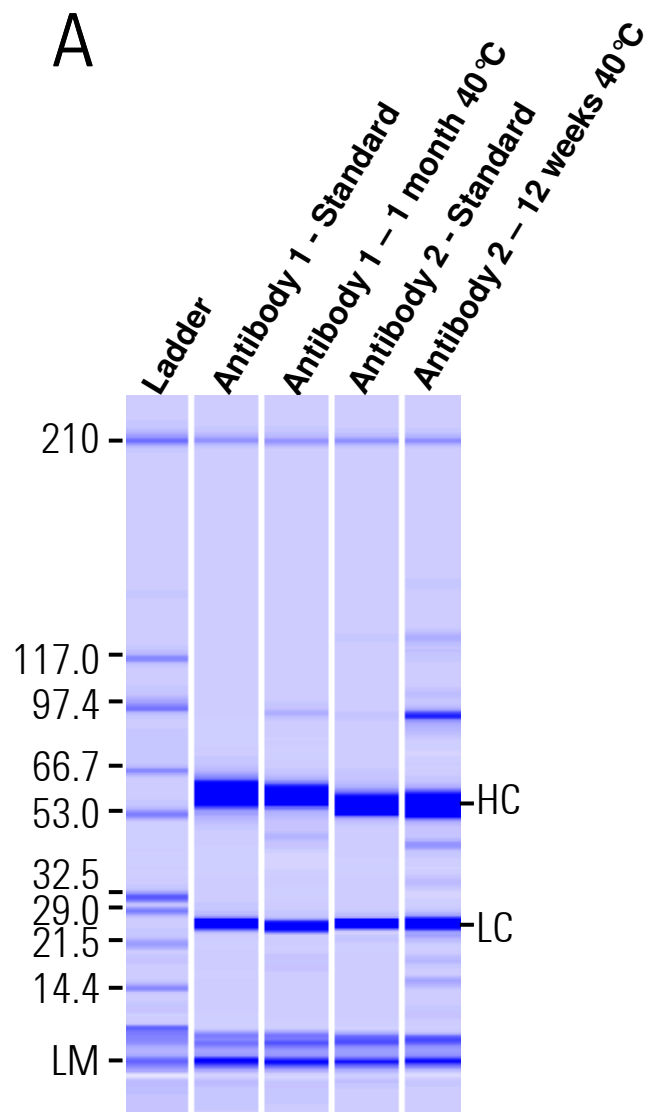
Antibody analysis under reducing and non-reducing conditions



Absolute Quantitation of IgG samples

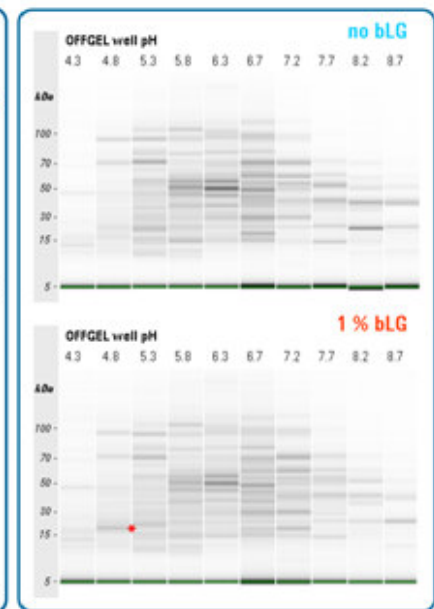
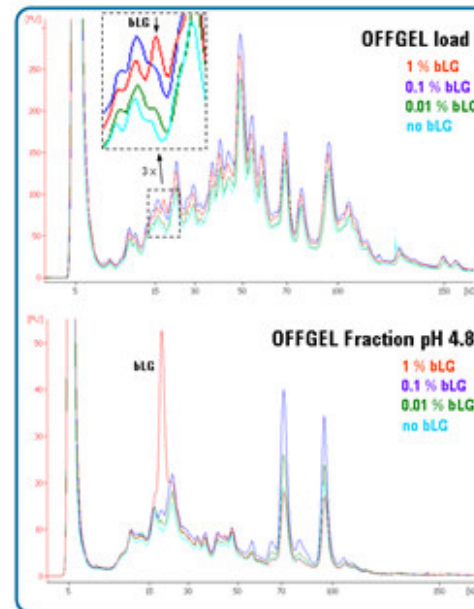
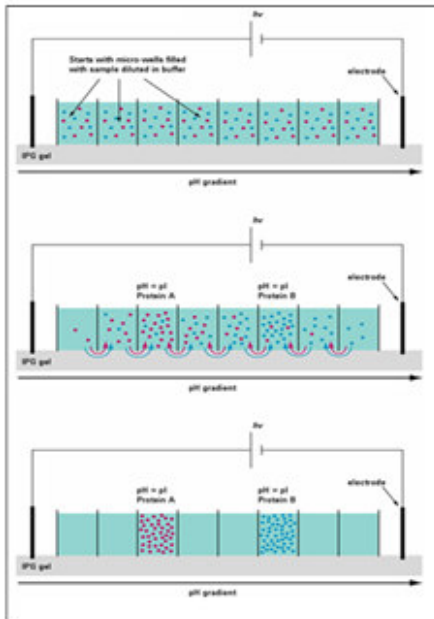
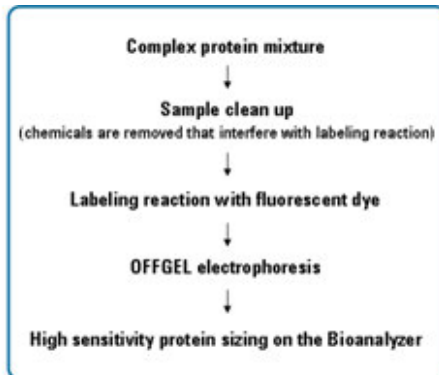


Analysis of Antibody Stability – stress test



Combination of IEF with SDS-PAGE

Agilent 3100 OFFGel Fractionator + 2100 bioanalyzer



Description of the new HSP-250 Assay (Direct labeling reaction, silver stain sensitivity)

- Reach and beat traditional „silver stain sensitivity“
- Offer solid quantitation for a large dynamic range

Target Applications:

- Protein QA/QC
reliable quantitation of main compound besides minor impurities
- Protein detection at lowest concentrations in research

High Sensitivity Protein 250 Kit 5067-1575 content is:

- 10 Chips (100 samples)
- Labeling Kit (Dye and Reagents)
- 2100 Separation Kit (Gel, Marker, Ladder, Buffer)
- User Documentation (Quick Start Guide & Labeling Protocol)

Extended experimental workflow

Sample

5-90 min

Transfer to suitable buffer
(precipitation, ultrafiltration, buffer exchange spin columns)

40 min

Labeling with dye
(N-hydroxy-succinimidyl ester chemistry)

labeling kit

5 min

Sample preparation for 2100
(SDS denaturation, dilution if desired)

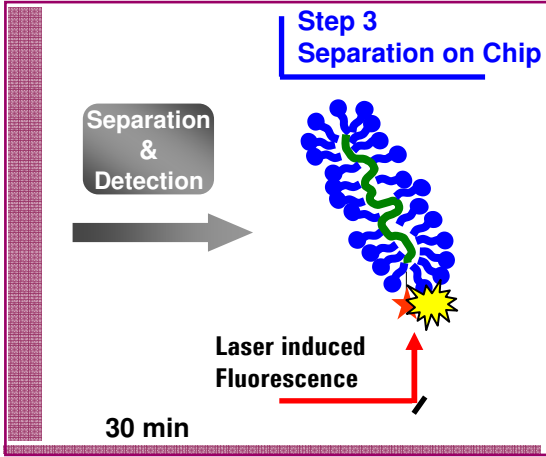
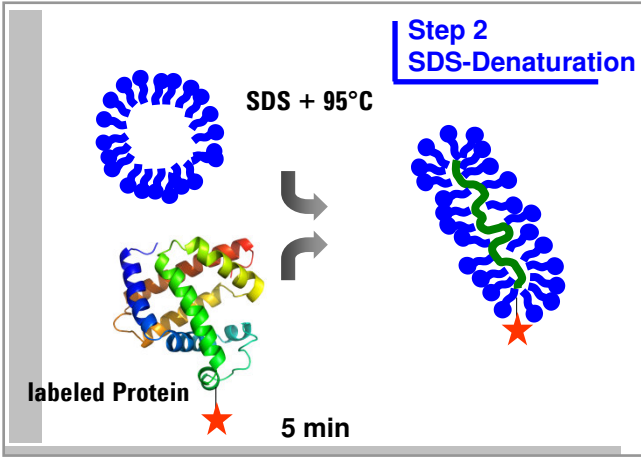
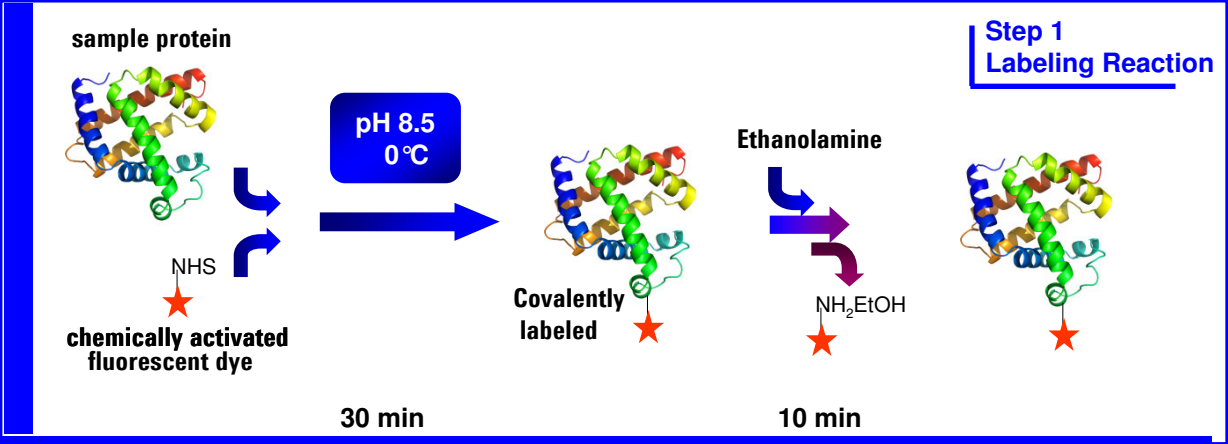
separation kit

35 min

Analysis on 2100

Data

Principle of High Sensitivity 250 Protein Staining

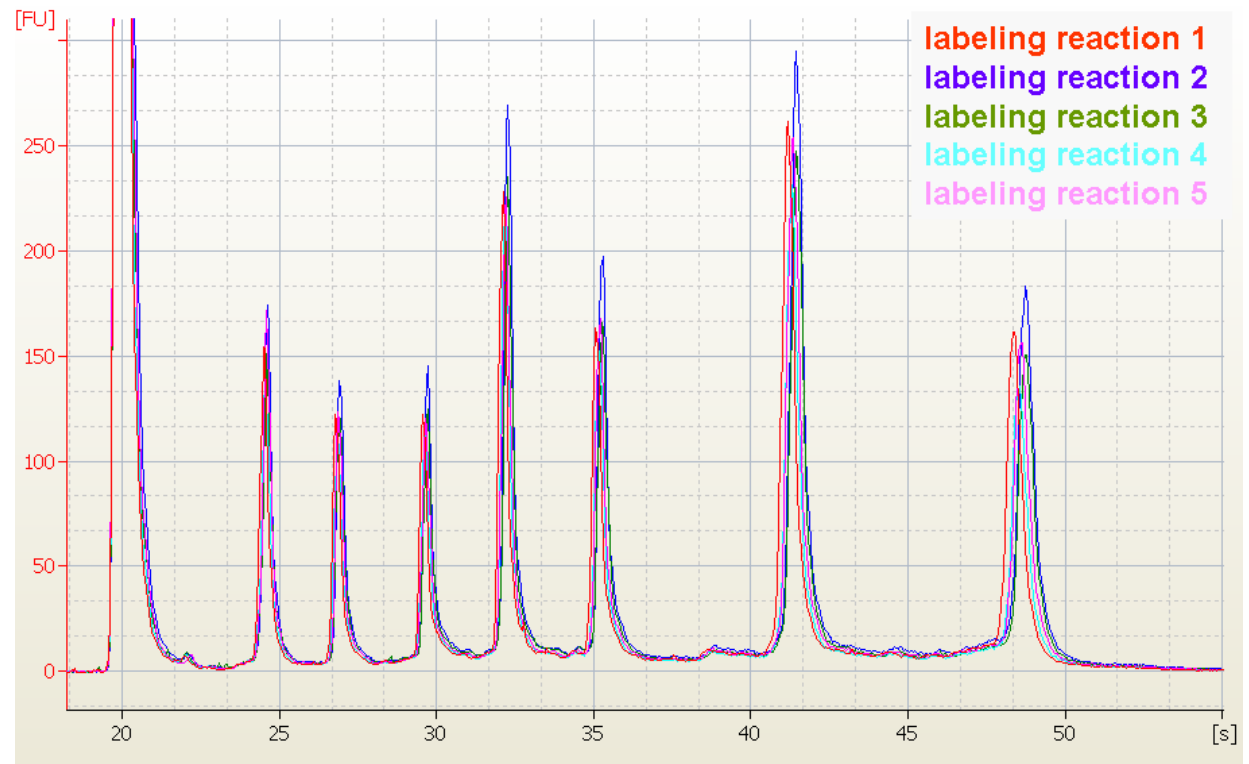


Reproducibility of Labeling Reaction: Ladder

Rugged Labeling reaction:

Reproducible reaction provides comparable signal intensities.
Homogenous labeling without extra bandbroadening

No deviation in peak width is indicating a constant number of dye per protein molecule and proves a stable protocol

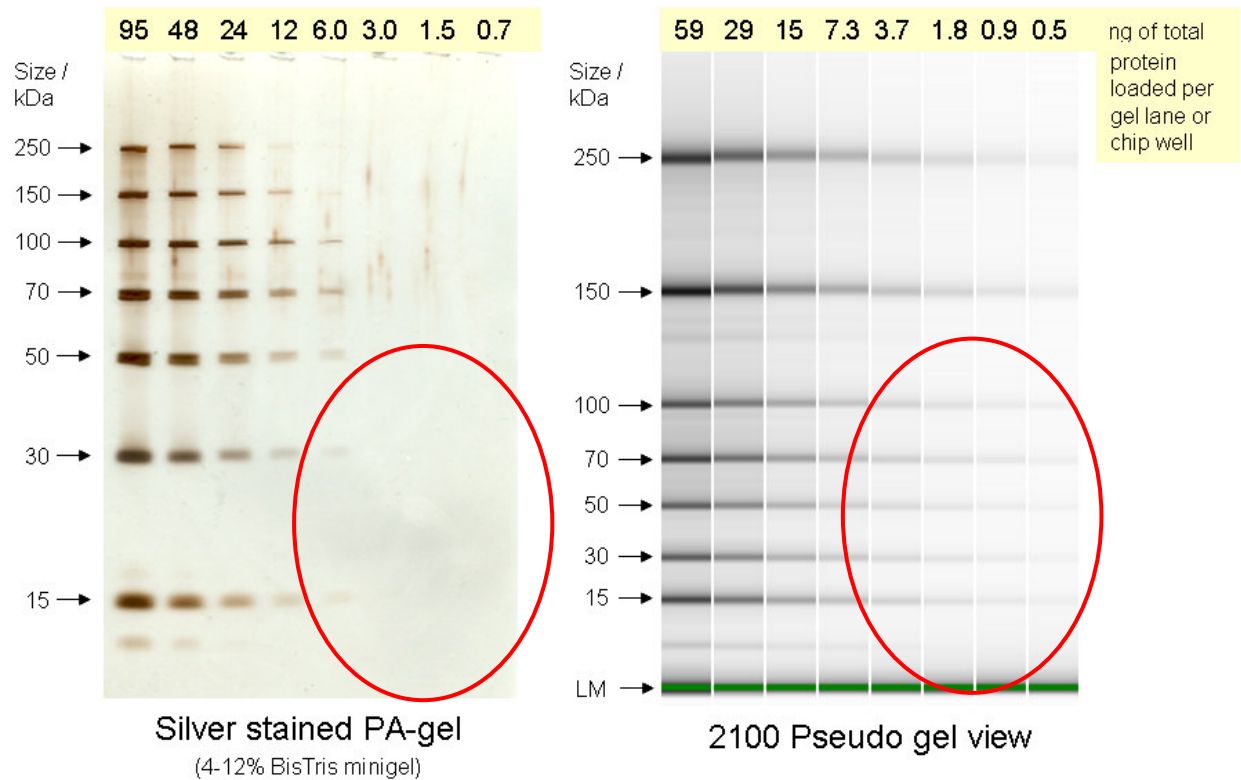


Sensitivity: Silver Staining vs. Bioanalyzer

Highest sensitivity:
Labeled proteins can be measured down to pg/ μ L concentrations loaded on Chip

Direct comparison of samples run on SDS-PAGE with Silver staining and on 2100 Bioanalyzer.

Concentrations are given per lane (as total concentration of 7 different proteins)



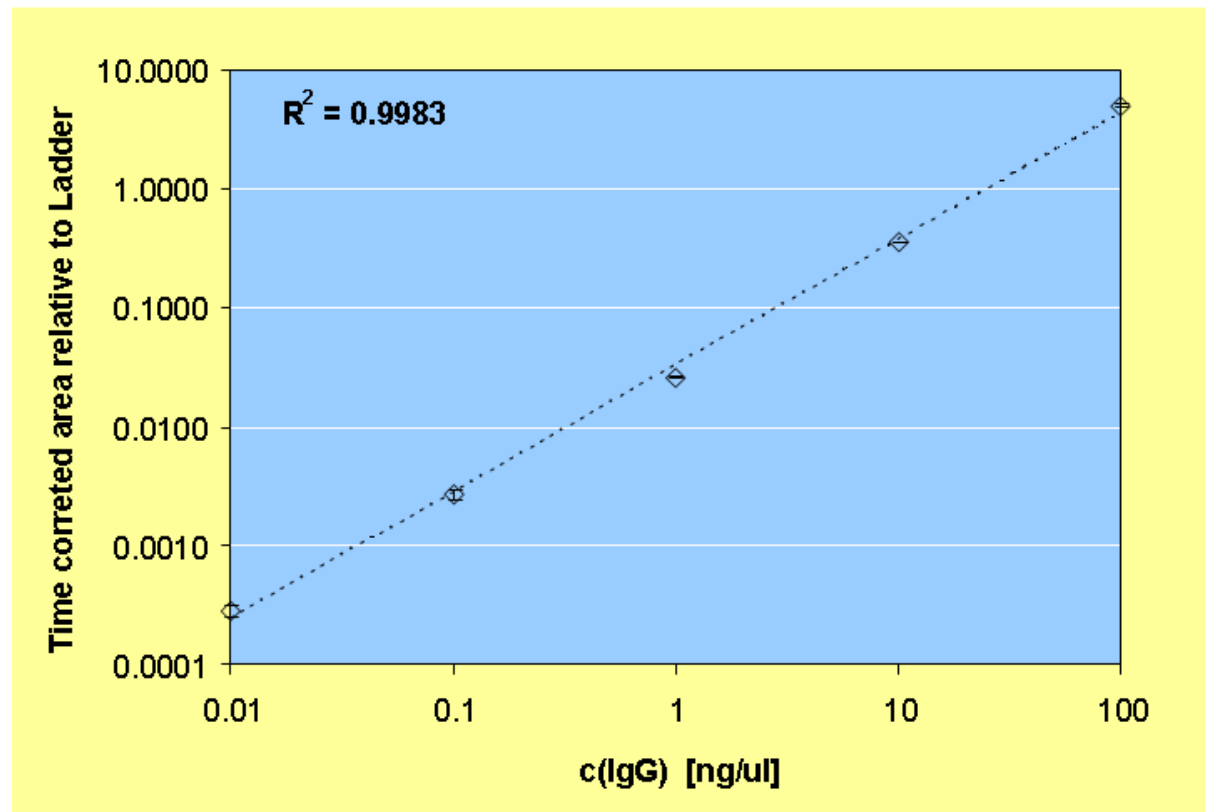
Linear Dynamic Range Test: IgG

Linear dynamic range:

Quantification of labeled IgG from 10 pg/ μ L to 100 ng/ μ L

averages \pm SD of 7 measurements (7 chips, 1 chip lots, 4 instruments)

4 orders of linear dynamic range allows to quantify an 0.05% impurity besides the main peak in a single run



2100 Bioanalyzer Compliance



2100 expert software

- One version for all assays
- Declaration of system validation



2100 expert security pack

- 21 CFR part 11 compliance
- Electronic records
- Electronic signatures
- Audit trails

2100 bioanalyzer

- IQ and OQ/PV services
- Declaration of conformity

Chips and reagents

- Declaration of conformity

Agilent Web pages with 2100 content

www.agilent.com/chem/labonachip

Agilent | Lab-on-a-Chip Products - Microsoft Internet Explorer provided by Agilent Technologies, Inc.

File Edit View Favorites Tools Help

Address <http://www.chem.agilent.com/Scripts/PCol.asp?Page=50>

Links Google Agilent Agilent Lab-on-a-Chip Products Frontier ePal Service Life Sciences Marketing Welcome ELSA

Agilent Technologies Our measure is your success.

Select a Country or Area **Contact Us**

Products & Services Technical Support Industries Buy About Agilent Search Register Login

Lab-on-a-Chip Products

Cutting edge Lab-on-a-Chip Products

Agilent Technologies is the leader in commercial microfluidic Lab-on-a-Chip technology. This technology utilizes a network of channels and wells that are etched onto glass or polymer chips to build mini-labs. Pressure or electrokinetic forces move pico liter volumes in finely controlled manner through the channels. Lab-on-a-Chip enables sample handling, mixing, dilution, electrophoresis and chromatographic separation, staining and detection on single integrated systems. The main advantages of Lab-on-a-Chip are ease-of-use, speed of analysis, low sample and reagent consumption and high reproducibility due to standardization and automation.

Buy

Request a quote
Where to buy
Store Home

Related Information

- Literature Library
- Applications
- Manuals
- more...
- Technical Support
- Frequently Asked Questions
- Technical Document Search
- RNA Integrity Database
- Downloads & Utilities
- 2100 Bioanalyzer User Forum
- Parts Information
- more...
- Education & Events
- Classroom Training Courses
- Customized On-Site Training
- e-Seminars
- more...

2100 Bioanalyzer

- 2100 Bioanalyzer
- 2100 Bioanalyzer Series II Kits
- 2100 Expert Software
- 2100 Expert Security Pack
- Compliance

1200 HPLC-Chip System

DNA Solutions

Protein Solutions

Cell Solutions

Services

Announcements

Are You Still Up-to-date?
Upgrade your Agilent 2100 Expert Software - and Save!

Order Now & Get 30% Discount!

Product Announcement
New Small RNA Assay for analysis of miRNA, siRNA and other oligonucleotides within 6-150 nt range

Special Offer
Save 30% on New Revision B.02.05 of 2100 Expert Software

Upcoming Events
2100 Bioanalyzer User Meetings in four US cities, Sept-Oct 2007

Other News
Issue 20 of 'Pharmaceutical Analysis' now available!

Open Genomics Home - Microsoft Internet Explorer provided by Agilent Technologies, Inc.

File Edit View Favorites Tools Help

Address <http://www.opengnomics.com/>

Links Google Agilent Agilent Lab-on-a-Chip Products Frontier ePal Service Life Sciences Marketing Welcome ELSA

The Latest in Microarray-Based Genomics Research

Presentations:

- Transcriptional Regulation of Embryonic Genomes
- Dr. Richard Young, Whitehead Institute for Biomedical Research, Massachusetts Institute of Technology
- Watch the Presentation Now
- Review the Presentation

Interviews:

- Hidden Problems - Uncovering Microdeletions Behind Uncovering Developmental Abnormalities
- An Interview with Dr. Cinzia Zuffanti, University of Pavia, Italy
- Listen to This Interview Now
- Review all Interviews

Research Reviews:

- Identifying New Combination Drug Classes for Hypertension and Diabetes
- A review of Molecular characterization of new selective neurotensin/proliferator-activated receptor (gammal) modulators with angiotensin receptor blocking activity
- Read More
- Using aCGH to Expand Your Cytogenetic Research
- A review of High-resolution global profiling of genomic alterations with long oligonucleotide microarray
- Read More

Publications Search

Enter keyword Search

Recent Publications

- Whole genome CHIP-chip analysis of Drosophila Twist and Giant suggests integration of diverse patterning processes in the Drosophila embryo
- Read More
- Direct and sensitive mRNA profiling from low-input total RNA
- Read More
- Amplified, lost, and fused genes in 11q23-25 simonin in acute myeloid leukemia, an array-CGH study
- Read More

News & Noteworthy

Agilent Technologies Offers First FFPE Kit for Oligo CGH Microarrays

Read More

Explore the power of CGH Analytics for array CGH data capture and interpretation

Watch the latest educational tutorial now

Visit the new bioanalyzer 2100 RNA Integrity Database to screen and validate your results based upon a multitude of annotated RNA profiles submitted by your peers

Access the RNA Integrity Database now

See our seamless workflow in action, from optimal bench set-up to array loading, hybridization, washing, and scanning

Watch an Agilent array experiment now

Current Promotions

Limited Time Offer on Human mRNA Microarrays

Get a 25% discount on human mRNA arrays and corresponding labeling/hybridization kits

Learn More

Introductory Pricing on the New Oligo aCGH FFPE Labeling Kit plus aCGH Array

Save 25% off your total order.

- Custom
- Custom

Subscribe to the Open Genomics Newsletter

Stay current with the latest in microarray-based genomics research.

Learn More

www.opengnomics.com

2100 kits for Protein applications

Protein Kits – Coomassie stain sensitivity

Number	Kit	Max # of samples
5067-1517	Agilent Protein 230 Kit	250
5067-1518	Agilent Protein 230 Reagents	
5067-1515	Agilent Protein 80 Kit	250
5067-1516	Agilent Protein 80 Reagents	

Protein Kits – Silver stain Sensitivity

Number	Kit	Max # of samples
5067-1575	High sensitivity Protein 250 Kit	100
5067-1576	High sensitivity Protein 250 Reagents	
5067-1577	High sensitivity Protein 250 Labeling Reagents	
5067-1578	High sensitivity Protein 250 Ladder	

2100 kits for Cell Assay and DNA applications

Cell Fluorescence Kits

Number	Kit	Max # of samples
5067-1519	Agilent Cell Kit	150
5067-1520	Cell Checkout Kit	

DNA Kits

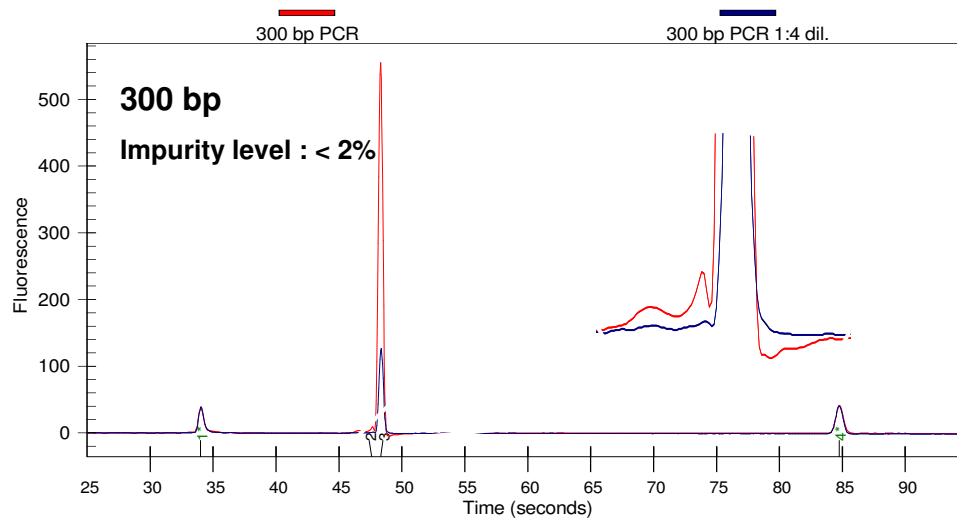
Number	Kit	Max # of samples
5067-1504	Agilent DNA 1000 Kit	300
5067-1505	Agilent DNA 1000 Reagents	
5067-1506	Agilent DNA 7500 Kit	300
5067-1507	Agilent DNA 7500 Reagents	
5067-1508	Agilent DNA 12000 Kit	300
5067-1509	Agilent DNA 12000 Reagents	

2100 kits for RNA applications

RNA Kits

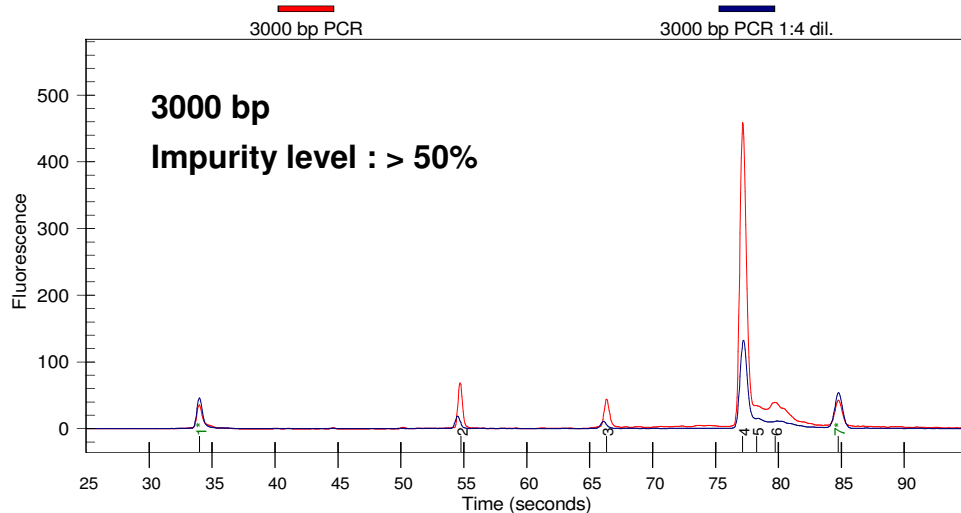
Number	Kit	Max # of samples
5067-1511	Agilent RNA 6000 Nano Kit	300
5067-1512	Agilent RNA 6000 Nano Reagents	
5067-1529	Agilent RNA 6000 Nano Ladder	
5067-1513	Agilent RNA 6000 Pico Kit	275
5067-1514	Agilent RNA 6000 Nano Reagents	
5067-1535	Agilent RNA 6000 Nano Ladder	
5067-1548	Agilent Small RNA Kit	275
5067-1549	Agilent Small RNA Reagents	
5067-1550	Agilent Small RNA Ladder	

Determination of PCR Product Impurity



Quantitative data from Agilent 2100 bioanalyzer

<u>Sample</u>	<u>c (DNA)</u>	<u>main peak</u>
300 bp PCR	41.4 ng/ul	40.7 ng/ul
300 bp PCR 1:4	9.6 ng/ul	9.6 ng/ul



Quantitative data from Agilent 2100 bioanalyzer

<u>Sample</u>	<u>c (DNA)</u>	<u>main peak</u>
3000 bp PCR	61.9 ng/ul	40.7 ng/ul
3000 bp PCR 1:4	14.8 ng/ul	9.8 ng/ul