



Live cell imaging in Target Validation and Lead Optimization

S. Prechtl

AD-HTS / High-Content Analysis

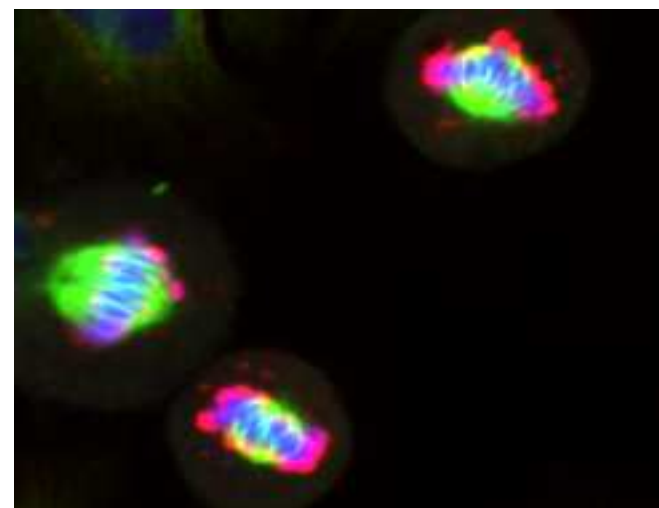
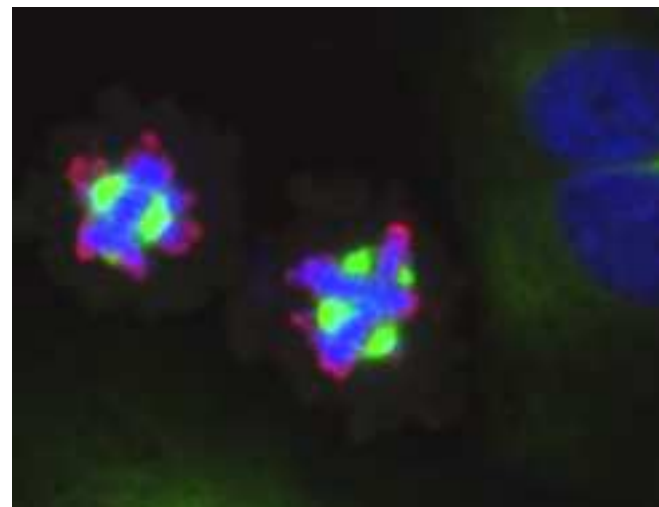
Assays and Cellular Targets (ACT) 2006
October 30 - November 1, 2006 - Green Valley Ranch Resorts - Las Vegas, NV

cell systems in modern drug research

- ▶ highly recommended scientific tools
 - ▶ examination of patho-physiological mechanism
 - ▶ investigation of compound interactions
- ⇒ within the cellular context

High-Content Analysis

- ▶ automated high-resolution imaging & image analysis
 - ⇒ exploits the detailed insight into cell system
- ▶ analysis of structural and metabolic changes
 - ⇒ functional high-quality information about targets and potential drug candidates
- ▶ quantification of complex drug actions at the level of the individual cell
 - ⇒ providing statistically significant results of greater reliability





HCA assays are applicable throughout nearly all stages of the drug discovery process

- ▶ target validation, lead discovery, lead optimization, secondary screening, *in vitro* toxicology studies, preclinical studies

HCA assays have the potential to accelerate drug discovery process by opening current bottlenecks

- ▶ performing higher-throughput cell-based assays providing high-quality data
- ▶ time-saving multiplexed / multilinear designed assays
- ▶ identifying genes or proteins involved in diseases
- ▶ qualifying drugability and therapeutic benefits of selected compounds
- ▶ identifying possible side-effects within cellular context

Target validation

Assay Development

HTS

Hit To Lead

Lead Optimization

High-Content Analysis

equipment consists of fully automated confocal HT imaging system, MT epi-fluorescence imager, motorized fluorescence microscope



OPERA
Evotec Technologies



Discovery 1
Molecular Devices

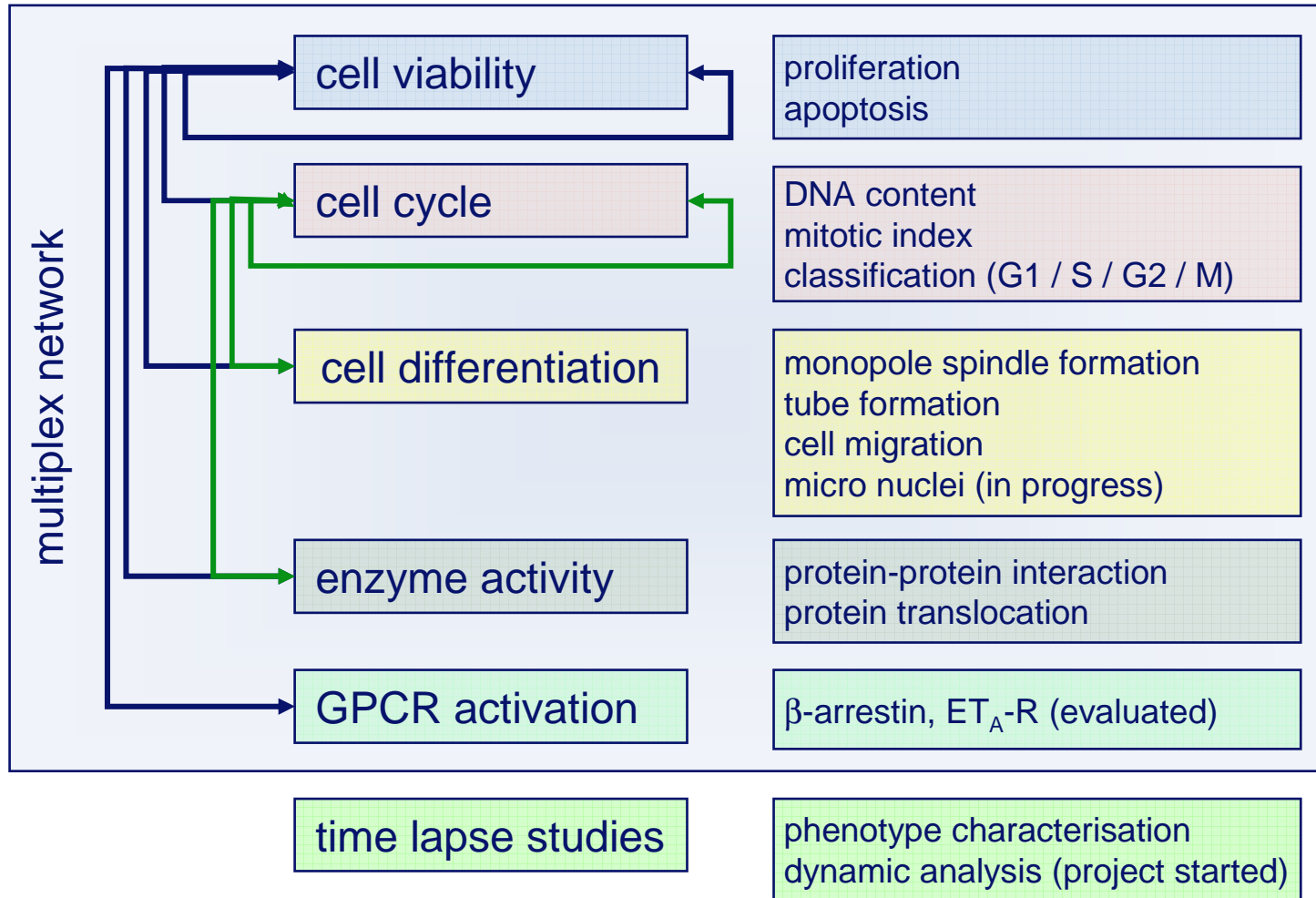


Axiovert 200
ZEISS



application panel

cell systems



- HeLa
- CHO
- DU145
- PC3
- LNCap
- MCF7
- T47D
- BT474
- SK-OV-3
- U-2 OS
- HUVEC
- MVEC
- PAEC
- S49
- Jurkat
- prim. MΦ

Target
validation

Assay
Development

HTS

Hit To Lead

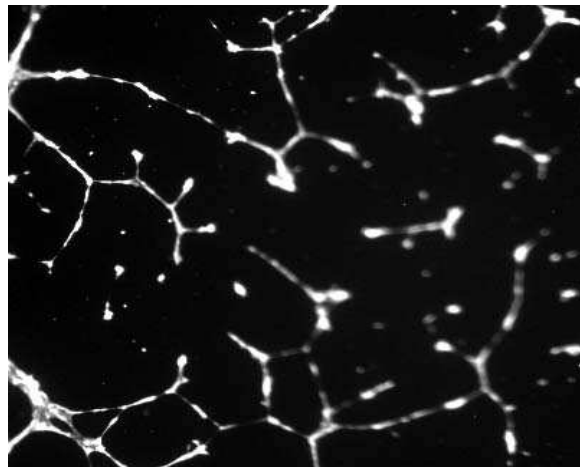
Lead
Optimization

High-Content Analysis

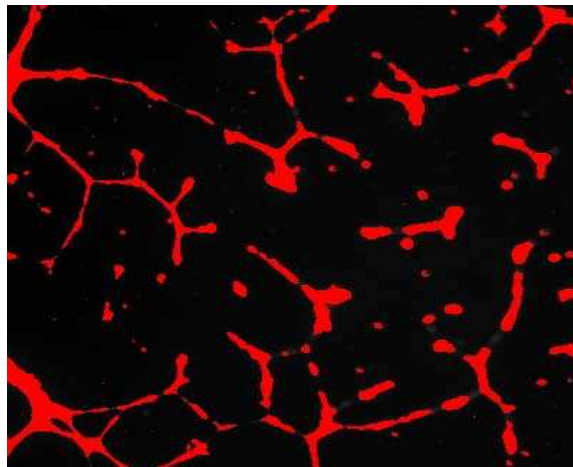
tube formation

- cellular model to quantify the differentiation of endothelial cells into capillary-like structures -> compounds modulating angiogenesis
- => automated *in vitro* tube formation assay enabling the screening of potential angiogenesis stimulators and inhibitors
- quantification of the extent of tube formation by measuring the total tube length

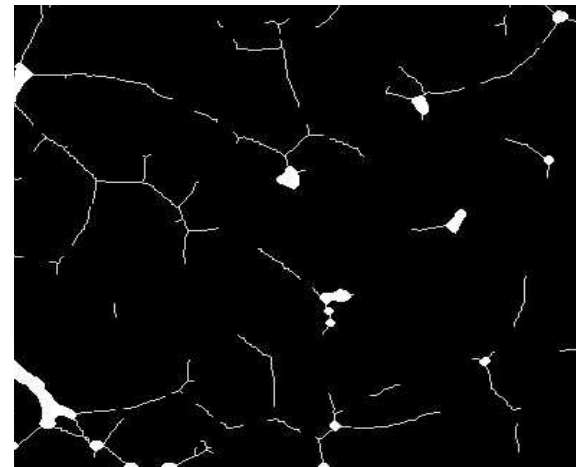
tube image



thresholding



analysis image



Target validation

Assay Development

HTS

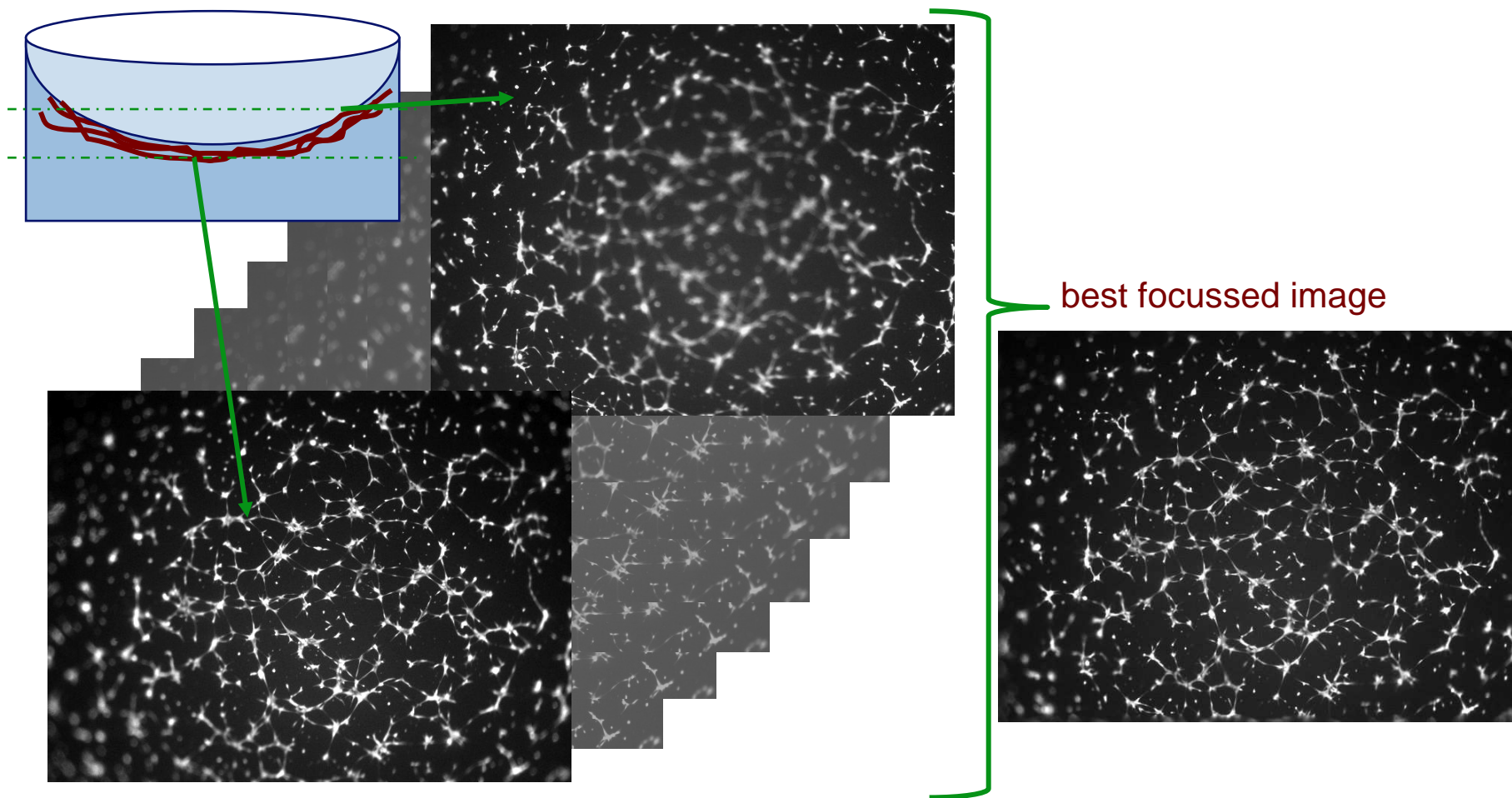
Hit To Lead

Lead Optimization

High-Content Analysis

“best focus” algorithm

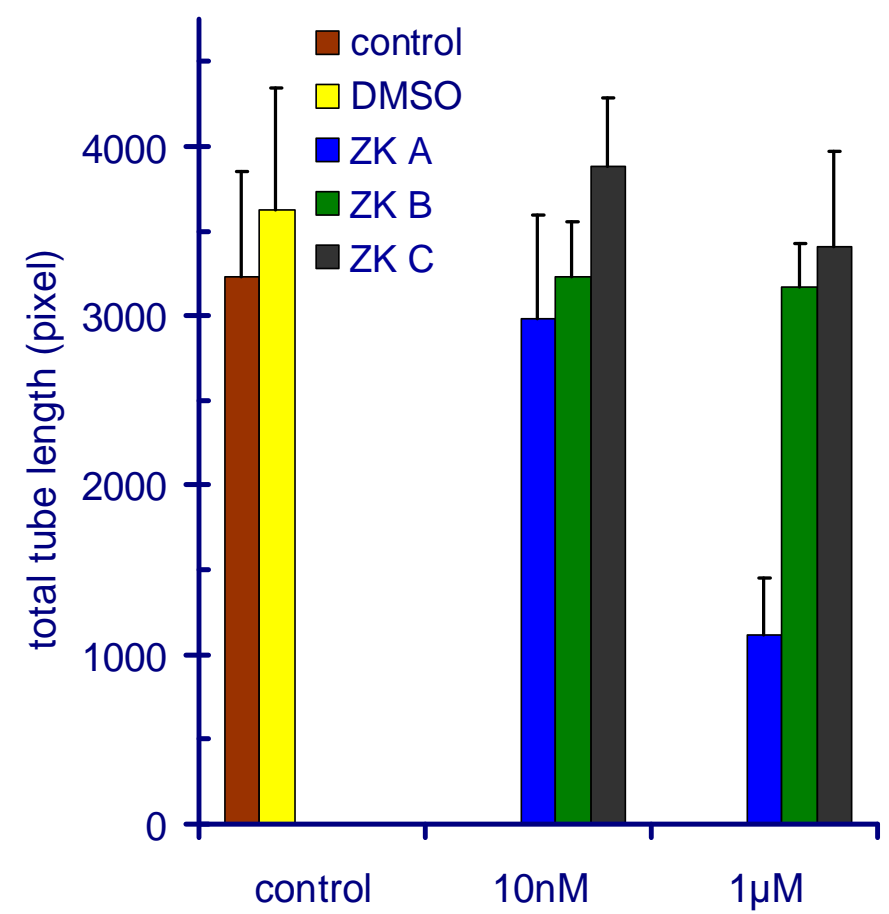
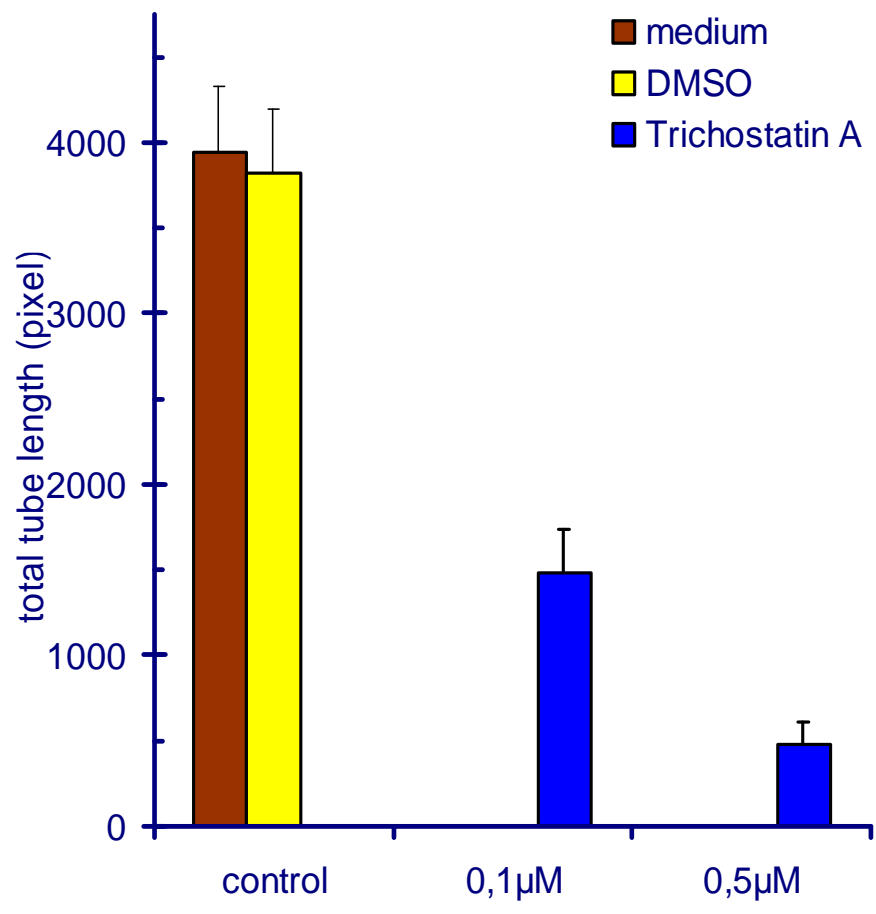
tube formation





MVEC

tube formation



Target validation

Assay Development

HTS

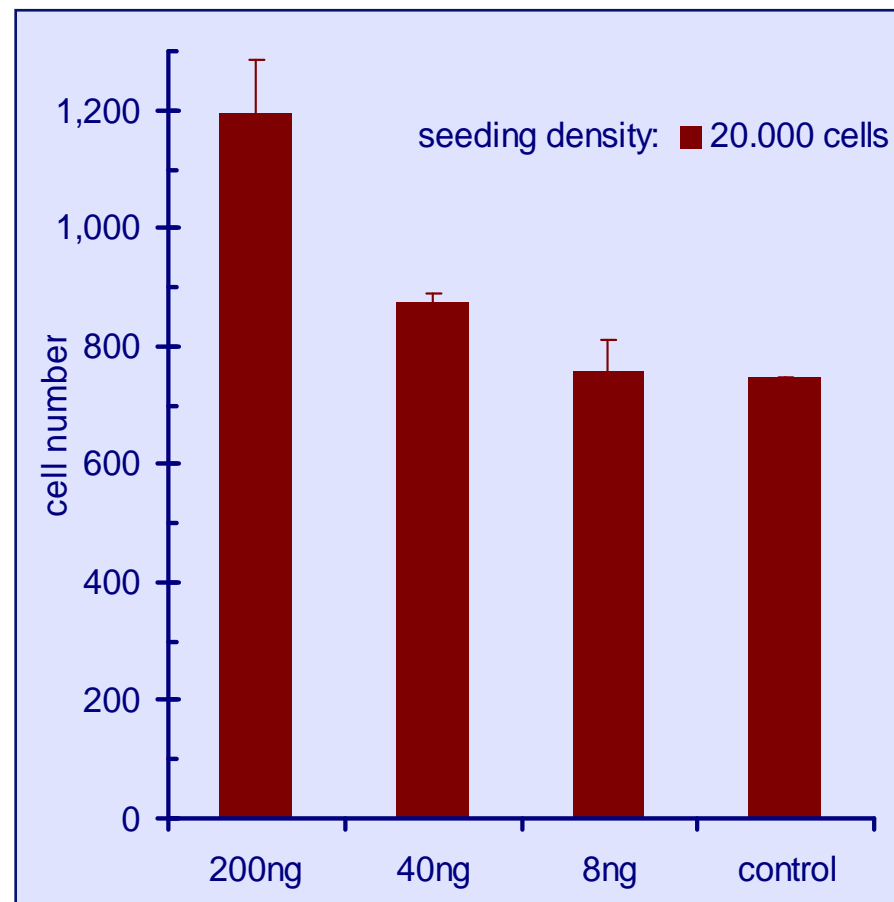
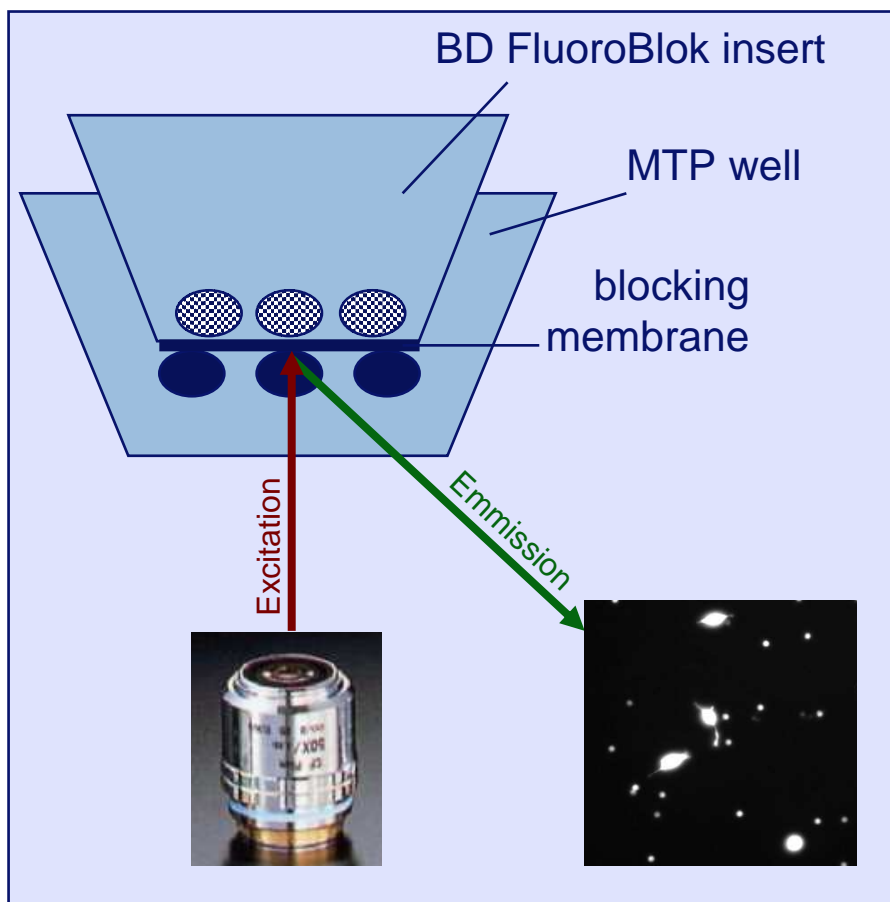
Hit To Lead

Lead Optimization

High-Content Analysis

migration

MVEC



Target validation

Assay Development

HTS

Hit To Lead

Lead Optimization

High-Content Analysis

DNA content

DNA content

Source image:

Approximate min width: $\mu\text{m} = 6$ pixels

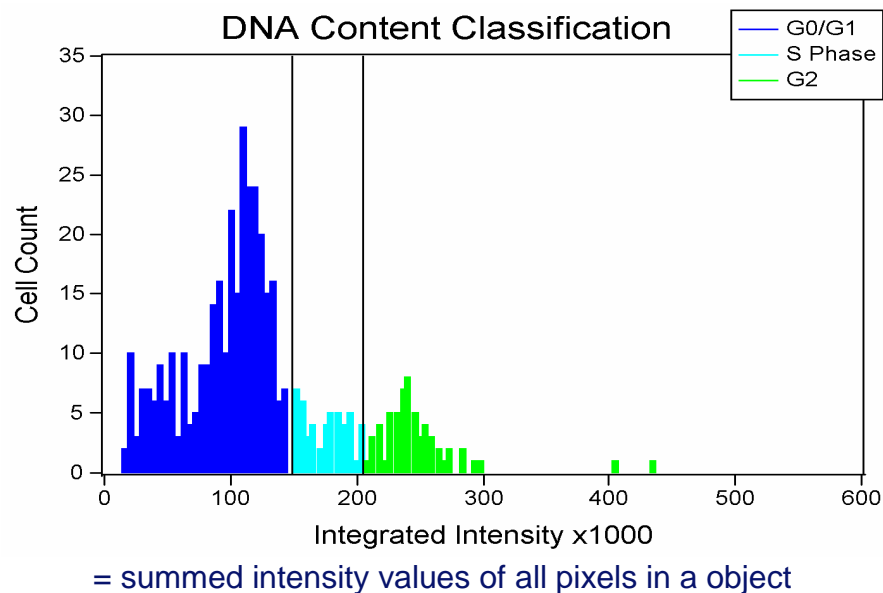
Approximate max width: $\mu\text{m} = 24$ pixels

Intensity above local background: graylevels

Background subtraction: at graylevels

Classification by integrated intensity (x1000)

G0/G1 (2N) S phase G2 (4N)



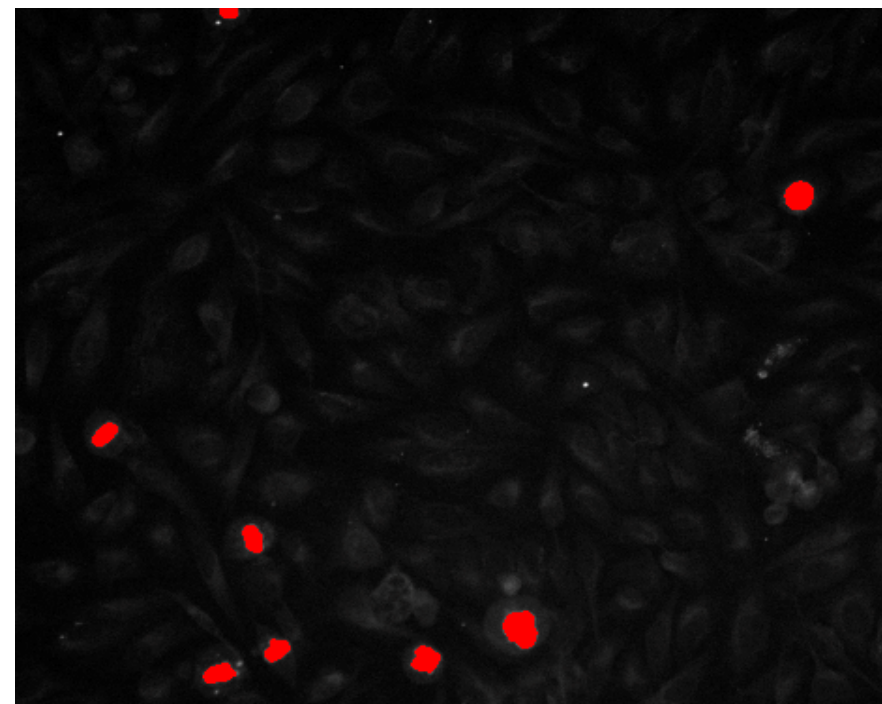
mitotic index

Mitotic classification

Mitotic-specific staining DNA average intensity

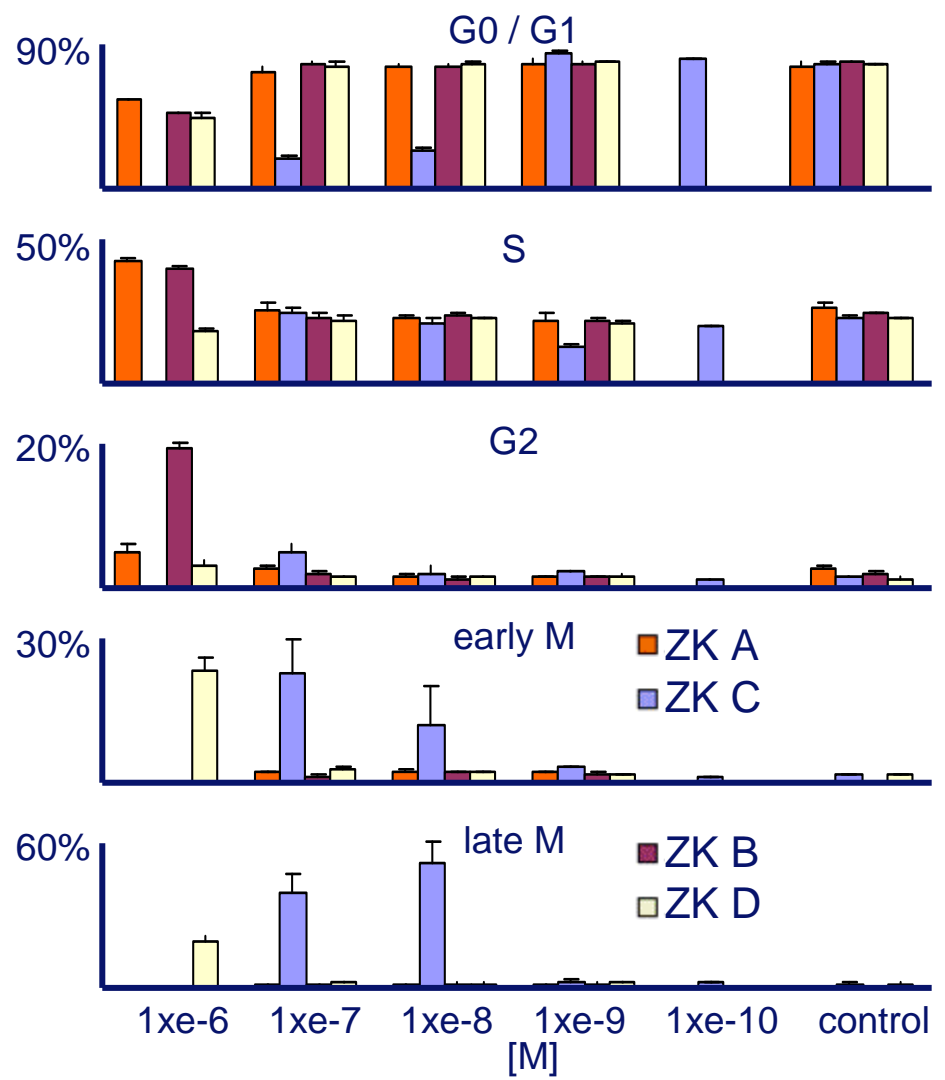
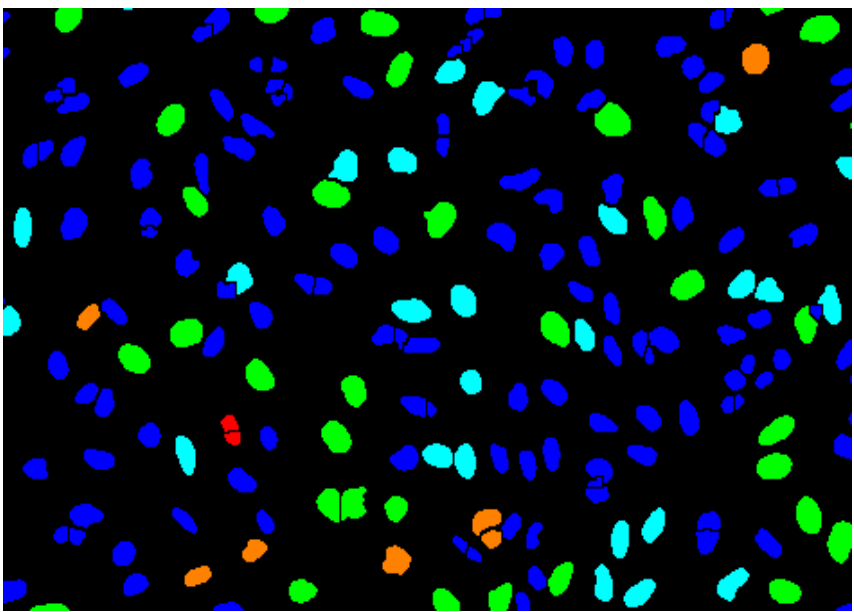
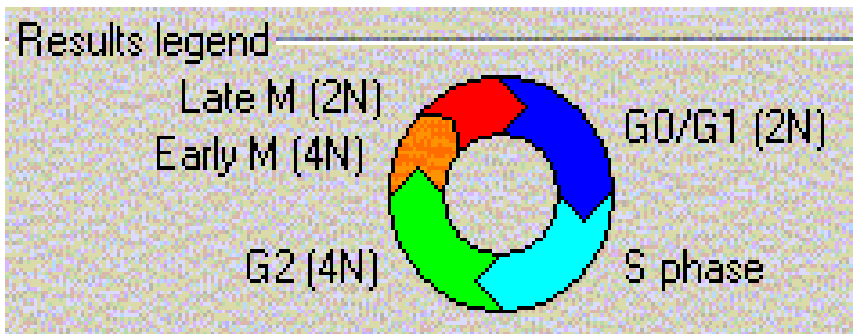
Source image:

Intensity above local background: graylevels





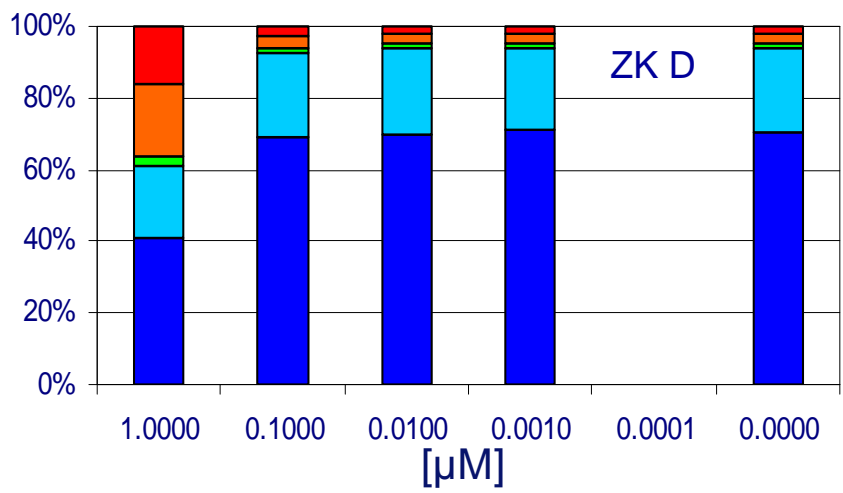
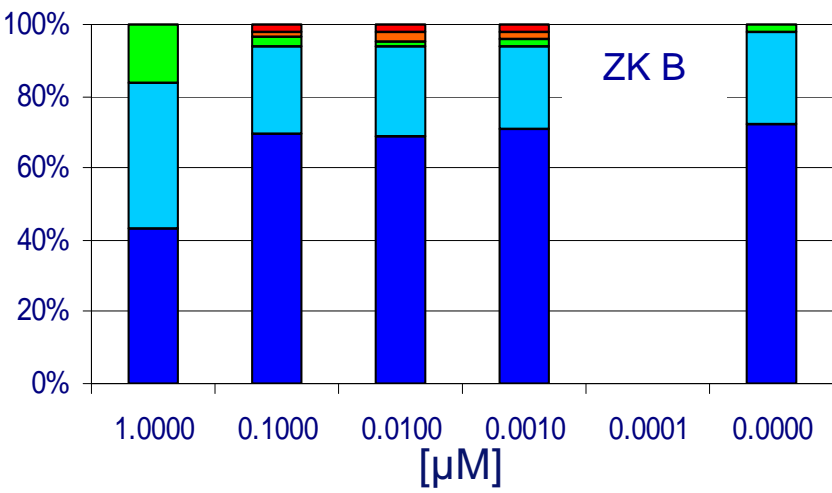
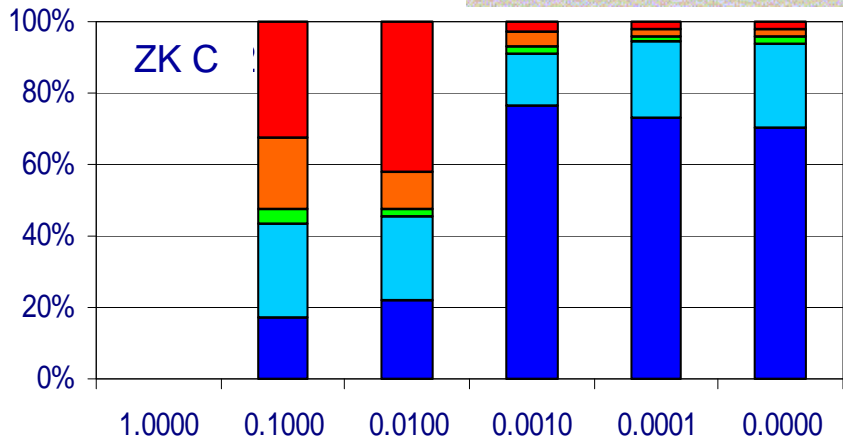
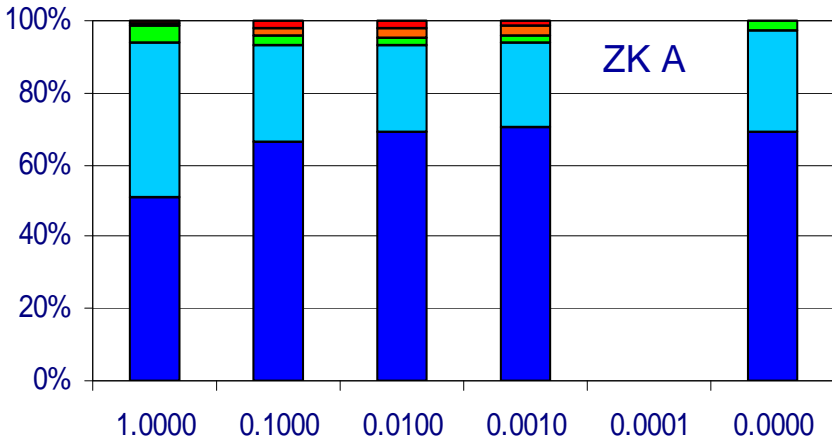
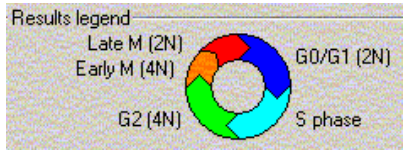
cell cycle classification





High-Content Analysis

cell cycle classification



Target validation

Assay Development

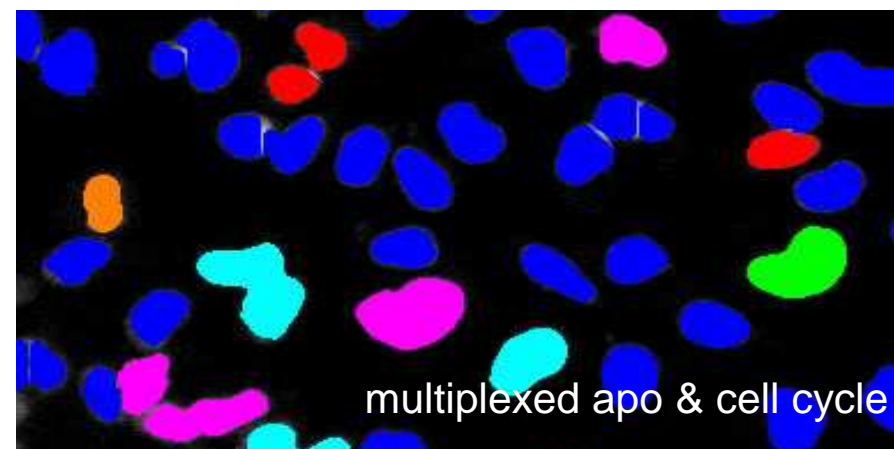
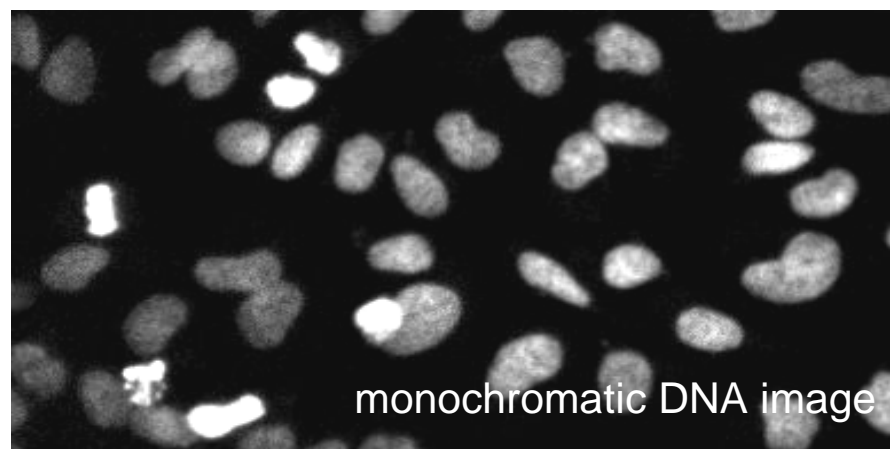
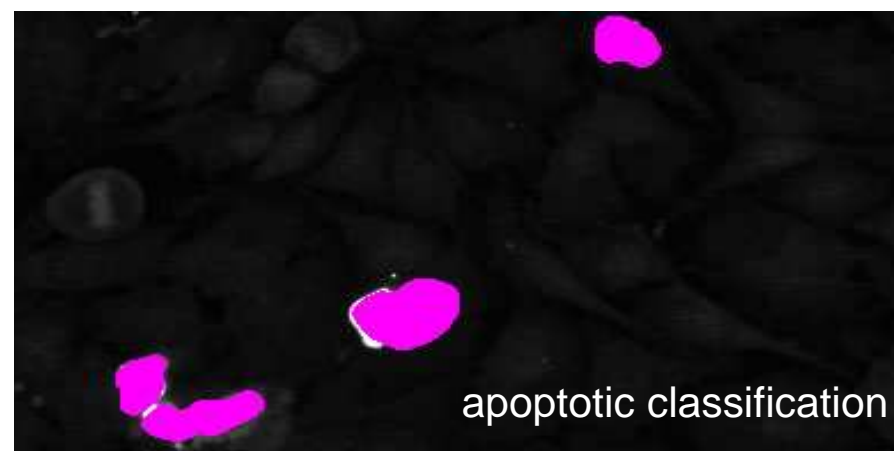
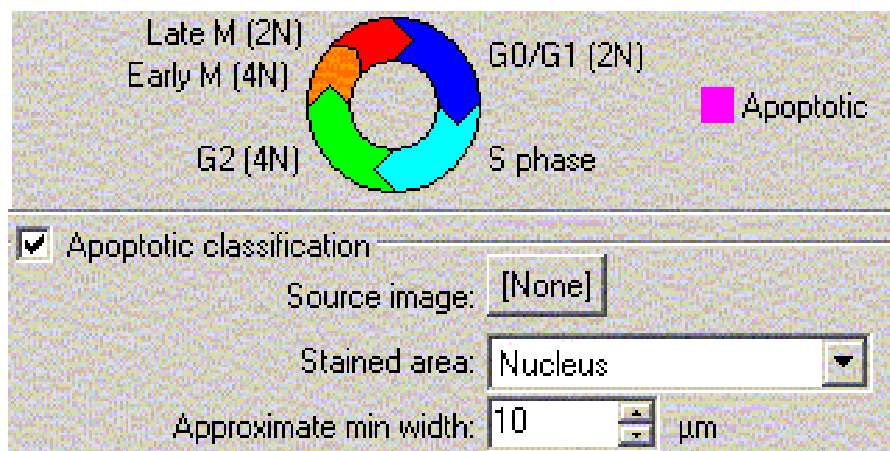
HTS

Hit To Lead

Lead Optimization

High-Content Analysis

apoptotic classification



Target validation

Assay Development

HTS

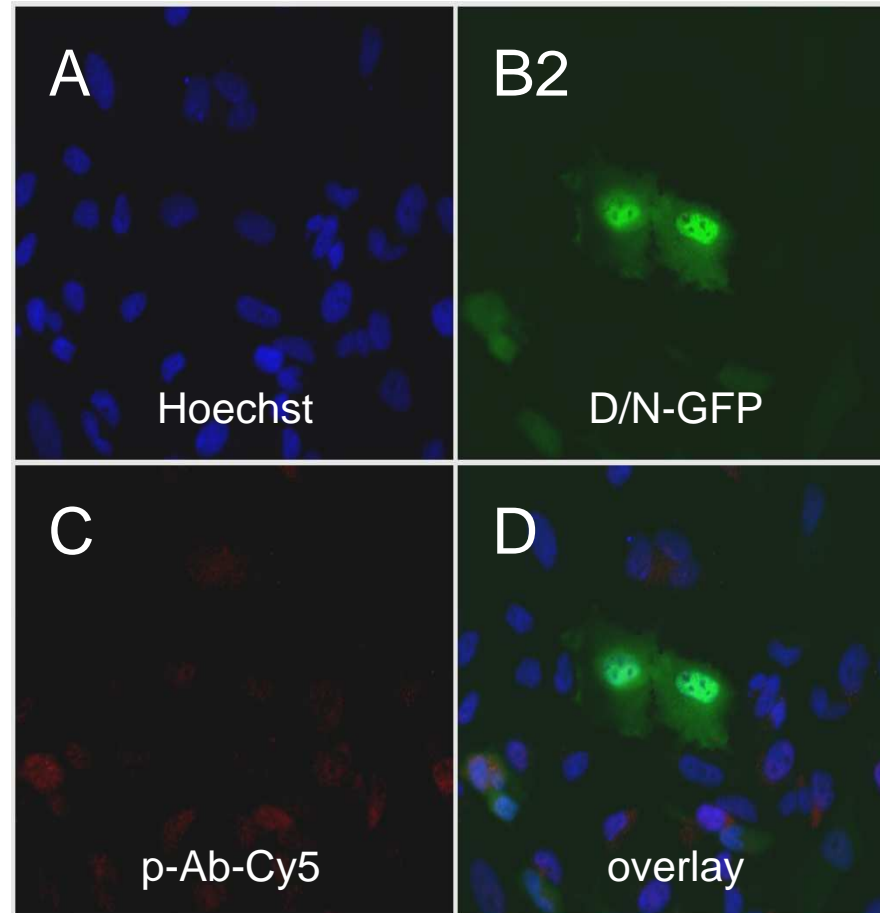
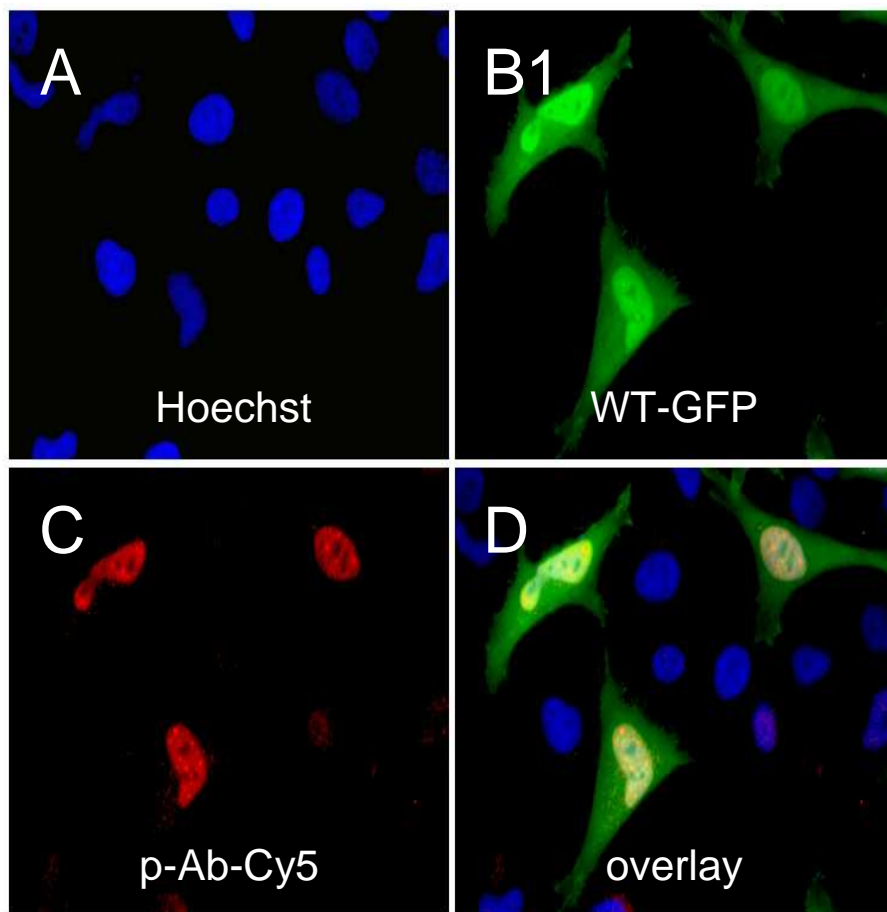
Hit To Lead

Lead Optimization

High-Content Analysis

protein-protein interaction

kinase vs. substrate



Target validation

Assay Development

HTS

Hit To Lead

Lead Optimization

High-Content Analysis

protein-protein interaction

kinase vs. substrate

transient transfection of GFP-tagged kinase (WT or D/N) and co-staining of phosphorylated known nuclear substrate

▶ **HOECHST dye (A)**

- visualisation of total number of nuclei

▶ **GFP expression (B)**

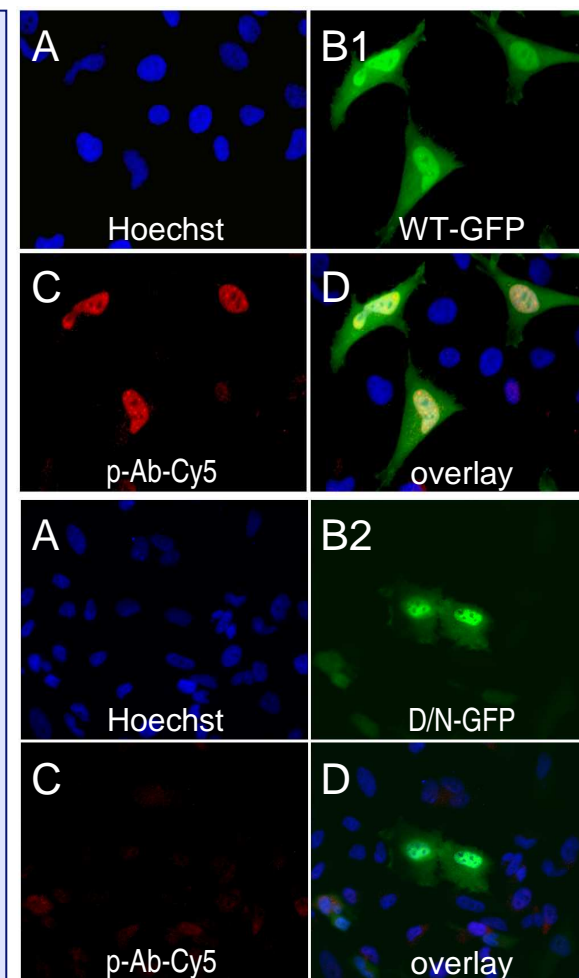
- identification of transfected cells
- distribution pattern of GFP-kinase
- B1: WT // B2: D/N

▶ **phosphorylated kinase substrate (C)**

- identification of substrate that has been phosphorylated at specific phosphorylation site

▶ **superimposition (D)**

- documentation of kinase dependent phosphorylation of substrate at specific site: only in WT-GFP positive nuclei

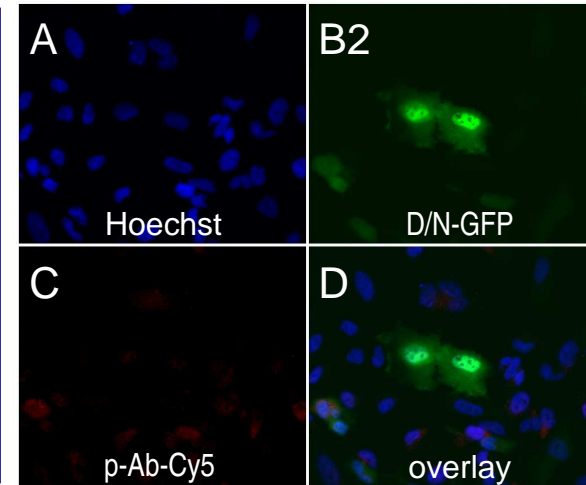
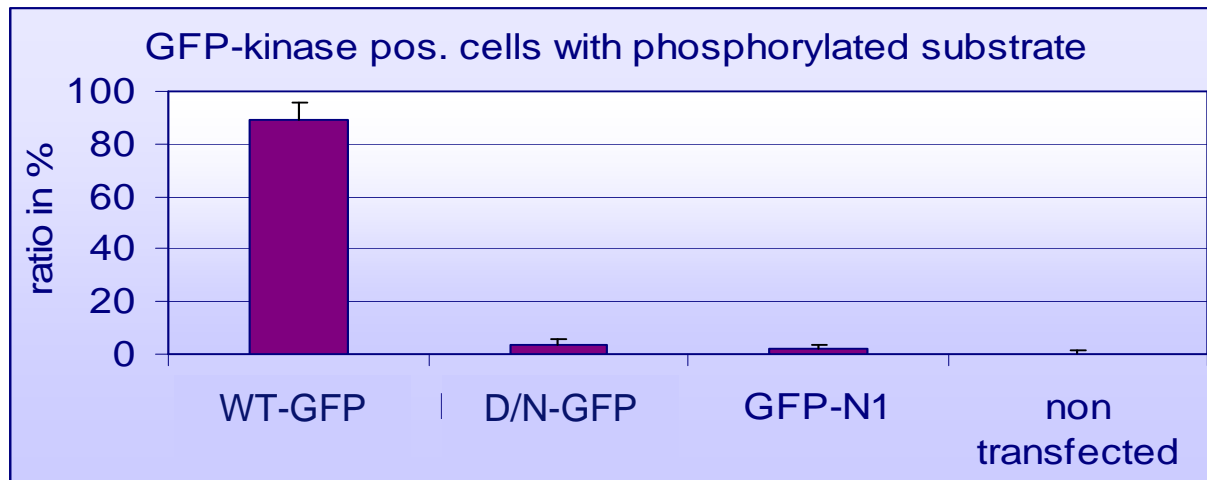
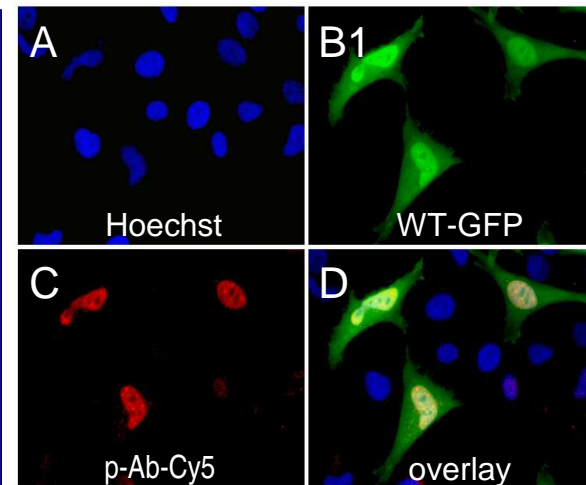
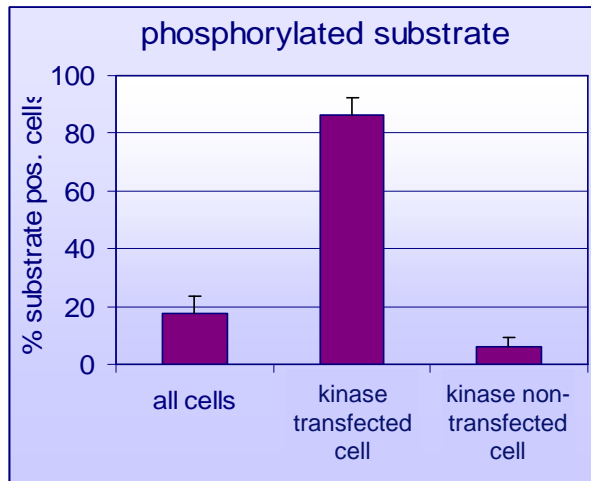
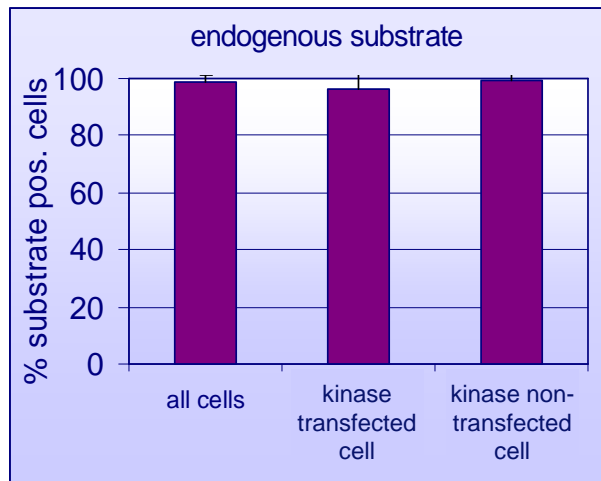




High-Content Analysis

protein-protein interaction

kinase vs. substrate



Target validation

Assay Development

HTS

Hit To Lead

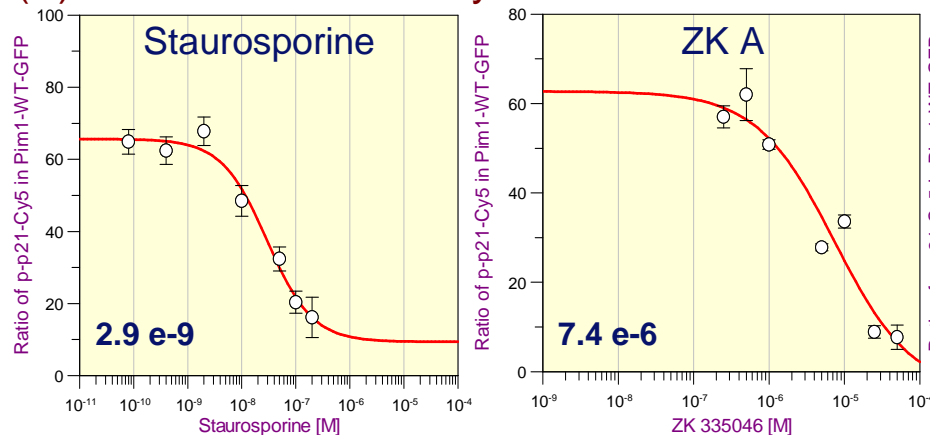
Lead Optimization

High-Content Analysis

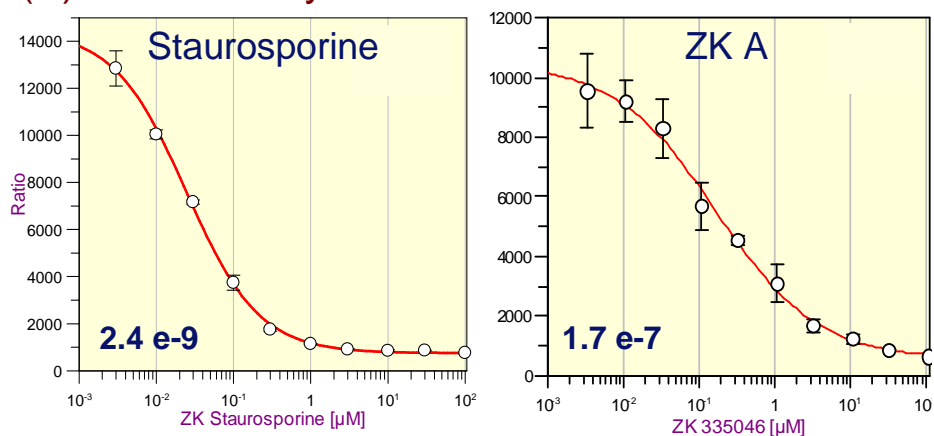
protein-protein interaction

kinase vs. substrate

(A) functional HCA assay



(B) HTRF assay



- functional HCA assay & biochemical HTRF assay revealed the same rank order for IC_{50} values of selected compounds
- IC_{50} values determined by HCA assay differ in 1 order of magnitude
- results from HCA assay contain
 - compound specificity
 - tox & site effects
 - cell permeability
 - MDR effects

⇒ results from HCA assay are of greater value

Target validation

Assay Development

HTS

Hit To Lead

Lead Optimization

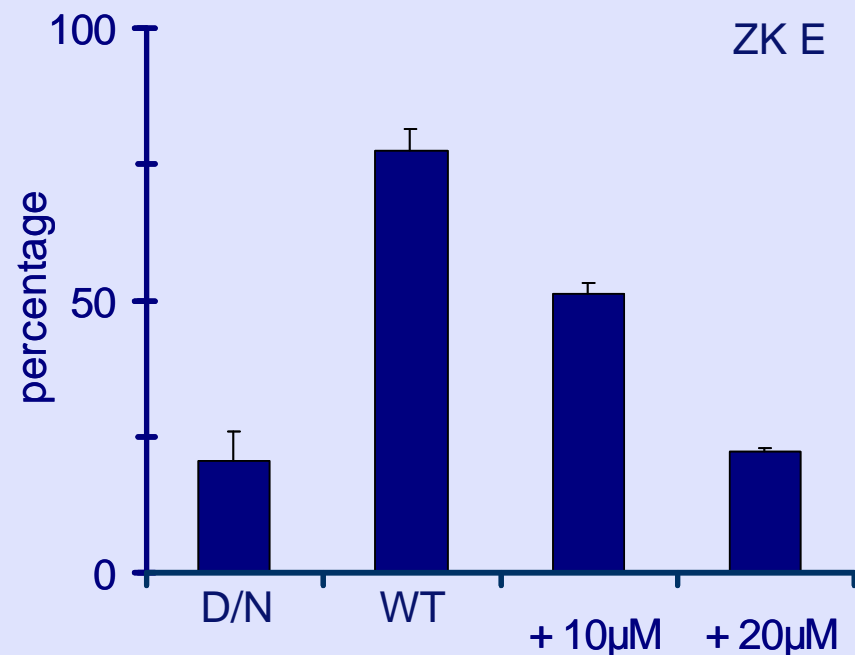
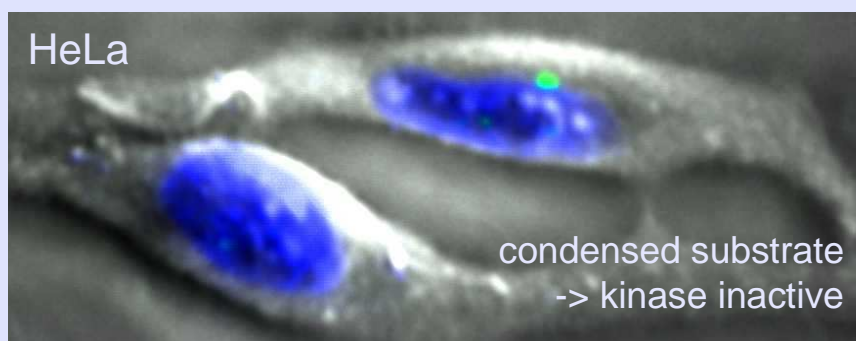
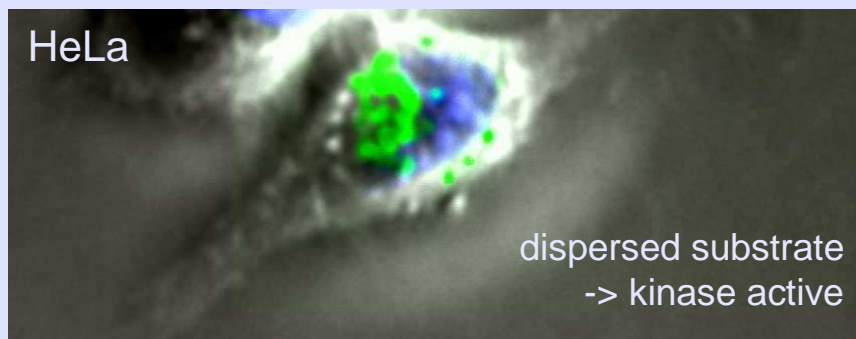
High-Content Analysis

protein-protein interaction

kinase vs. substrate

phosphorylation of substrate by kinase leads to a specific morphological alteration

- identification of kinase inhibitors
- quantification of potency



Target validation

Assay Development

HTS

Hit To Lead

Lead Optimization

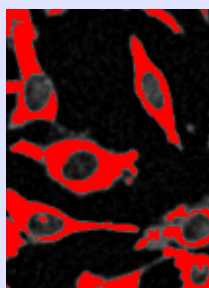
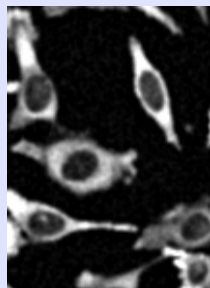
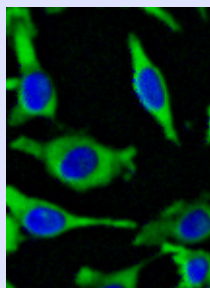
High-Content Analysis

protein translocation

protein activation often correlates with dislocation from origin to point of interaction

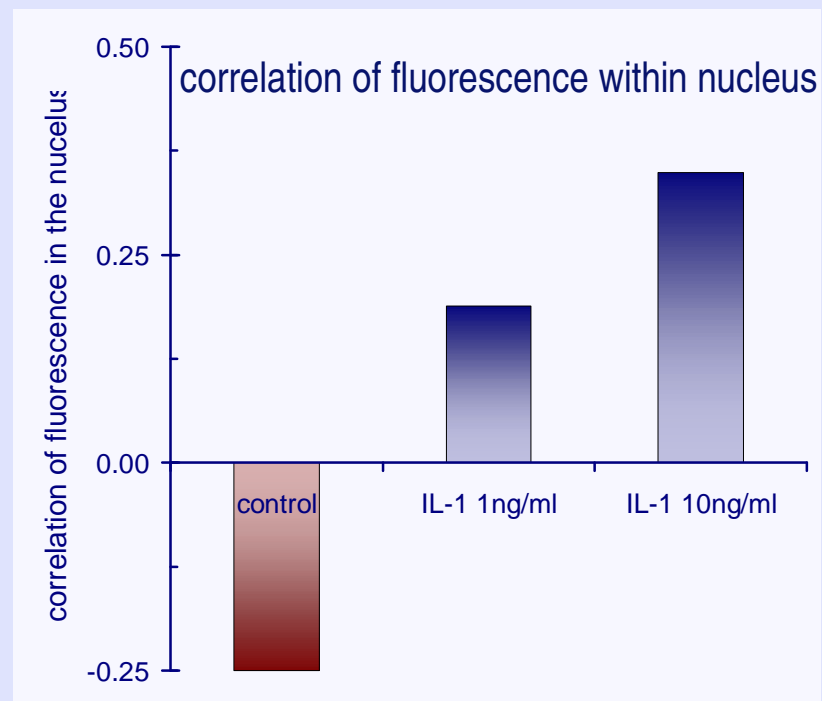
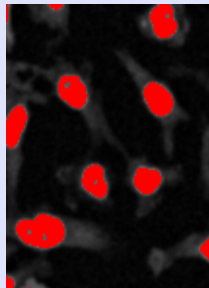
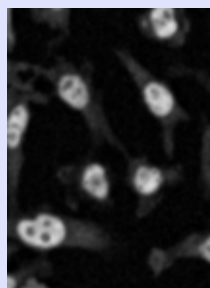
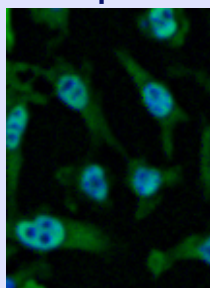
- phenotypical parameter which can be utilized by HCA
- identification of inhibitors / inducers

control



HeLa

compound



Target
validation

Assay
Development

HTS

Hit To Lead

Lead
Optimization

High-Content Analysis

receptor internalization

GPCR activation

Transfluor

- technology which can be used to screen for ligands or drugs regulating GPCR activity
- GPCR desensitization occurs very early in a common pathway and is coupled with binding to β -arrestin
- can be monitored by cycling from membrane to cytoplasm and spot formation

Endothelin A receptor internalization (ET_A-R EVOTEC Technologies)

- after agonistic stimulation, ET_A-R-GFP translocates from membrane to cytoplasm
- accumulation of ET_A-R-GFP in endosomes can be quantified by HCA

⇒ functional receptor internalization assays can be utilized for primary HTS

Target validation

Assay Development

HTS

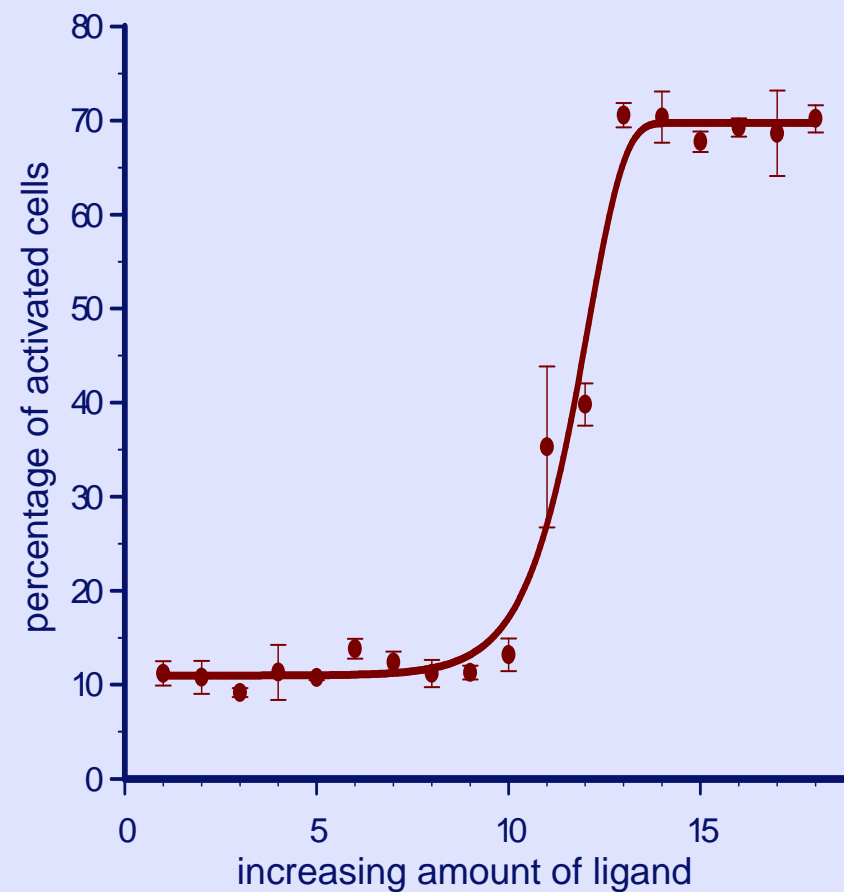
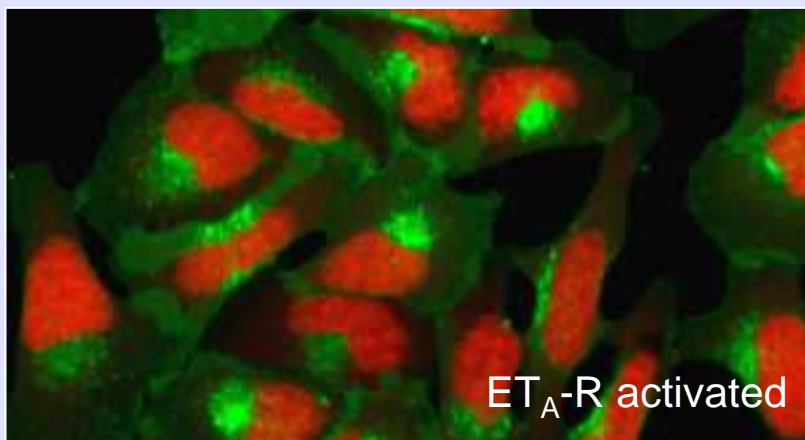
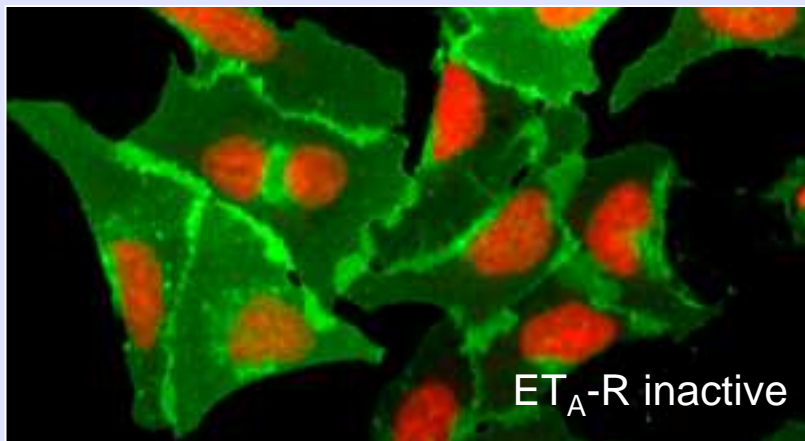
Hit To Lead

Lead Optimization

High-Content Analysis

receptor internalization

OPERA



EVOTEC Technologies

Target
validation

Assay
Development

HTS

Hit To Lead

Lead
Optimization

High-Content Analysis

time lapse studies

utilization of dynamic HCA approaches during preclinical drug research

- enlighten complex subcellular processes and pathways
- establish a profound knowledge data base
- improve specificity of compounds, efficiency of drugs and compliance

live-cell assays

- increasingly requested during target validation & lead optimization
- identification of important check points e.g. during cell cycle progression
- improve the understanding of cellular mechanisms
- qualify compound specificity in selected cell systems
- document the uniqueness of drug candidates

⇒ require improved data management, data storage, data handling & automated data analysis solutions

Target validation

Assay Development

HTS

Hit To Lead

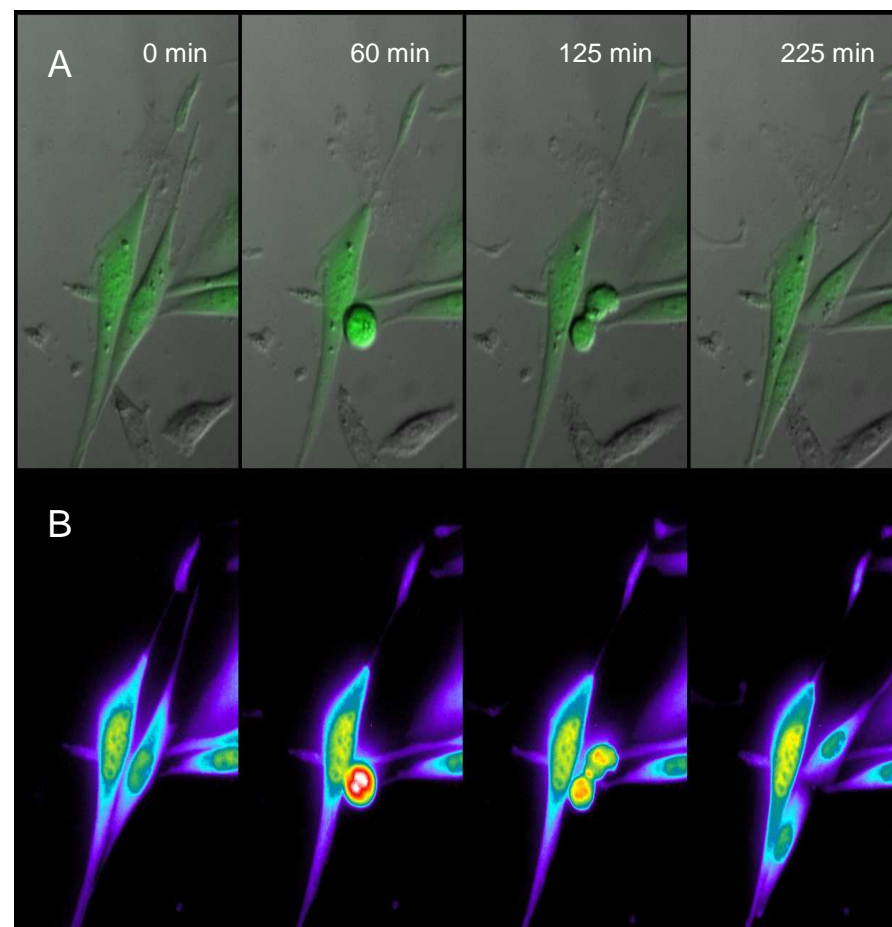
Lead Optimization

High-Content Analysis

cell cycle kinase

GFP-labeled kinase

- ▶ intracellular localization during cell cycle progression
 - superimposition (A) of DIC and GFP channel
 - intensity profile of GFP fluorescence (B)
- ▶ interphase (0 min)
 - predominant nuclear localization
- ▶ mitosis
 - localisation in specific subcellular regions
- ▶ metaphase (60 min)
 - bipolar localisation
- ▶ cytokinesis (125 min)
 - area of the reassembling nuclei and region of the contractile ring



Target
validation

Assay
Development

HTS

Hit To Lead

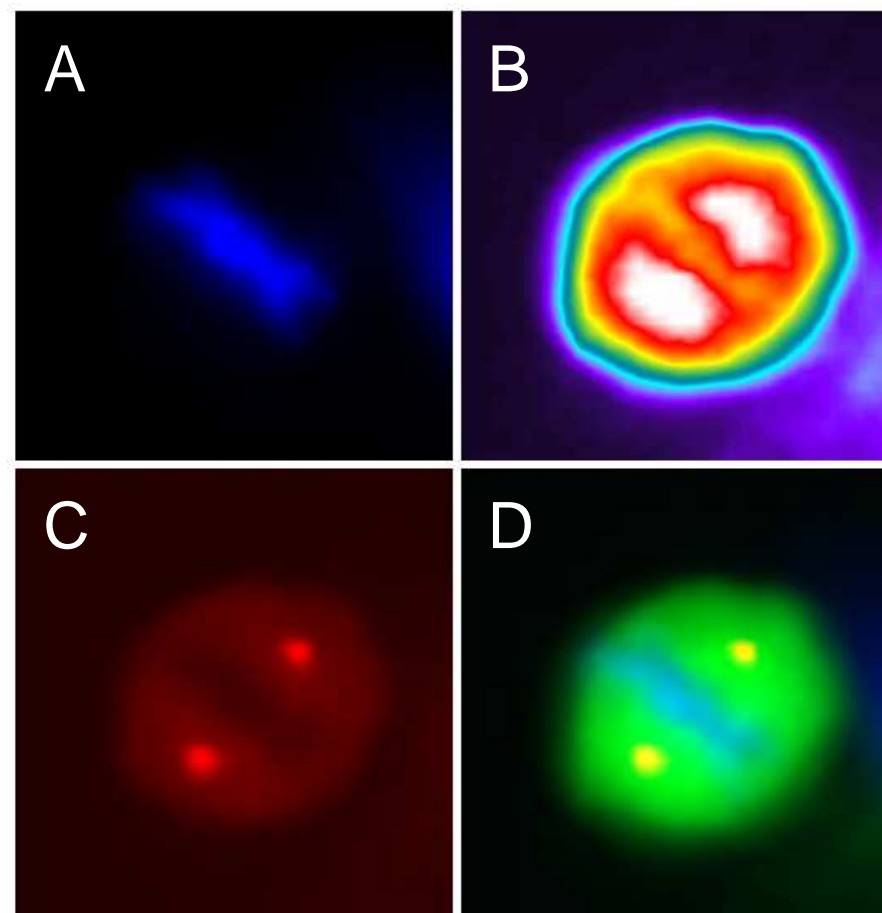
Lead
Optimization

High-Content Analysis

cell cycle kinase

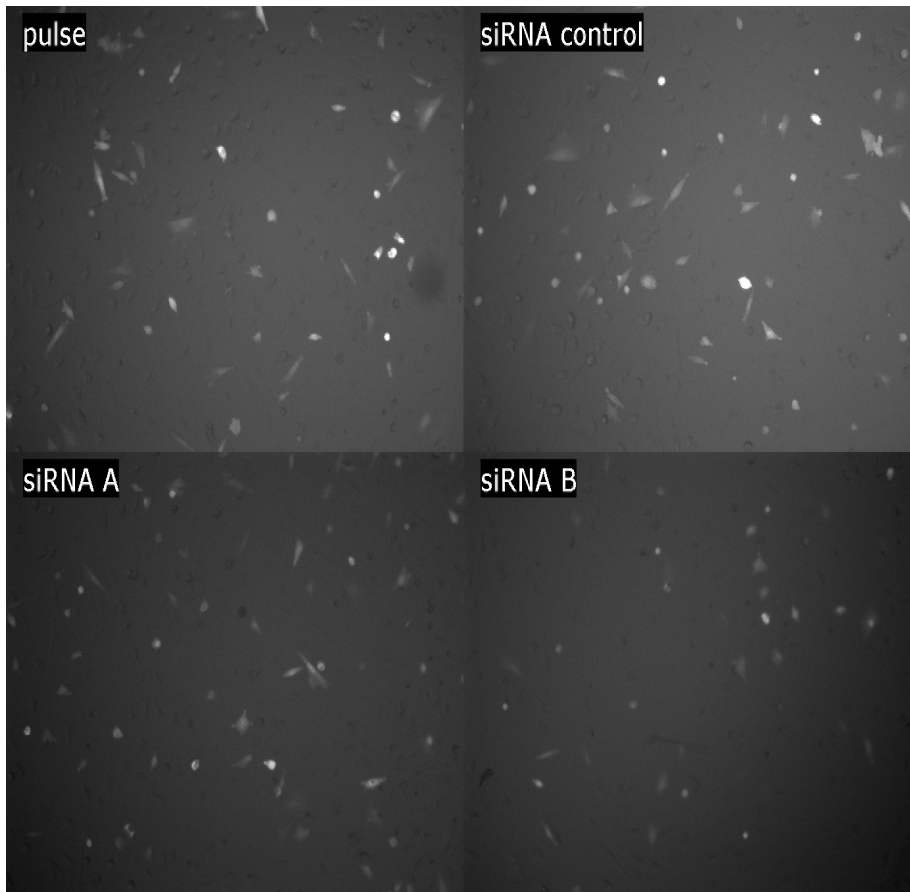
GFP-labeled kinase

- ▶ intracellular localization of GFP-tagged kinase during metaphase
- ▶ **chromosomes (A)**
 - condensed localisation in equatorial plate
- ▶ **GFP-tagged kinase (B)**
 - intensity profile of GFP expression
- ▶ **γ -tubulin (C)**
 - bipolar localisation of spindle poles
- ▶ **superimposition (D)**
 - indication of a predominant localisation of GFP-kinase fusion protein around spindle poles during metaphase

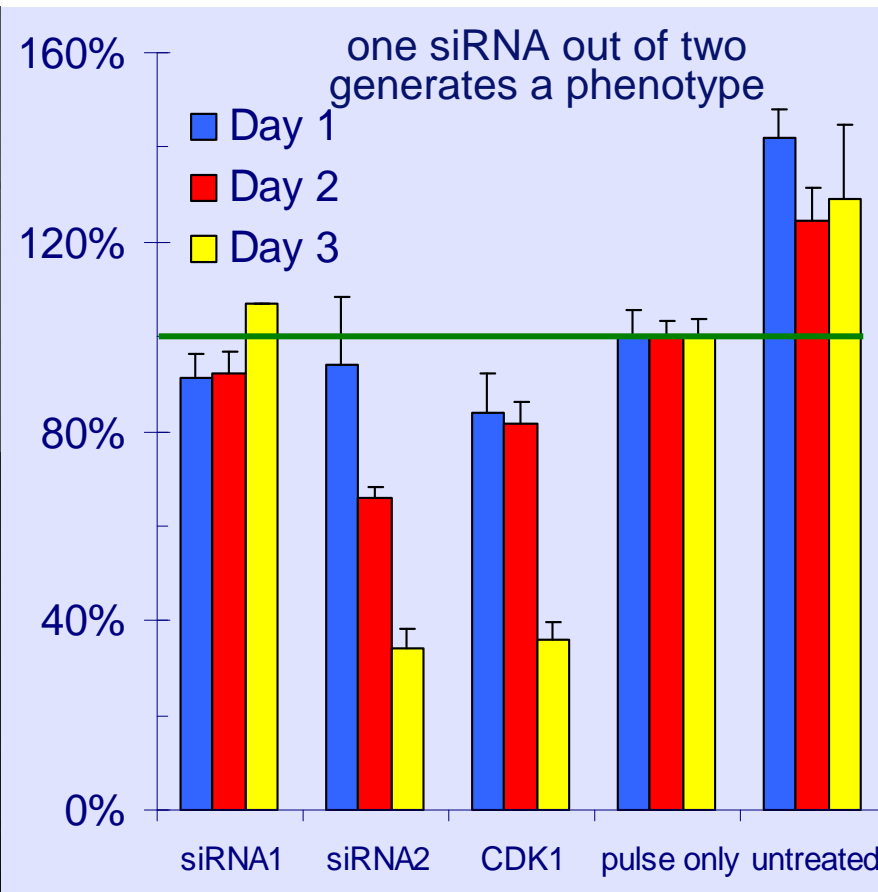




siRNA knockdown of a GFP-kinase in CHO cells



siRNAi k.d. of endogenous kinase in MCF7 cells



in collaboration with Claudia Merz

Target validation

Assay Development

HTS

Hit To Lead

Lead Optimization

High-Content Analysis

time lapse studies

cell cycle kinase

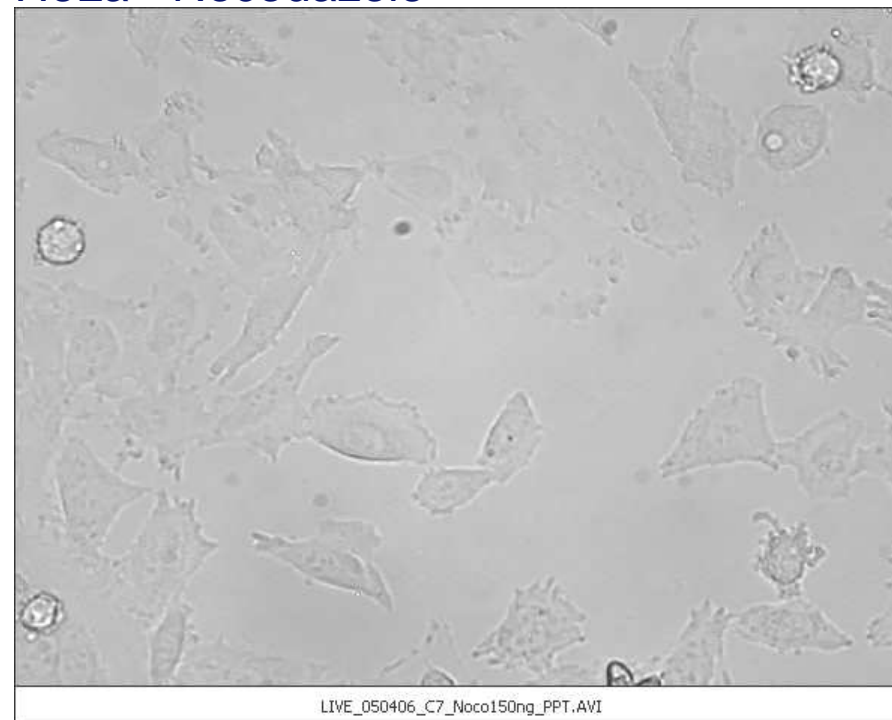
inhibition of mitotic spindle check point protein leads to G2/M arrest

- documented by Nomarski contrast method, 24h monitoring, images taken every 5min

HeLa - ZK D



HeLa - Nocodazole



Target validation

Assay Development

HTS

Hit To Lead

Lead Optimization

High-Content Analysis

time lapse studies

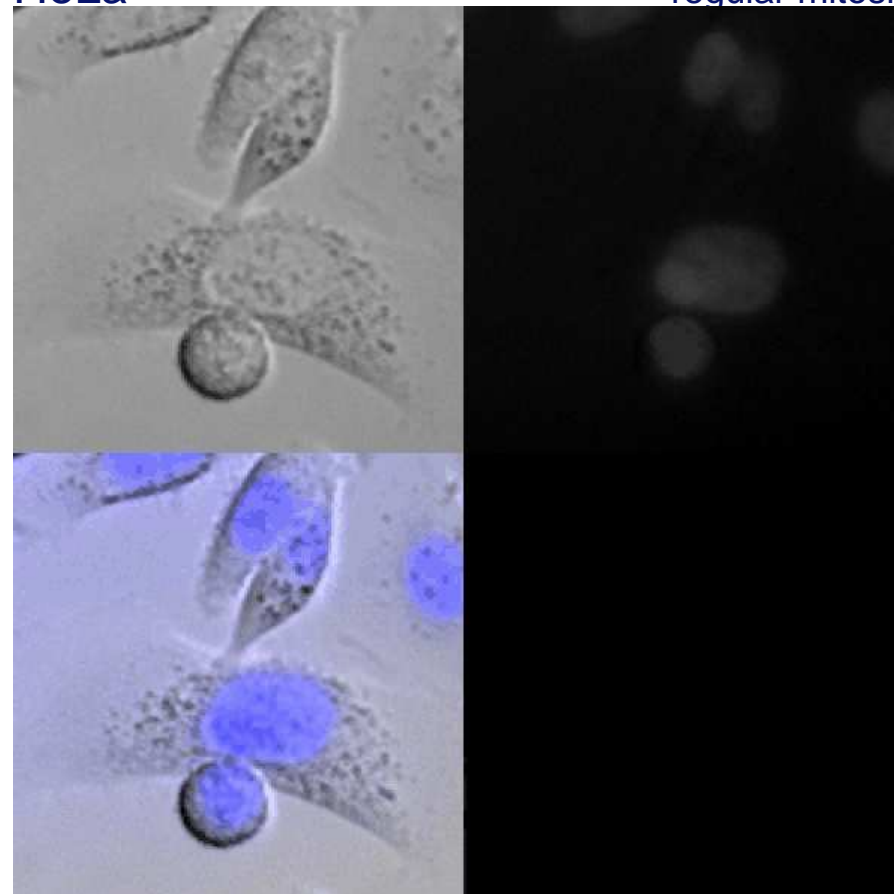
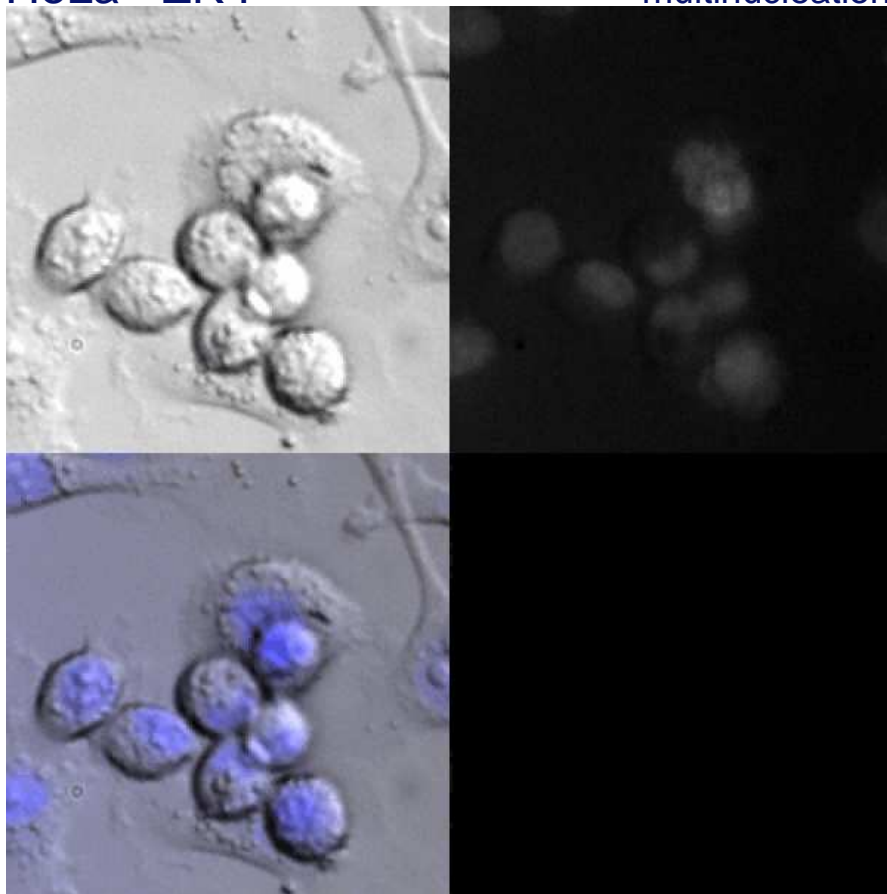
cell cycle kinase

HeLa - ZK F

multinucleation

HeLa

regular mitosis



Target
validation

Assay
Development

HTS

Hit To Lead

Lead
Optimization

High-Content Analysis

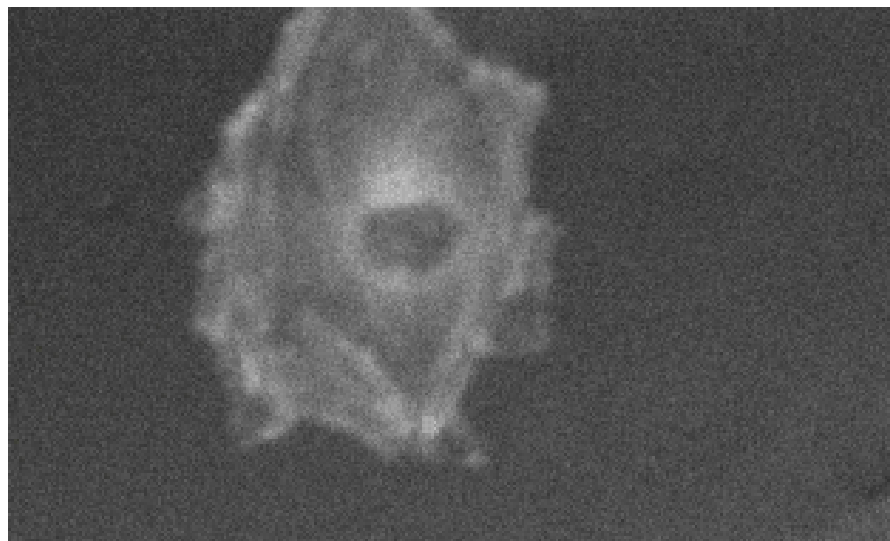
time lapse studies

Discovery 1

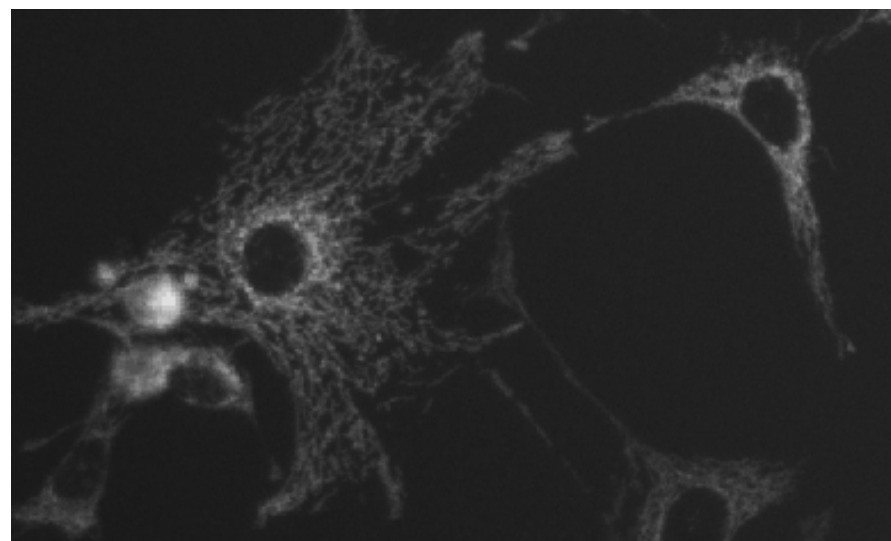
stably transfected cells expressing AFP tagged cell compartment or target protein

- phenotypical characterisation ⇒ improved black box screening
- functional pathway analysis ⇒ quantification of compound specificity
- ⇒ accounting the dynamic and spatial context

PtK2-PhiYellow-Actin



HeLa-TurboGreen-Mitochondria



Target validation

Assay Development

HTS

Hit To Lead

Lead Optimization

High-Content Analysis

3D sectioning

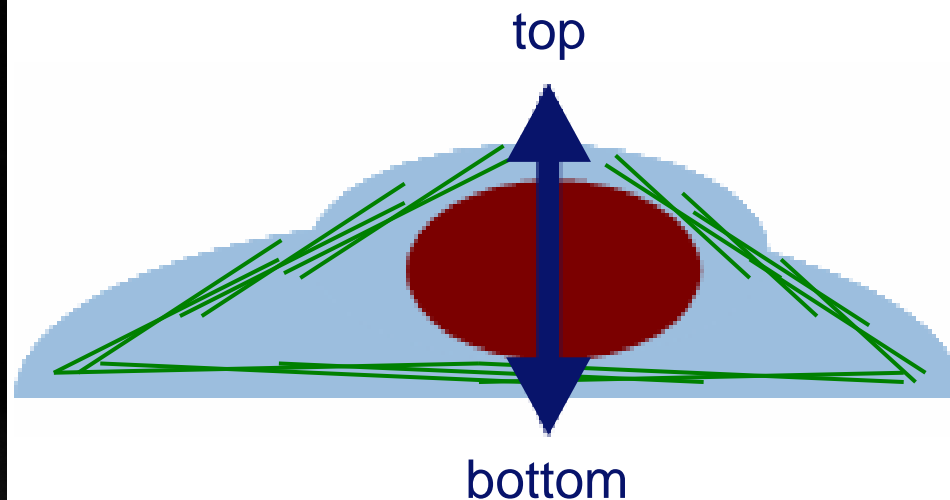
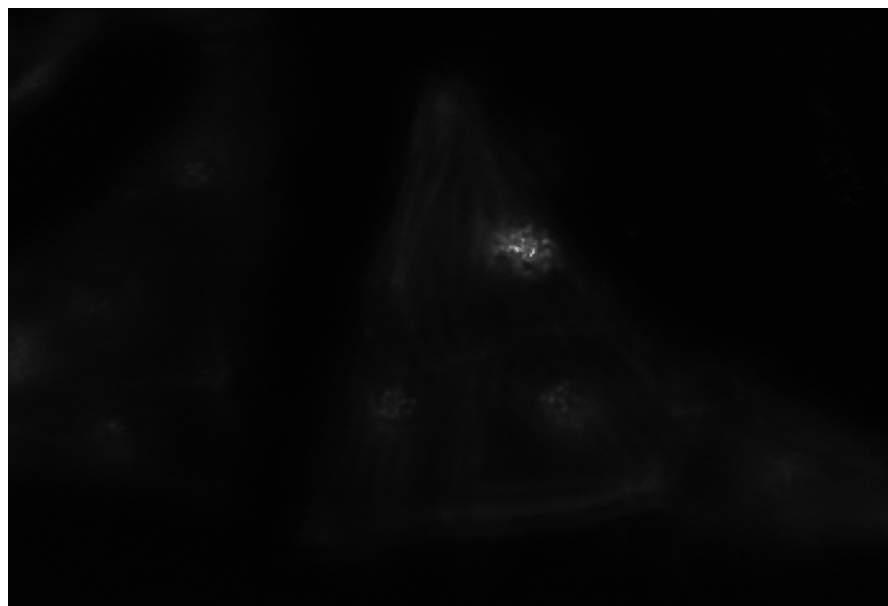
OPERA

confocal imaging enables the dissection of a cell into a stack of discrete planes

- cells can be analysed in a three-dimensional context
- 3D sectioning combined with time-lapse generates a powerful novel HCA tool

PtK2-PhiYellow-Actin

40x



Target validation

Assay Development

HTS

Hit To Lead

Lead Optimization

High-Content Analysis

supported projects

applications

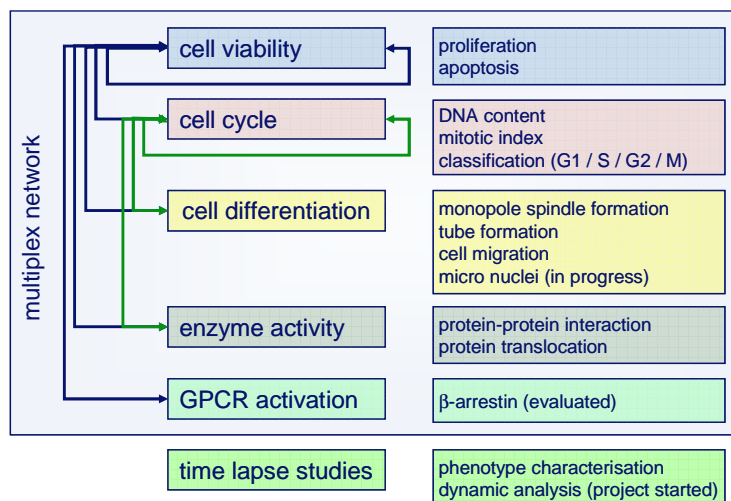
HT-RNAi project

- phenotypical screening: proliferation, apoptosis, DNA content, mitotic index, cell cycle classification, monopole spindle formation, etc.

Oncology

- siRNA / cDNA transfection: proliferation, apoptosis, cell cycle classification, protein-protein interaction, protein translocation, time lapse studies

application panel



cell systems

HeLa
CHO
DU145
PC3
LNCap
MCF7
T47D
BT474
SK-OV-3
U-2 OS
HUVEC
MVEC
PAEC
S49
Jurkat
prim. M Φ



High-Content Analysis

supported projects

applications

Gynecology & Andrology

- proliferation, migration, primary macrophage classification

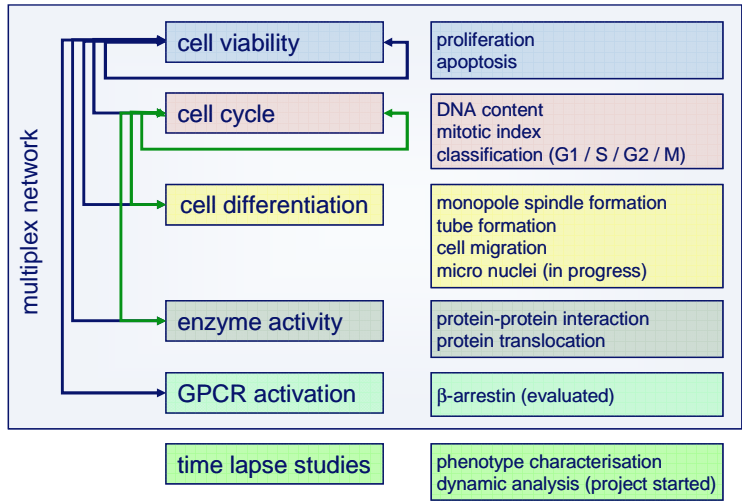
Inflammation

- proliferation, apoptosis

Oncology

- proliferation, apoptosis, tube formation, mitotic index, DNA content, cell cycle classification, protein-protein interaction, time lapse studies

application panel



cell systems

- HeLa
- CHO
- DU145
- PC3
- LNCap
- MCF7
- T47D
- BT474
- SK-OV-3
- U-2 OS
- HUVEC
- MVEC
- PAEC
- S49
- Jurkat
- prim. MΦ



supported projects

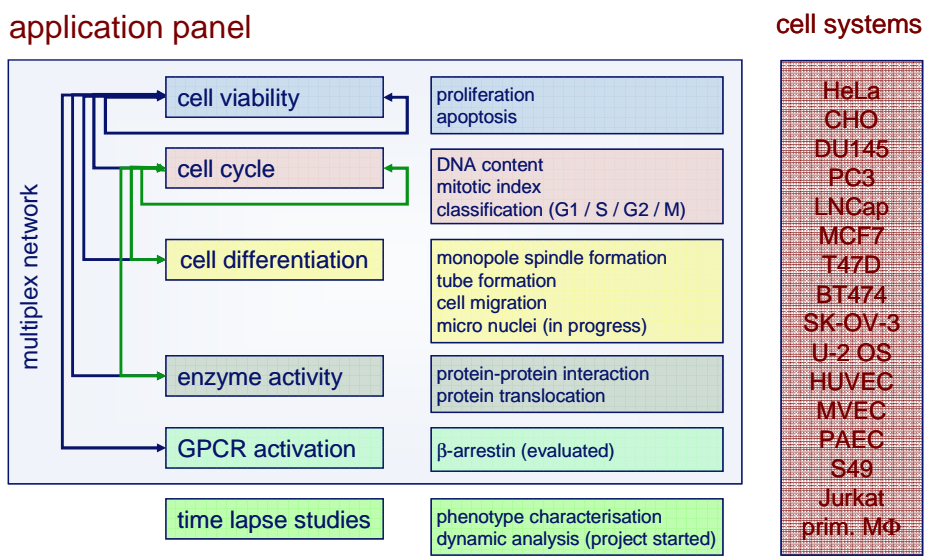
applications

Inflammation

- proliferation, apoptosis

Oncology

- proliferation, apoptosis, mitotic index, DNA content, cell cycle classification, protein-protein interaction, time lapse studies



Target validation

Assay Development

HTS

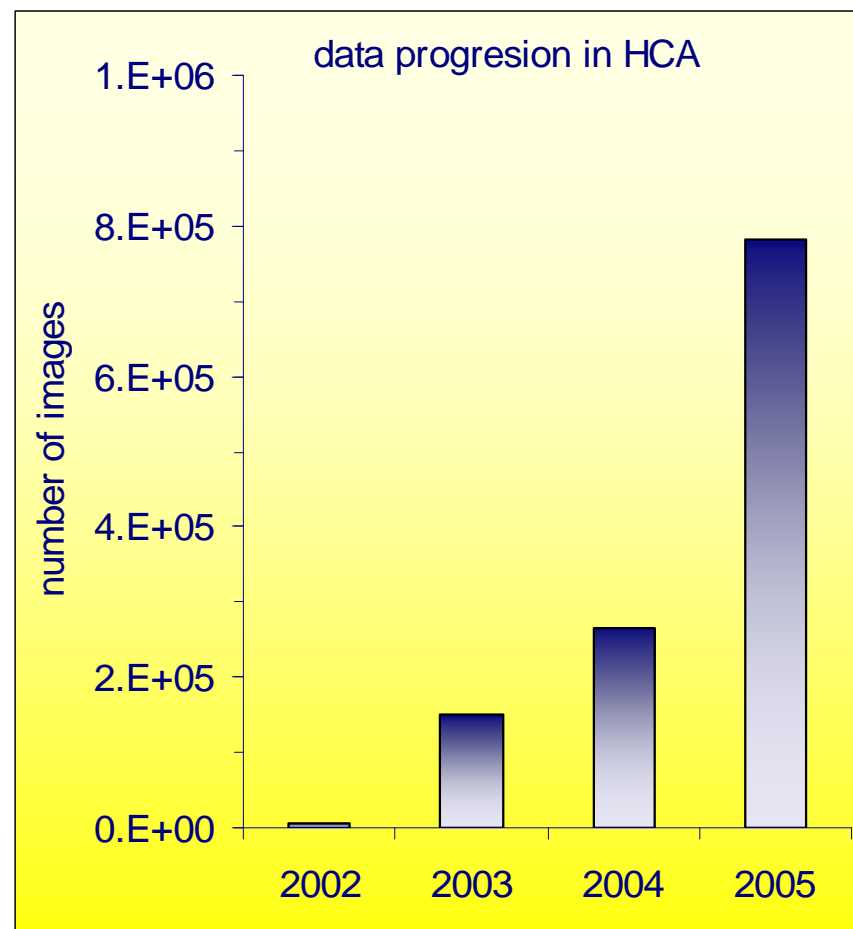
Hit To Lead

Lead Optimization

High-Content Analysis

HCA at Schering AG in 2006

- ▶ broad panel of HCA applications established that meet CRBA needs
 - statistically secured and relevant data confirmed HCA as a reliable technology
- ▶ highly requested technology
- ▶ Target Validation
 - RNAi k.d. & transient transfection studies for all major target families
- ▶ LD / LO support
 - multiplexed & functional assays established -> target specific mechanistic profiling
- ▶ contributions to primary HTS possible
 - target oriented assays, improved black box



acknowledgements

- Philip Denner
- Alexander Dimerski
- Sebastian Schäfer
- Janine Schmalowsky

