



## ***Cell-based kinase assays in HTS***

*Potential and limitations for primary and secondary screening*

***Benjamin Bader***

*HTRF-symposium Avignon 25. April 2013*

- Introduction
- General considerations cell-based kinase assays
- EFC-assays for Tyr-Kinases
- TR-FRET systems: HTRF and Lanthascreen assays for the PI3K / Akt / mTOR pathway
- Summary & conclusions

# Kinases as drug targets



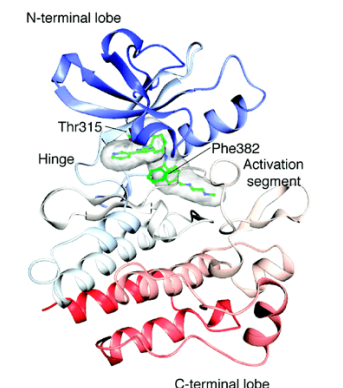
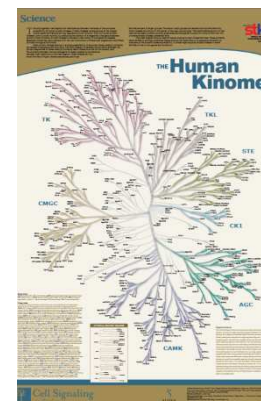
## Registered kinase inhibitors

Compound	Kinase target	Cancer target	Company
Imatinib (Glivec, Gleevec, STI571)	ABL 1-2, PDGFR, KIT	CML, Ph+ B-ALL, MML, CEL, GIST	Novartis
Gefitinib (Iressa, ZD1839)	EGFR	NSCLC	AstraZeneca
Erlotinib (Tarceva, OSI-774)	EGFR	NSCLC, pancreatic cancer	OSI, Genentech Inc, Roche
Lapatinib (Tykerb, GW2016)	EGFR, ERBB2	Breast cancer	Glaxo SmithKline
Dasatinib (Sprycel, BM-354825)	ABL1-2, PDGFR, KIT, SRC	CML	Bristol Myers
Nilotinib (Tasigna, AMN107)	ABL1-2, PDGFR, KIT	CML	Novartis
Sunitinib (Sutent, SU11248)	VEGFR1-3, KIT, PDGFR, RET, CSF1R, FLT3	RCC, GIST	Pfizer
Sorafenib (Nexavar, Bay 43-9006)	VEGFR2, PDGFR, KIT, FLT3, BRAF	RCC	Onyx and Bayer Pharmaceuticals
Pazopanib (Votrient, GW-786034)	VEGFR1-3, PDGFR, KIT,	RCC	GlaxoSmithKline
Everolimus (Afinitor, Rad001)	mTOR	RCC	Novartis
Temsirolimus (Torisel, CCI-779)	mTOR	RCC	Wyeth

(Fabbro 2012)

## recent additions

Vandetanib	VEGFR, EGFR, RET	thyroid	AstraZeneca
Vemurafenib	B-RafV600E	melanoma	Roche/Plexxikon
Regorafenib	multiKinase	colorectal	Bayer
Critozinib	ALK, ROS1	NSCLC	Pfizer
Bosutinib	BCR/Abl	CLL	Pfizer
Ruxolitinib	JAK1/2	Myelofibrose	Incyte/Novartis



cAbl and Gleevec (Noble 2004)

- human kinome: 518 protein kinases + 20 lipid kinases (Manning 2002)
- currently, ~150 kinase targeted drugs are in clinical development (Fabbro 2012)
- most registered kinase drugs target Tyrosine kinases, with more Ser/Thr kinase targeted drugs in the pipeline
- most kinase drugs target the ATP-pocket
- main indication: oncology



# Kinase Inhibitors in the BAYER Development Pipeline

Phase I (11)	Phase II (8)	Phase III (14)
Cancer / CDK Inhibitor ★	Cancer / PI3K Inhibitor ★	Thyroid Cancer / Sorafenib ★
Cancer / Mesothelin-ADC	Cancer / Regorafenib* ★	Breast Cancer / Sorafenib ★
Cancer / PSMA BiTE Antibody	Cancer / MEK-Inhibitor ★	Adjuvant HCC / Sorafenib ★
Cancer / PI3K $\alpha/\beta$ Inhibitor ★	Cancer / Radium-223 Dichloride	Adjuvant RCC / Sorafenib ★
Anemia / HIF-PH	Additional Indications / Sorafenib ★	Major Adverse Cardiac Events / Rivaroxaban
Heart Failure / Partial Adenosine A1 Agonist	CHF / MR Antagonist	Hemophilia / peg rFVIII**
Heart Failure / Vasopressin Receptor Antag.	PH / Riociguat (sGC Stimulator)	Hemophilia / rFVIIa**
Heart Failure / sGC Stimulator	Gram-neg. Pneumonia / Amikacin inhale	Myopic CNV / Aflibercept
Bronchiectasis / Neutrophil Elastase Inhibitor		DME / Aflibercept
Sympt. Uterine Fibroids / S-PRAnt		Contraception / LCS 16
Endometriosis / BAY 1026153		VV Atrophy / Vaginorm
		Submental fat removal / ATX-101
		Lung Infection / Cipro Inhale
		Skin and Lung Infections / Tedizolid

\*Regorafenib is a Bayer compound developed solely by Bayer. In 2011, Bayer entered into an agreement with Onyx Pharmaceuticals, Inc. under which Onyx will receive a royalty on any future global net sales of regorafenib in oncology

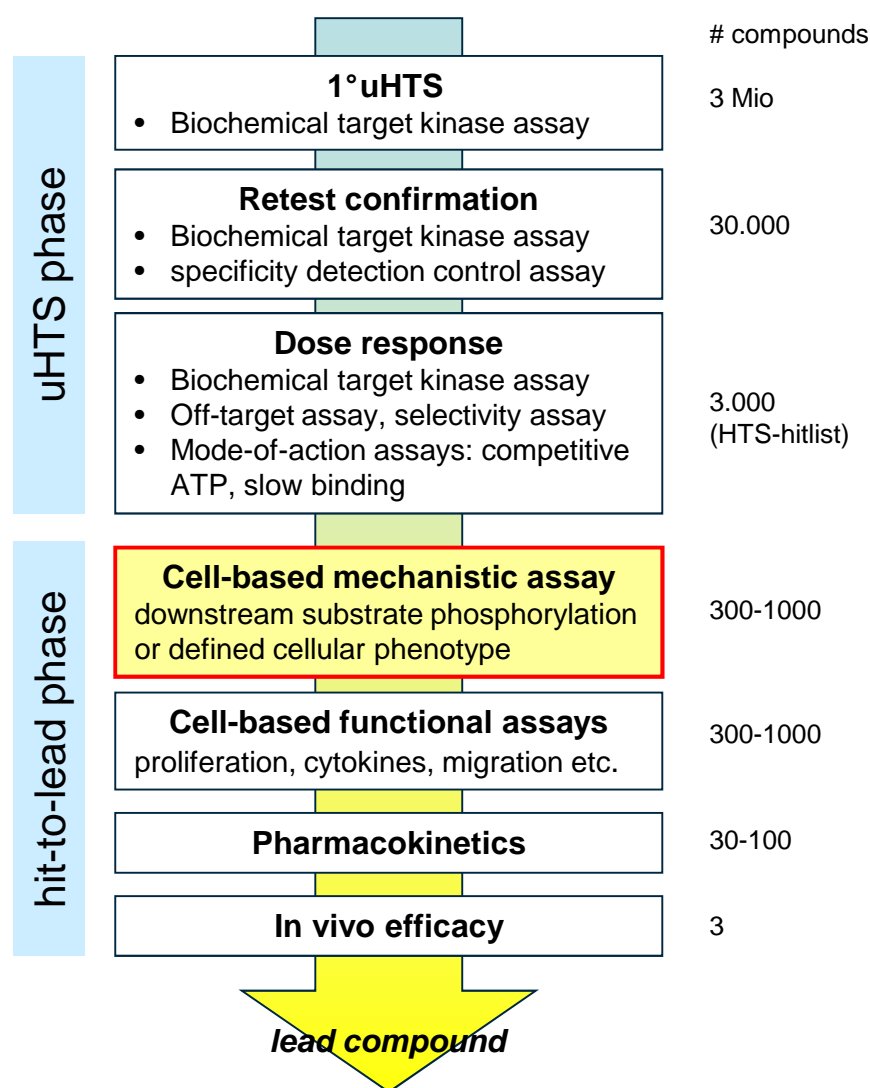
\*\* Combined Phase II/III

Status February 2013

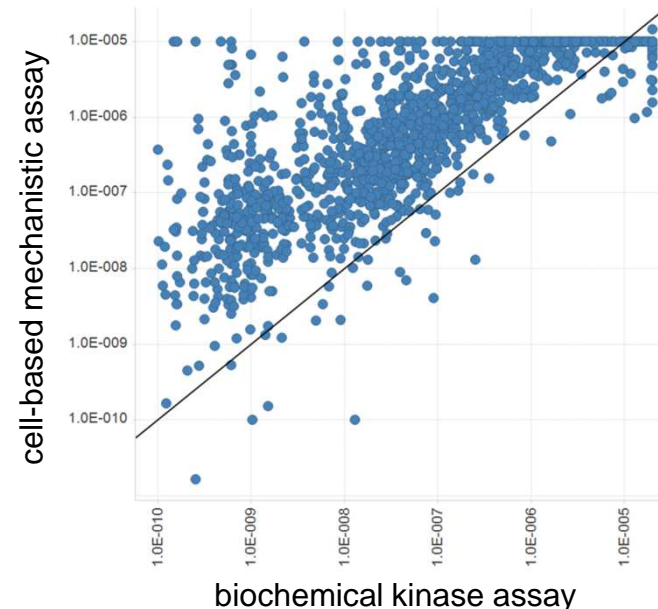
Selection of major Pharma pipeline projects in clinical Phase I to III

- █ Oncology
- █ Cardiovascular Diseases
- █ Women's Healthcare
- █ Hematology
- █ Ophthalmology
- █ Others

# Features of a typical early kinase drug discovery project



correlation mechanistic – biochemical



- specific cell-based kinase assays are essential for hit profiling during hit-to-lead phase
- integration of cell-based kinase assays as early as possible, best case even before hitlist delivery



# Requirements for efficient assay support in HTS and Hit-to-lead phase @ BAYER



## ***Key objectives for every assay***

(except HCA staining procedures)

- assay-ready plates
- 1536 well format, at least 384 well
- homogenous addition only
- endpoint assays
- frozen cells
- short
- robust

automated preparation of  
assay-ready plates



cells and reagent  
dispensing into 1536 well  
plates → addition-only



endpoint detection  
with imaging-based  
multimode-readers



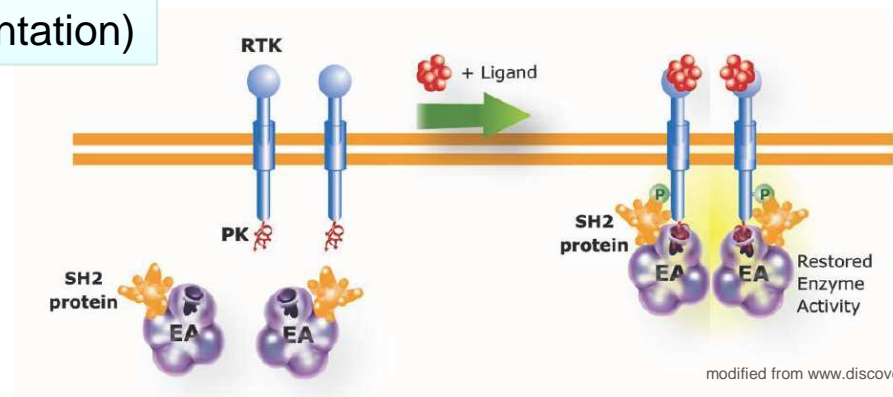
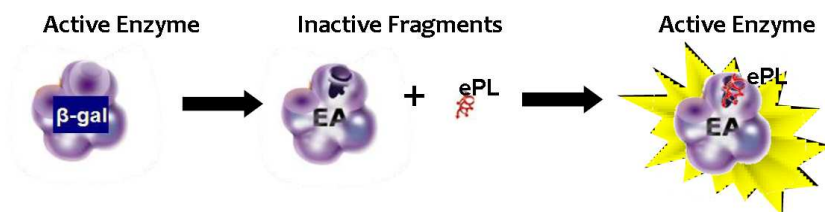
# General options for cell-based kinase assays



Assay technology	Provider	comments	critical tools
Western Blot	various	heterogenous low throughput	phospho-antibody
Incell Western	various	high content imaging	phospho-antibody
ELISA-type	various	heterogenous	antibody pair
Luminex	Millipore	heterogenous	antibody pair
Surefire / ALPHAscreen	PerkinElmer	homogenous	antibody pair
HTRF	Cisbio	homogenous	antibody pair
Lanthascreen	LifeSciences	homogenous	phospho-antibody recombinant cell-line
EFC - Enzyme Fragment Complementation	DiscoverRx	homogenous	antibody-free recombinant cell-line
Survival assay	Advanced Cellular Dynamics / Carna	Tyr-kinase activity restores survival of Ba/F3 cells	antibody-free recombinant cell-line
Pathway reporter assays	various	downstream gene activation	antibody-free recombinant cell-line

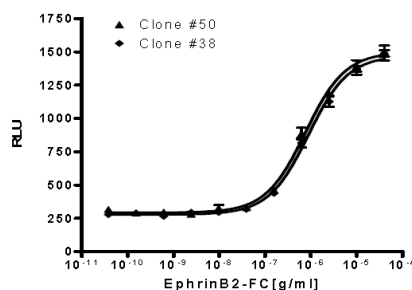
# EFC-assays for Tyr-Kinases (DiscoverRx)

## Technology EFC (enzyme fragment complementation)

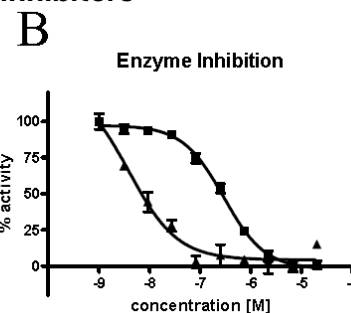
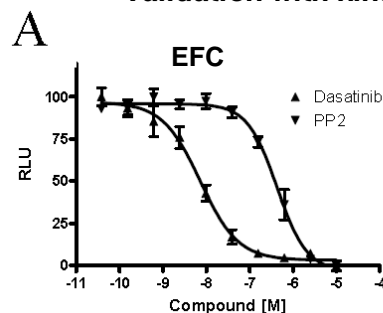


## EphB4 cellular kinase activity assayed using an enzymatic protein interaction system (Wehrman et al. 2013 ADT)

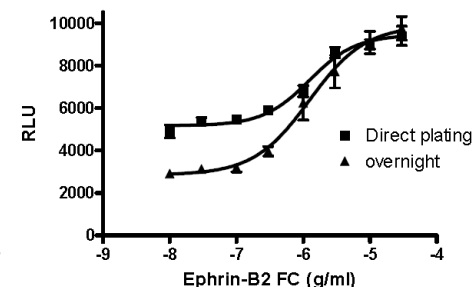
### Agonist stimulation



### Validation with kinase inhibitors



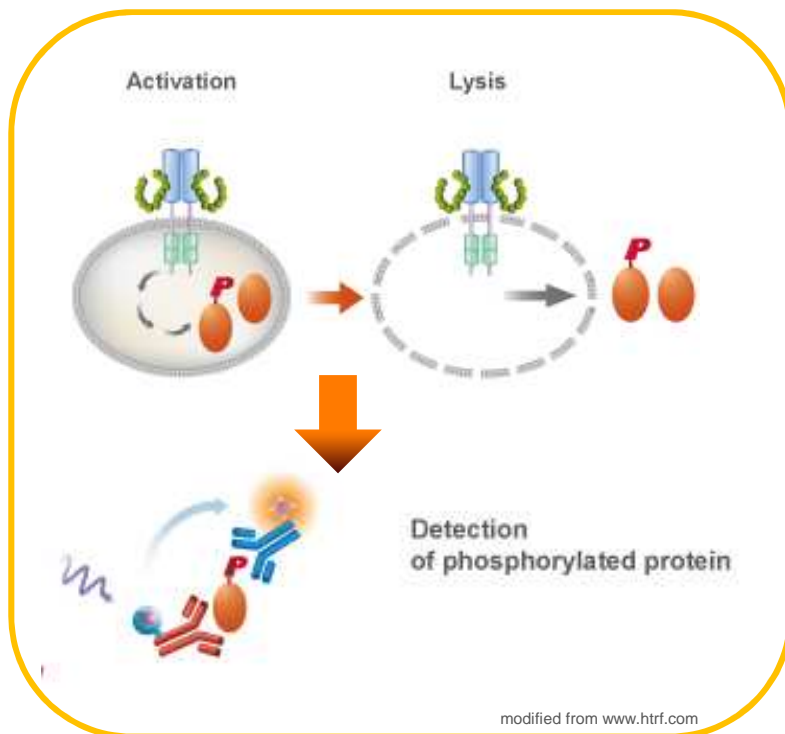
### Frozen cell and direct assay format



- EFC-kinase technology is suitable for miniaturized frozen cell assays
- limitations: largely restricted to Tyrosine kinases, requires recombinant cells

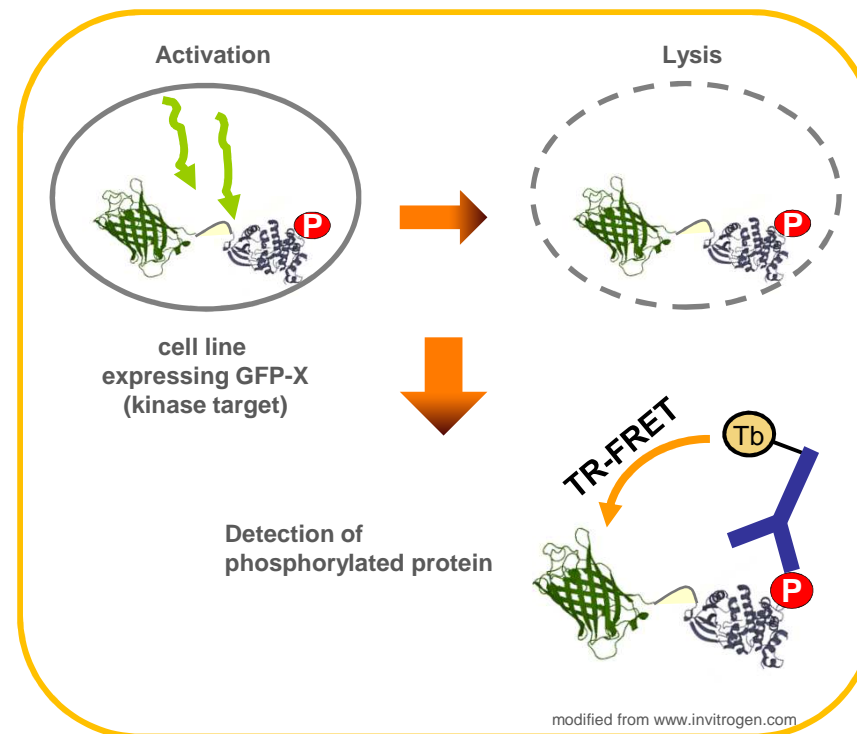


# TR-FRET systems: HTRF and Lanthascreen



## **HTRF**

- endogenous phospho-protein
- any cell line expressing protein
- antibody pair required
- HTRF-detection at 620/665 nm

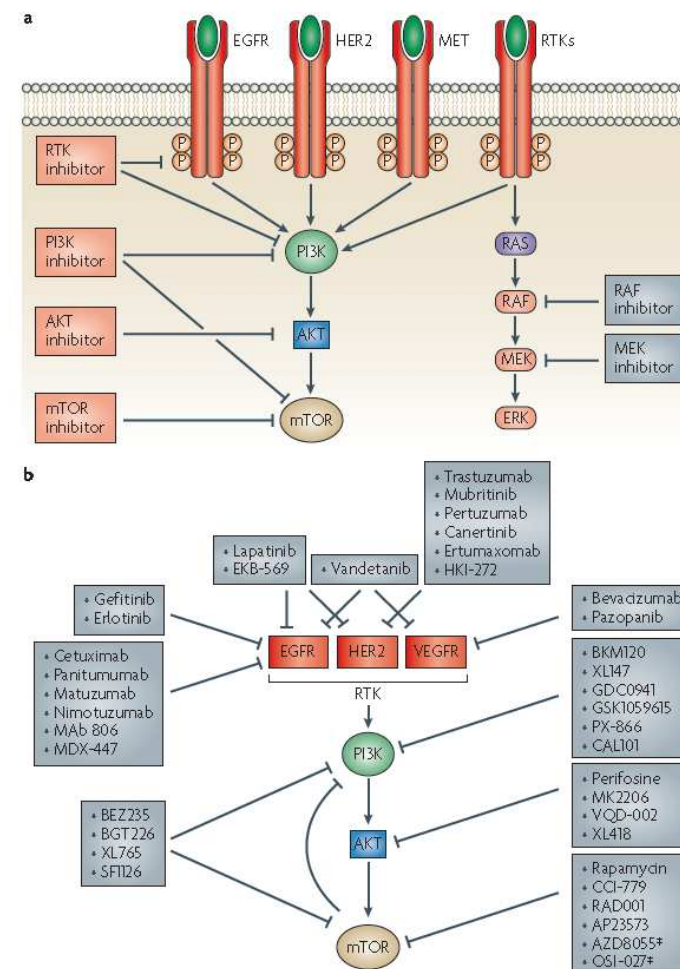
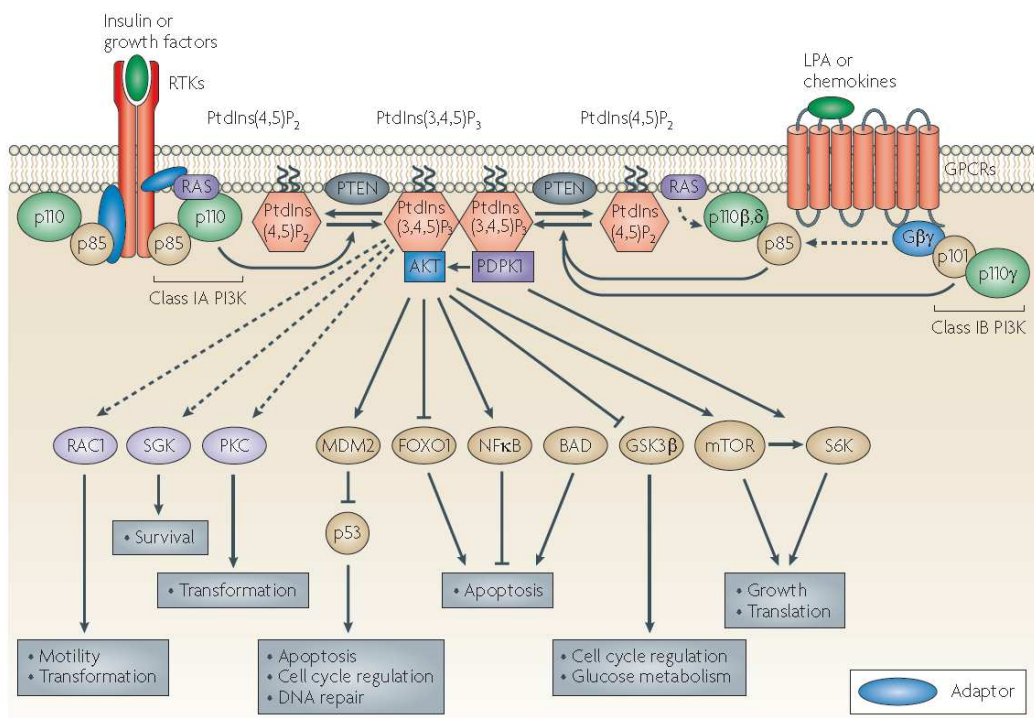


## **Lanthascreen**

- overexpressed GFP-phospho-protein
- stable cell line or BacMam transient
- only phospho antibody required
- Lanthascreen detection at 490/520

***HTRF and Lanthascreen assays for  
the PI3K / Akt / mTOR pathway***

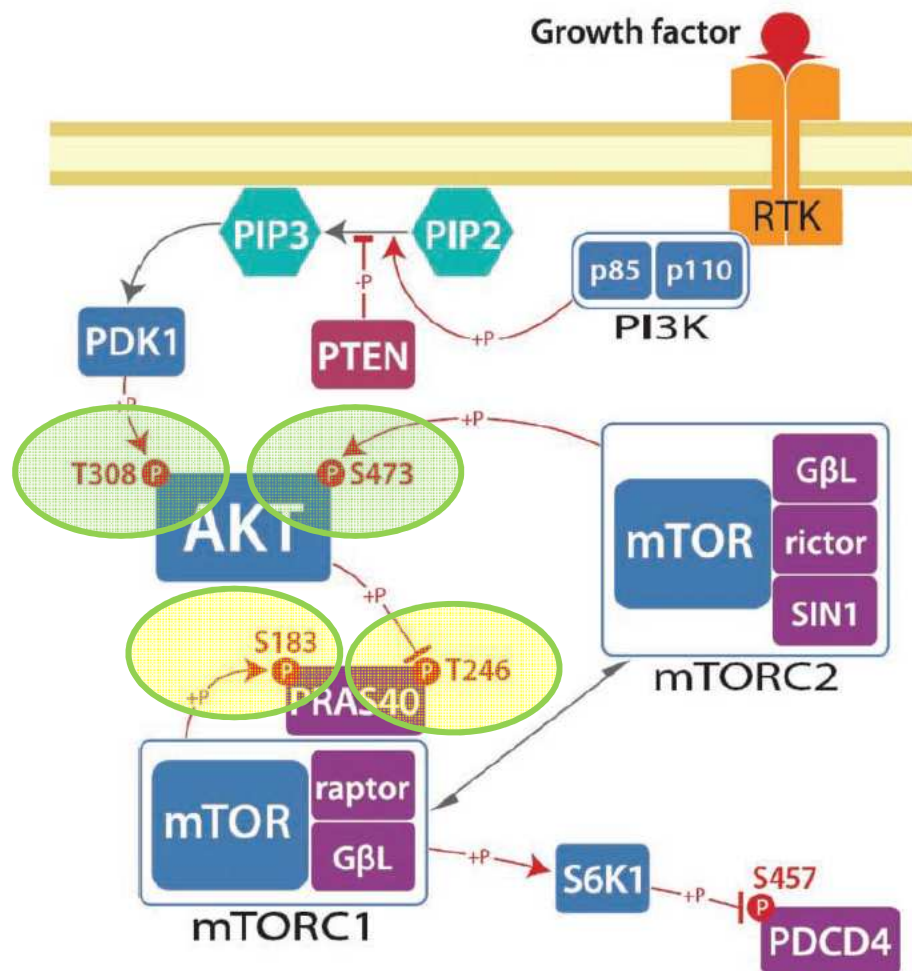
# Introducing the PI3K / Akt / mTOR pathway



- PI3K / Akt pathway is central to cancer formation
- Chemotherapeutic approaches against multiple targets are in the pipelines

Liu 2009 Nat. Rev Drug Discovery

# Assaying the PI3K / Akt / mTOR pathway



(modified from Carlson 2009)

## *Lanthascreen study goals:*

- Evaluate BacMam Lanthascreen technology
- GFP-Akt @ two sites (pT308 and pS473)
- GFP-PRAS40 @ two sites (S183 and T246)
- test different cell backgrounds
- validate with reference inhibitors

## *HTRF study goals:*

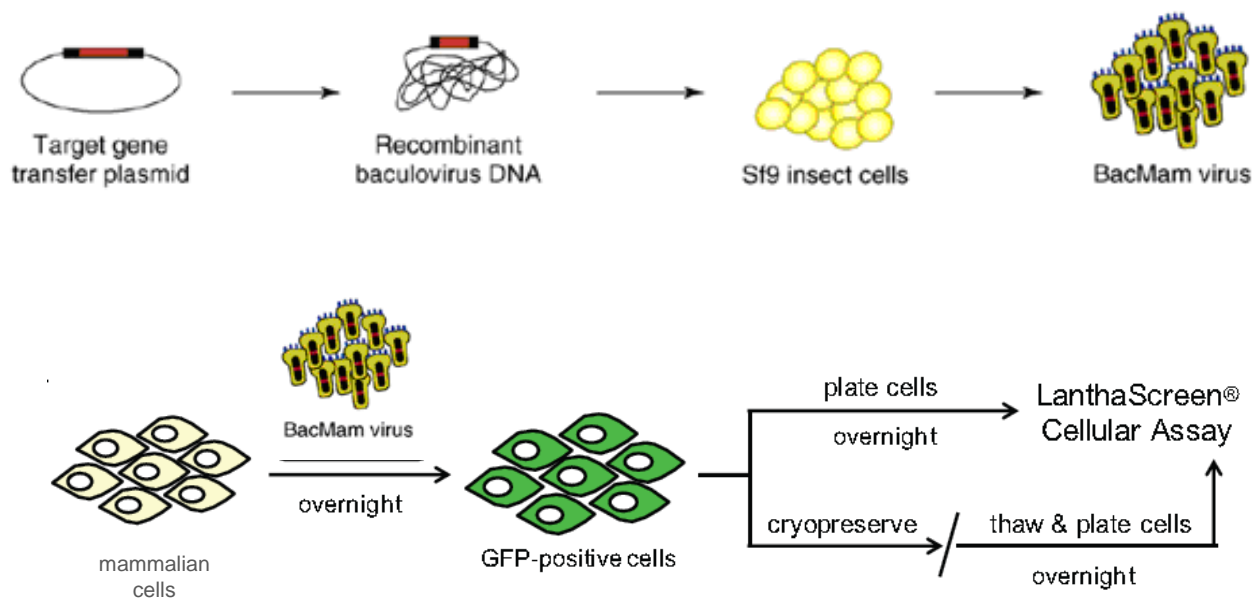
- set up cell-based mTOR kinase activity assay for HTS
- identify optimal cancer cell background
- validate with reference inhibitors

→ validate HTS-compatibility with both formats using a miniaturized 384 well focussed screen

# Lanthascreen Assay Development



*BacMam technology combined with  
Lanthascreen Cellular assay*



(modified from Kost 2007, Carlson 2010)



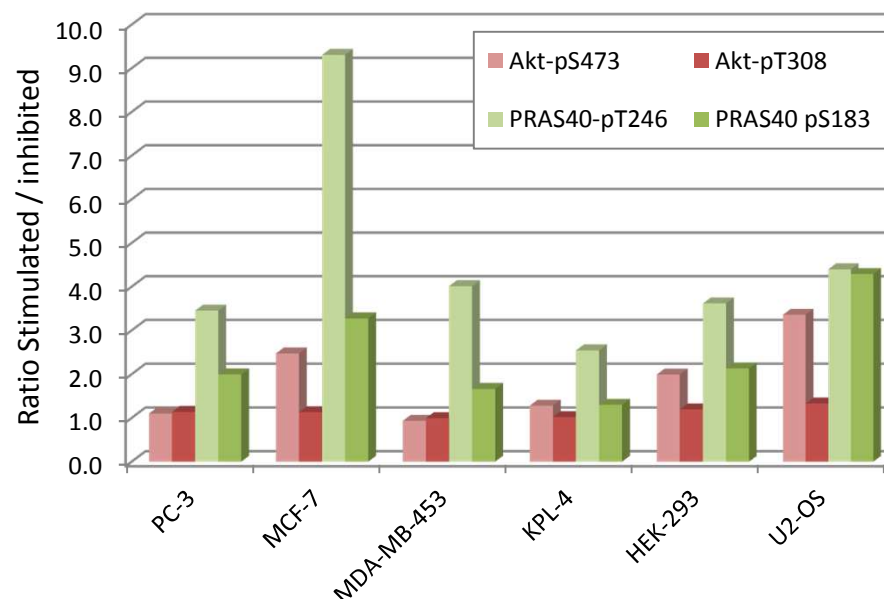
# Lanthascreen – cellular background



## Lanthascreen study goals:

- Evaluate BacMam Lanthascreen technology
- GFP-Akt @ two sites (pT308 and pS473)
- GFP-PRAS40 @ two sites (S183 and T246)
- test different cell backgrounds
- validate with reference inhibitors

Cell line	Type	PI3K Pathway mutation
PC-3	Human prostate cancer	PTEN negative
MCF-7	Human breast cancer	PI-3-Kinase mutation E545K
MDA-MB-453		PI-3-Kinase mutation H1047R
KPL-4		
HEK-293	Human embryonal kidney	WT
U2-OS	Human osteosarcoma	WT

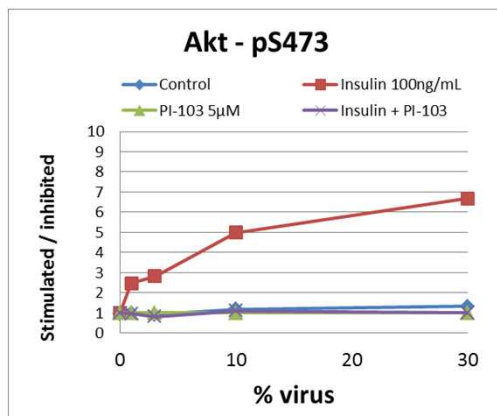


→ MCF7 PRAS40-pT246 selected for further optimization work

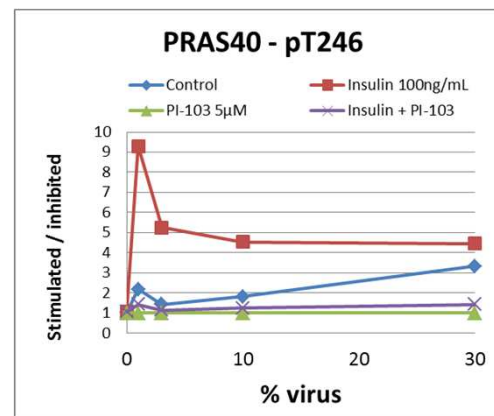
# Lanthascreen – virus efficiency



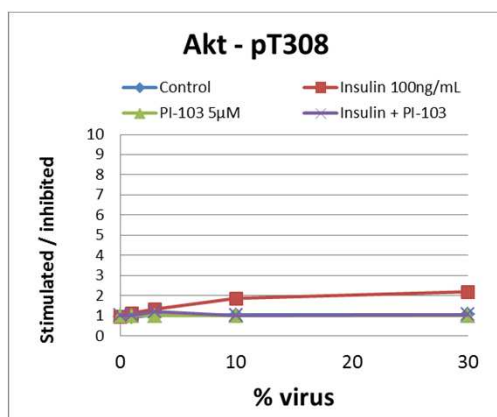
30% virus  
 $S/B_{\max} = 7$



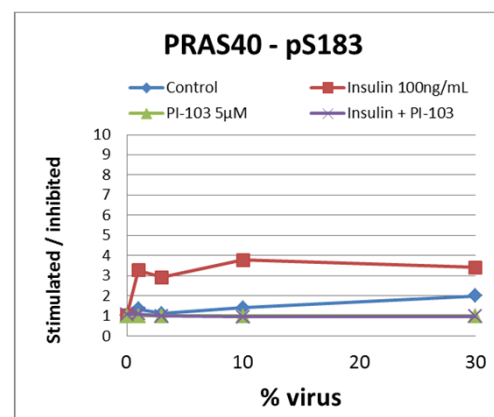
1% virus  
 $S/B_{\max} = 9$



10% virus  
 $S/B_{\max} = 2$



1% virus  
 $S/B_{\max} = 3$



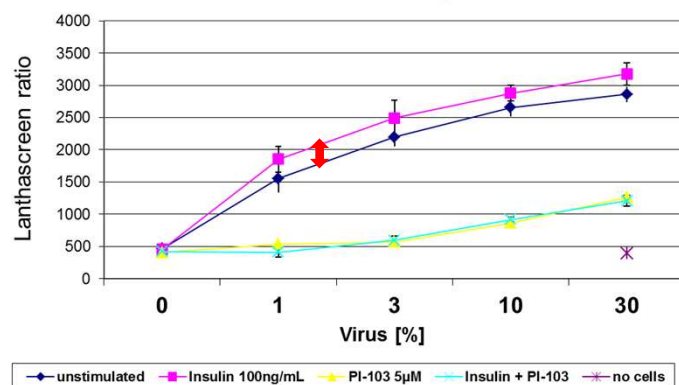
- different virus efficiency: PRAS40 with highly efficient expression ( $S/B_{\max}$  reached at 1%)
- differences between phosphosites → antibody quality (?)

# Lanthascreen – cellular background



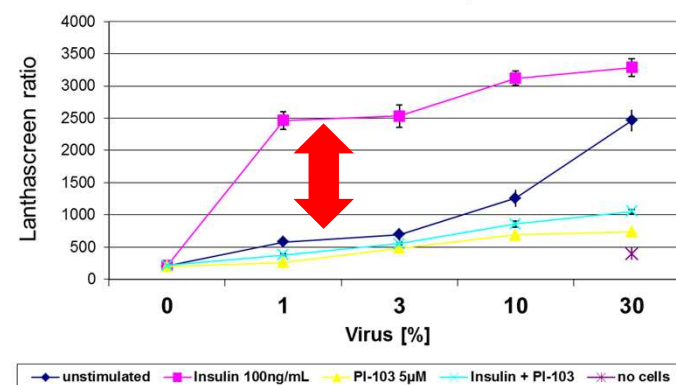
PTEN negative

PC3 cells - PRAS40 pThr246

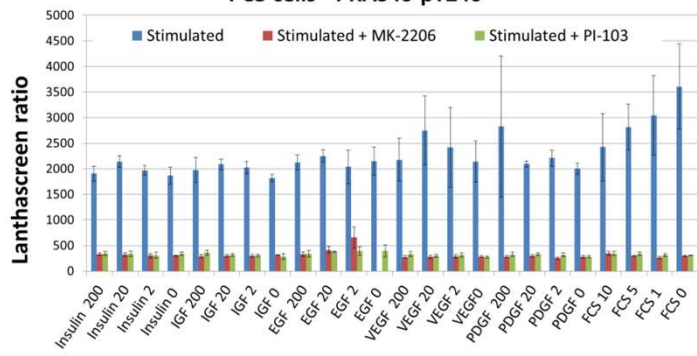


PI3K E545K

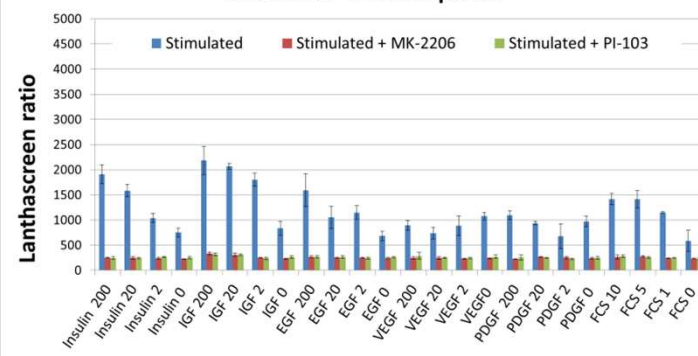
MCF7 cells - PRAS40 pThr246



PC3 cells - PRAS40-pT246



MCF7 cells - PRAS40-pT246

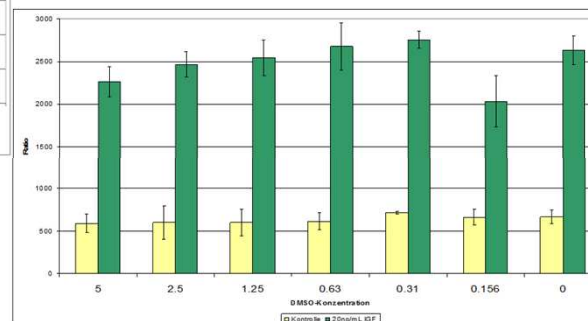
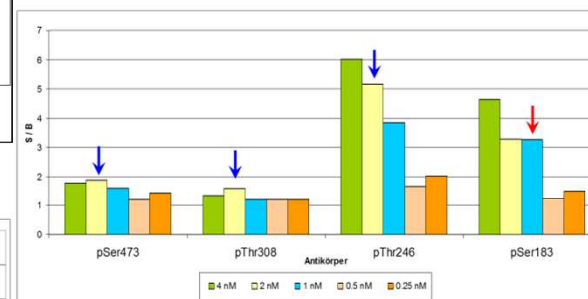
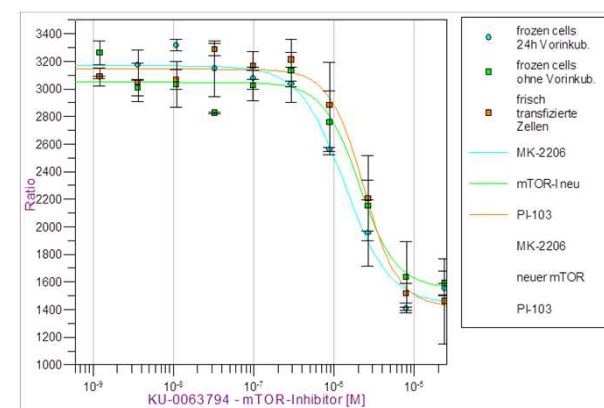
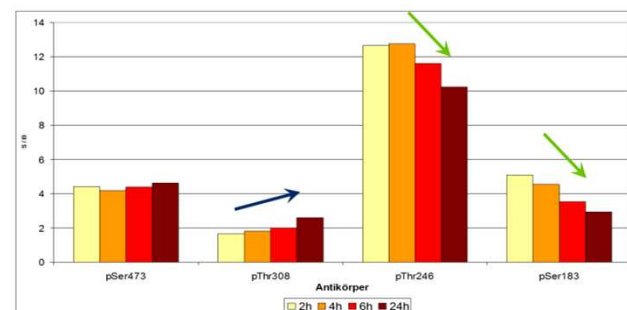
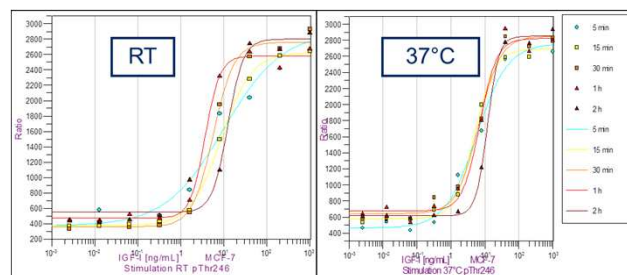


- PC3-cells - in contrast to MCF7 - have fully stimulated pathway  
cave → different pathway mutations have different impact on basal activation

# Lanthascreen - optimization



- stimulation time
- stimulation temperature
- antibody concentration
- time cell lysis → read
- DMSO sensitivity
- Volume 10 µl → 5 µl
- frozen cells



# Lanthascreen - optimized assay protocol

day 1:

seed cells into of MCF7 cells T-flask (frozen or continuous culture)

day 2:

transduce MCF7 cells with 1% PRAS40 virus in T-flask

day 3:

harvest transduced cells and prepare cell suspension in medium + 1% FCS

dispense 3  $\mu$ l (5000 cells/well) into assay-ready MTP containing 50 nl cpd.



2 h @ 37°C

add 1  $\mu$ l lysis/detection mix



2 h @ RT

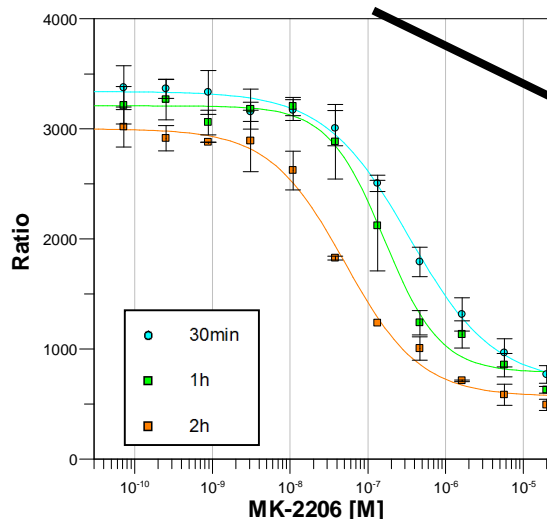
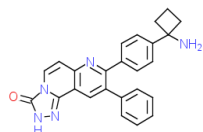
read TR-FRET



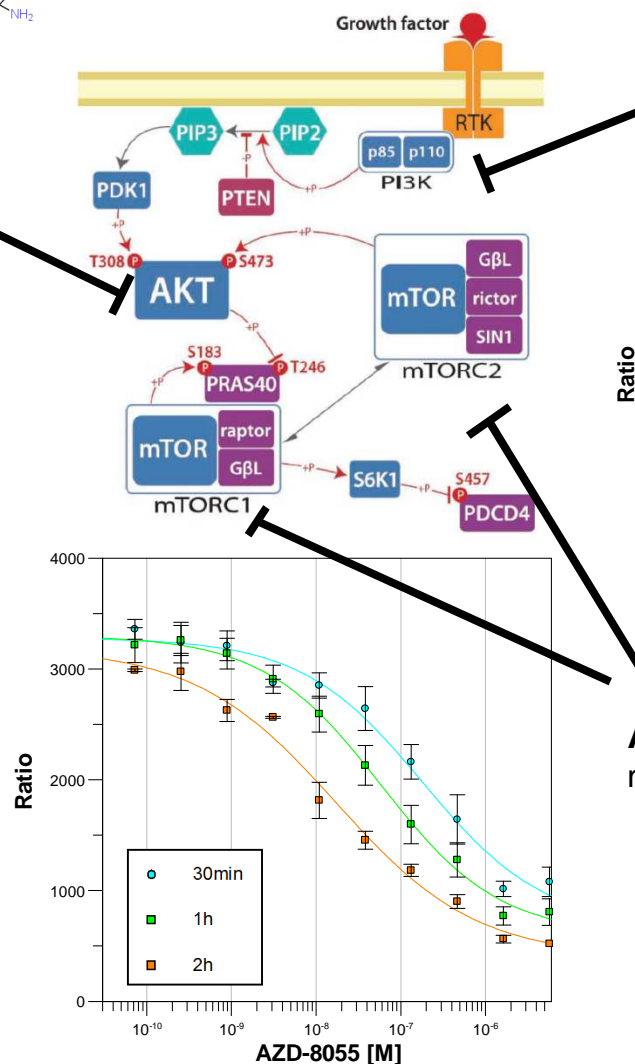
# Lanthascreen - inhibitor validation PRAS40-pT246



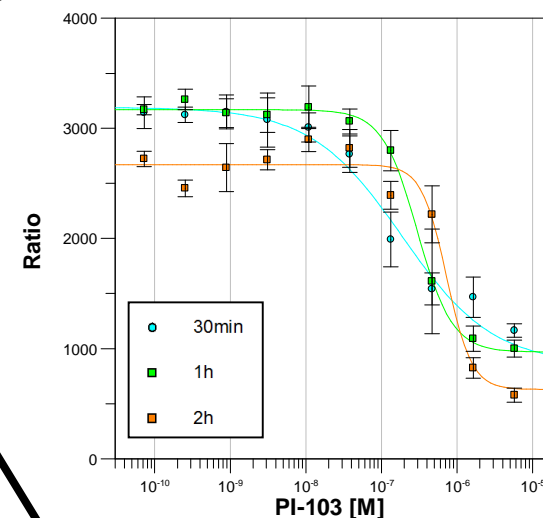
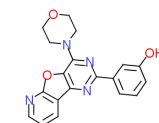
**MK-2206**  
allosteric Akt inhibitor



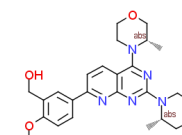
- potency of pathway inhibitors depends on incubation time
- Akt and mTOR inhibitors improved IC50 after 2 h



**PI-103**  
PI3K kinase inhibitor



**AZD-8055**  
mTOR kinase inhibitor

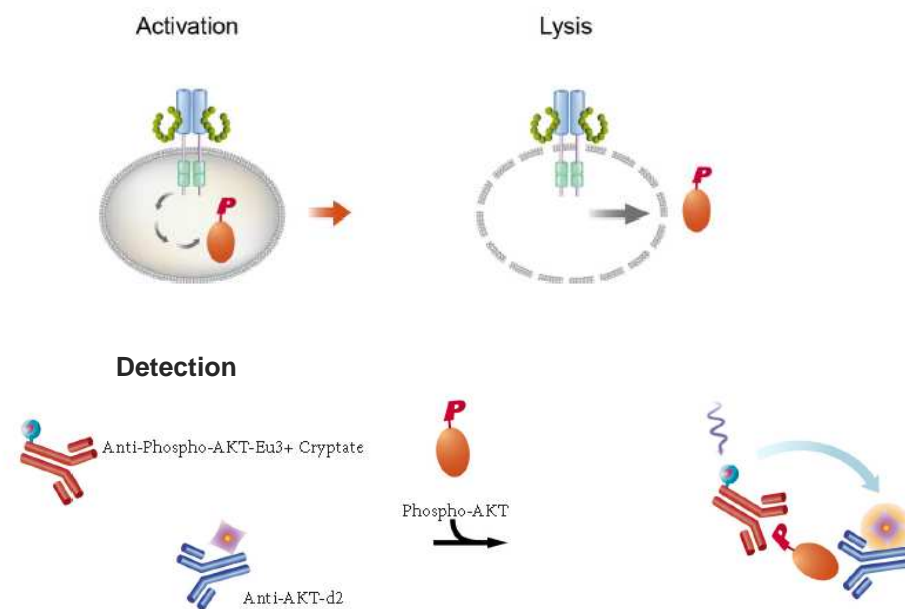


# HTRF – assay

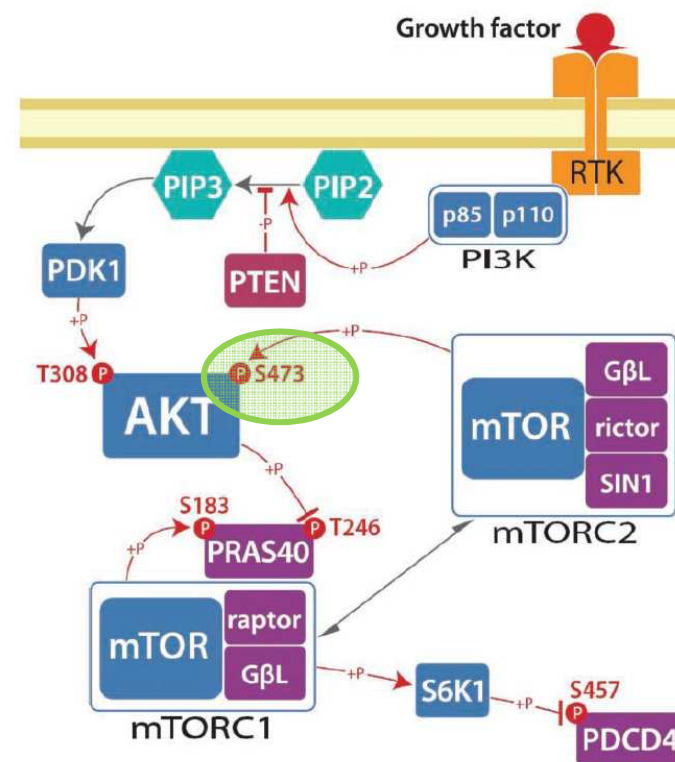


## *HTRF study goals:*

- set up cell-based mTOR kinase activity assay for HTS using pAkt Ser473 readout
- identify optimal cancer cell background
- validate with reference inhibitors



(from Cisbio product insert)

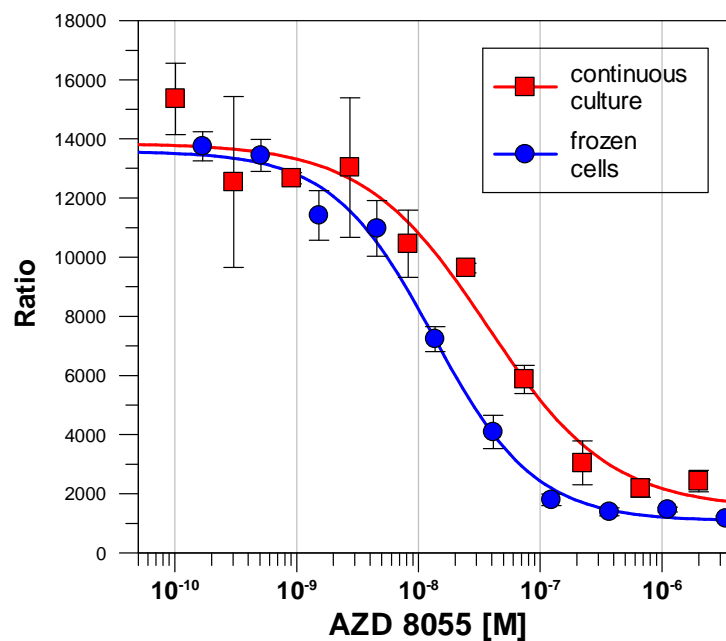
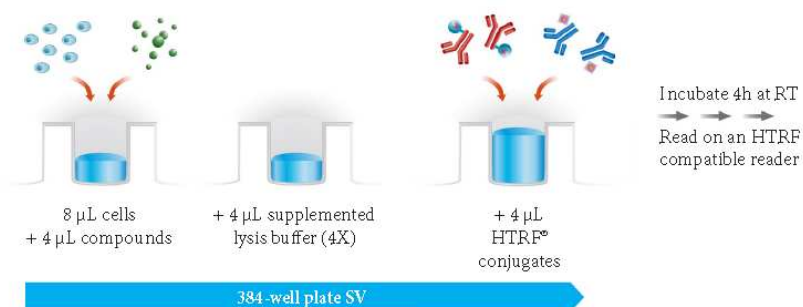


(modified from Carlson 2009)

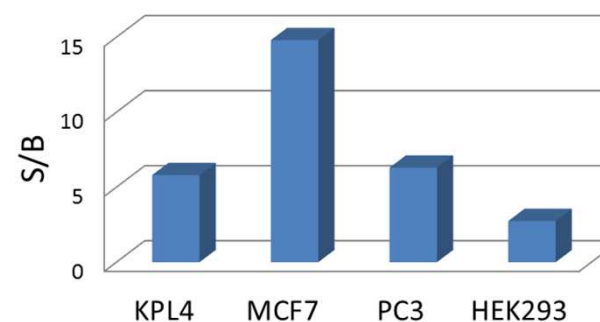
# HTRF – cell lines



One-plate assay protocol



HTRF - pAkt-S473



- MCF7 cells strongest S/B
- frozen cell assay
- suspension cell format
- miniaturized to 1536 well

# HTRF - optimized assay protocol



day 1:

thaw frozen MCF7 cells and prepare cell suspension in medium + 1% FCS

dispense 3  $\mu$ l (4000 cells/well) into assay-ready MTP containing 50 nl cpd.



30 min @ 37°C

add 1  $\mu$ l lysis buffer



1 h @ RT shake

add 4  $\mu$ l antibody detection mix

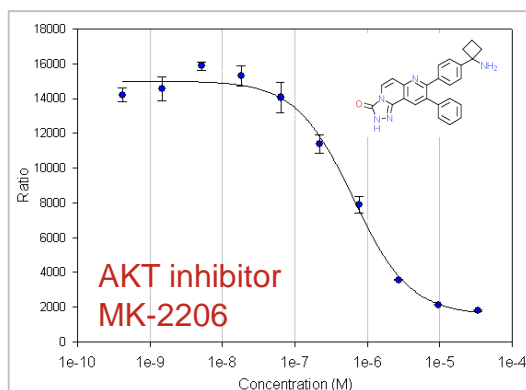


20 h @ RT

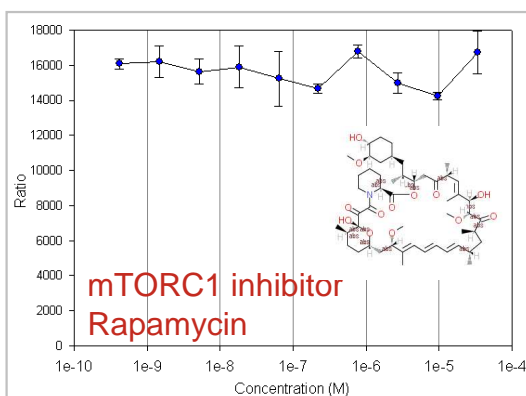
day 2:

read TR-FRET

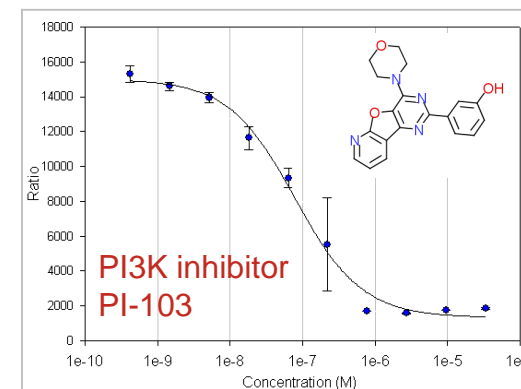
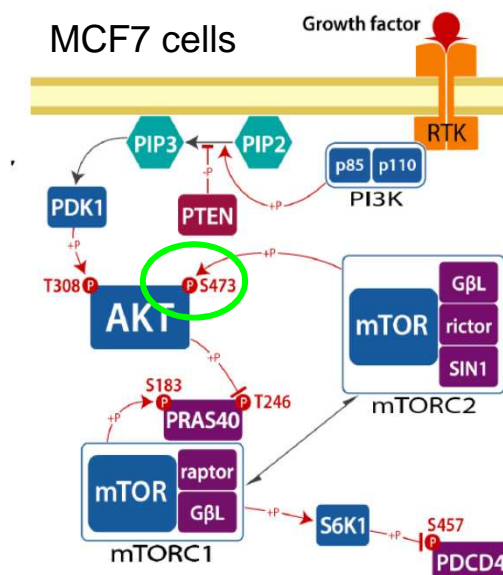
# HTRF – pathway inhibitor validation



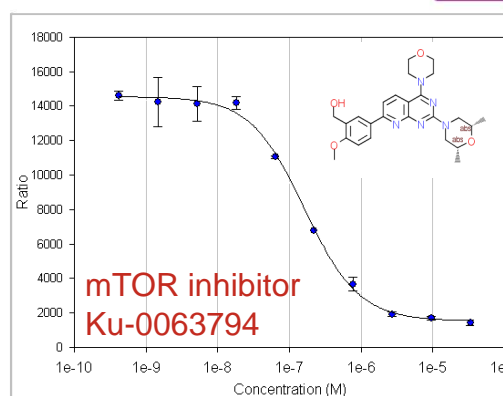
IC50 = 635nM



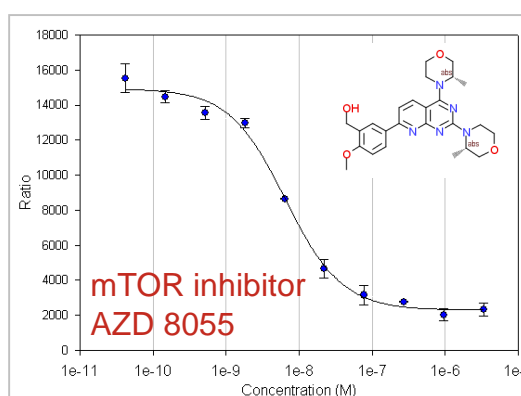
no effect



IC50 = 79nM



IC50 = 163nM



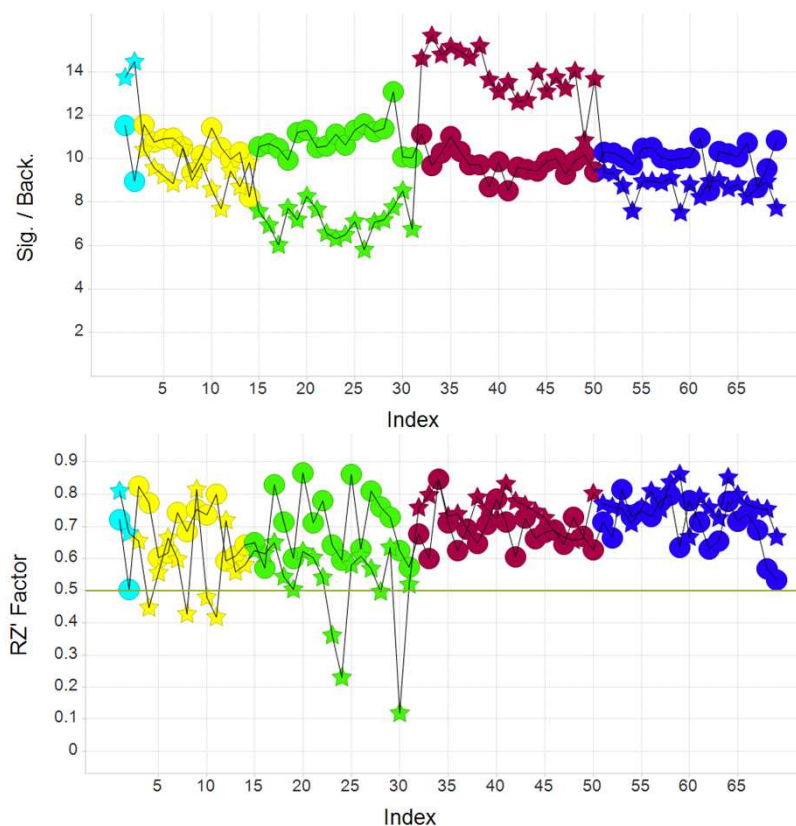
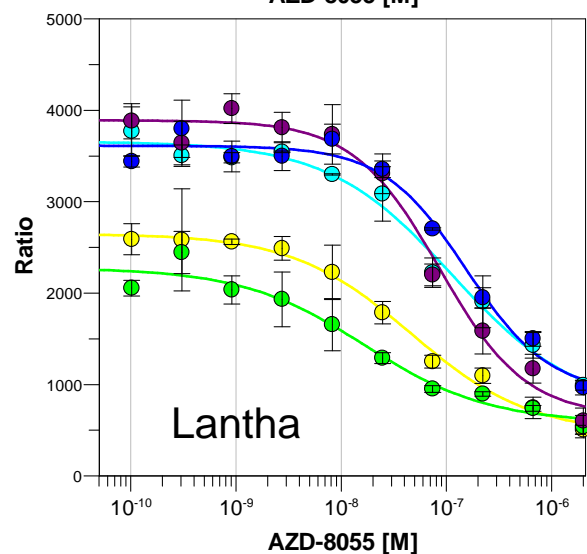
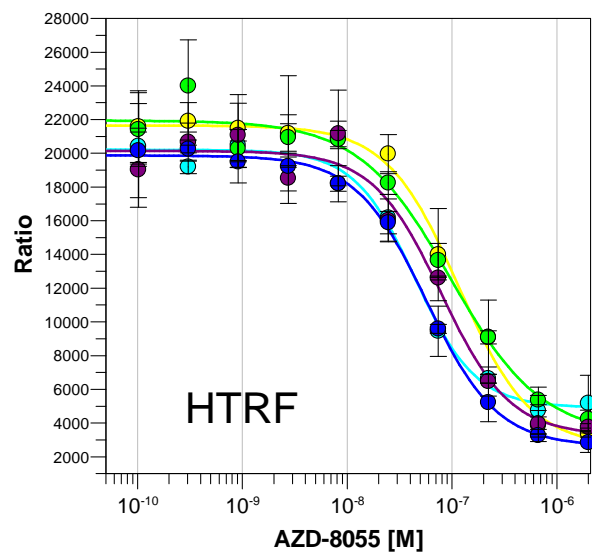
IC50 = 7nM



### ***Focussed medium throughput screen:***

- BacMAM Lantha PRAS40-pT246
  - HTRF Akt-pS473
- 24.300 compounds, kinase targeted library
  - 384 well single format
  - 69 MTPs
  - 5 parallel assay runs (Lantha and HTRF)

# Focussed medium throughput screen



shape by assay:

● HTRF

★ Lanthascreen

color by run

● Run 1

● Run 2

● Run 3

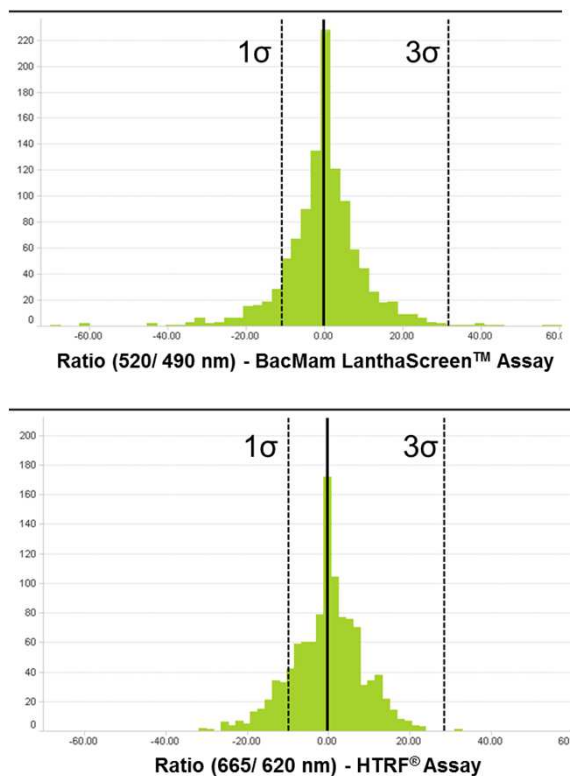
● Run 4

● Run 5

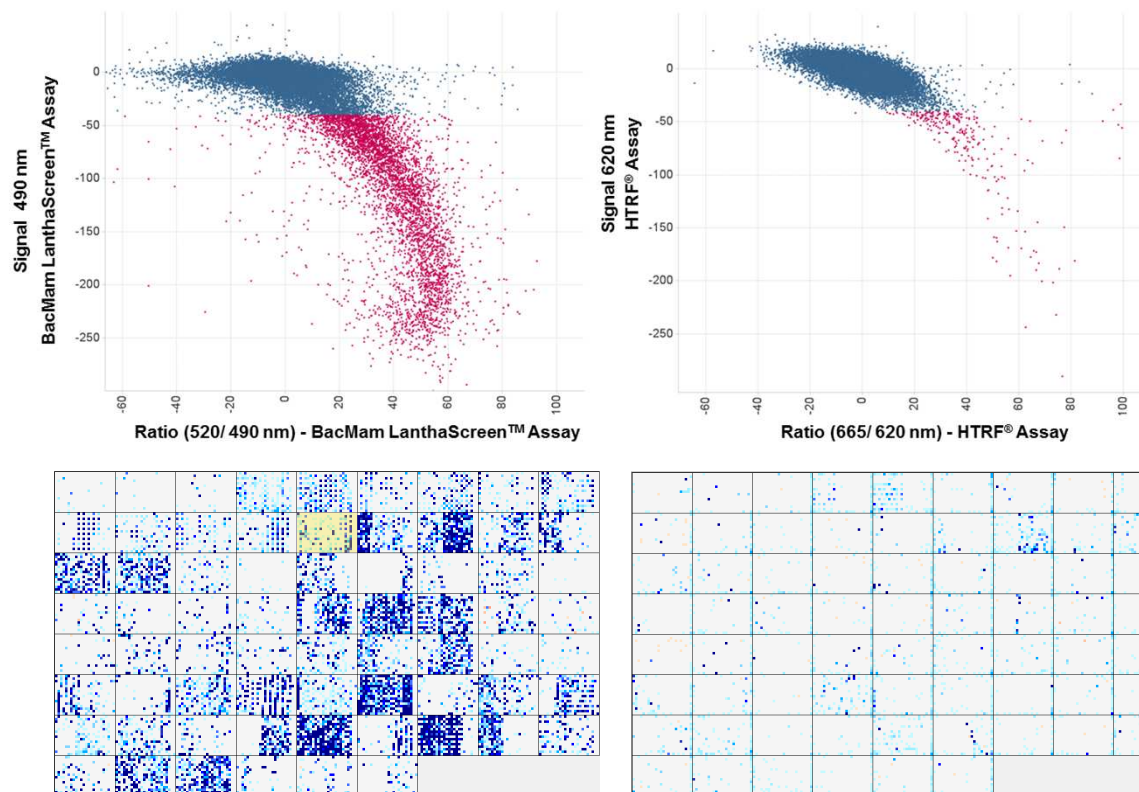
- Similar sensitivity of both assays against mTORi
- HTRF: stable and robust assay performance
- Lanthascreen: day-to-day variability in S/B

# Focussed medium throughput screen

Distribution neutral controls:

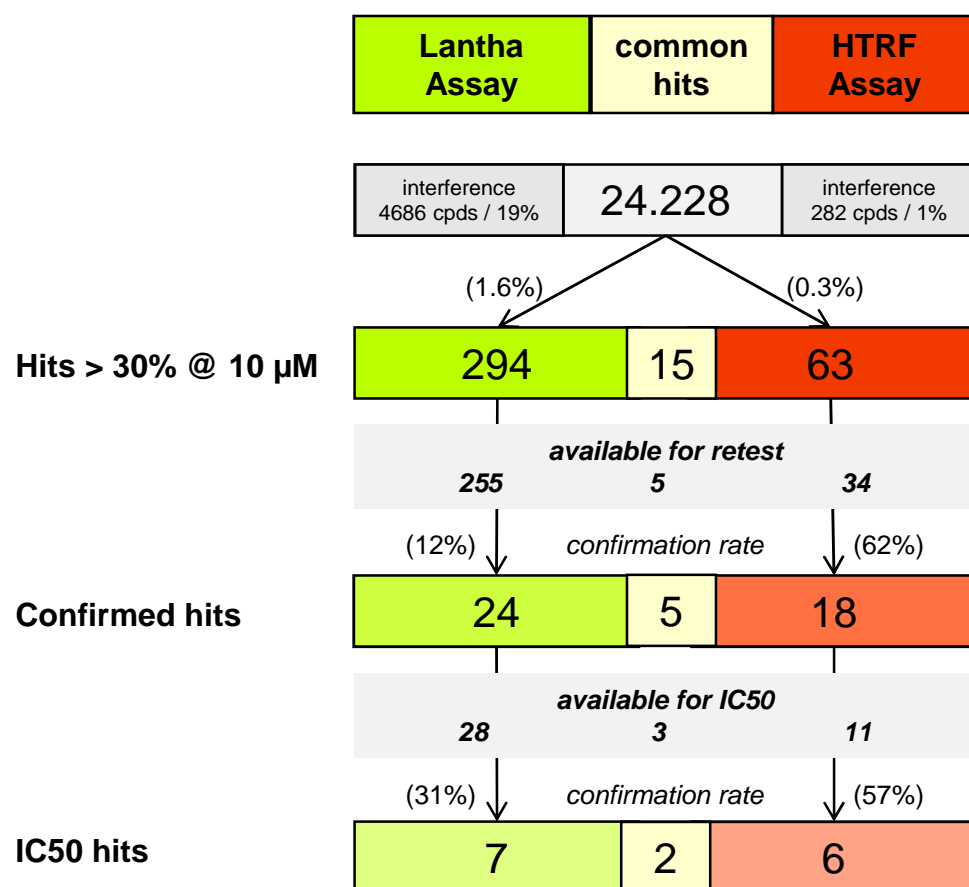


Distribution compounds / donor interference:



- Neutral controls: similar distribution for both assays
- Compounds: strong interference in BacMAM donor channel

# Focussed medium throughput screen



	Lantha	HTRF
Interference	high (19%)	low (1%)
Hit rate	high (1.6%)	low (0.3%)
Confirmation rate	low (12%)	high (62%)

# Summary & Conclusions



- Kinases remain interesting target class in pharmaceutical research
- Specific and efficient cell-based kinase assays are essential in pharma research projects
- Homogenous assay systems fit best to Bayers lead discovery platform
- Positive experience at Bayer with EFC-technology, HTRF and Lanthascreen
- EFC-technology interesting for Tyr-Kinase uHTS
- Lanthascreen technology positioned for secondary testing
- HTRF positioned for uHTS and secondary testing



# Acknowledgements



Anja Kretzschmar

Sabrina Reimann

Katrin Nowak-Reppel

Dagmar Zeggert-Springer

Karsten Parczyk

Monika Gross (Beuth University of Applied Science)



data published in master thesis of Anja Kretzschmar, November 2011



Science For A Better Life

Thank you!