



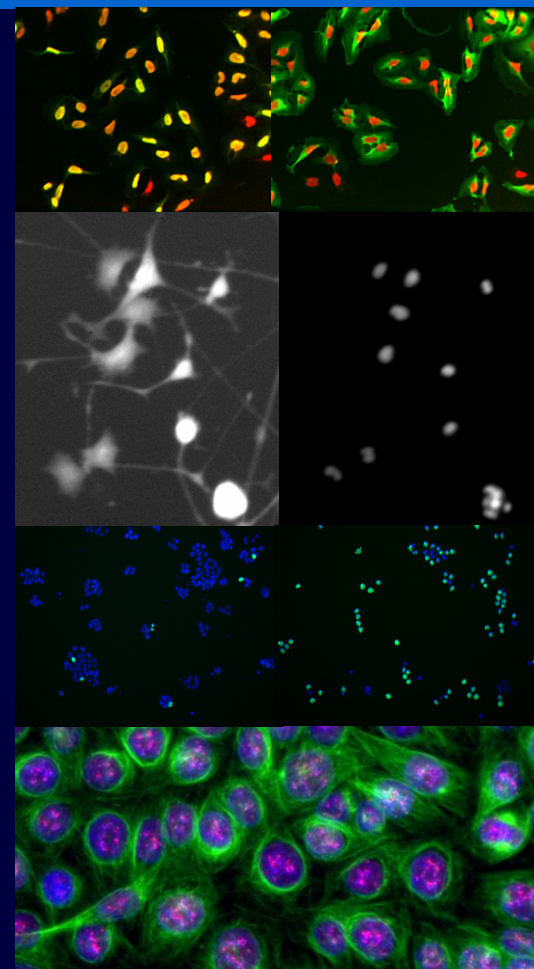
Breakthrough science. Breakthrough medicine.SM

Techniques in High-Content Screening and Assay Development Workshop

Developing a Robust Automated Image-Based Assay

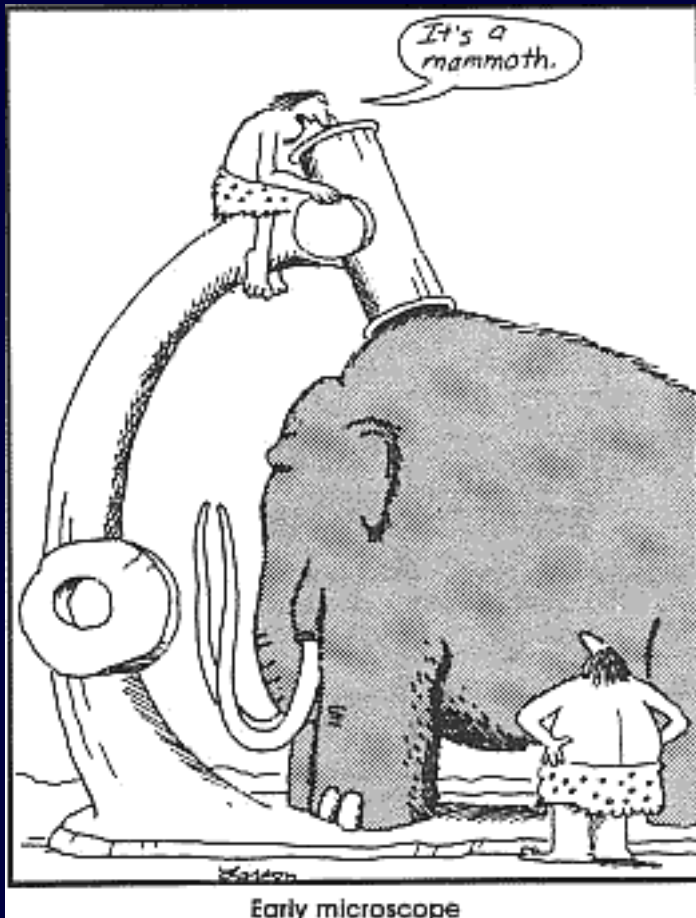
Doug Bowman,
Sr. Mgr Imaging Sciences
Millennium Pharmaceuticals

Drug Discovery Technology
August 7, 2006



The Big Picture

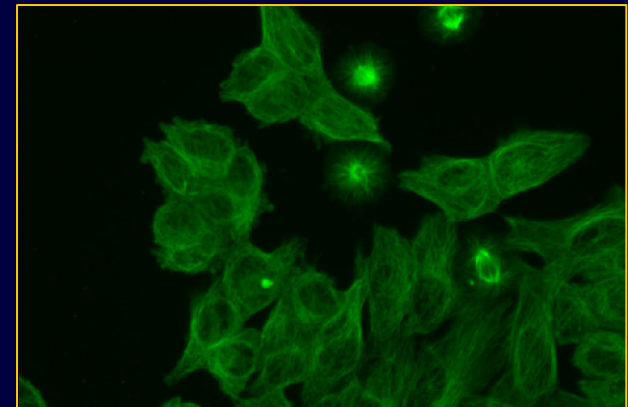
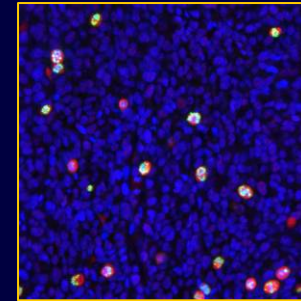
The Far Side *by Gary Larson*



- Cell-based assay
- Immunofluorescence
- Technologies / Instrumentation
- Screening Assay
- People

HCS Assays using Immunofluorescence

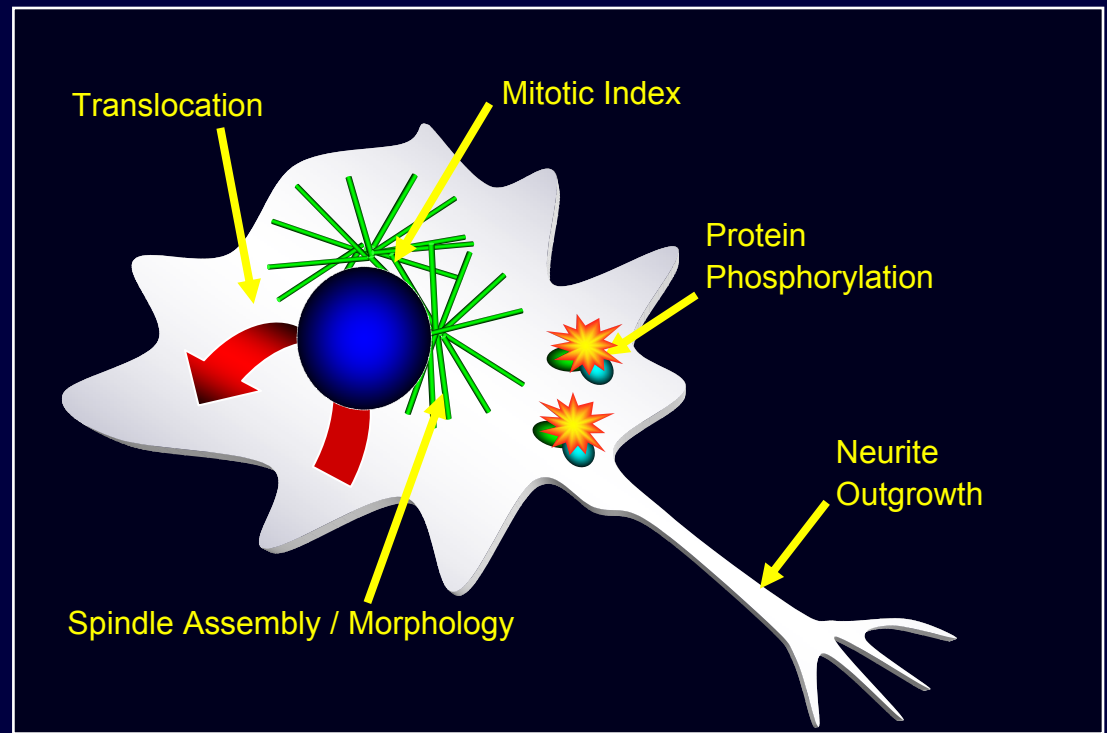
- IF Staining
 - Fix cells to substrate
 - Permeabilize membranes
 - Block
 - Add primary
 - Add secondary
- Image Acquisition
- Image Analysis



Cellular “High Content” Imaging

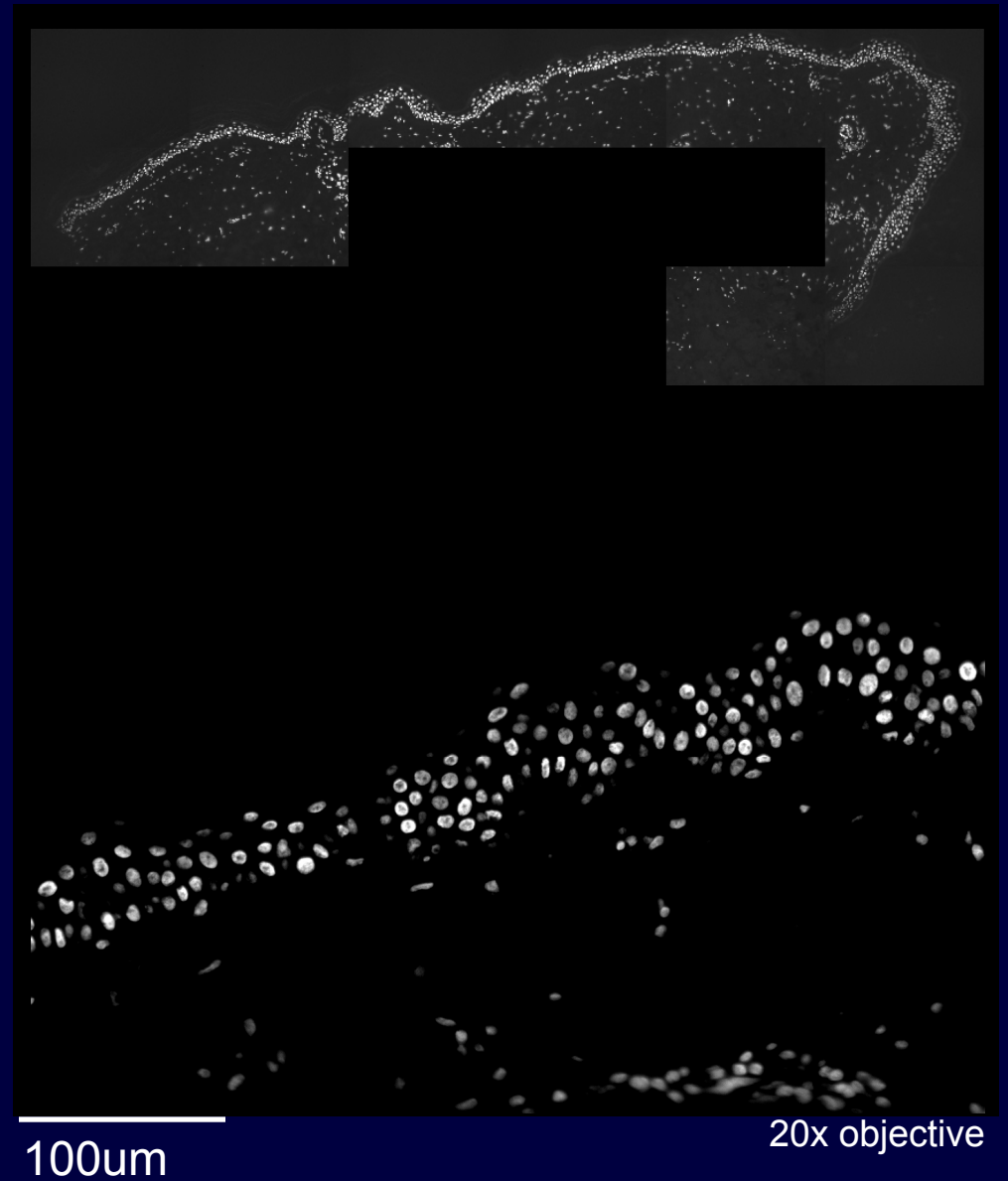
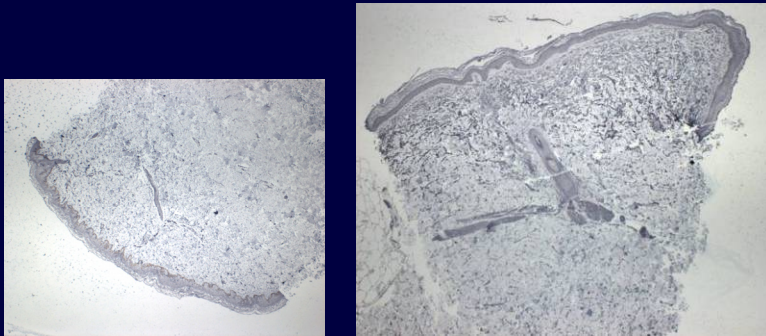
- Provide physiological context for assaying drug activity via the quantitative analysis of cellular and sub-cellular events.
- Analysis also allows for the classification of cellular phenotypes based on morphological parameters.

“HIGH CONTENT”



Tissue...

- High content also!
 - Many of same assays
- Similar issues
- High resolution scanning



Technologies

- Instrumentation

- Liquid handling robotics
- Automated microscope-based systems capable of acquiring high resolution images

Discovery-1



Opera



Automated
Microscope



InCell1000



ArrayScan



- Software algorithms



MetaXpress



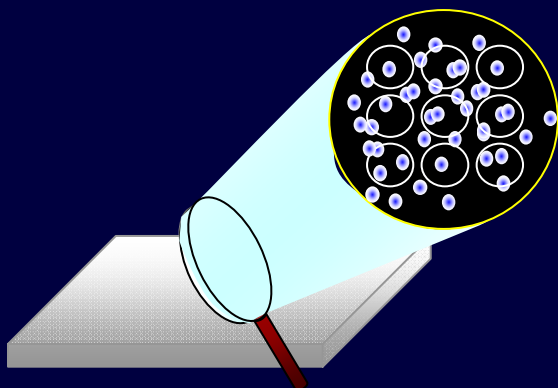
ProGuru



Acapella

Image-based assays

- Large volume of data



= 2λ / image * 9 sites / well * 96 wells / plate

= 1728 images / plate

= 1GB+ / plate

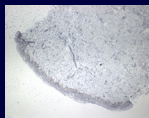
3 plates / hr (Opera) 1 plate / hr (Discovery-1)



= 2λ / image * 5 sites / sample * 3 mice / timepoint

= 30 images / timepoint

= 400+MB / 5 timepoints

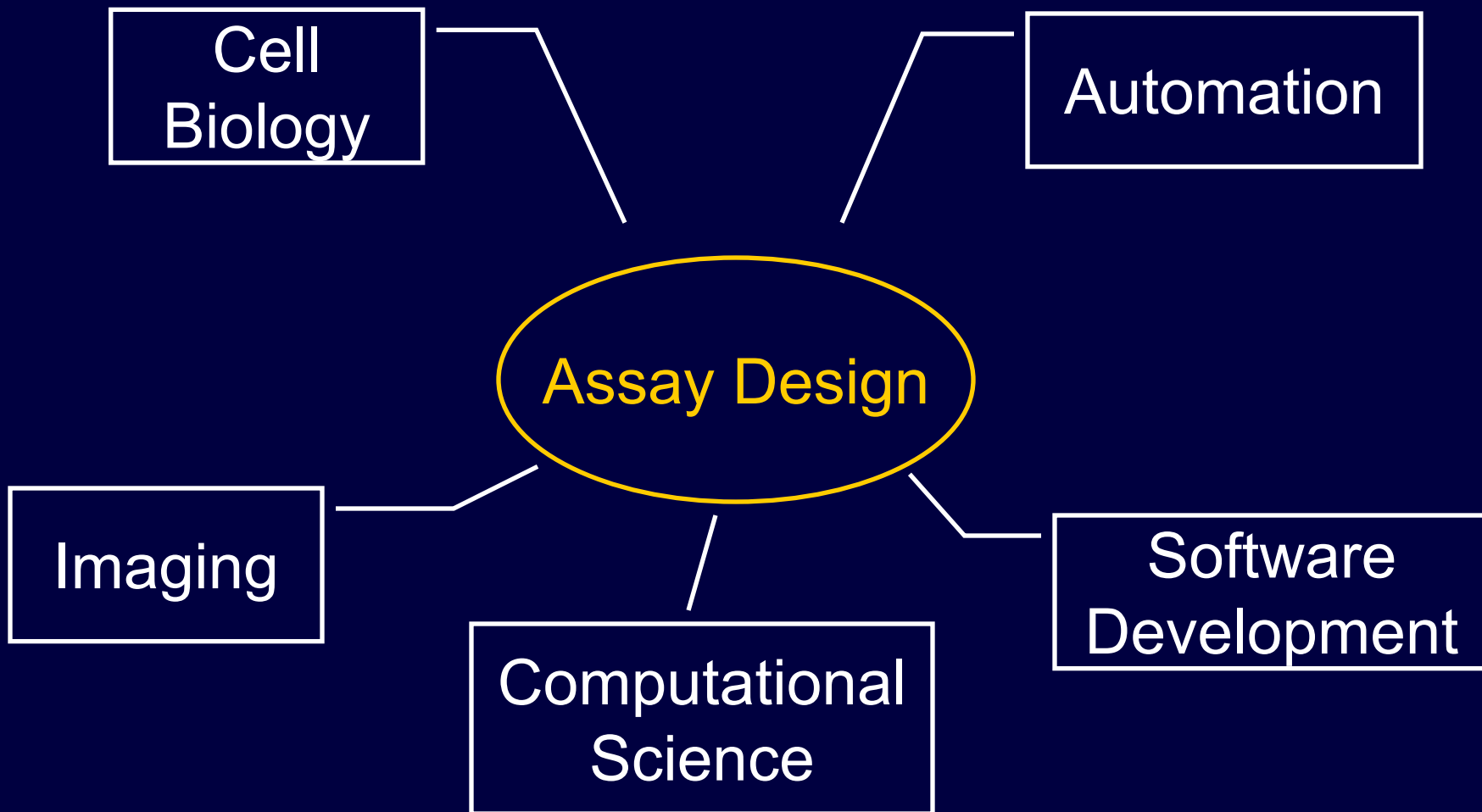


= 3λ / image * 15 sites / membrane

= 30 images / sample

= 120MB / sample

Working Group



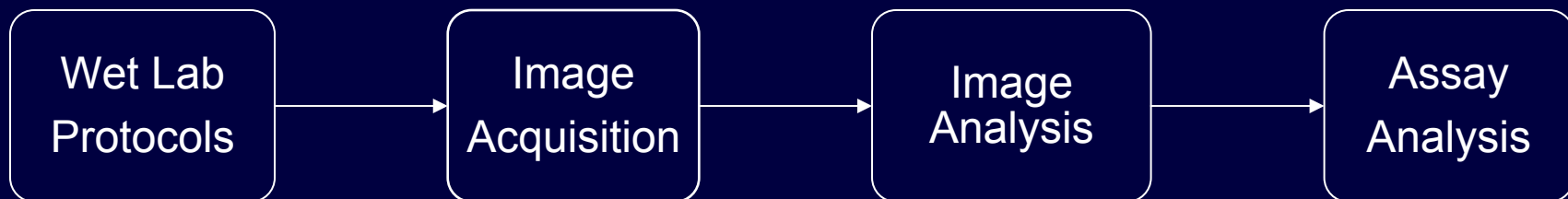
Developing a Robust Automated Image-Based Assay: Guidelines

- Use automation where possible
- Visualize images
- Develop robust process controls
- Use data-driven approach

A data-driven approach to assay design

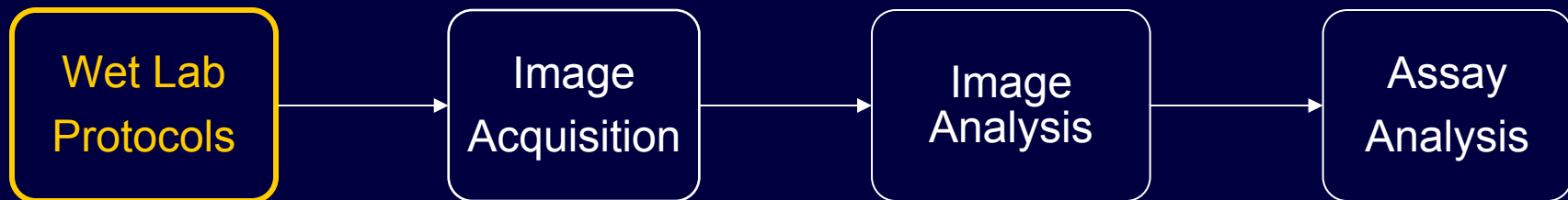
- Dissect components of the assay
- Using control experiments
 - Identify drivers of variability
 - Minimize them experimentally or account for them statistically
 - Optimize parameter selection (robust)
- Assess performance of assays in project-relevant context

HCS Assay Components



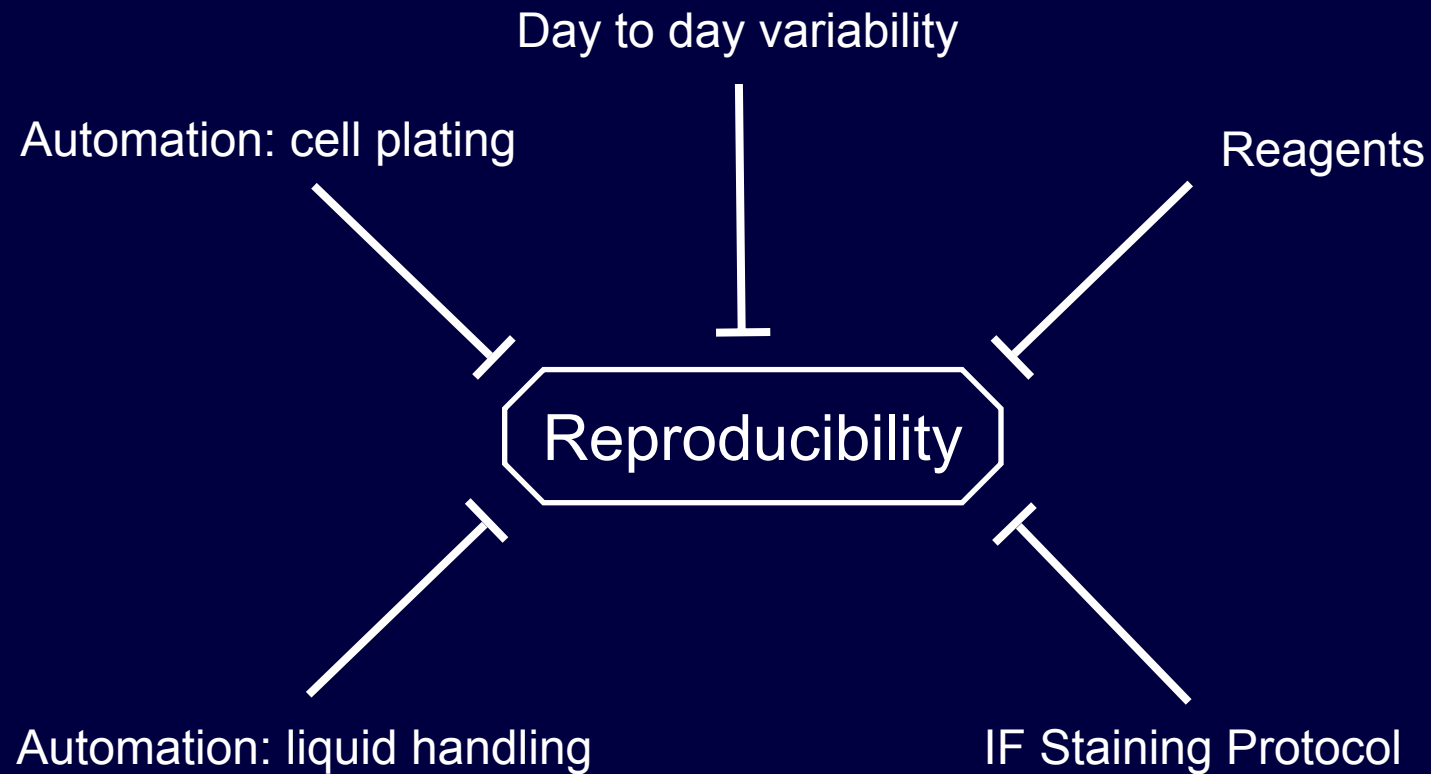
- Dissect components
- Using control experiments
 - Identify drivers of variability
 - Minimize them experimentally or account for them statistically

HCS Assay Components



- Dissect components
- Using control experiments
 - Identify drivers of variability
 - Minimize them experimentally or account for them statistically

Wet Lab Protocols



Wet Lab Protocols:

Testing liquid handling using Artel system

- Accuracy and Reproducibility
 - Compound addition / dilution
 - Cell Plating
 - Antibody addition / protocols



Artel MVS

		<u>CVs</u>
100ul	manual (12 channel Multipipette)	4.43%
100ul	Multidrop	1.67%
50ul	Platamate	0.75%
50ul	Tecan	0.80%
10ul	Precision 2000	1.34%

Older Multidrop tube cassettes found inaccurate
– *relative inaccuracy* ~25% and *CVs* ~2%

New cassettes purchased and validated –
relative inaccuracy ~2% and *CVs* ~0.7%

Wet Lab Protocols:

IF Staining Protocol

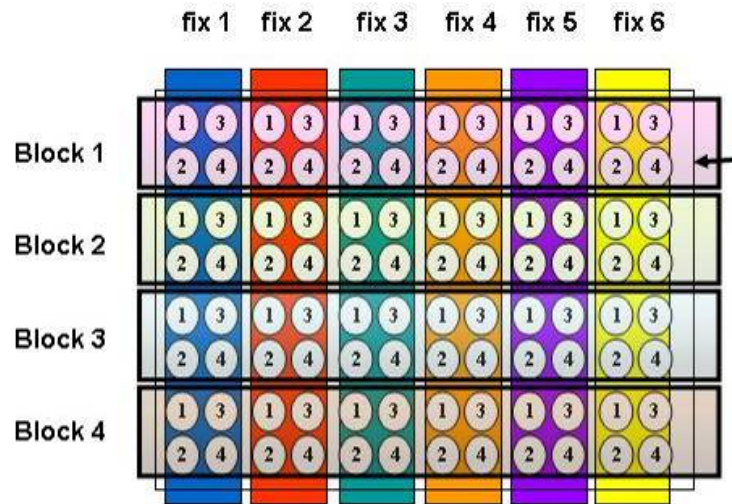
Facilitate a systematic comparison of a large number of assay parameters to maximize data quality while minimizing the time required to find optimal parameters.

- fix
- perm
- block
- primary ab dilution
- secondary ab dilution
- buffers
- wash buffers
- number of washes
- time/temp of incubation....etc

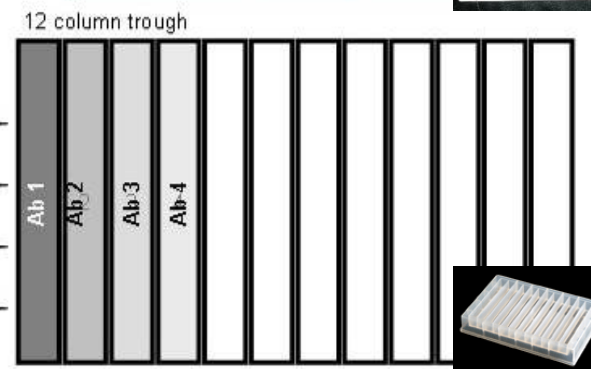
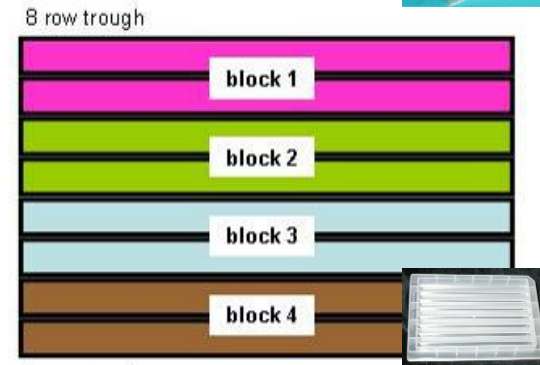
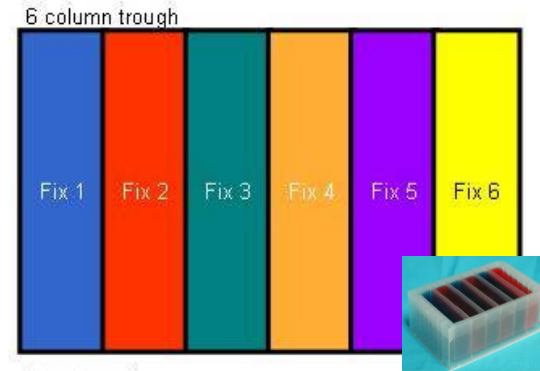
Antibody Optimization

Matrix of 6 fixatives - 4 blocks - 4 ab dilutions to assay

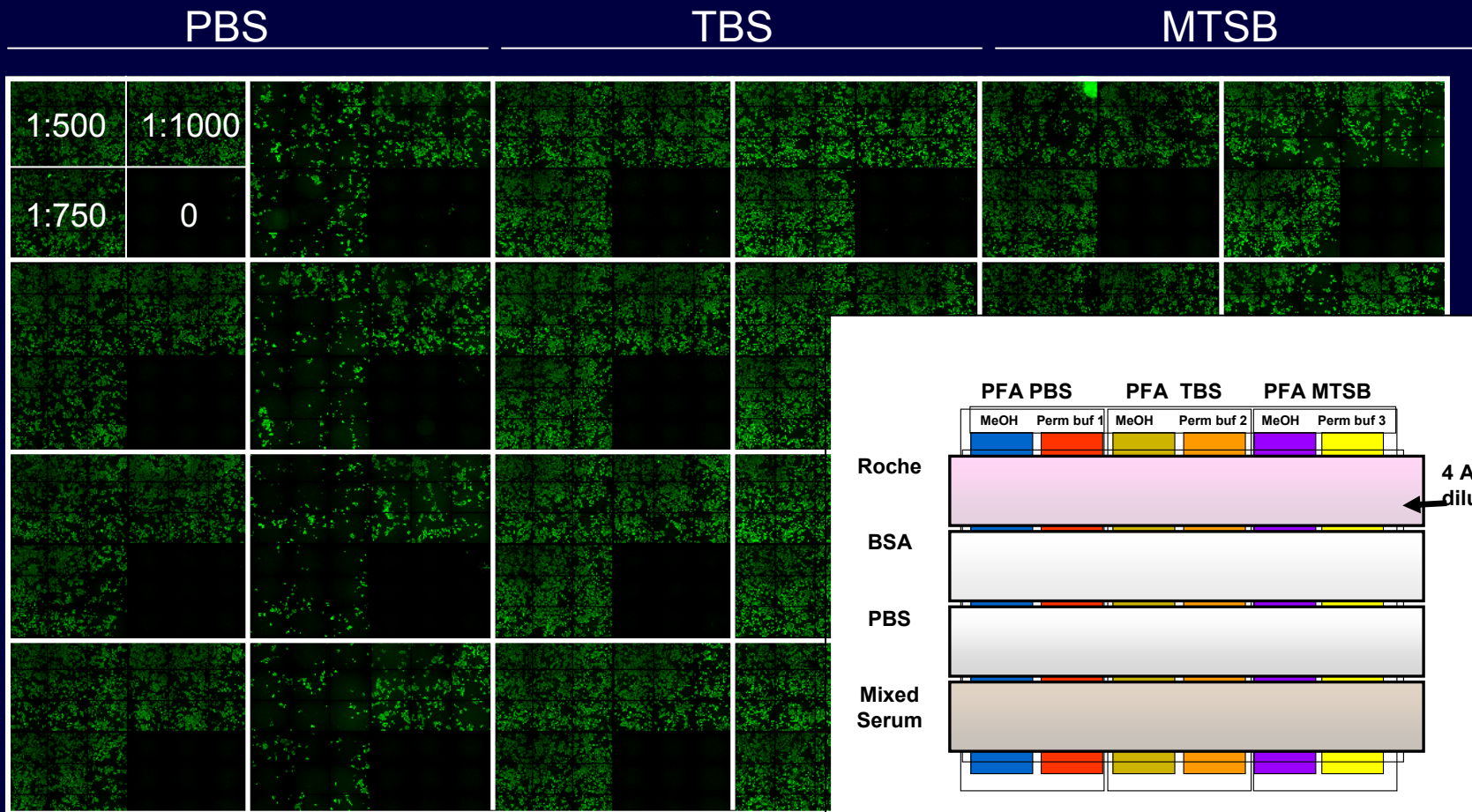
96 individual conditions



4 Ab dilutions in quadrants

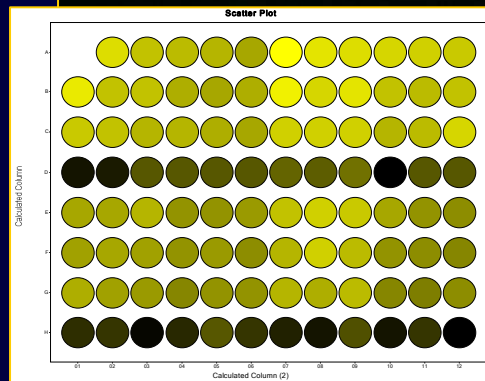
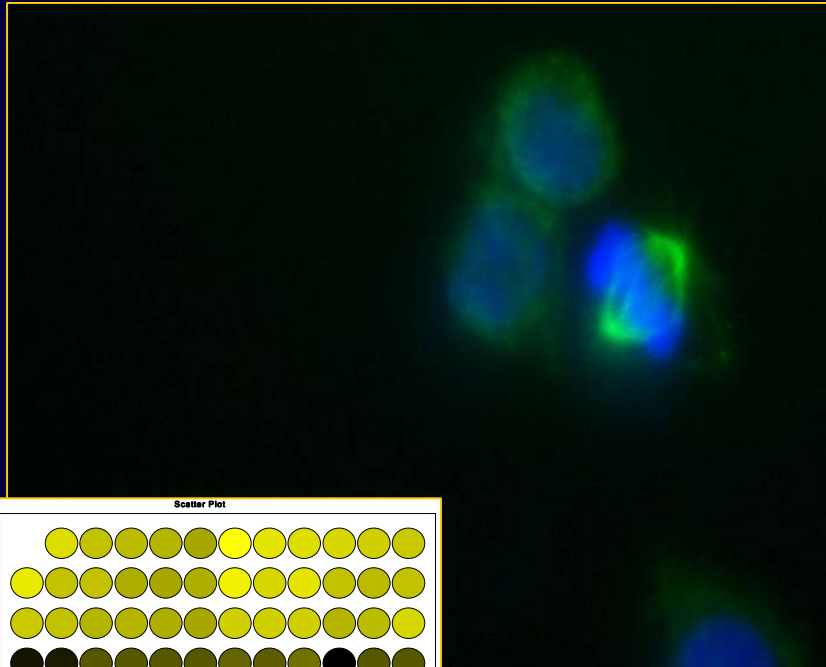


Wet Lab Protocols: Antibody Optimization



Wet Lab Protocols: Antibody Optimization

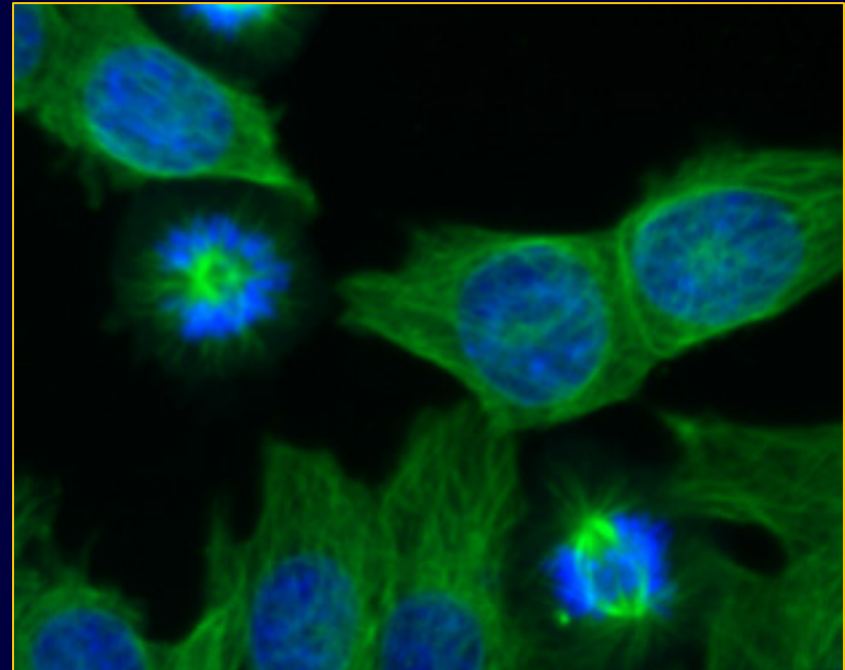
Before



S/N

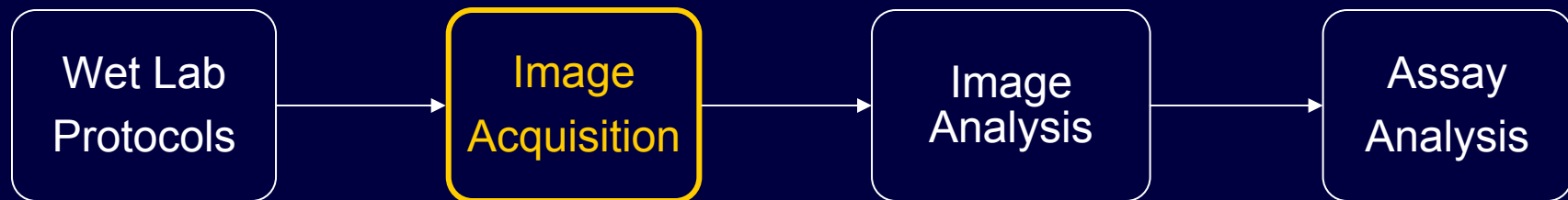
DNA / Tubulin

After



- Comparison of > 400 parameters (~2 weeks)
- Enhances analysis of spindle morphology
- Save \$ Antibody concentration

HCS Assay Components



- Dissect components
- Using control experiments
 - Identify drivers of variability
 - Minimize them experimentally or account for them statistically

Image Acquisition

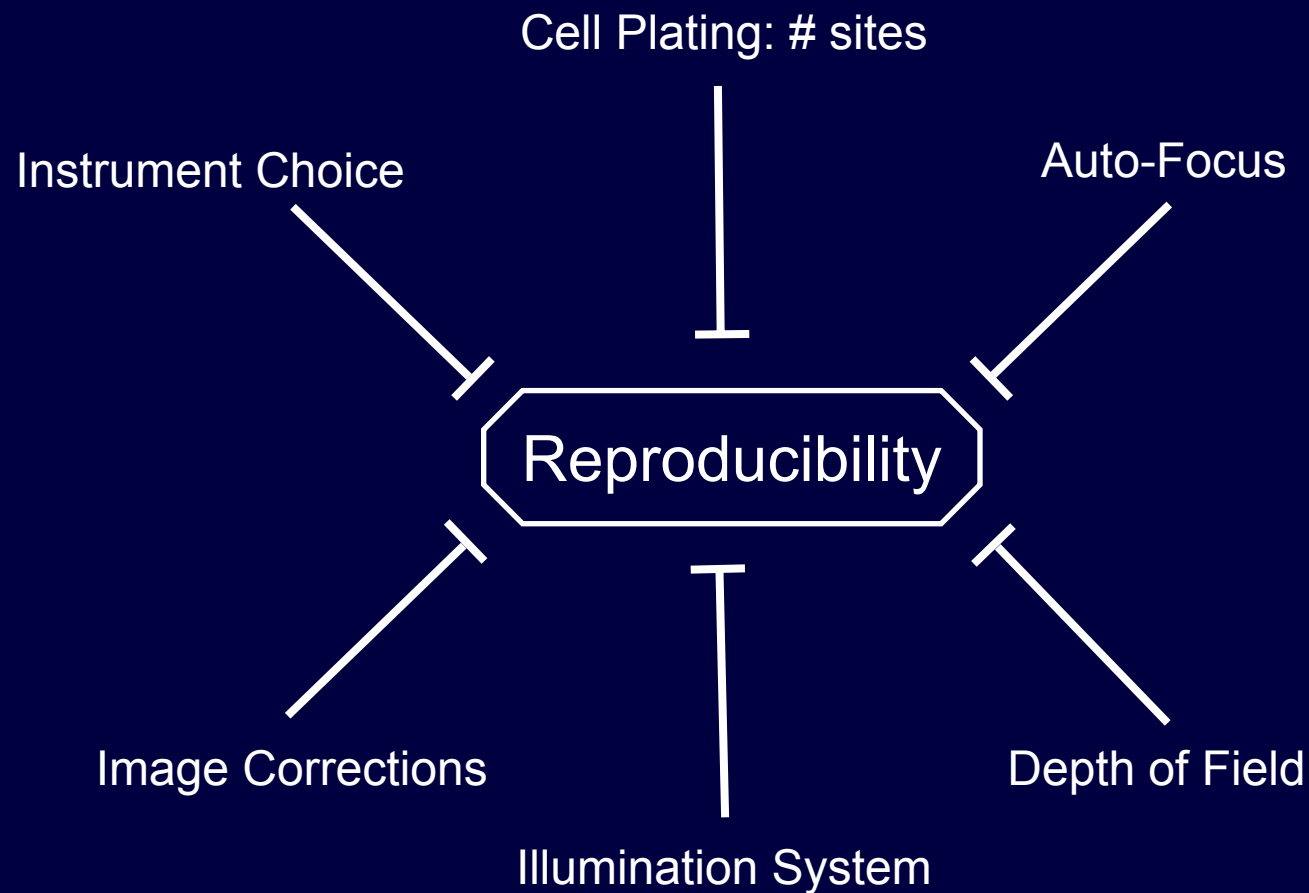


Image Acquisition

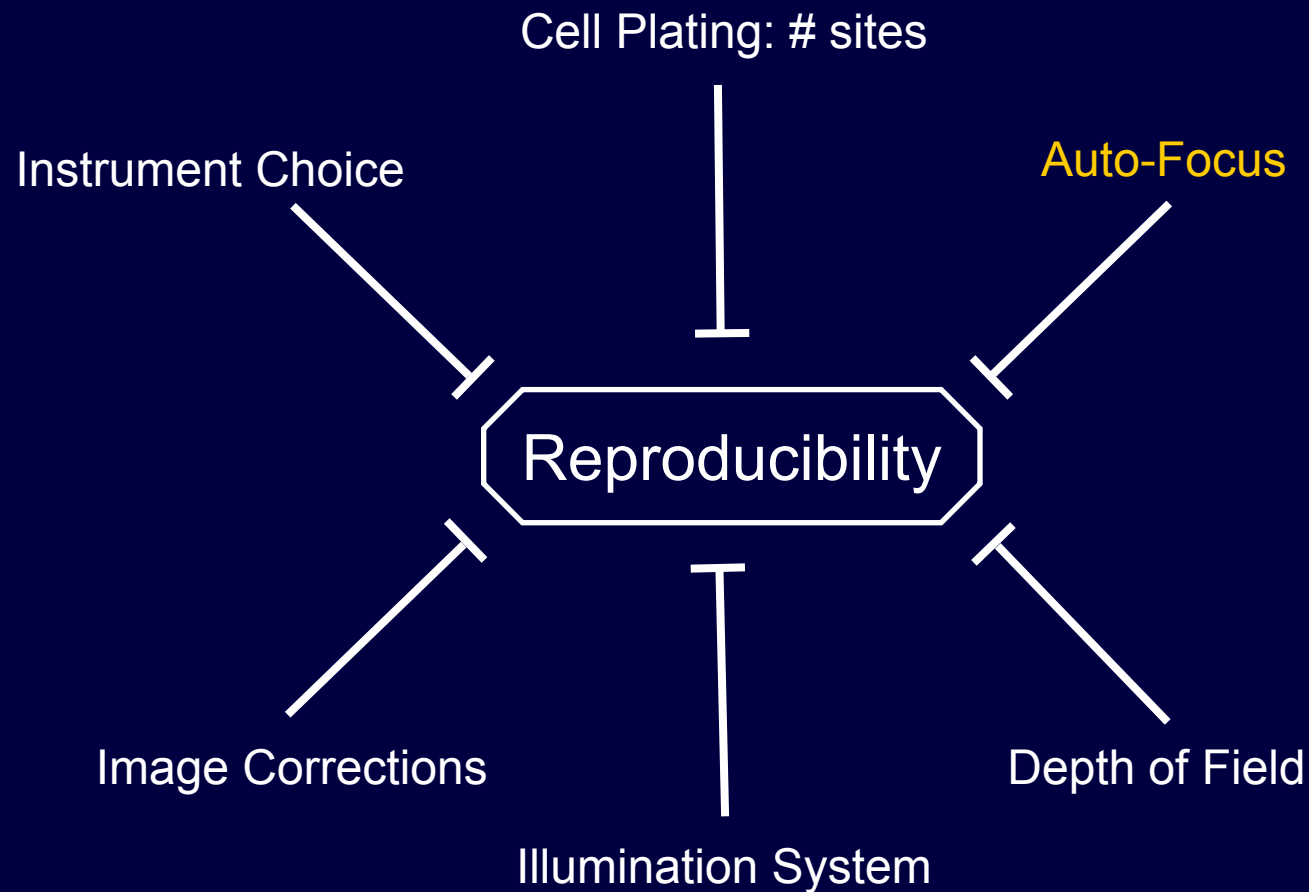


Image Acquisition: AutoFocus

- Image-based and Laser-based

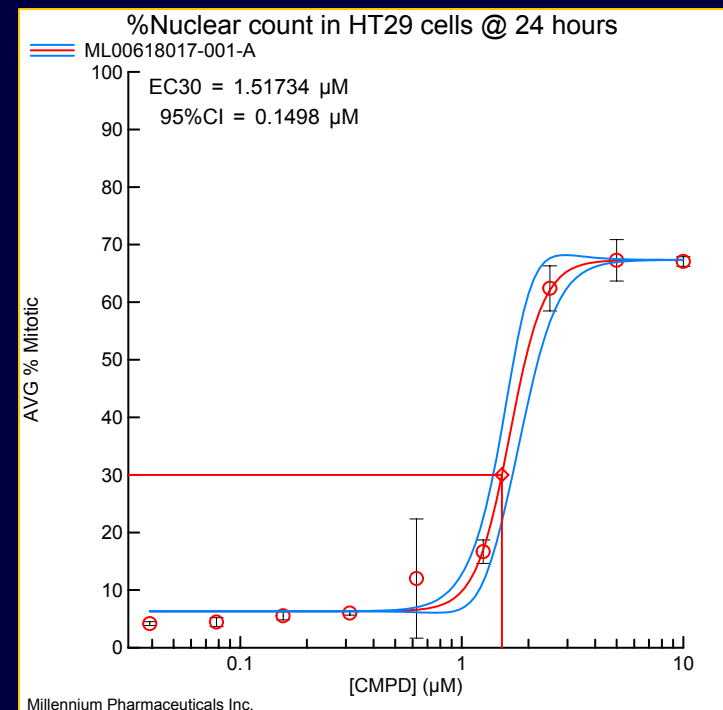
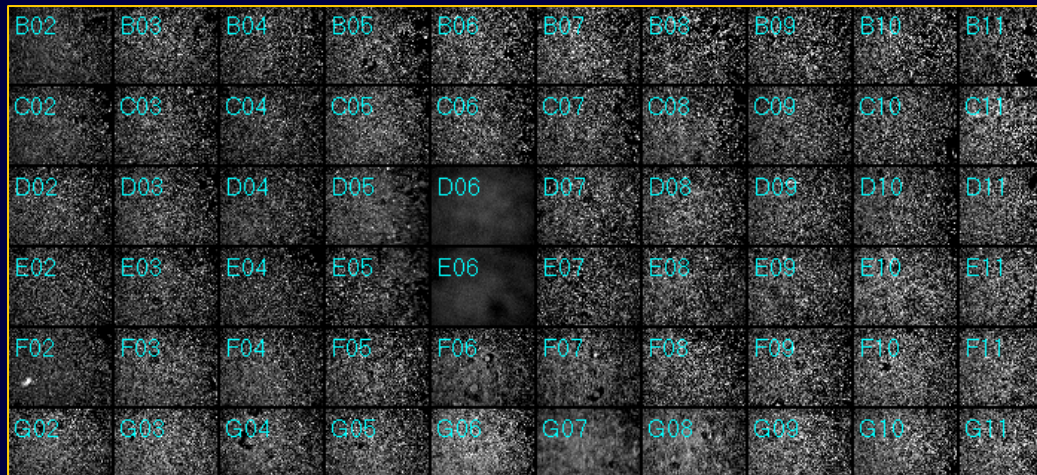
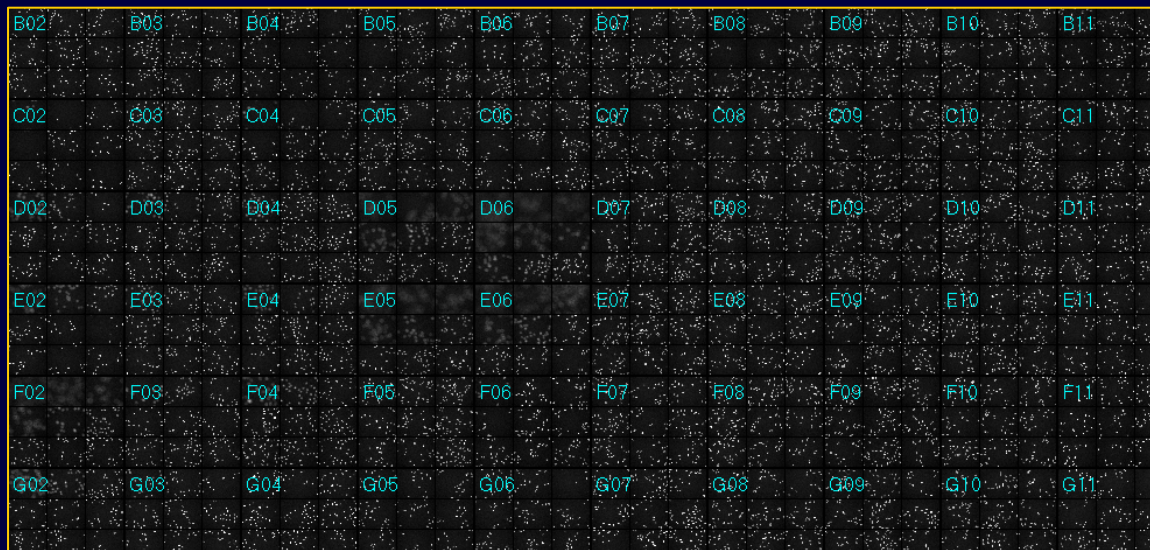


Image Acquisition: AutoFocus



60 wells, 9 sites / well

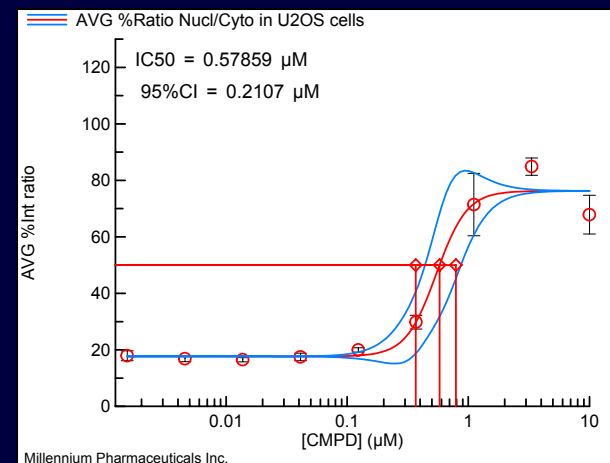


Image Acquisition:

AutoFocus

- Minimize experimentally
 - Instrument parameters
 - Plate selection
- Identify, Remove: Auto-focus Classifier
 - Classify several hundred images by hand
 - Use image processing measures to extract quantities that vary based on edges and intensity
 - Train a classifier to call out-of-focus wells

Image Acquisition: AutoFocus

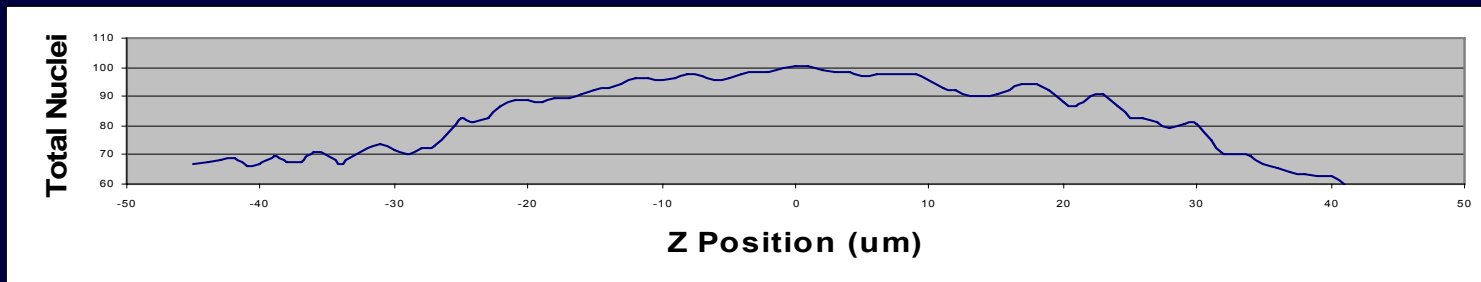
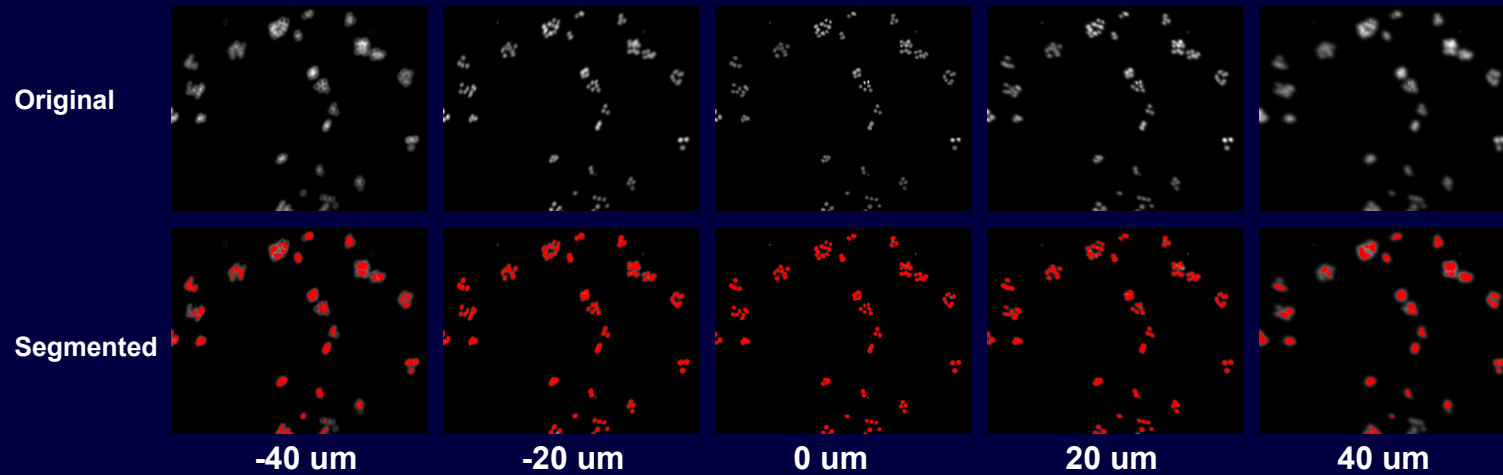


Image Acquisition:

AutoFocus

- BayesNet classifier learns differences between in-focus and out-of-focus images
- Algorithm maximizes separation of classes on training set by:
 - automatically assigning weights for each predictor
 - automatically selecting threshold

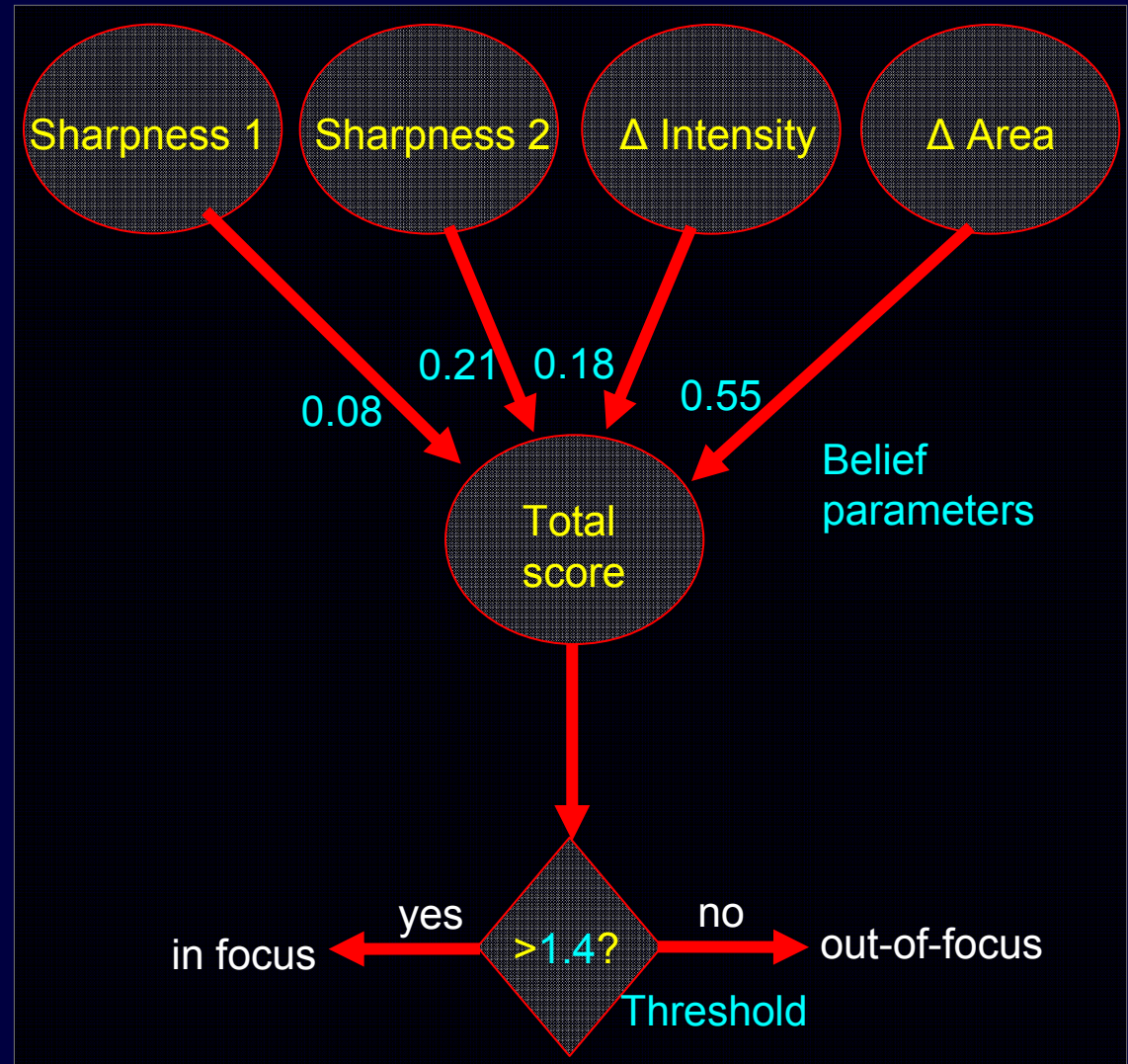


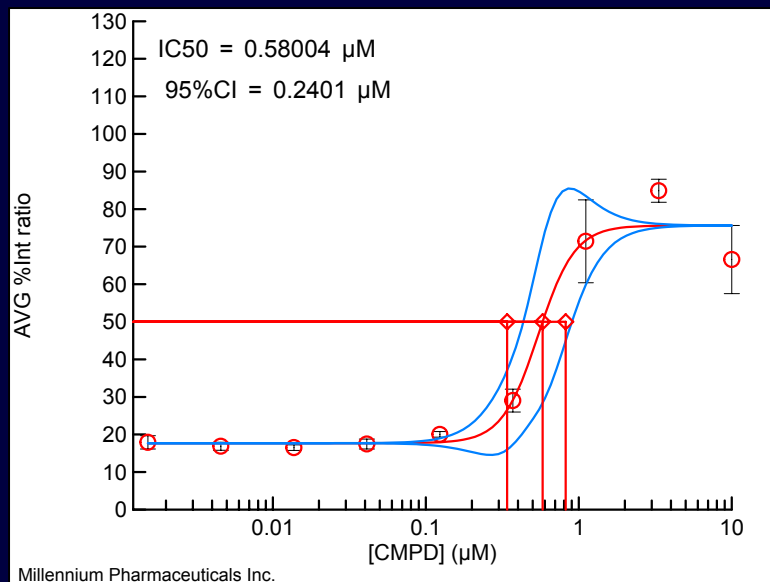
Image Acquisition:

AutoFocus

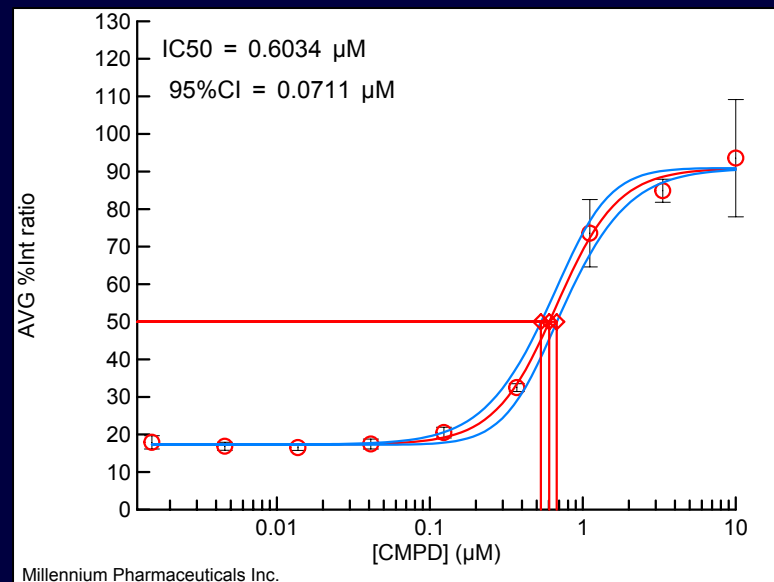
- Results of BayesNet assessed using cross-validation:
 - Algorithm is trained on 9/10ths of the data and tested on remaining 1/10th
 - Process is repeated 10 times
 - Error rates are averaged
- 93% to 98% of the wells were called correctly (for 3 different cell types).
- Nearly all classification errors occurred in 'marginal' category
- Zero-R (baseline classification error)= 32%

Image Acquisition: AutoFocus

Filtering out-of-focus images via machine learning



Original Data



Filtered Data

Image Acquisition

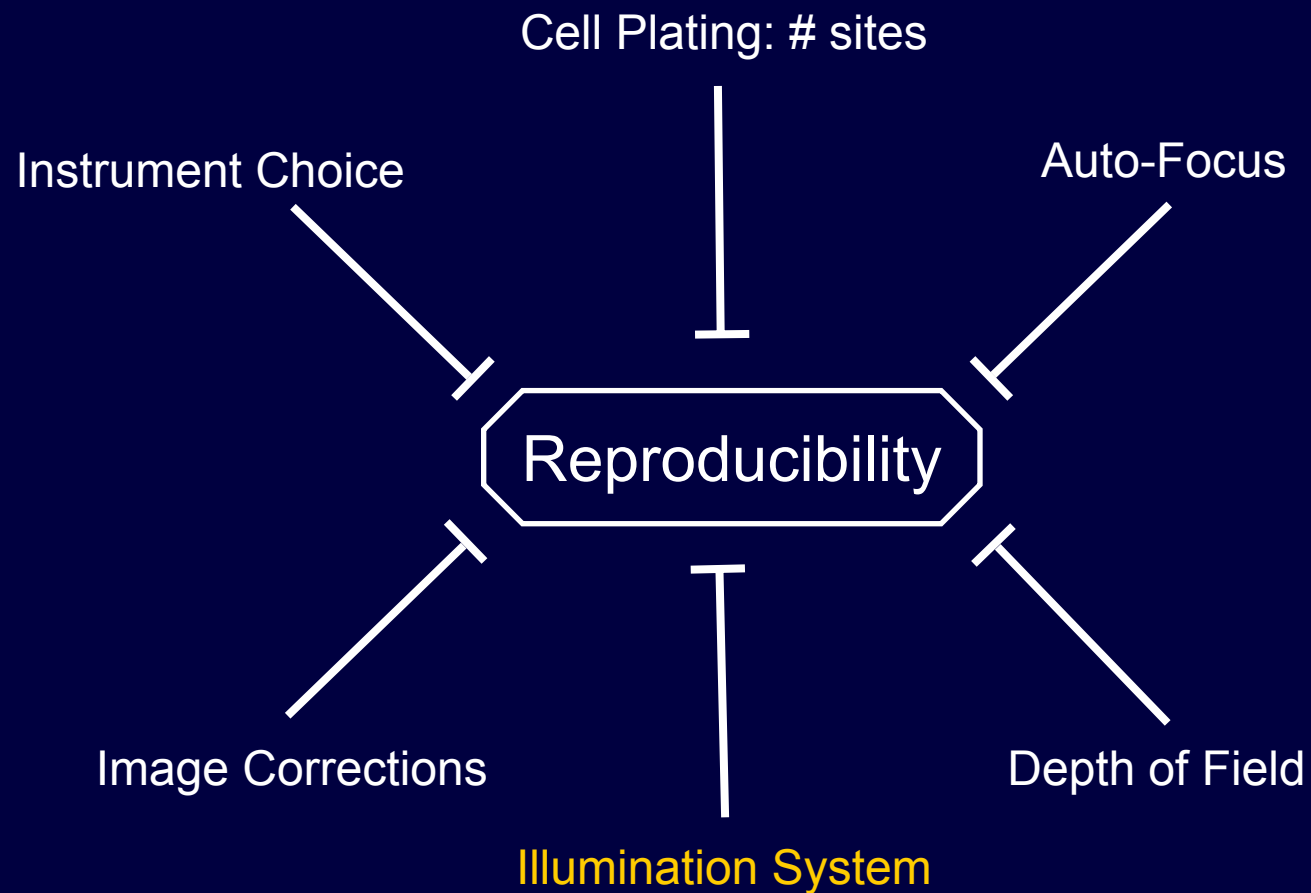
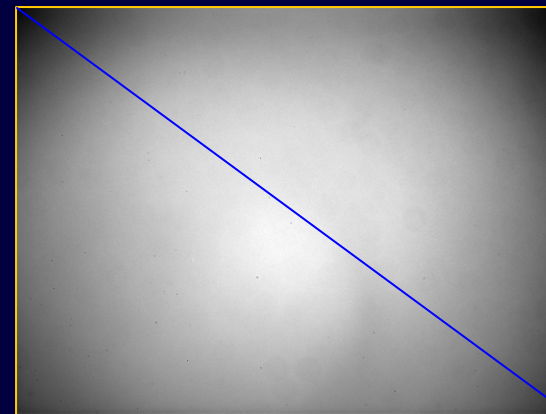
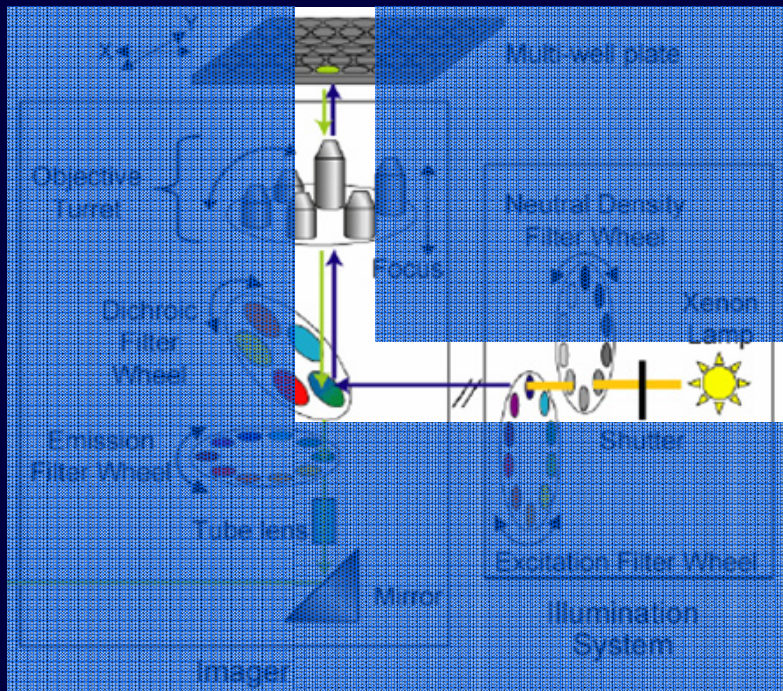


Image Acquisition:

Illumination system variation

- What is it?
 - Fluorescence illumination system



Uniform sample

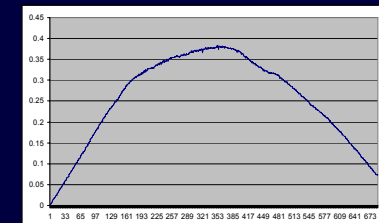


Image Acquisition:

Illumination system variation

- Is it a problem?
 - ↓ Cell counting assays
 - ↑↑ Cell intensity assays
 - ↑↑ Phenotyping

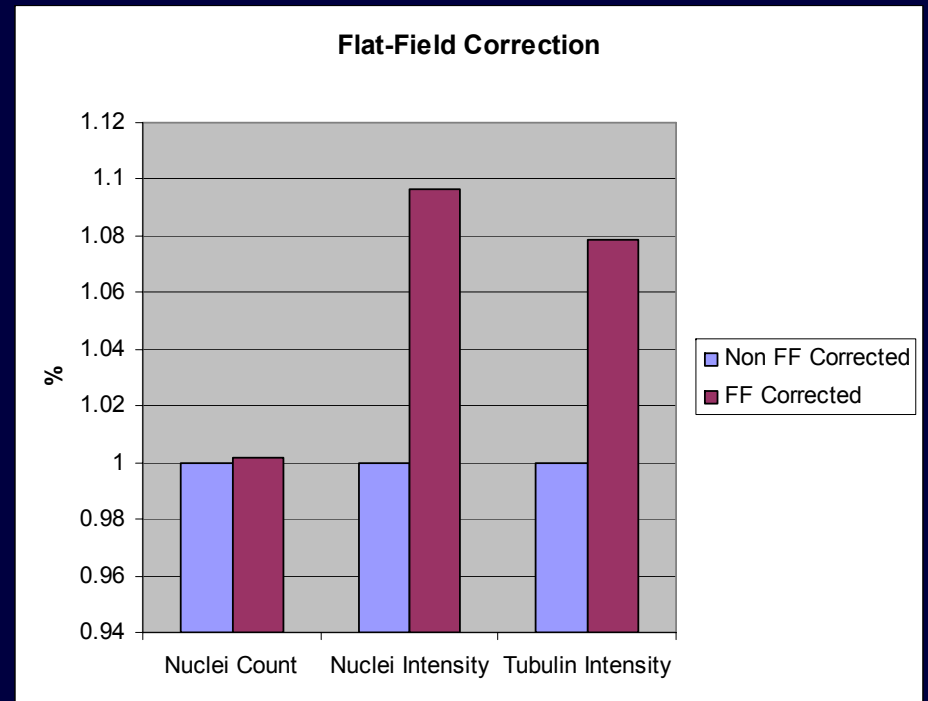


Image Acquisition

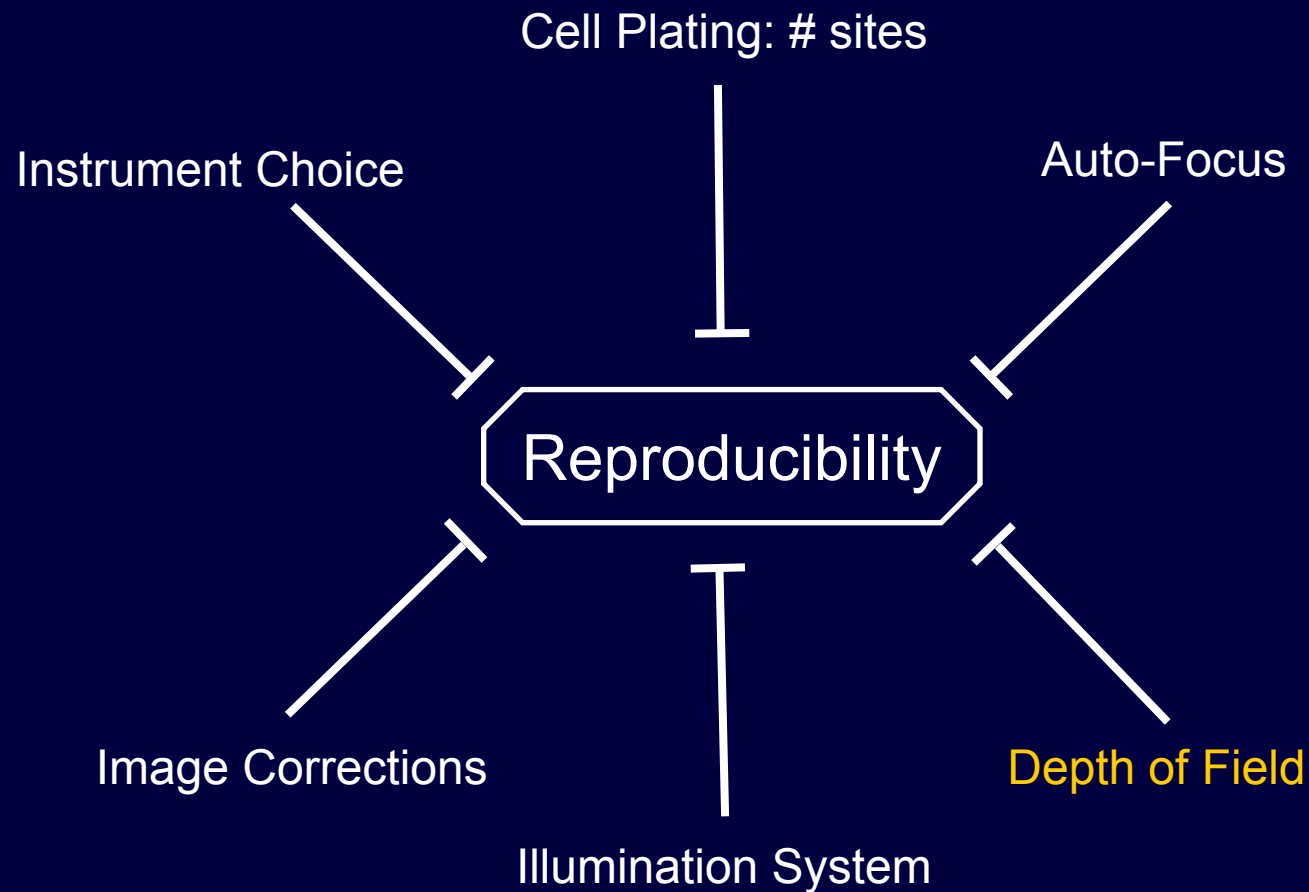
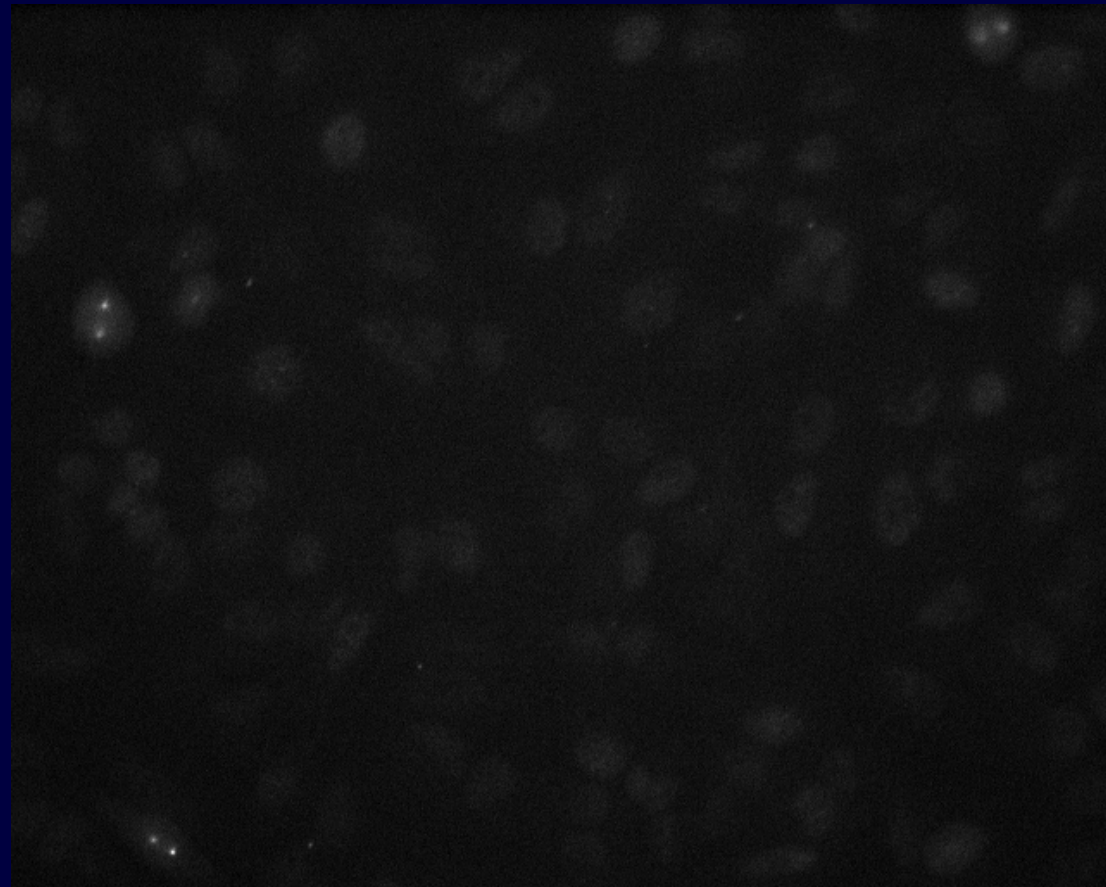


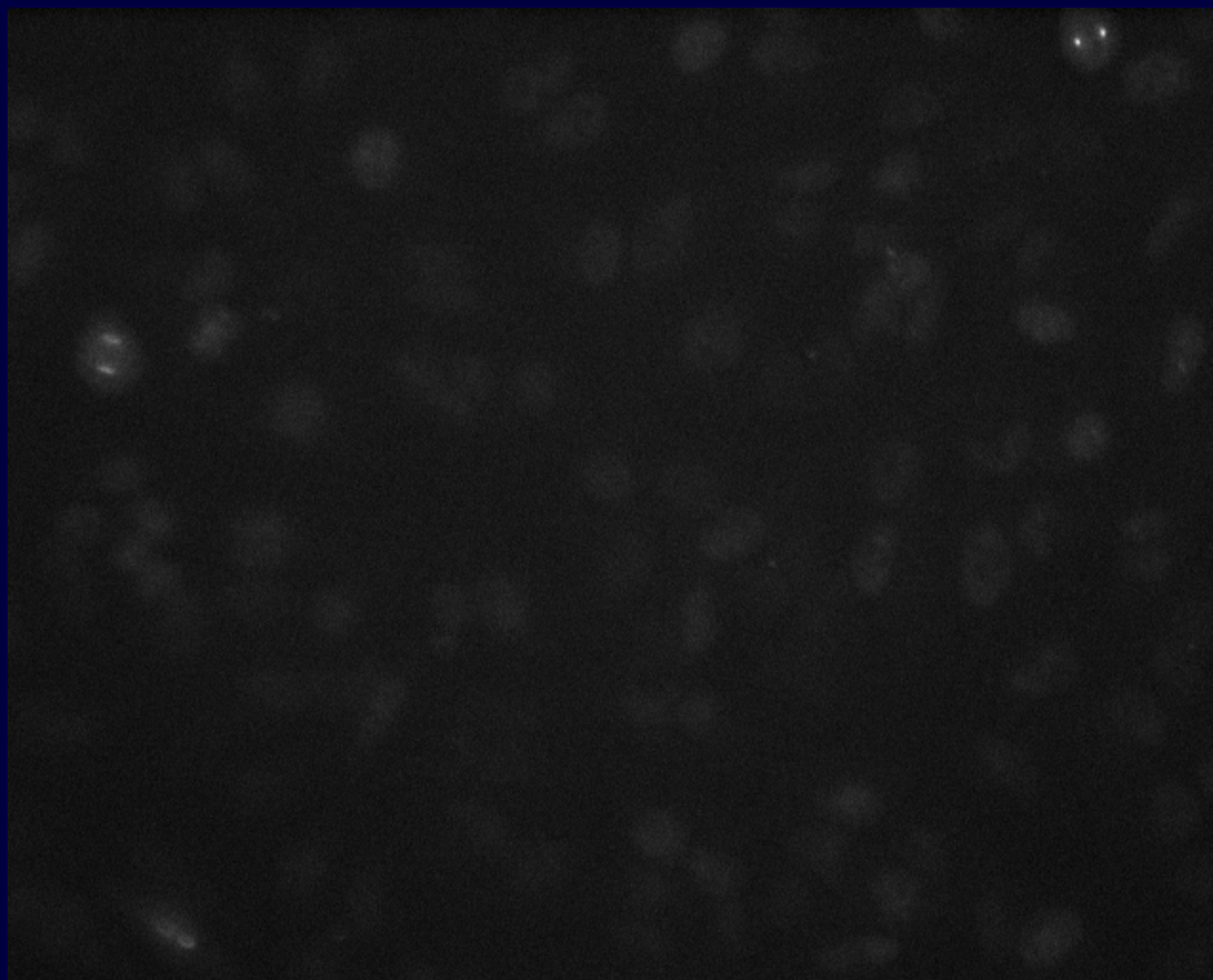
Image Acquisition:

Depth of Field



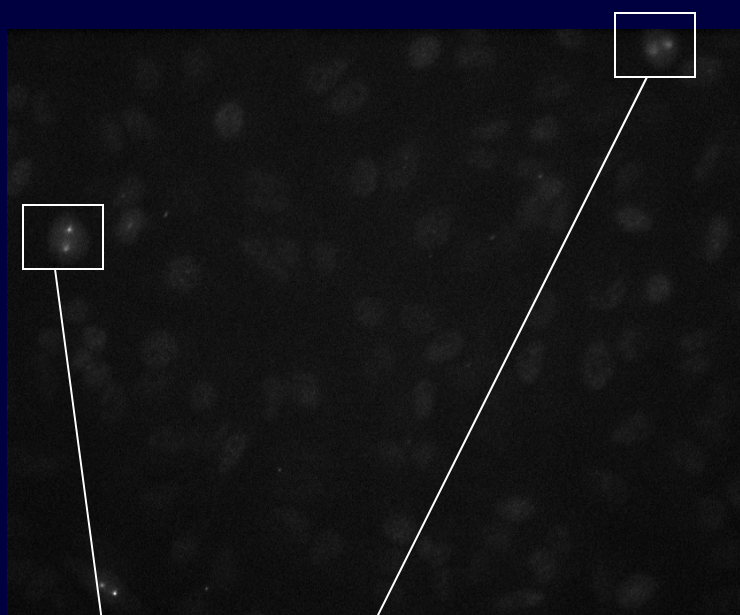
20x

Image Acquisition: Depth of Field

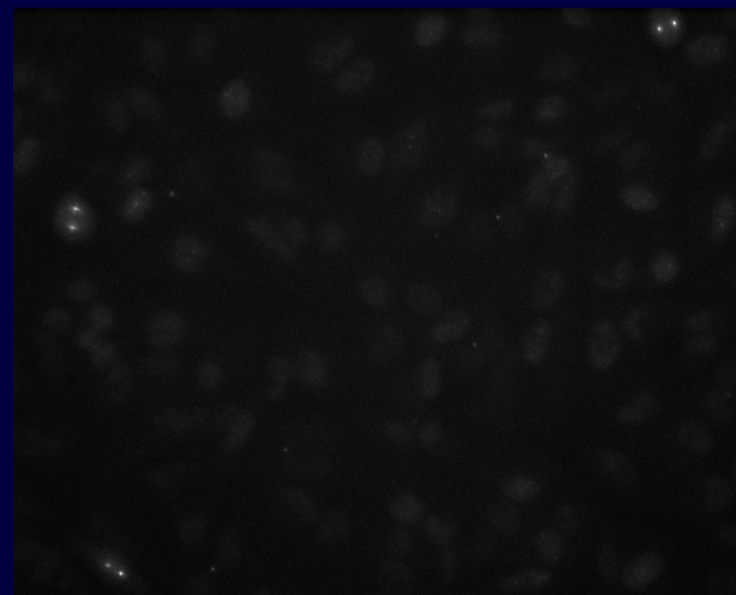


- 2.0 μm

Image Acquisition: Depth of Field



Single Plane



“Best Focus”



Z = -4.0um

Z = -2.0um

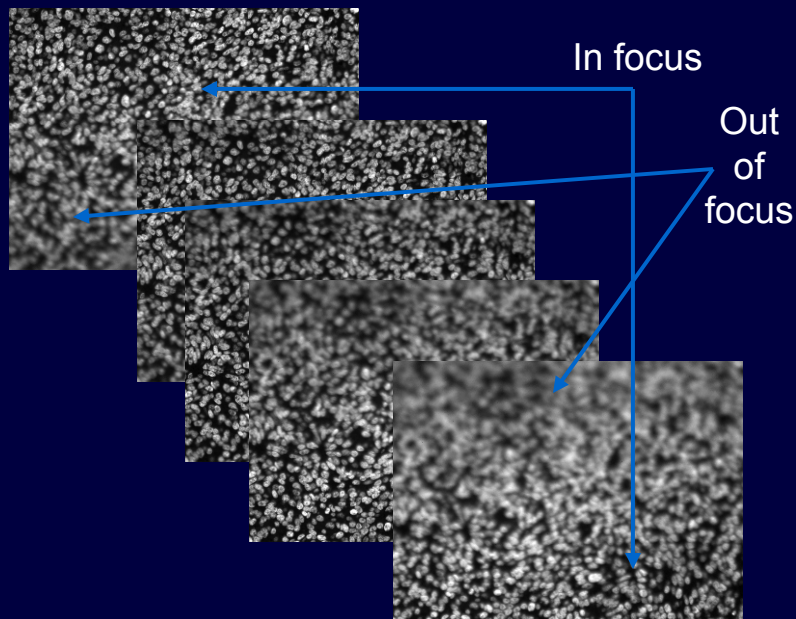
Z = 0.0um

Z = 2.0um

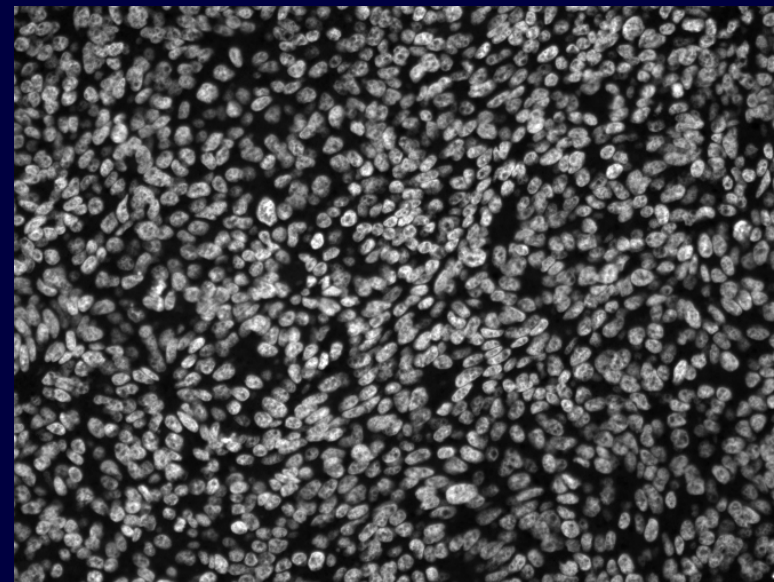
Z = 4.0um

Tissue Imaging

- Challenge: depth of focus of the objective is less than the flatness of the specimen. This results in portions of a single image plane being out of focus.

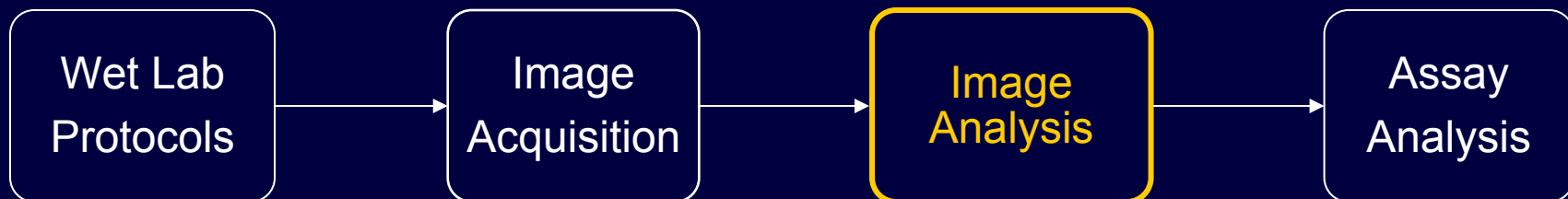


Multiple optical sections @ 2.0um intervals



"Best Focus" Image

HCS Assay Components



- Dissect components
- Using control experiments
 - Identify drivers of variability
 - Minimize them experimentally or account for them statistically

Image Analysis

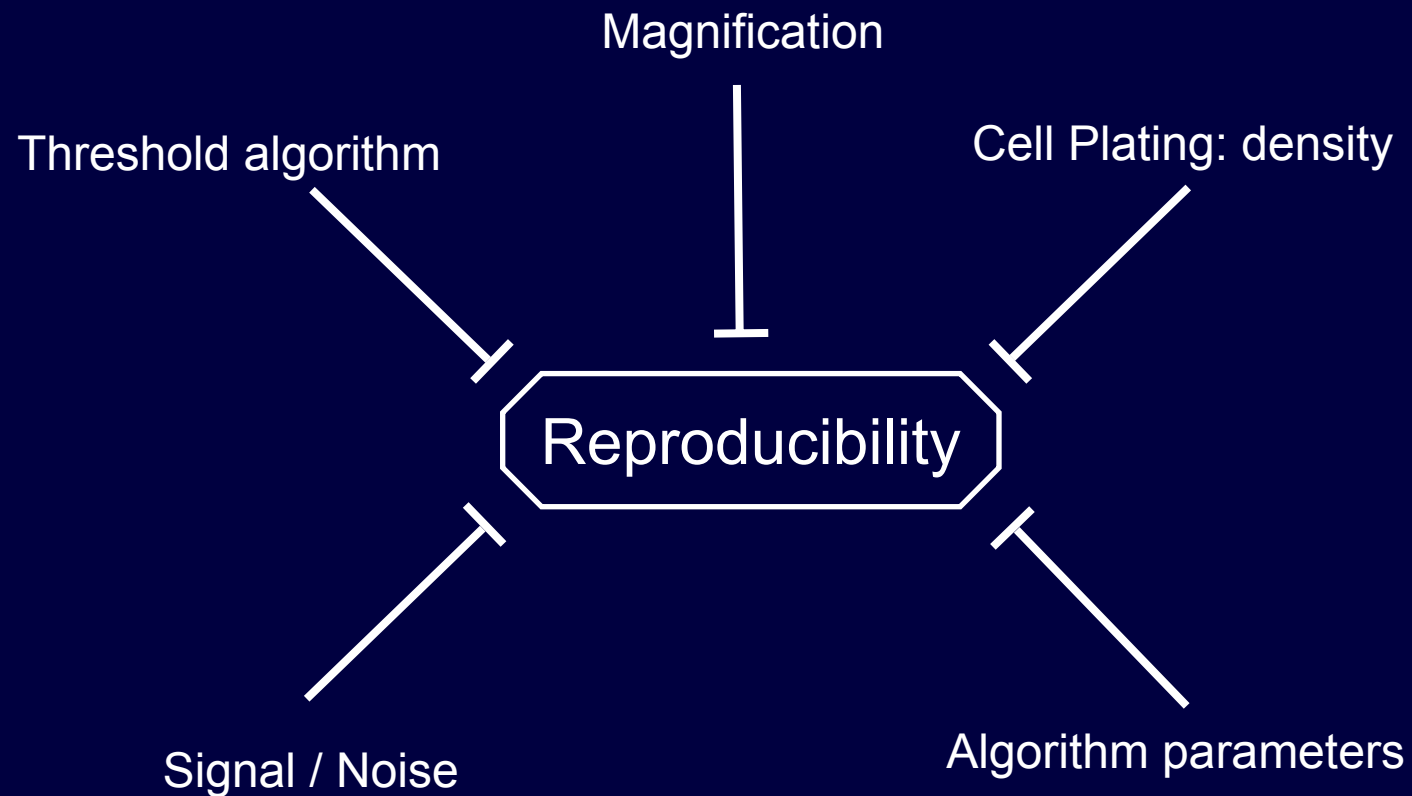


Image Analysis:

Cell Plating Density

- Assess
 - Optimal cell plating density to minimize segmentation errors
- DOE:
 - Serial dilution
 - DAPI
 - 6 replicates

	1	2	3	4	5	6	7	8	9	10	11	12
A	X	X	X	X	X	X	X	X	X	X	X	X
B	X	← Cell plating density →									X	
C	X	← Cell plating density →									X	
D	X	← Cell plating density →									X	
E	X	← Cell plating density →									X	
F	X	← Cell plating density →									X	
G	X	← Cell plating density →									X	
H	X	X	X	X	X	X	X	X	X	X	X	X

Image Analysis:

Cell Plating Density

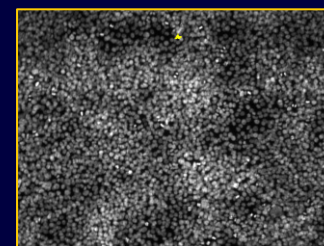
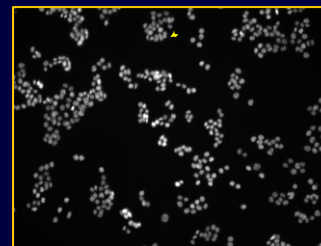
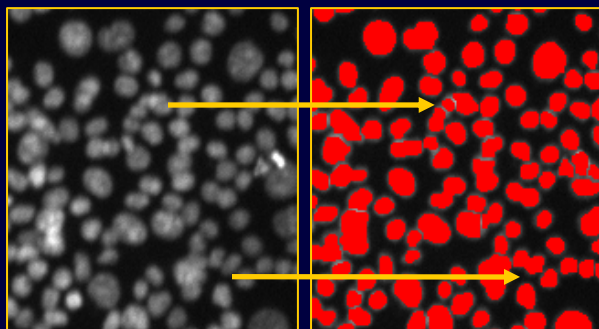
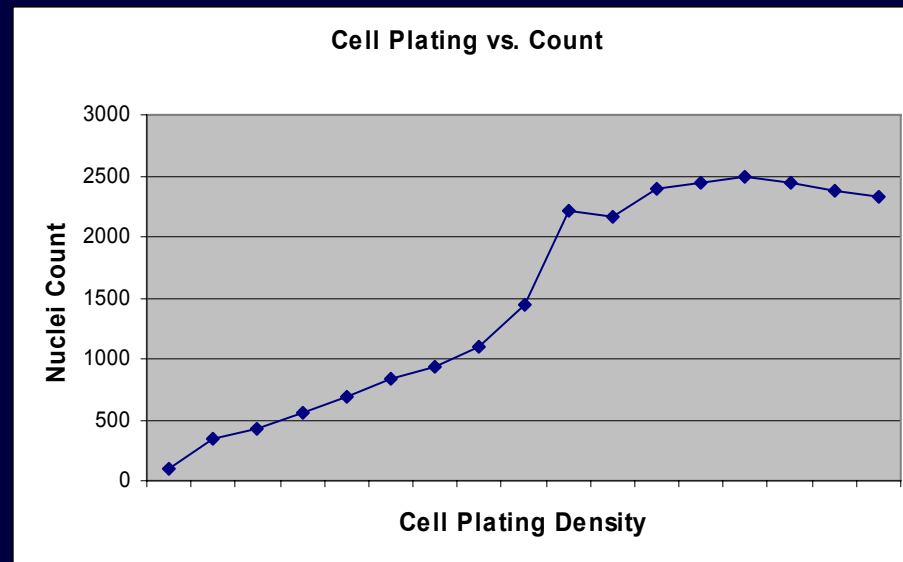


Image Analysis:

Threshold algorithm

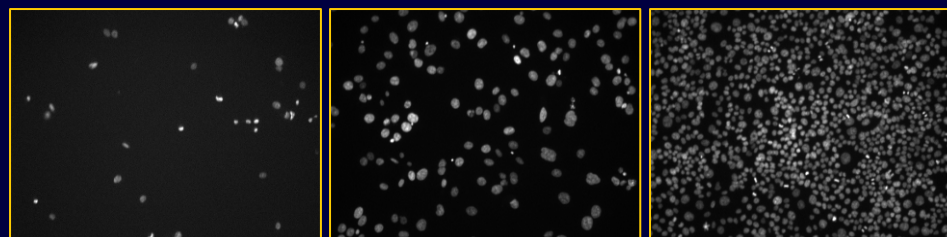
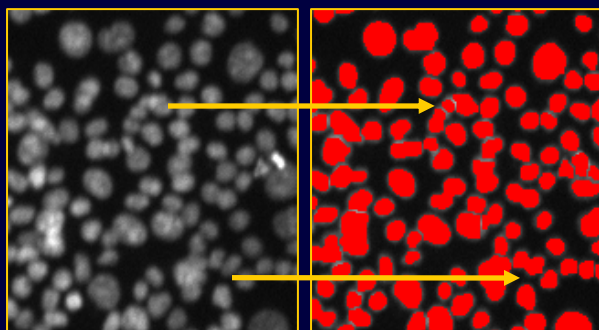
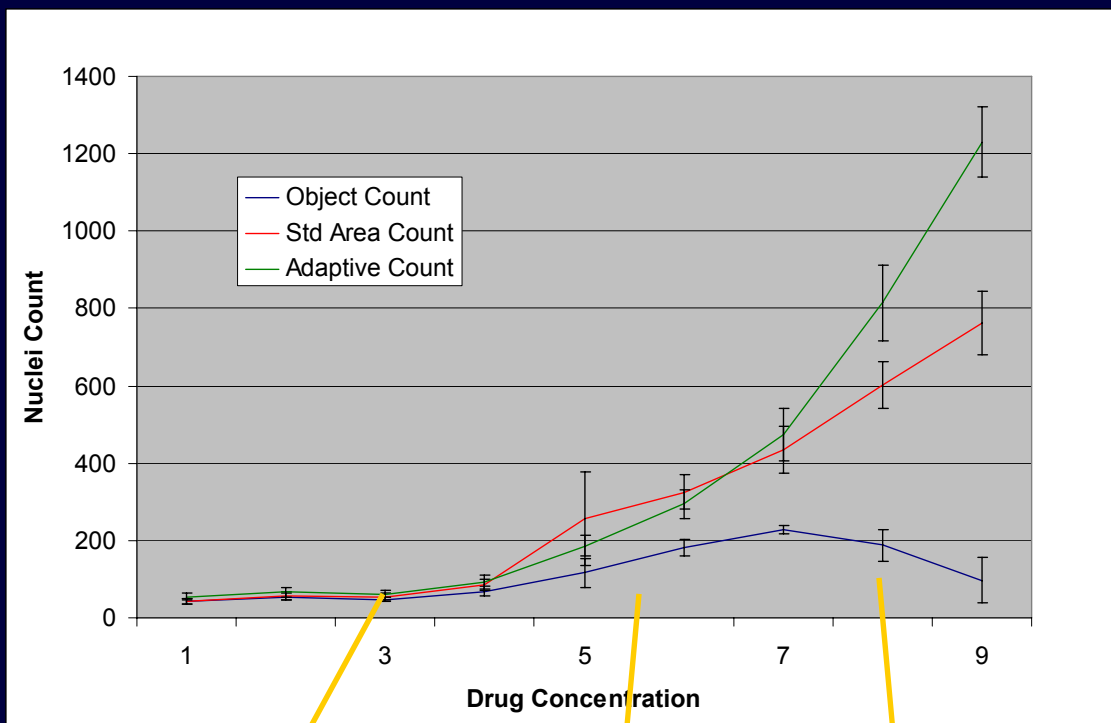


Image Analysis:

Signal / Noise

- Assess

- Optimal exposure times to minimize acquisition time and maximize Signal/Noise

- Image-based autofocus

- Intensity measurements

- Segmentation

- DOE:

- Uniformity Plate

- DAPI

- Multiple exposures

	1	2	3	4	5	6	7	8	9	10	11	12
A	X	X	X	X	X	X	X	X	X	X	X	X
B	X											X
C	X											X
D	X											X
E	X											X
F	X											X
G	X											X
H	X	X	X	X	X	X	X	X	X	X	X	X

Image Analysis: Signal / Noise

- Single Uniformity Plate, 9 sites/well

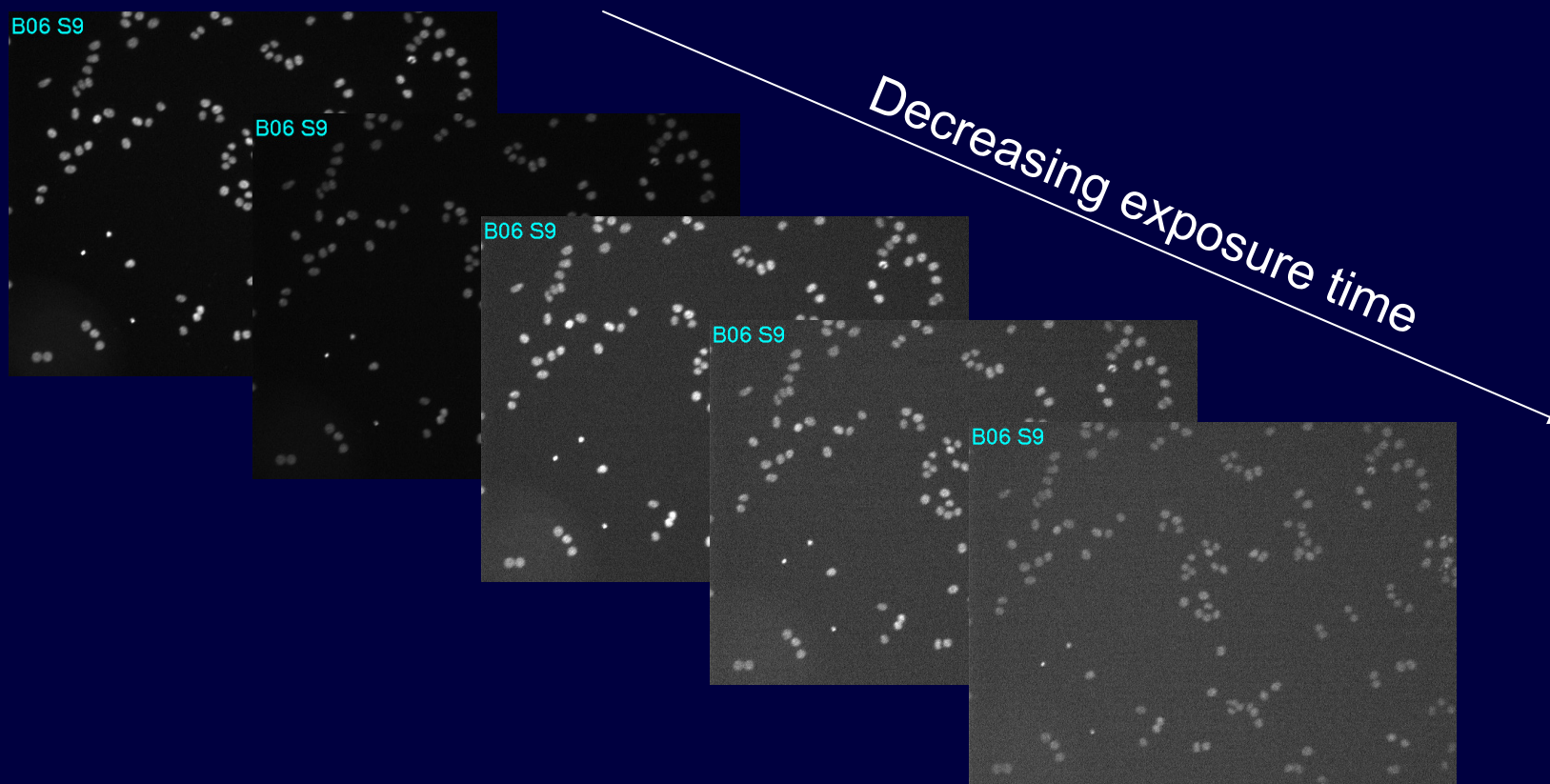


Image Analysis: Signal / Noise

- Effect on thresholding

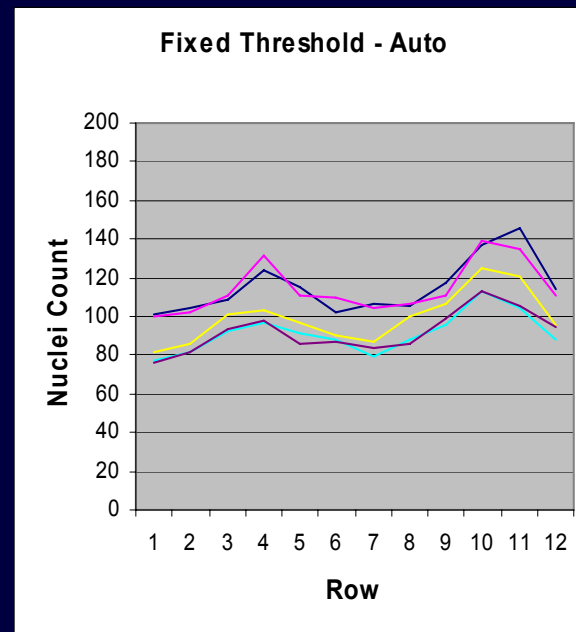
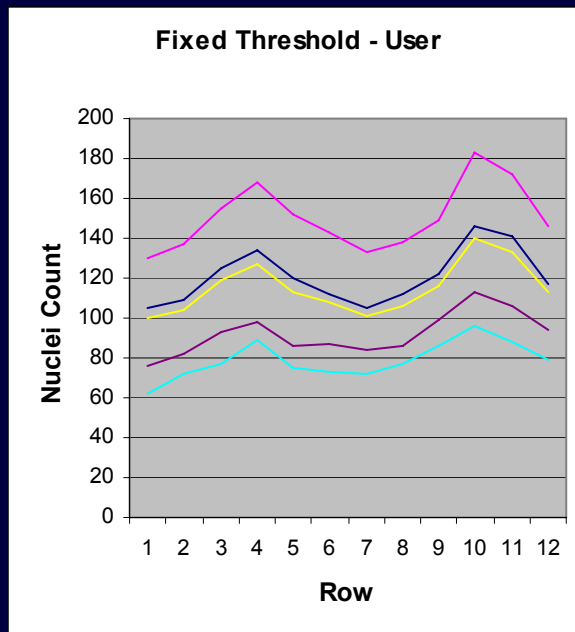
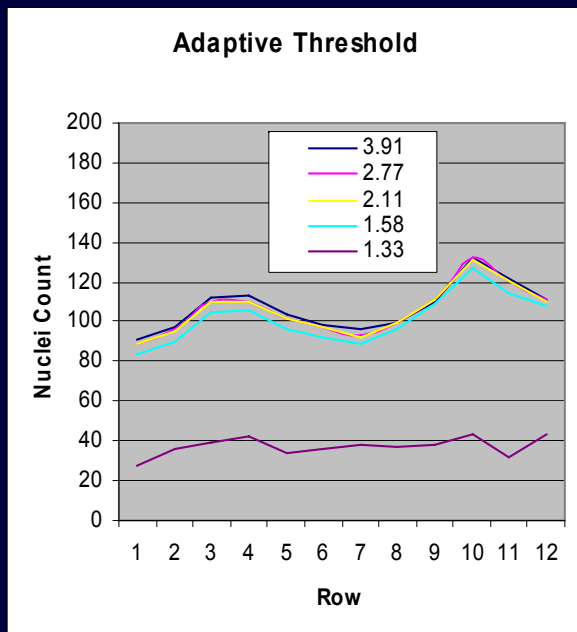
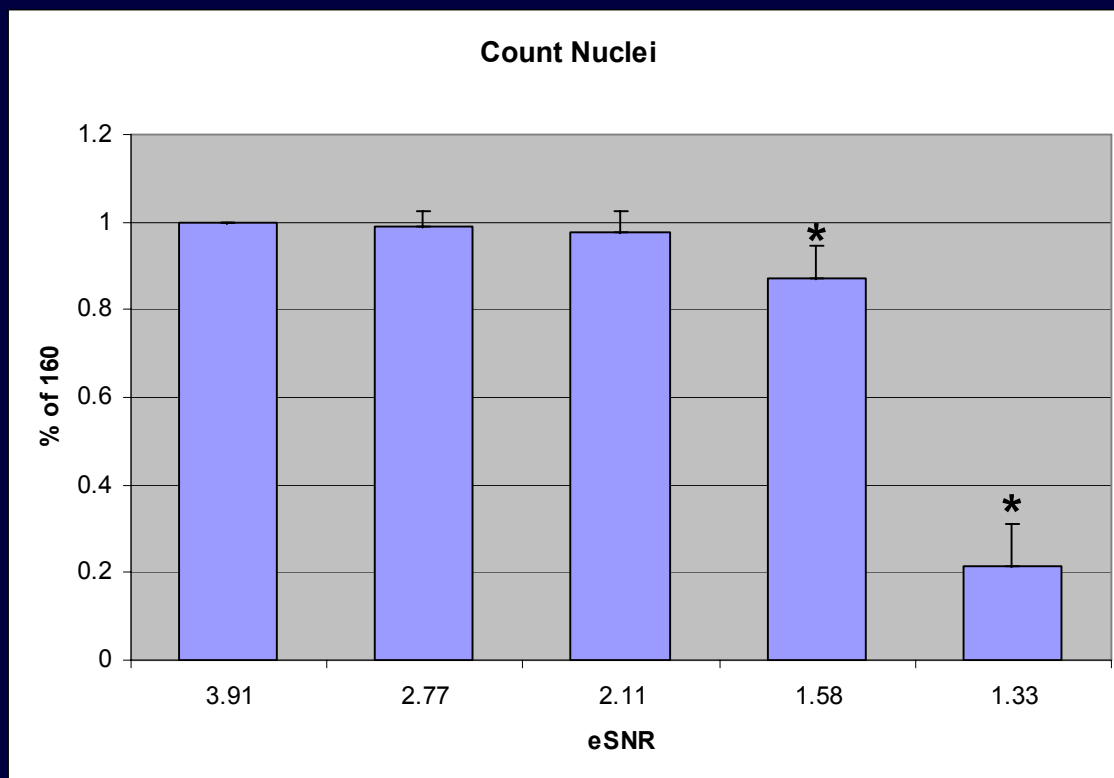


Image Analysis: Signal / Noise



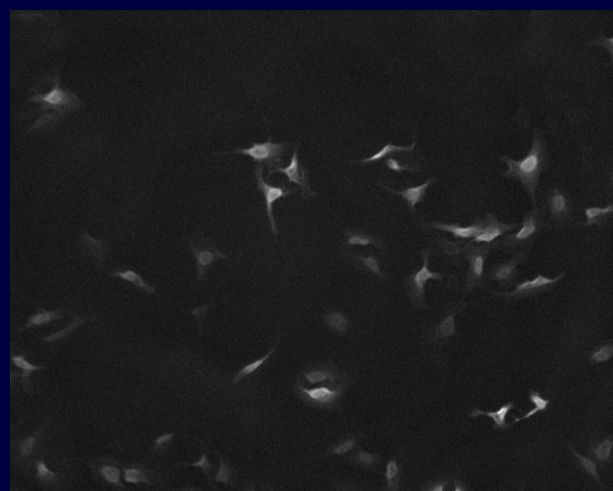
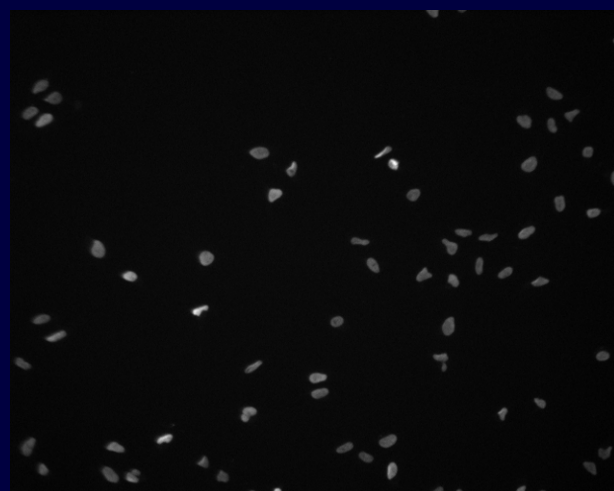
*: $p < 10^{-100}$,
Welch's two-tailed t-test
Error bars are standard deviations

Image Analysis:

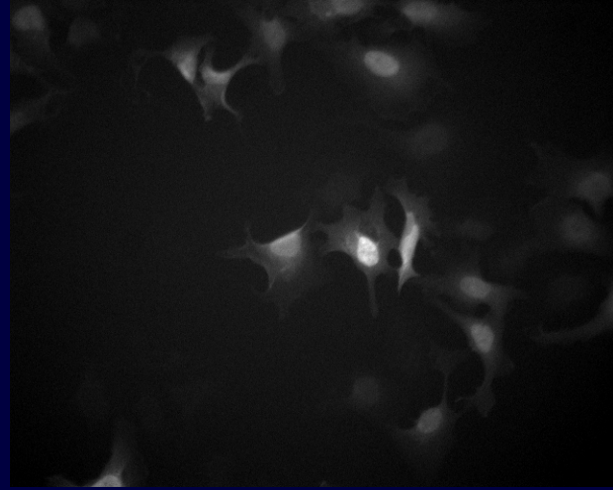
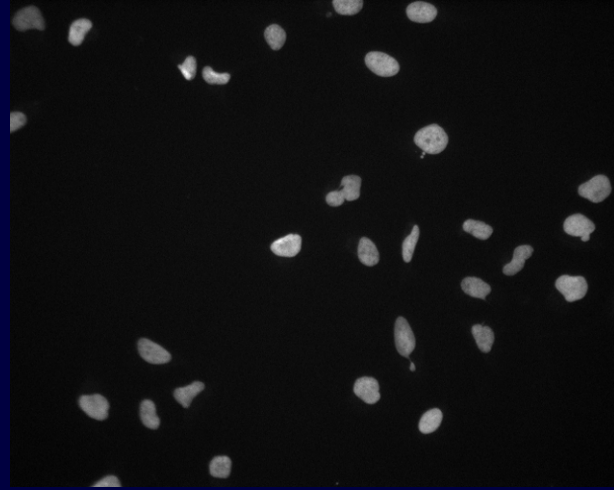
Magnification

- What objective to use?
 - Segmentation accuracy
 - Dose response curve re-sampling

Image Analysis:
Magnification



10x



20x

DNA

FKHR-GFP

Image Analysis:
Magnification

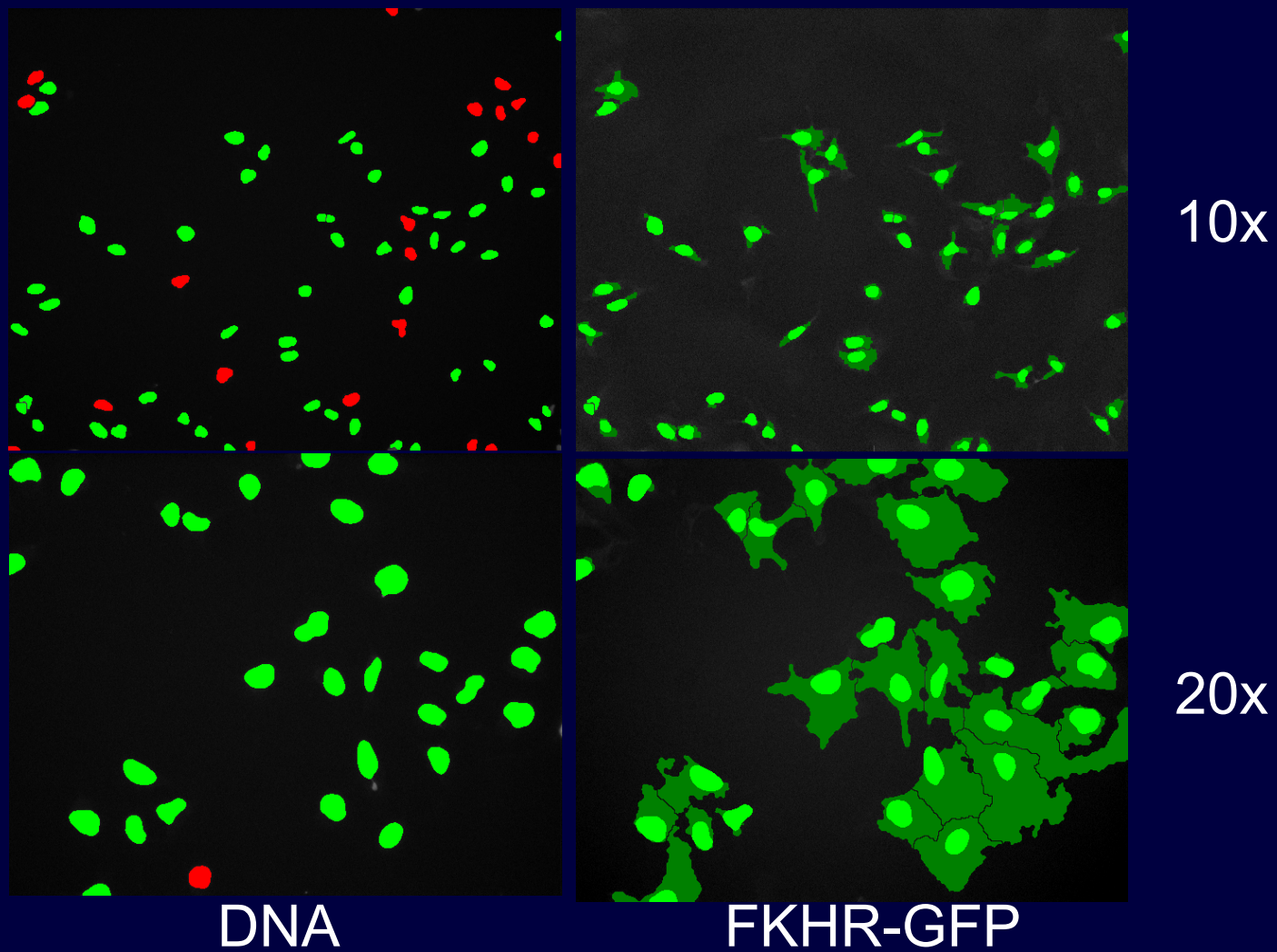
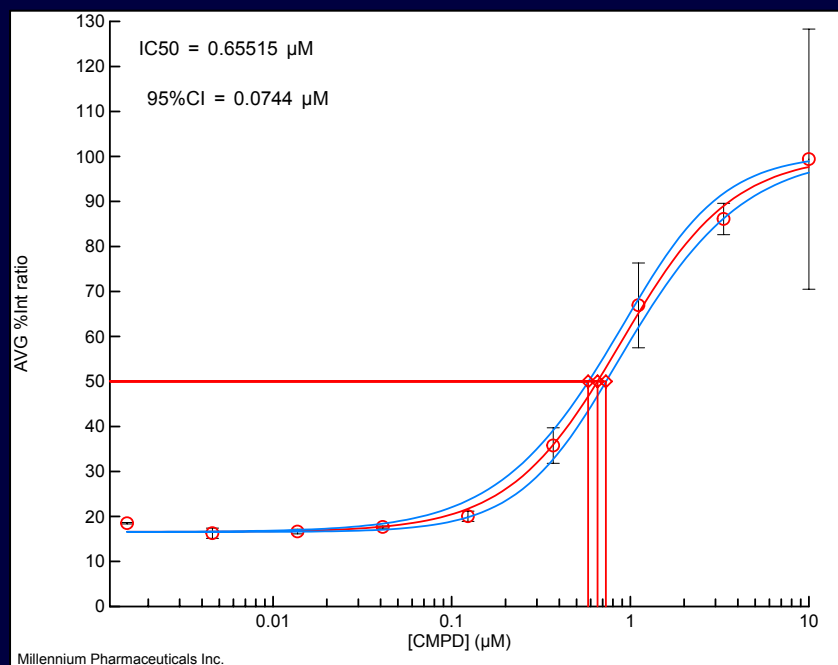
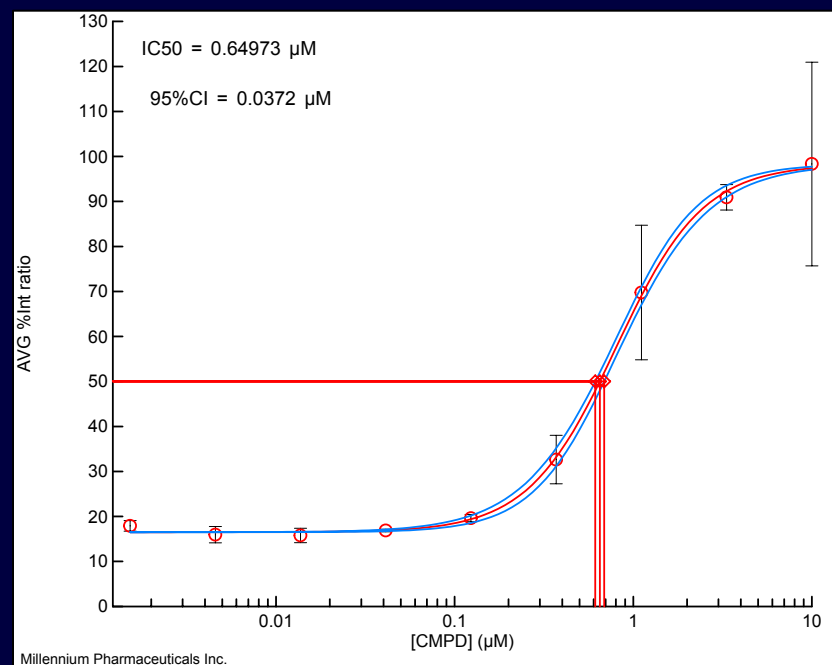


Image Analysis: Magnification



10x



20x

Image Analysis:

Algorithm parameters

- Assess
 - Effect of variability of user-defined parameters
 - Different software algorithms (vendors)
- DOE:
 - Dose Response
 - Multiple algorithm parameters

	1	2	3	4	5	6	7	8	9	10	11	12
A	X	X	X	X	X	X	X	X	X	X	X	X
B	X	← dose response →									C	X
C	X	← dose response →									C	X
D	X	← dose response →									C	X
E	X	← dose response →									C	X
F	X	← dose response →									C	X
G	X	← dose response →									C	X
H	X	X	X	X	X	X	X	X	X	X	X	X

Image Analysis: Algorithm parameters

- User variability?

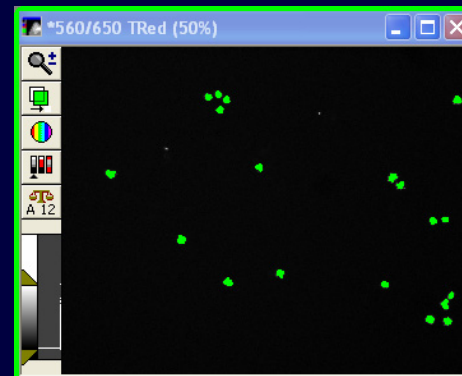
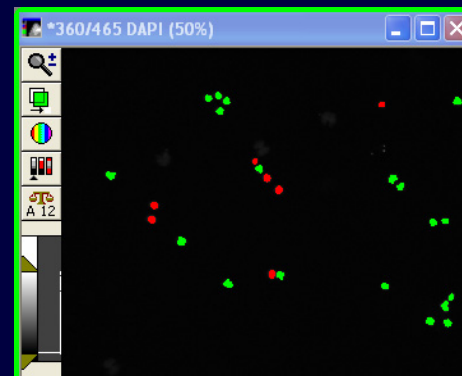
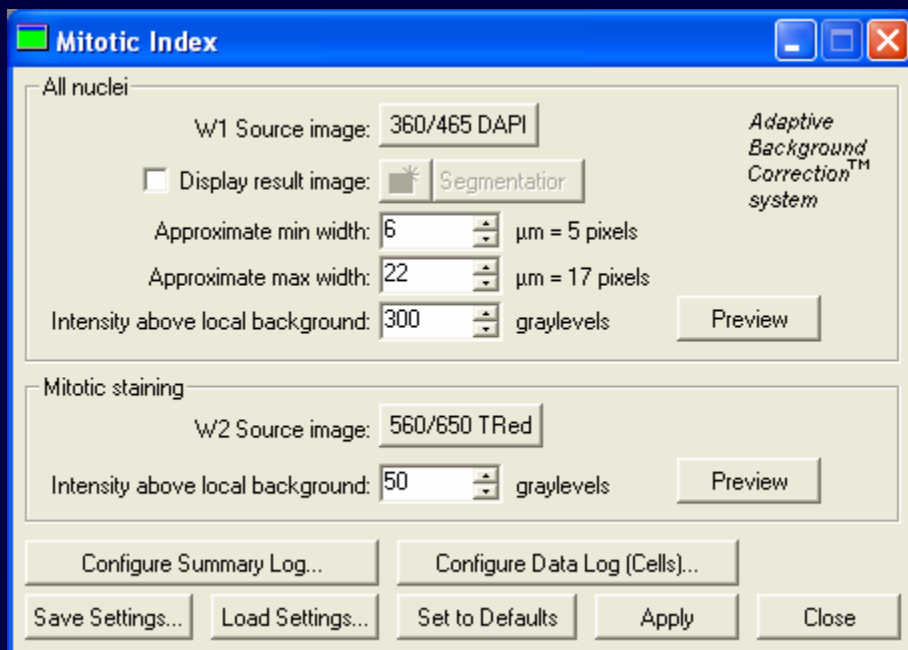
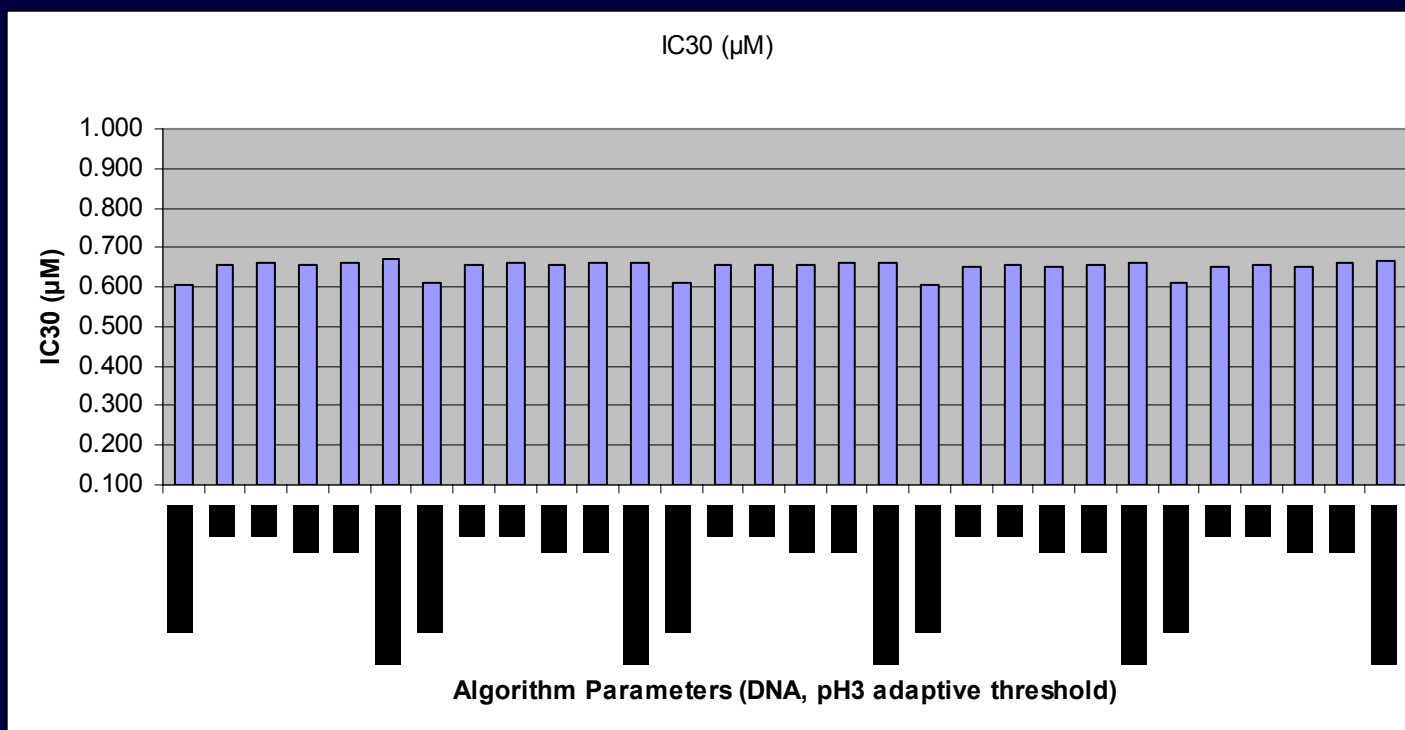
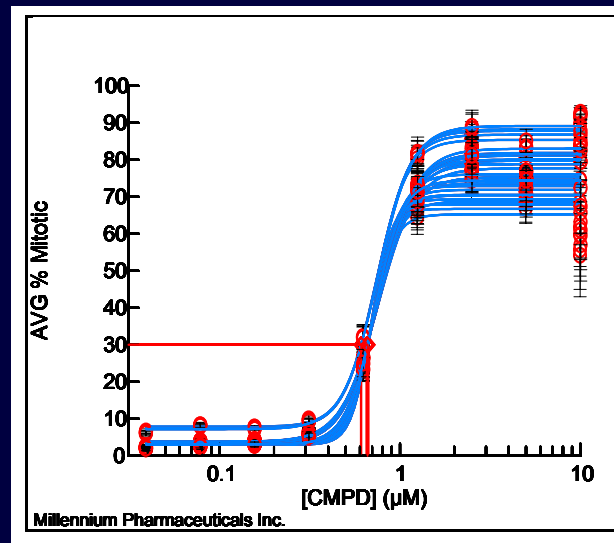
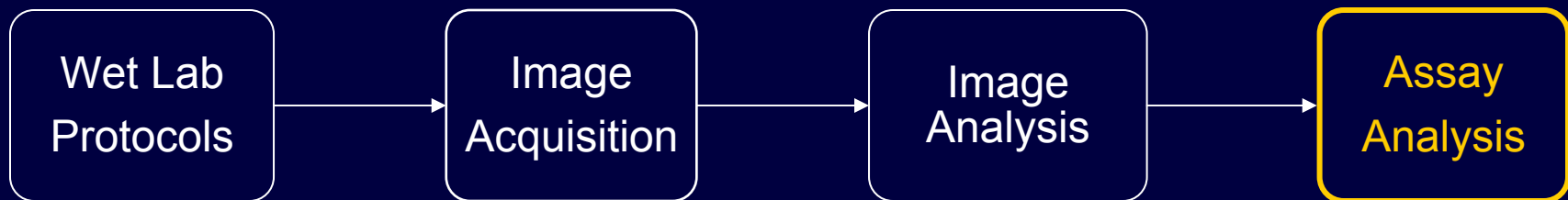


Image Analysis: Algorithm parameters

- Single Plate, 9 sites/well, dose-response
- 30 variations of DNA, pH3 threshold

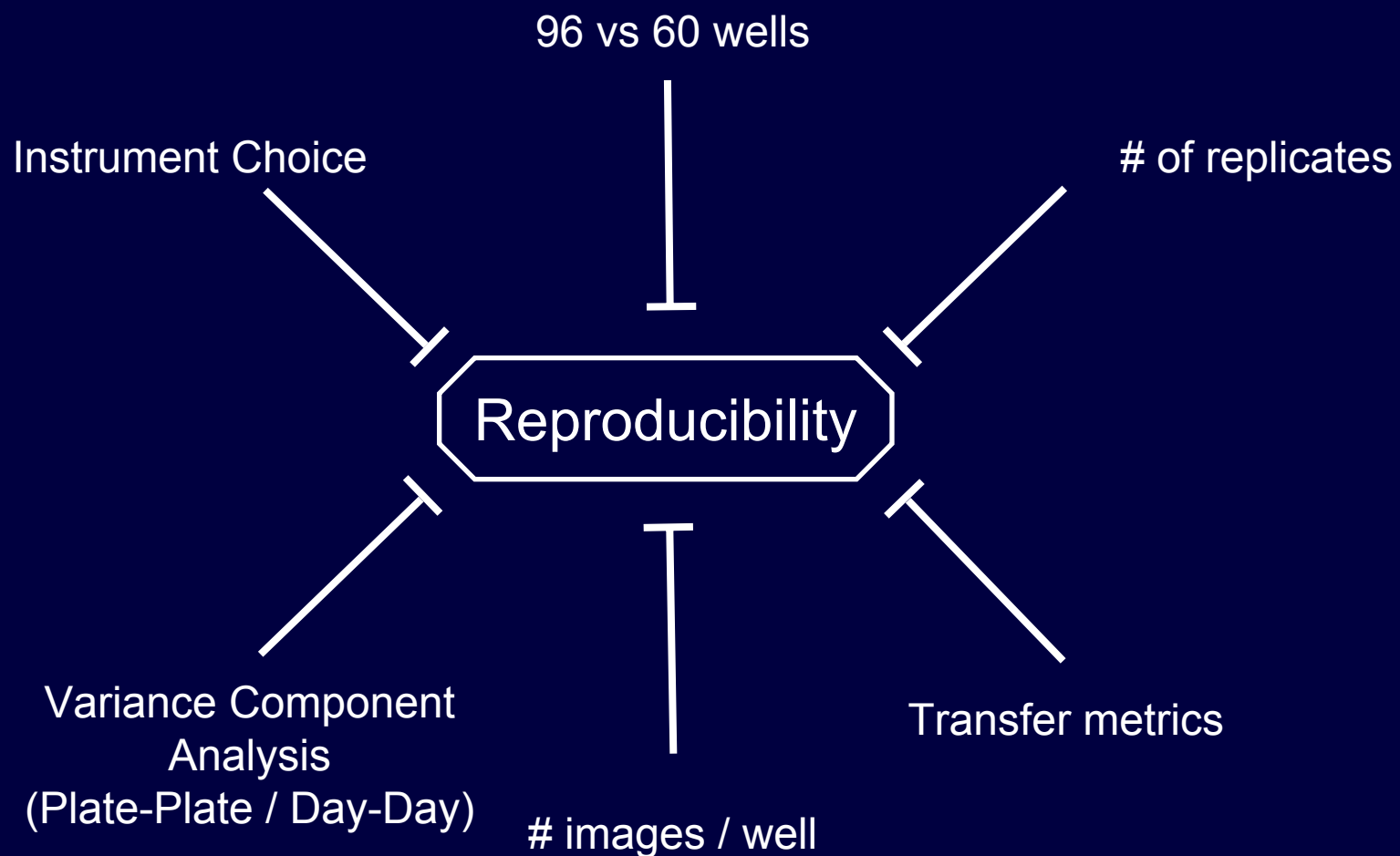


HCS Assay Components

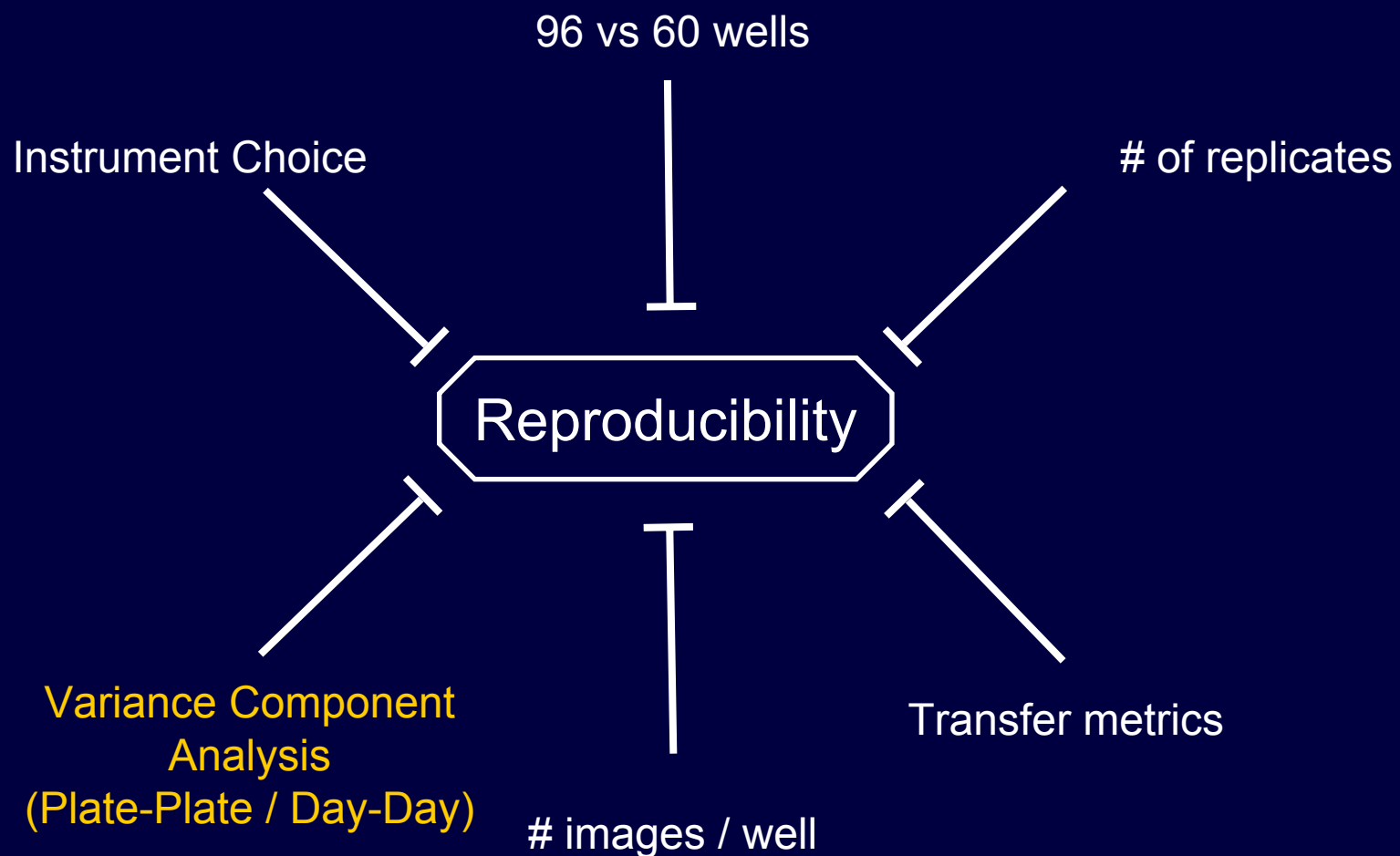


- Dissect components
- Using control experiments
 - Identify drivers of variability
 - Minimize them experimentally or account for them statistically

Assay Analysis



Assay Analysis

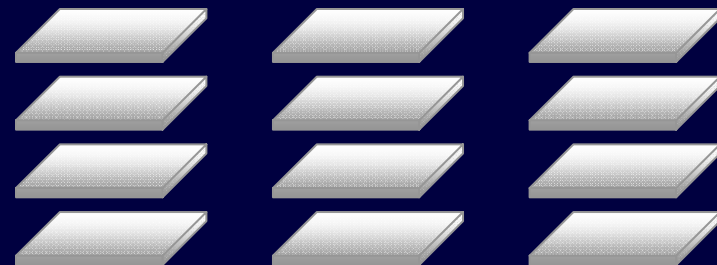


Assay Analysis:

Variance Component Analysis

- Assess variability

- Row / column
- Plate-plate
- Day-day



Day 1

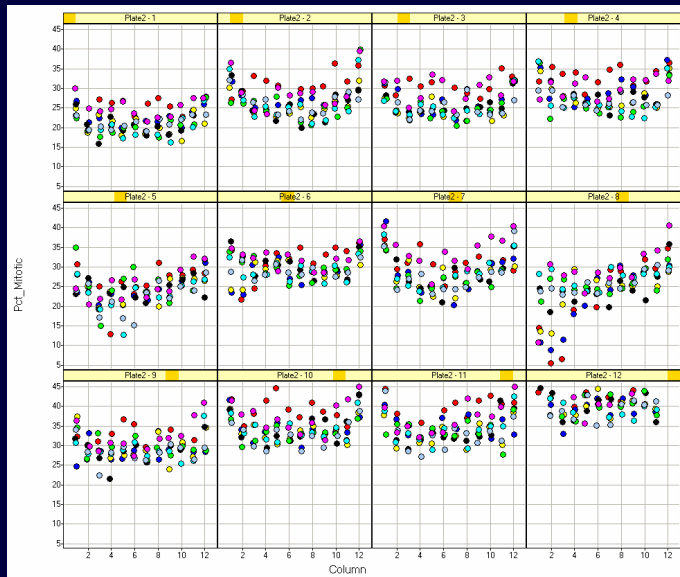
Day 2

Day 3

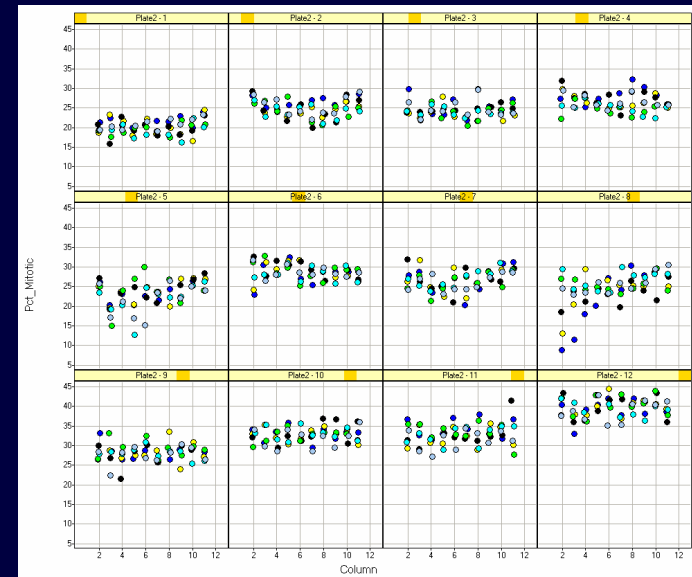
- DOE:

- Single drug concentration (Uniformity plates)
- 4 plates x 3 days

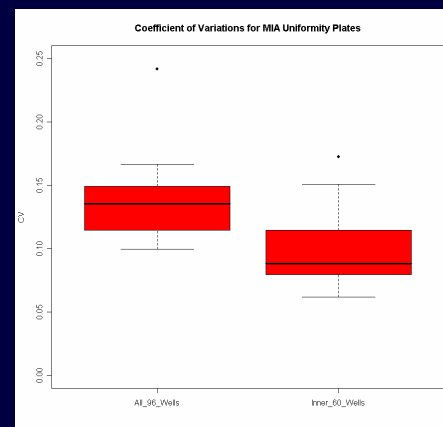
Assay Analysis: Uniformity Data



96 wells



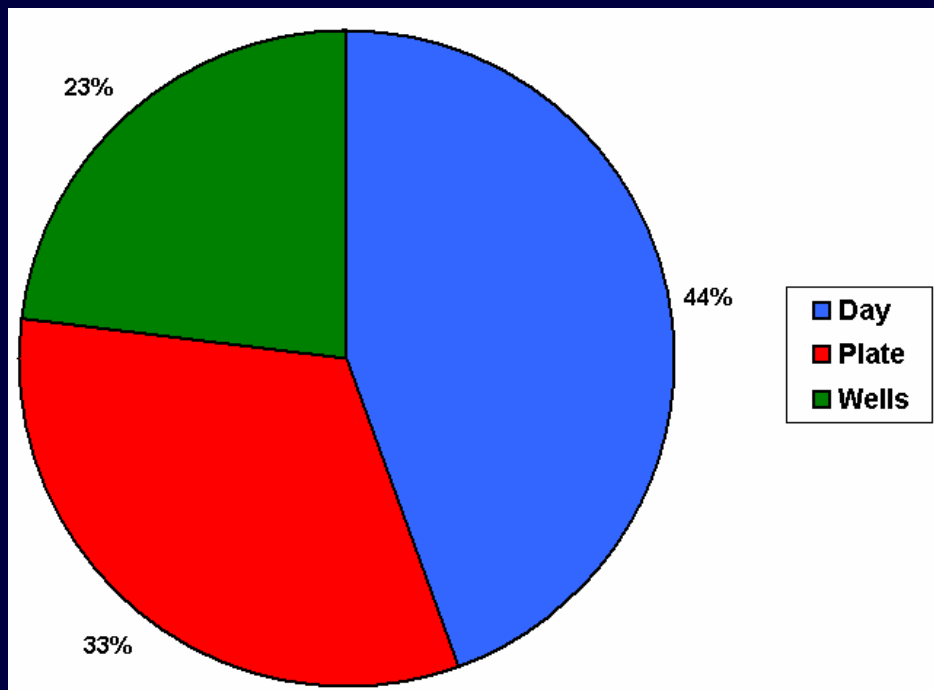
60 wells



- Inner 60 wells had median CV which was 35% lower than that for all 96 wells

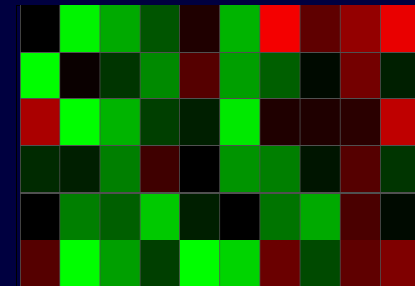
Assay Analysis: Uniformity Data

Variance Component Analysis



12 Plates

- Variability among days was largest component followed by plate
- Replicate dose response curves should be run on different plates on different days to account for largest amount of variability.



Single Plate

Total Mean: 27.7

Total Variance: 32.0 (std dev: 5.7)

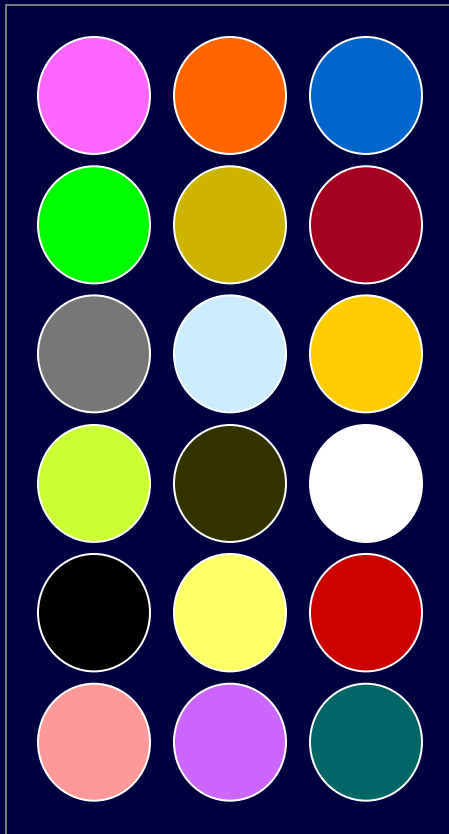
Coefficient of Variation: 20.4%

Assay Analysis: Re-sampling

- Determine
 - How many sites / well?
 - How many replicate wells?
 - Is the assay robust?
 - Does assay perform well across facilities?
- Sample
 - Uniformity Plates
 - Control Compound
 - Multiple Compounds

Bootstrapping

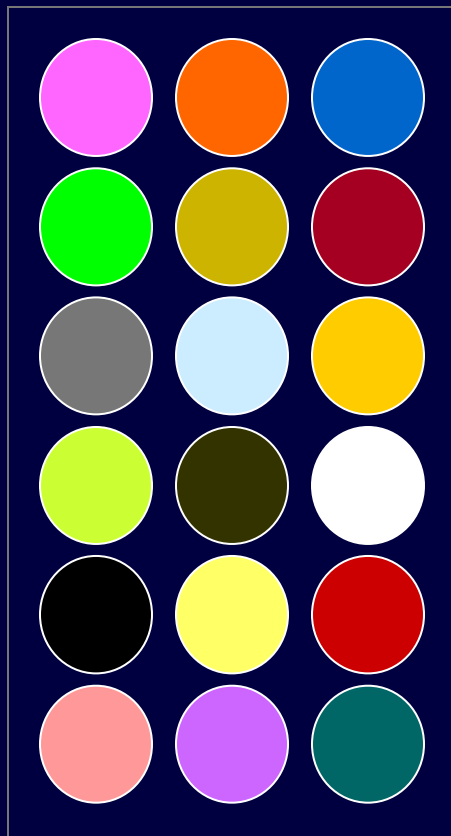
- Method for estimating the sampling distribution of an estimator by re-sampling with replacement from the original sample



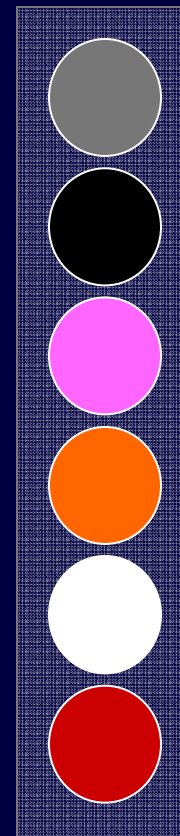
Population

Bootstrapping

- Method for estimating the sampling distribution of an estimator by re-sampling with replacement from the original sample



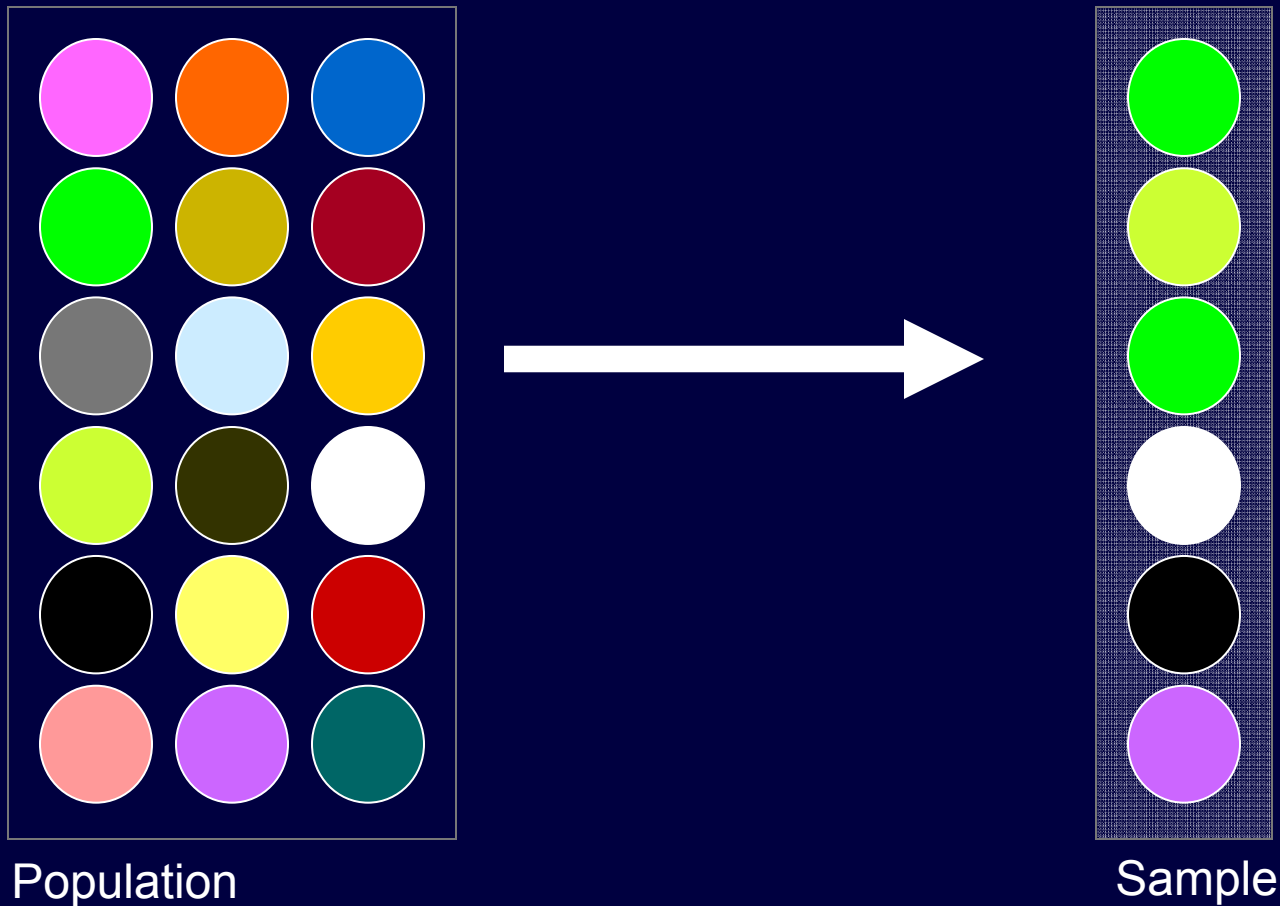
Population



Sample

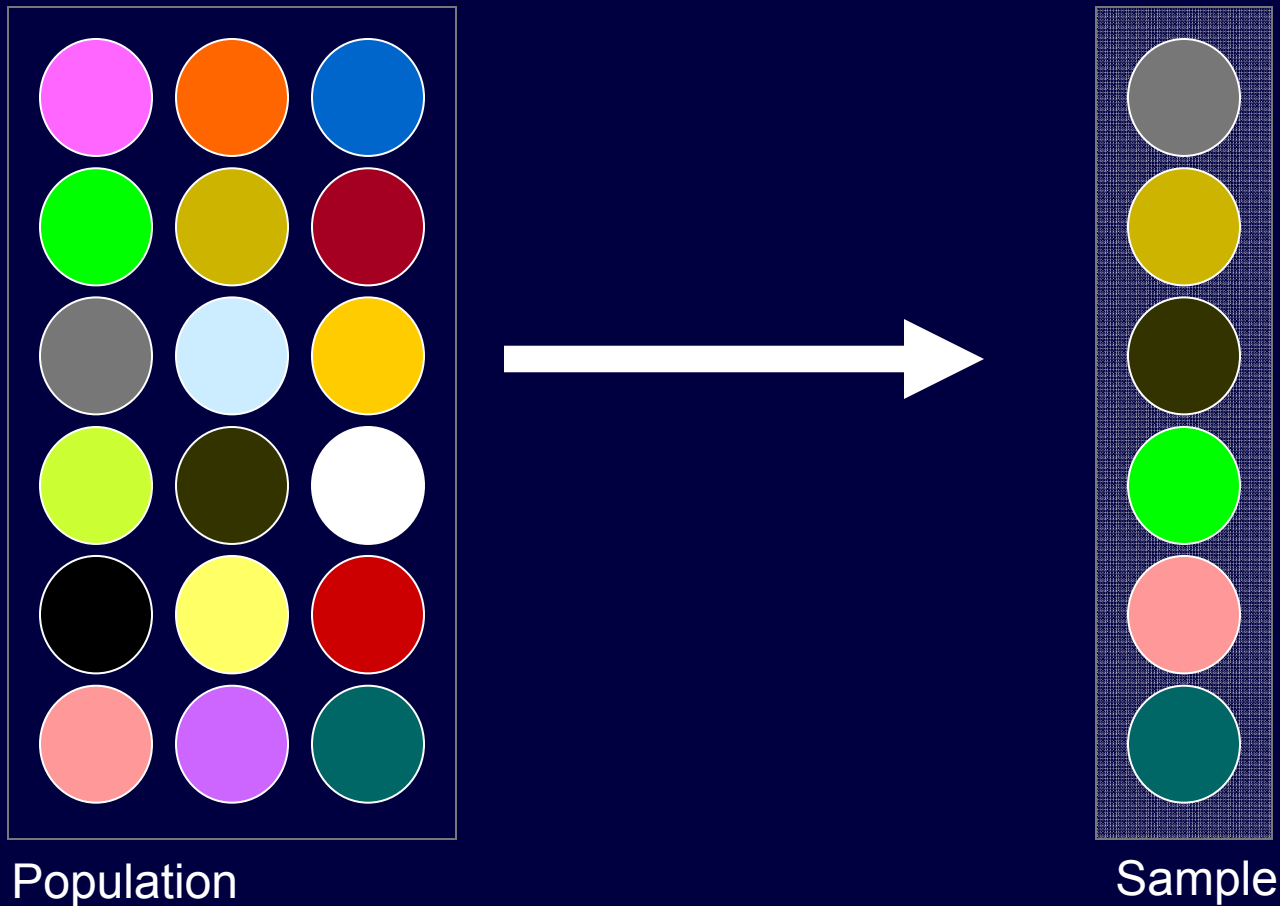
Bootstrapping

- Method for estimating the sampling distribution of an estimator by re-sampling with replacement from the original sample



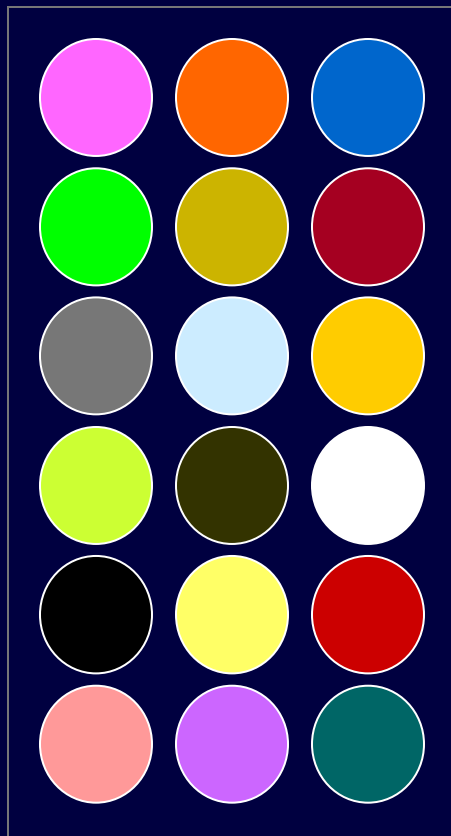
Bootstrapping

- Method for estimating the sampling distribution of an estimator by re-sampling with replacement from the original sample

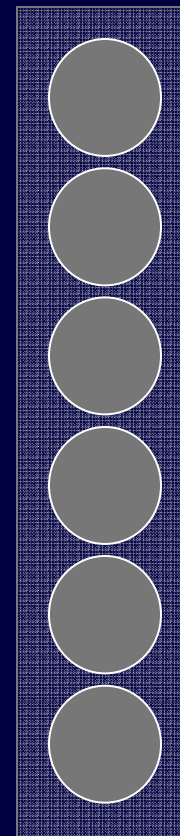


Bootstrapping

- Method for estimating the sampling distribution of an estimator by re-sampling with replacement from the original sample



Population



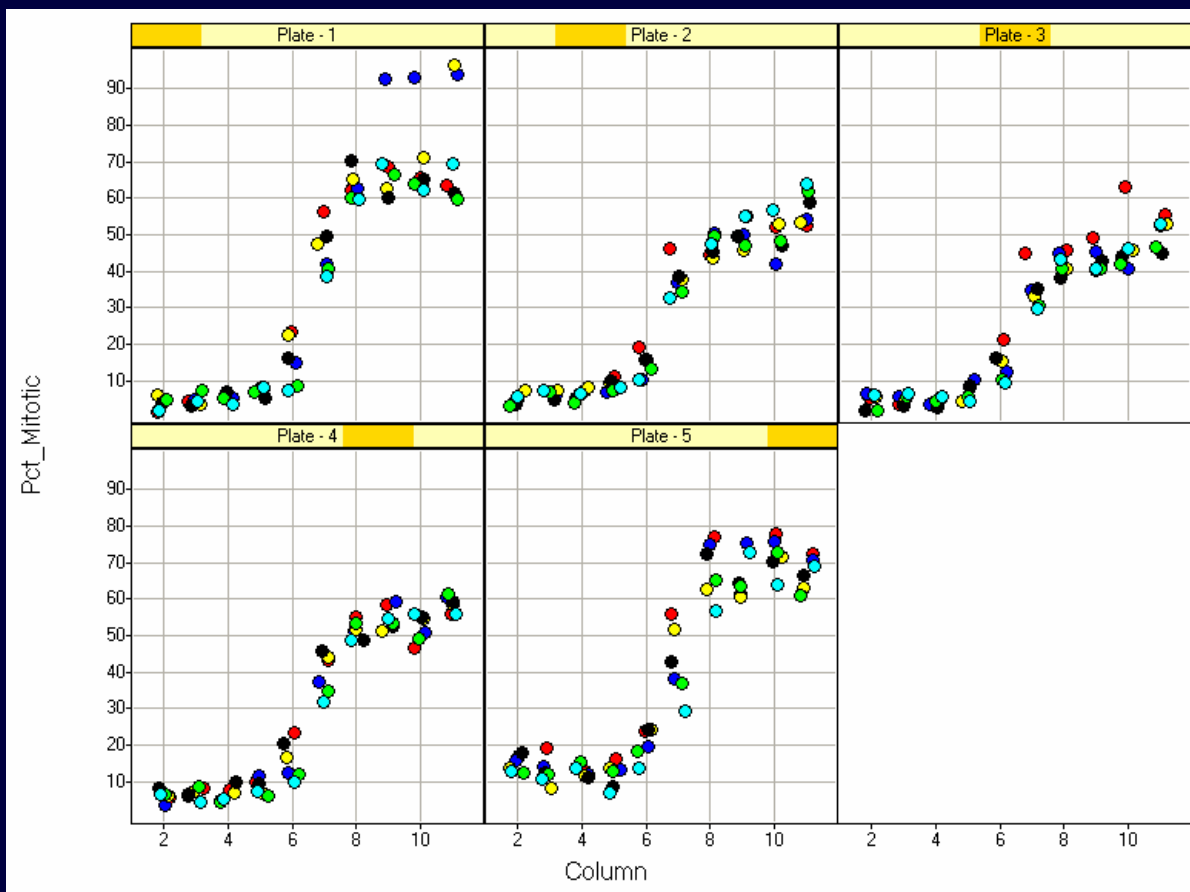
Sample

Assay Analysis: Control Compound

- Assess
 - Effect of number of replicate curves upon EC30 estimates
 - Effect of number of images per well upon EC30 estimates
- DOE:
 - 6 dose response curves
 - 5 plates

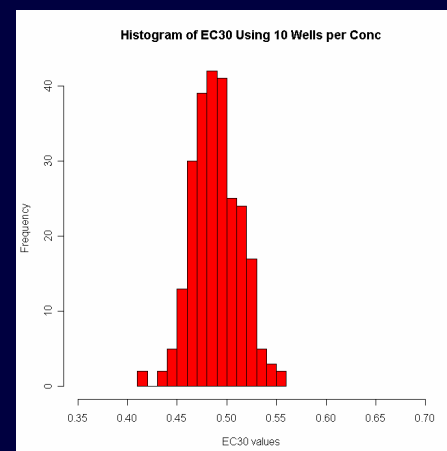
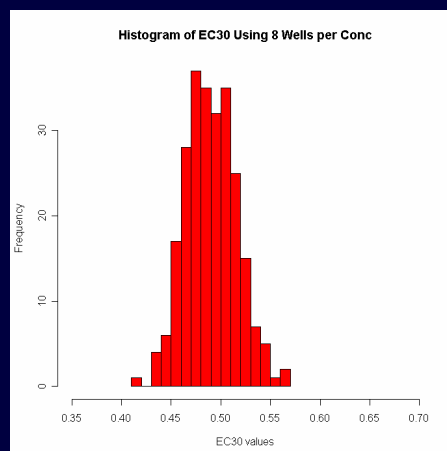
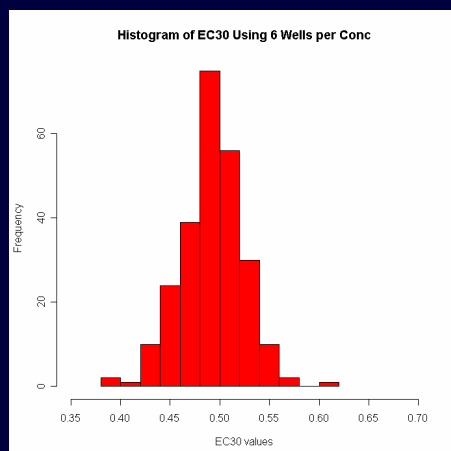
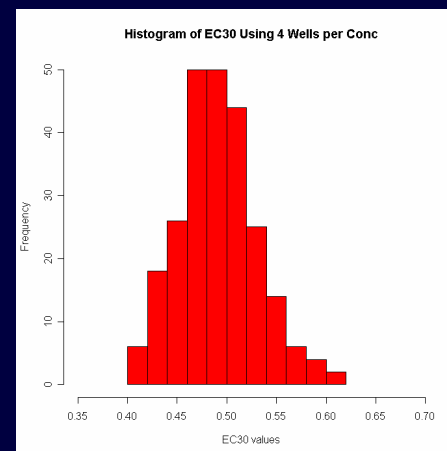
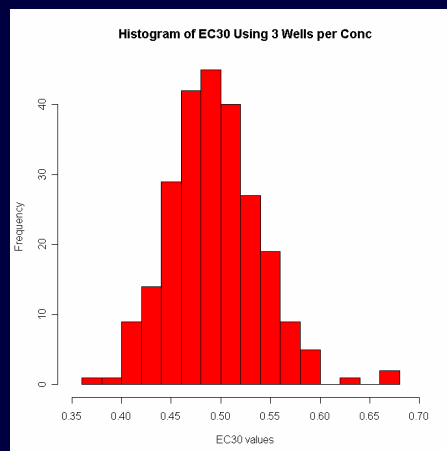
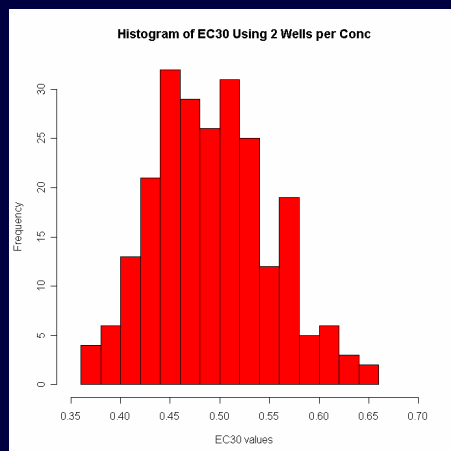
	1	2	3	4	5	6	7	8	9	10	11	12
A	X	X	X	X	X	X	X	X	X	X	X	X
B	X	← dose response →									C	X
C	X	← dose response →									C	X
D	X	← dose response →									C	X
E	X	← dose response →									C	X
F	X	← dose response →									C	X
G	X	← dose response →									C	X
H	X	X	X	X	X	X	X	X	X	X	X	X

Assay Analysis: Control Compound Data

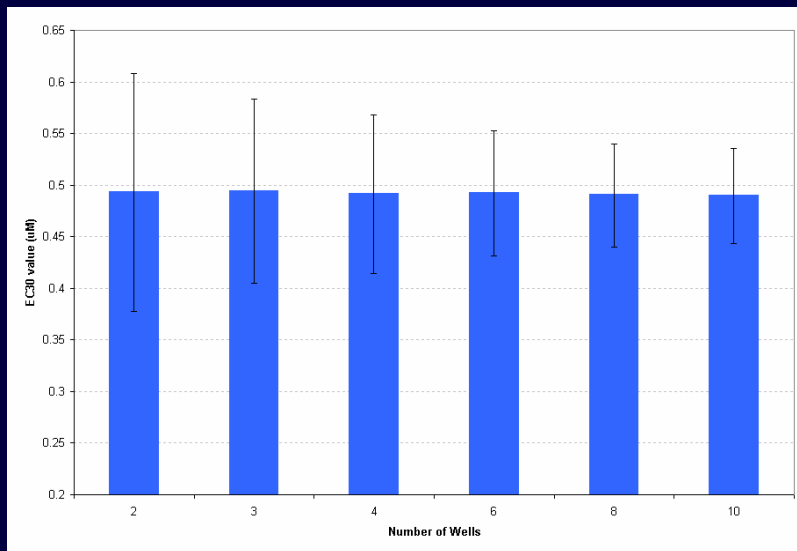


Re-sampling Results

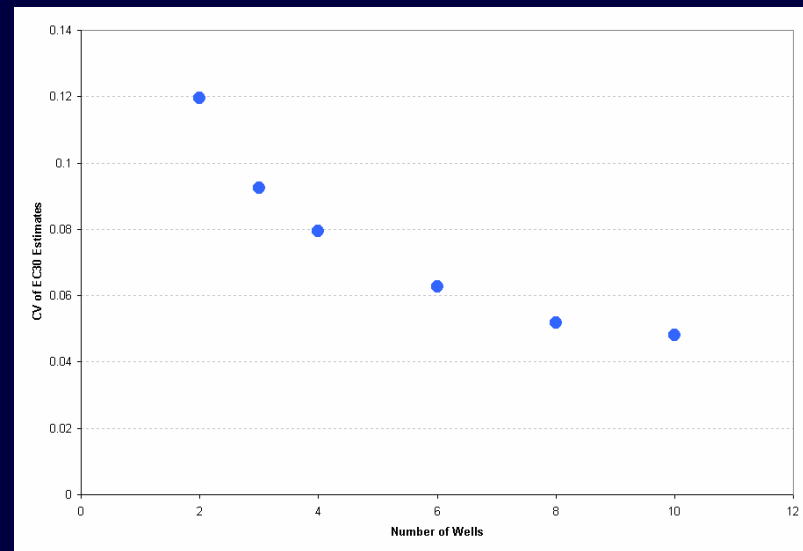
Distribution of EC30 values



Re-sampling Results: Determine # wells



Error bars are 95% CI

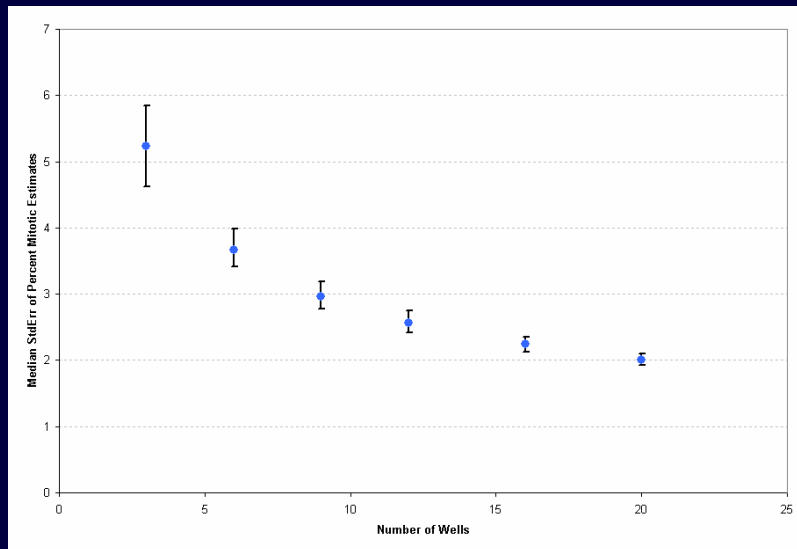


All CVs are less than 12%

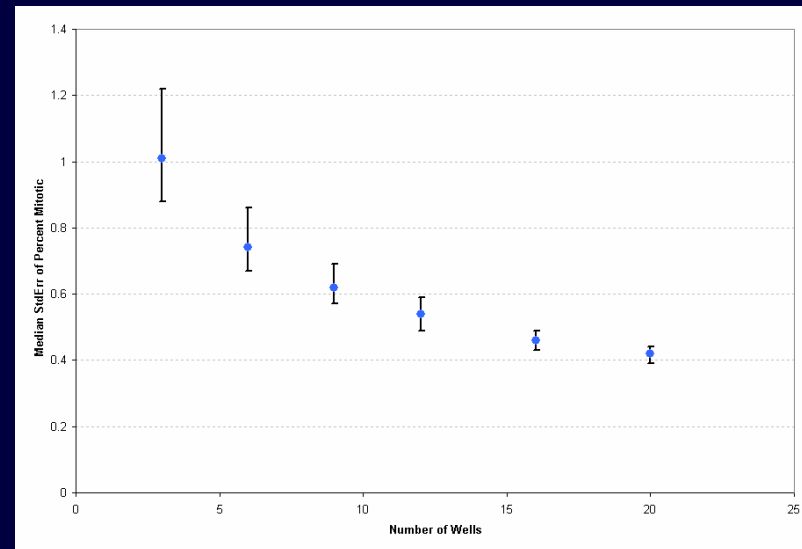
-Results incorporate differences among plates and there were four observations with high values.

Re-sampling Results: Determine # sites/well

10uM

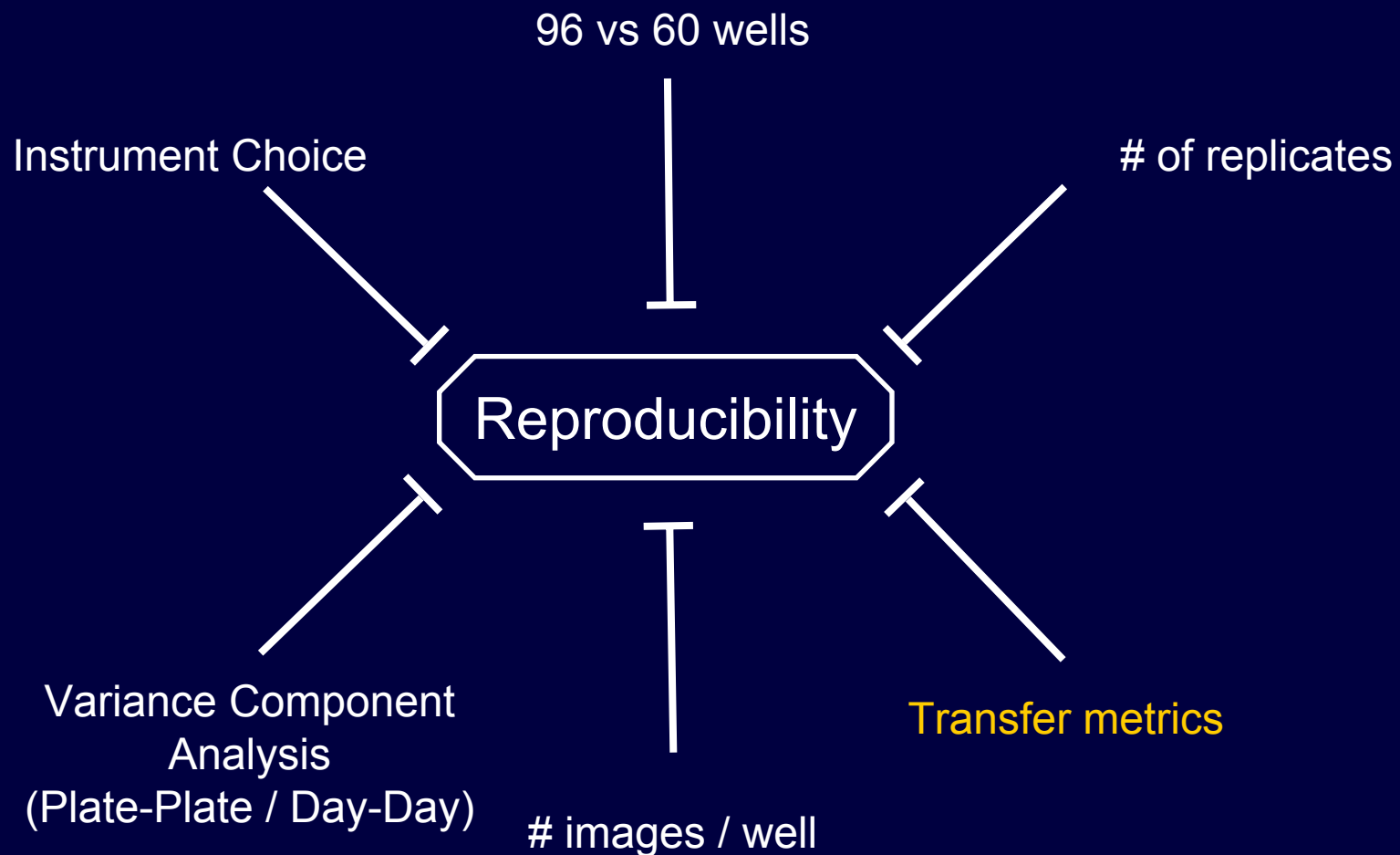


DMSO



Going from 6 images to 9 has almost as much impact on standard error as going from 9 to 16.

Assay Analysis



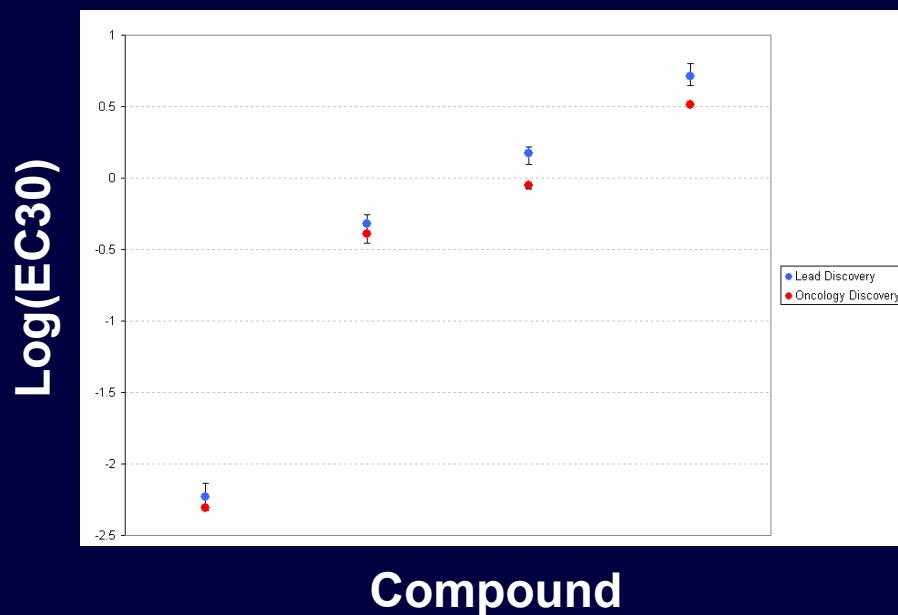
Assay Analysis:

Transfer

- Assess
 - Compare re-sampled EC30 estimates to estimates generated in Oncology Discovery
- DOE
 - 4 compounds
 - 3 replicates / compound
 - Each compound was run on 4 plates = 12 curves
 - All plates run in two facilities

Assay Analysis: Transfer

Four compound re-sampling results

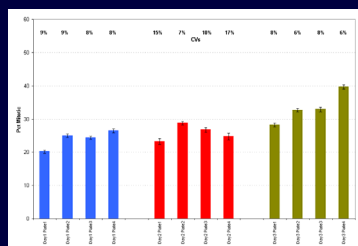


- Three wells from each concentration were sampled.
 - Fit dose response curve and calculated EC30.
 - Total of 250 curves
 - Examine distribution of EC30 values
- A and B were not significantly different, C and D were approximately 60% higher

Assay Analysis:

Instrument selection

- Uniformity Plates

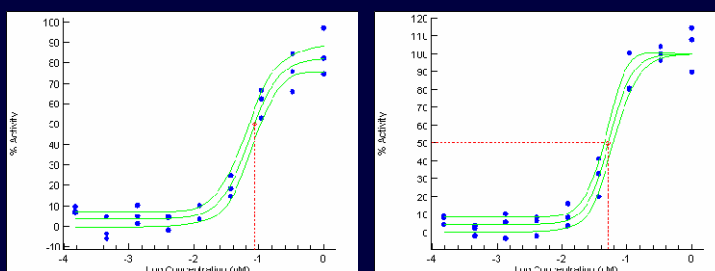


Discovery-1



Opera

- Single Compound



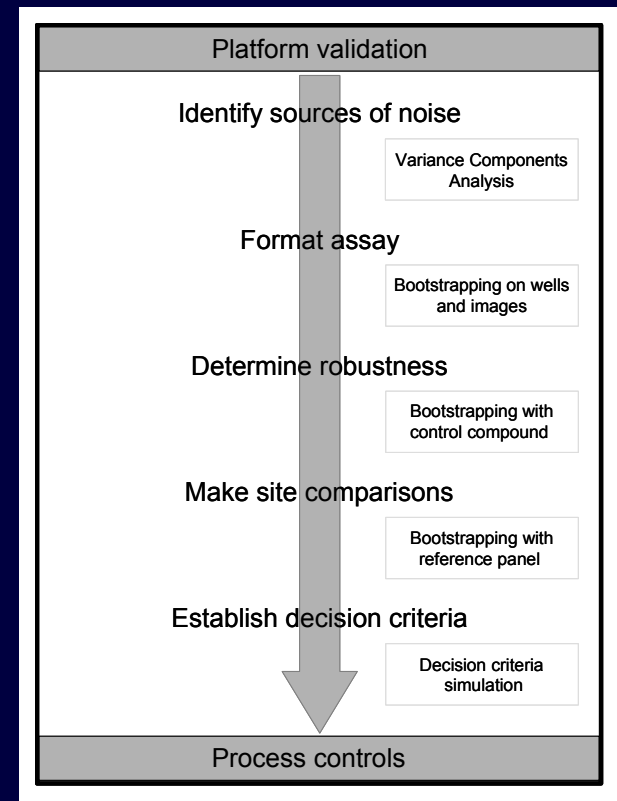
- Signal/Noise, acquisition time, CVs

- Z', EC50 CVs, confidence interval

* determination based on signal/noise, speed, robustness, availability

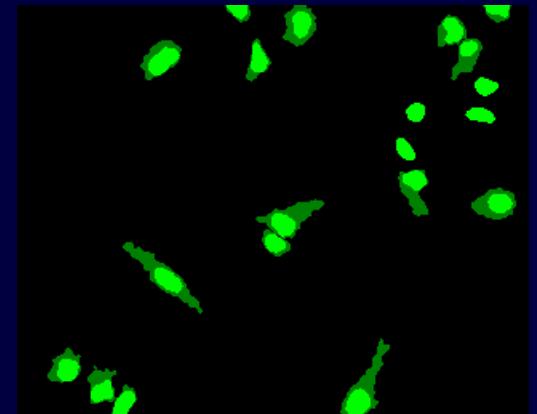
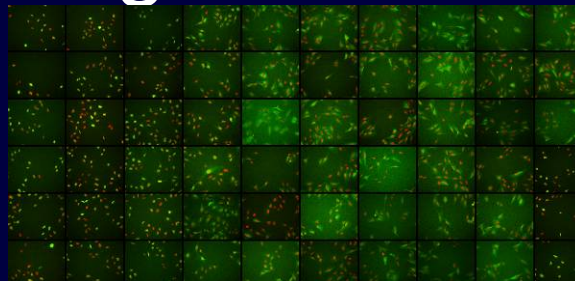
Overall Assay Evaluation

- ~2 months
- Confidence
 - Immunofluorescence staining protocol
 - Automation / liquid handling
 - Instrument / Instrument QC
 - Software Algorithm
 - Robust assay
 - Transfer from Discovery to Platform Group



Summary

- Look at the images



- Run uniformity plates to identify and minimize noise (biology, instrument, and algorithm)
- Use analysis modules throughout process

Process Deliverables

- Development Guidelines
 - QC metrics: instruments, algorithm
 - IF staining protocol optimization
 - Image analysis optimization
- Production Guidelines
 - Statistical analysis for transfer
 - Ongoing QC metrics

Acknowledgements

- Molecular & Cellular Oncology
 - Ben Amidon
 - Jeff Ecsedy
 - Natalie Roy D'Amore
 - Liz Carideo
 - Kara Hoar
 - Claudia Rabino
 - Michelle Tighe-Nestor,
 - Denise Driscoll
 - Deborah Wysong
- Cancer Pharmacology
 - Arijit Chakravarty
 - Katherine Galvin
- Lead Discovery
 - John Ringeling
 - John Donovan
- Computational Sciences
 - Sudeshna Das
 - Mike Pickard

Breakthrough science. Breakthrough medicine.SM

