



Bayer HealthCare
Bayer Schering Pharma

Cell-based Receptor Tyrosine Kinase Screening Reloaded

Development of a new cellular assay system for a target protein so far resistant against HTS formats

Benjamin Bader

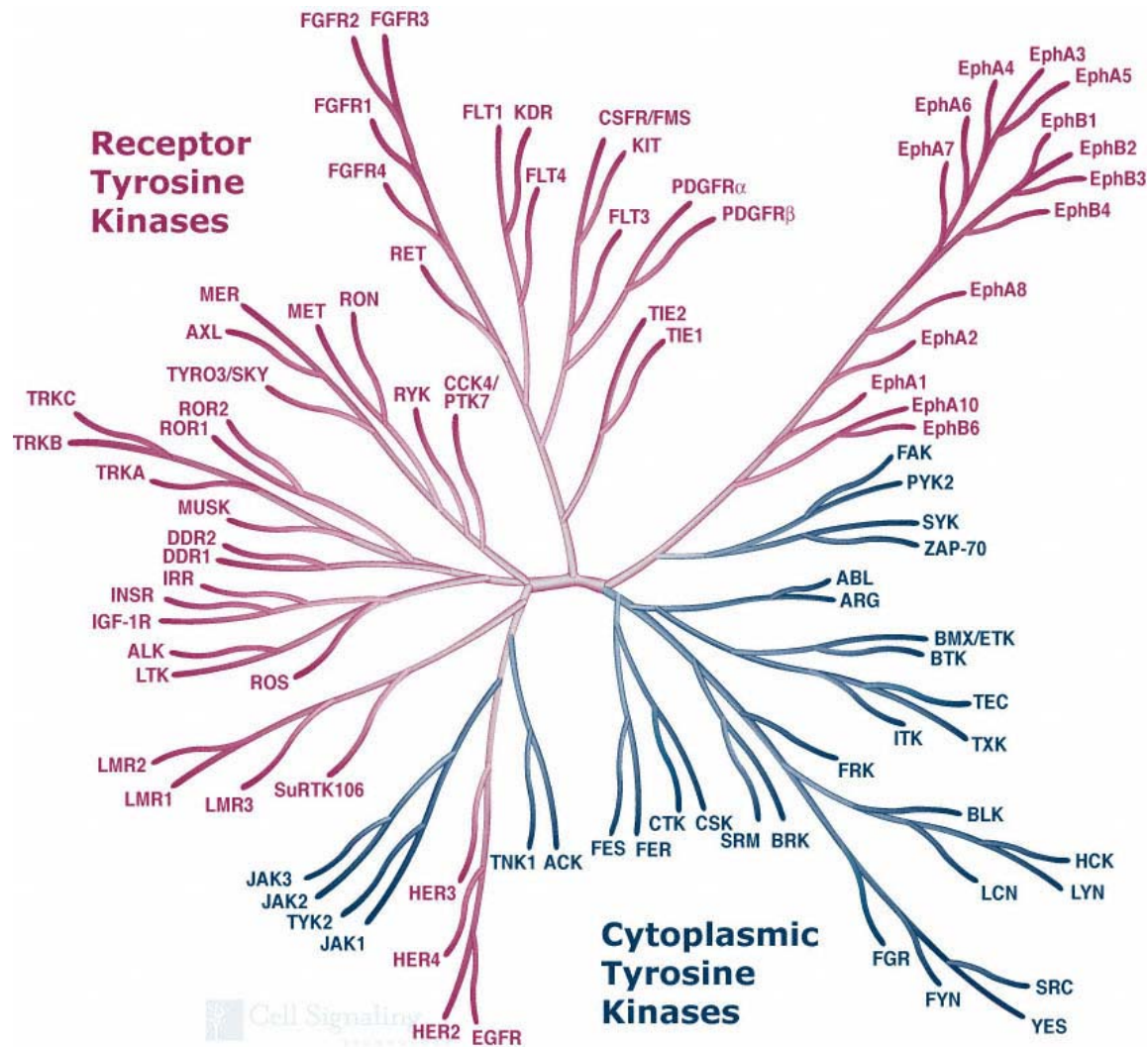
SBS-meeting Lille 2009

- Rationale for cell-based screening of RTKs
- Some project history
- The DiscoverRx workplan
- Assay development data
- Assay transfer, adaptation and optimization
- Summary and Outlook



Bayer HealthCare
Bayer Schering Pharma

Receptor Tyrosine Kinases



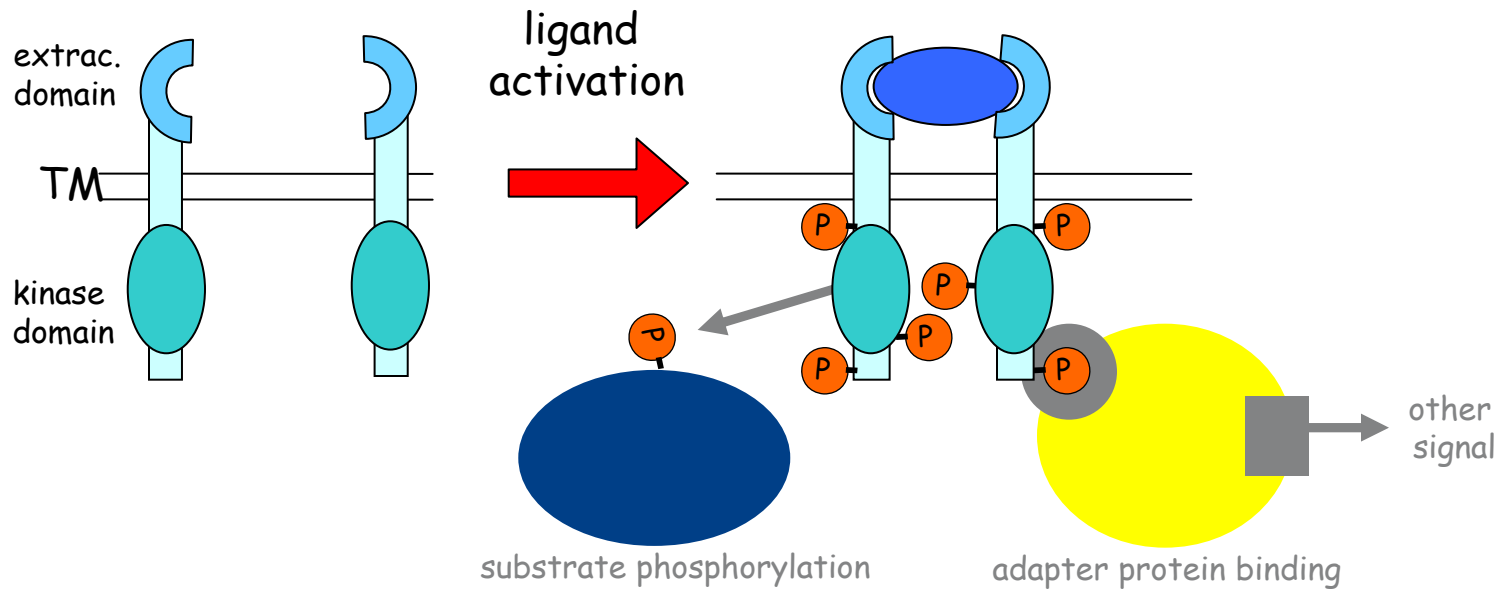
RTKs play important roles in fundamental cellular processes like

- proliferation,
- differentiation,
- migration,
- metabolism
- survival

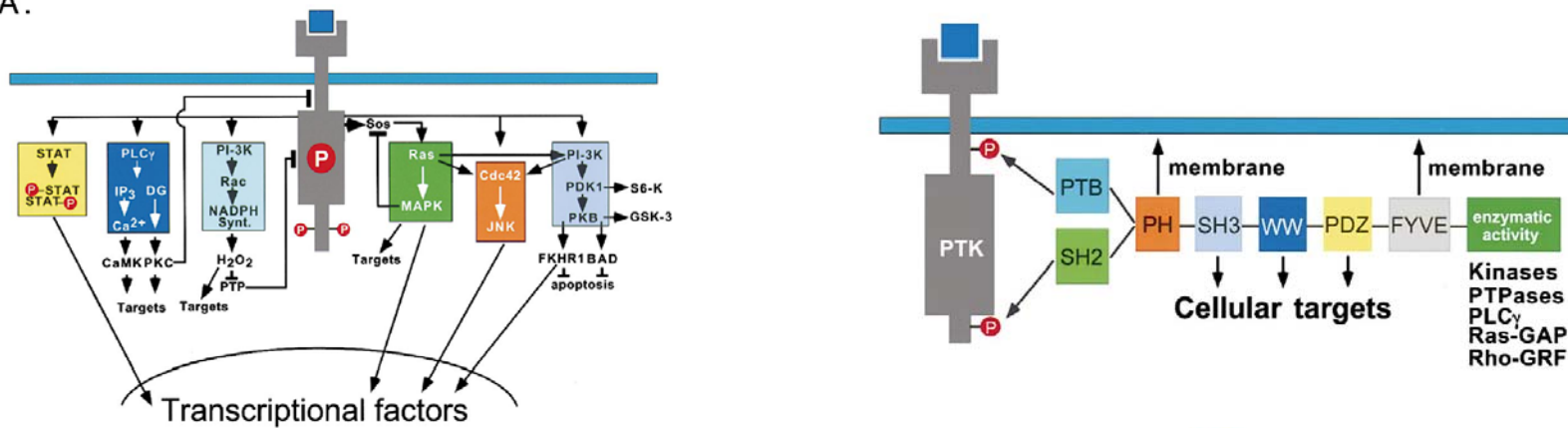


Bayer HealthCare
Bayer Schering Pharma

General scheme of activation and signal transduction of RTKs



A.



Adapted from: Schlessinger, 2000, Cell 103, 211-225

SBS 2009, Bader, page 4



Bayer HealthCare
Bayer Schering Pharma

Advantages of a cell-based RTK assay in comparison to enzymatic kinase assays

- **full length protein**, not only kinase domain
- physiological **membrane environment** of RTK
- **native substrate** compared to peptides in enzyme assays
- the RTK is present in its **inactive state** when compounds are added
- possibility to find **prodrugs** which can be activated inside the cell
- **inhibitors of ligand binding site** can be found



***Chance to find new inhibitors
with different binding modes***



Bayer HealthCare
Bayer Schering Pharma

Project history

RTK target protein has been screened using enzymatic assays



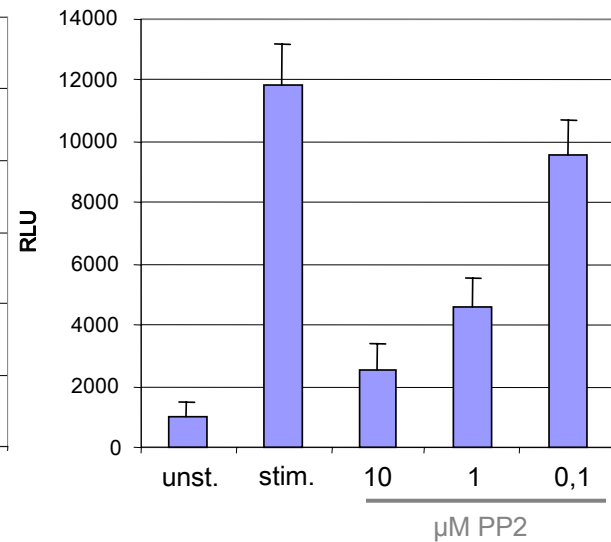
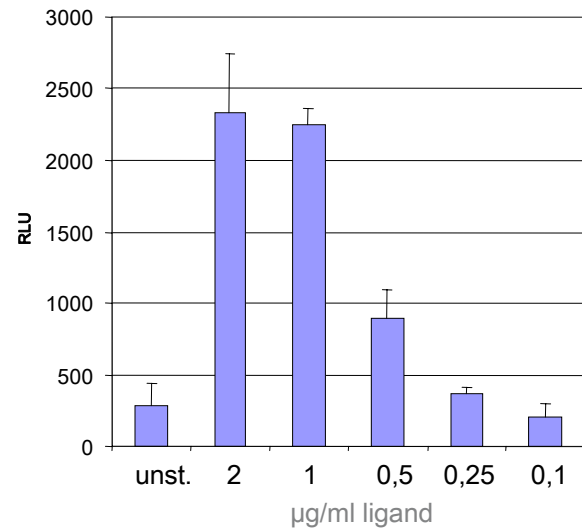
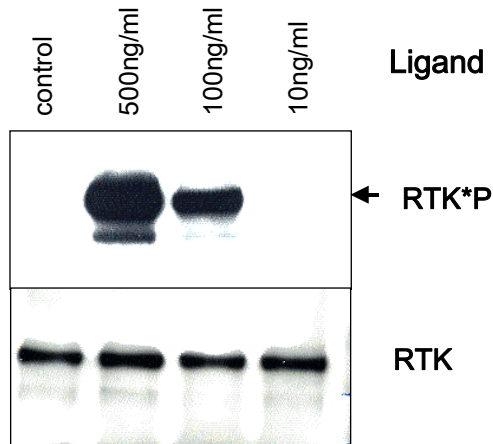
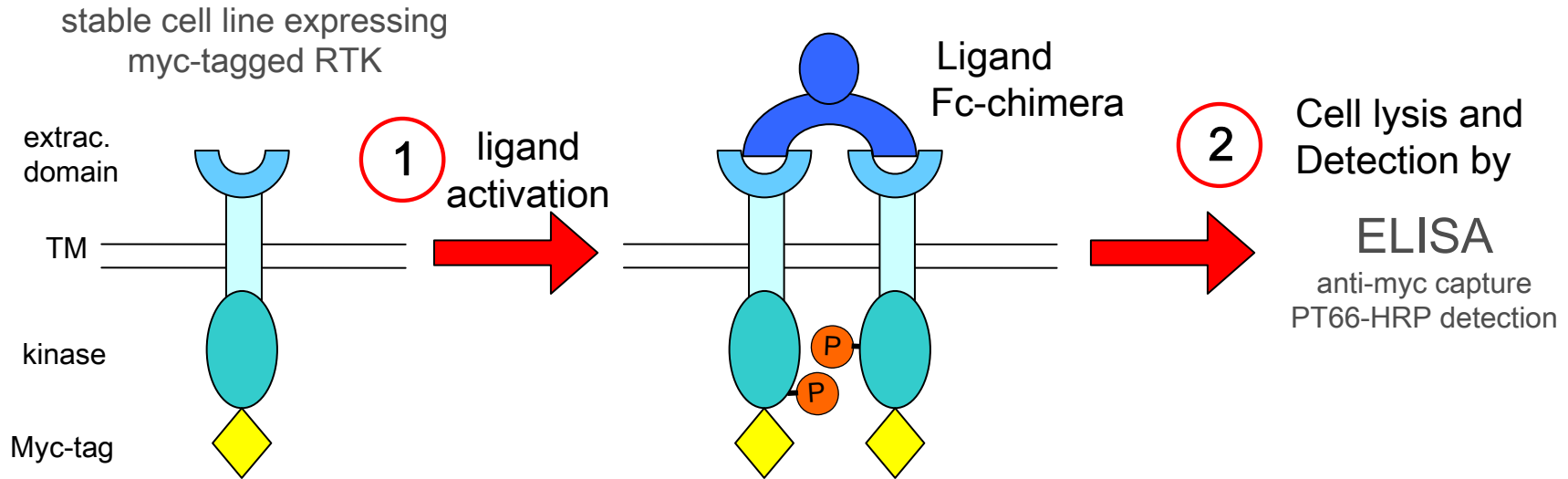
- considerable amount of IC50 hits with activity in cellular assays
- but: no lead structure identified

Cellular ELISA assay has been set-up as secondary assay



Bayer HealthCare
Bayer Schering Pharma

Cellular autophosphorylation ELISA in 384 well



Project history

RTK target protein has been screened using enzymatic assays



- considerable amount of IC50 hits with activity in cellular assays
- but: no lead structure identified

Cellular ELISA assay has been set-up as secondary assay



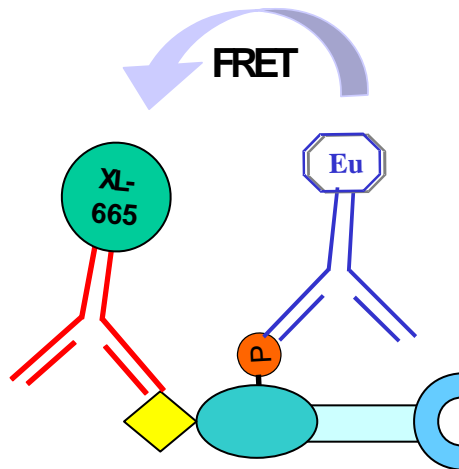
- worked well in 384 well
- but: too complicated for full HTS
- and: ELISA in 1536 well !?!

Homogenous formats were developed (HTRF)

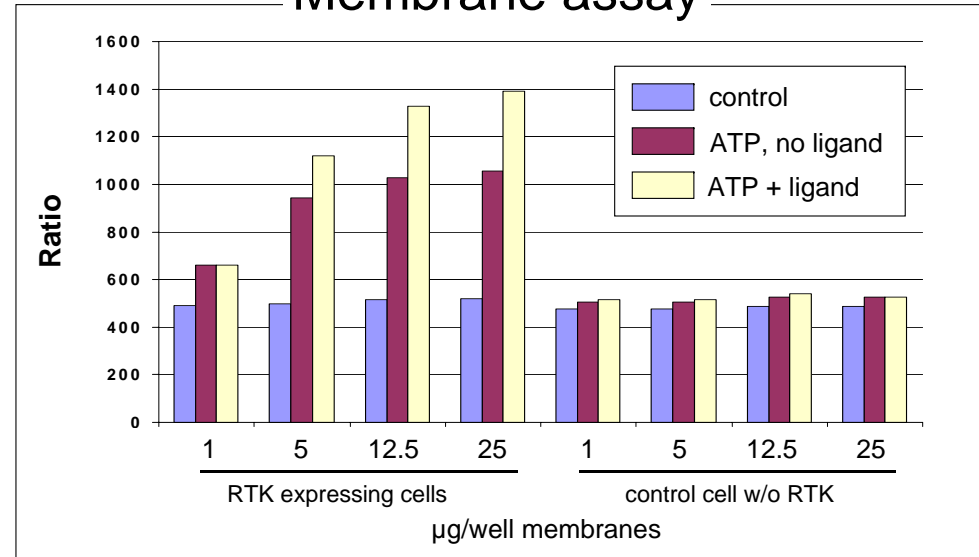


Bayer HealthCare
Bayer Schering Pharma

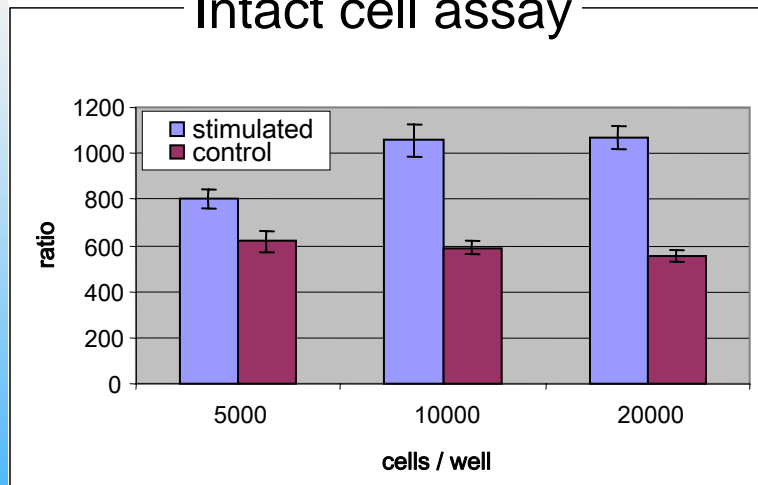
HTRF detection



Membrane assay



Intact cell assay



Membrane assay:

- Homogenous format possible
- Strong ligand independent signal

Intact cells assay:

- S/B too low for 384 well assay
- Too many cells needed
- protocol contains several wash steps



Bayer HealthCare
Bayer Schering Pharma

Project history

RTK target protein has been screened using enzymatic assays



- considerable amount of IC50 hits with activity in cellular assays
- but: no lead structure identified

Cellular ELISA assay has been set-up as secondary assay



- worked well in 384 well
- but: too complicated for full HTS
- and: ELISA in 1536 well !?!

Homogenous formats were developed (HTRF)



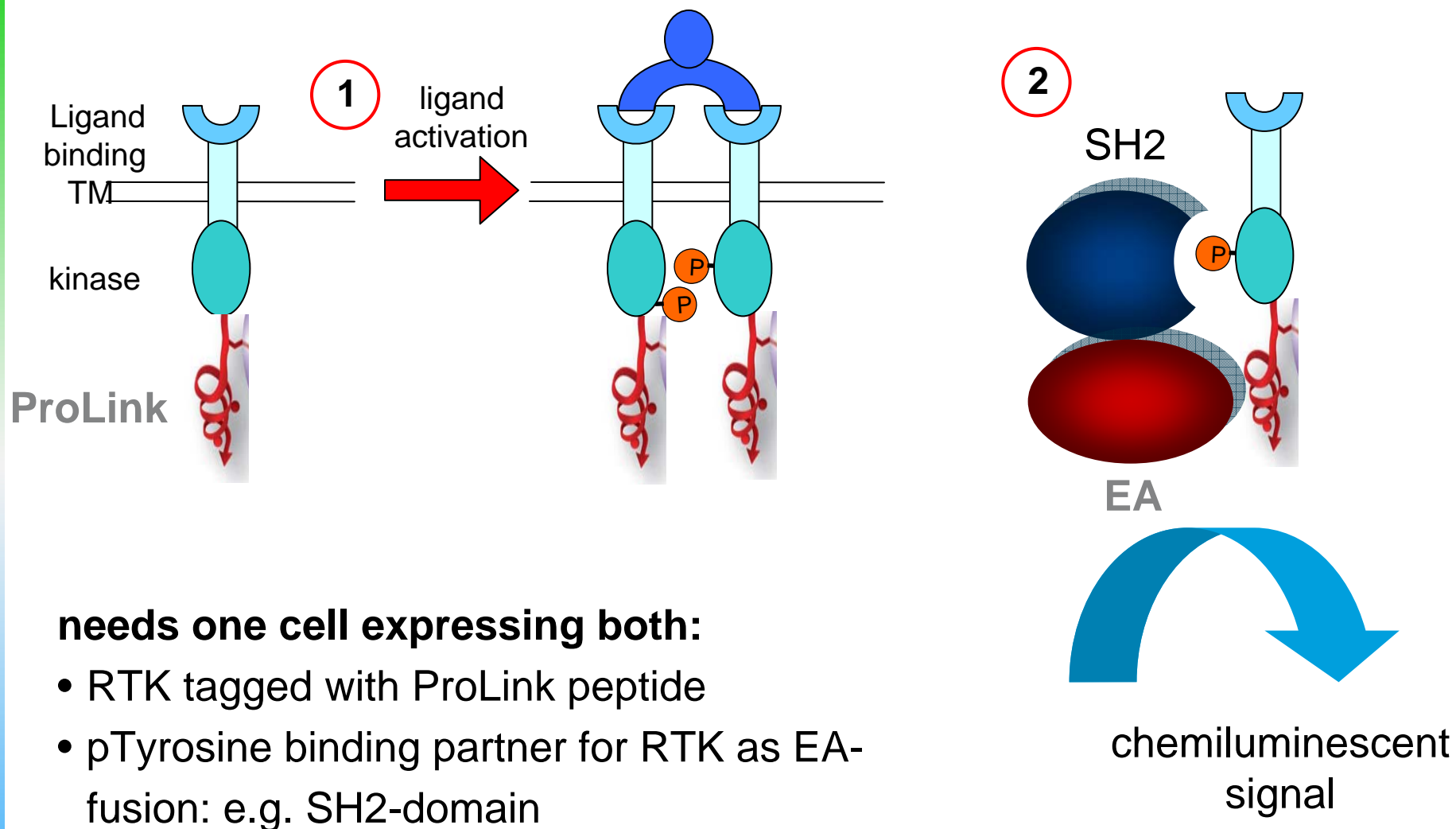
- Membranes feasible in 384 well
- but: ligand independent signal was not accepted by project team

New option: DiscoverRx



Bayer HealthCare
Bayer Schering Pharma

Concept for RTK activation assay (DiscoverRx):



needs one cell expressing both:

- RTK tagged with ProLink peptide
- pTyrosine binding partner for RTK as EA-fusion: e.g. SH2-domain



Bayer HealthCare
Bayer Schering Pharma

The DiscoverRx workplan

Aim: build a PathHunter protein-protein interaction assay for monitoring autophosphorylation of the target RTK through the interaction of an SH2 binding domain

Milestone 1

cloning of vectors
protein expression in CHO-K1



Milestone 2

retroviral co-infections of
RTK + SH2-domain (3 pairs)
2 week selection and functional test



Milestone 3

Show ligand-mediated auto-phosphorylation
of RTK in Western Blot



Milestone 4A-C

stable cell line generation of one RTK + SH2 pair:
clonal stability, functional tests, reference compounds,
DMSO-tolerance, HTS-compatibility (%CV, S/B,
cells/well), viability, Mycoplasma test



Milestone 5

reproduce data at BSP
Screening Berlin

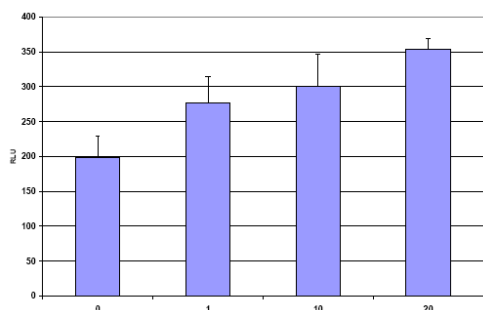


Bayer HealthCare
Bayer Schering Pharma

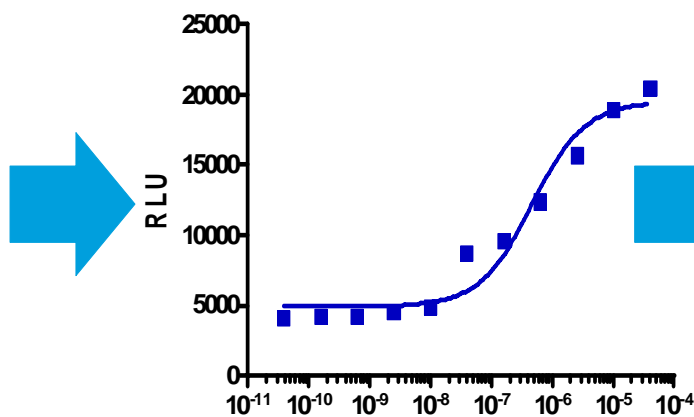
The data – Milestone 2

Aim: Find suitable RTK + SH2 pair and show functionality in the EFC-assay

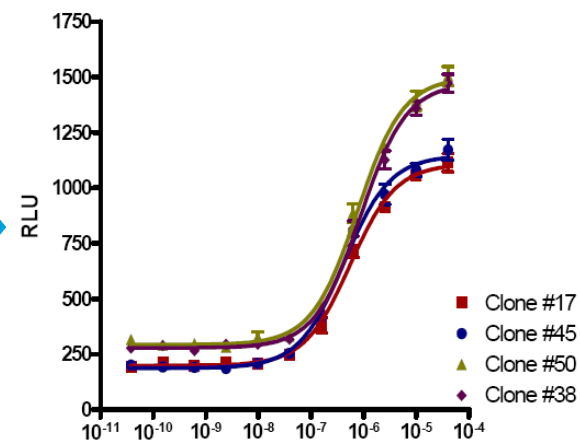
- First of all: this was the hardest part
- More than 3 different SH2-domains were tested
- Other cell backgrounds were tested



CHO RTK/Shc1
S/B = 1.7



U2OS RTK/Shc1
S/B = 4



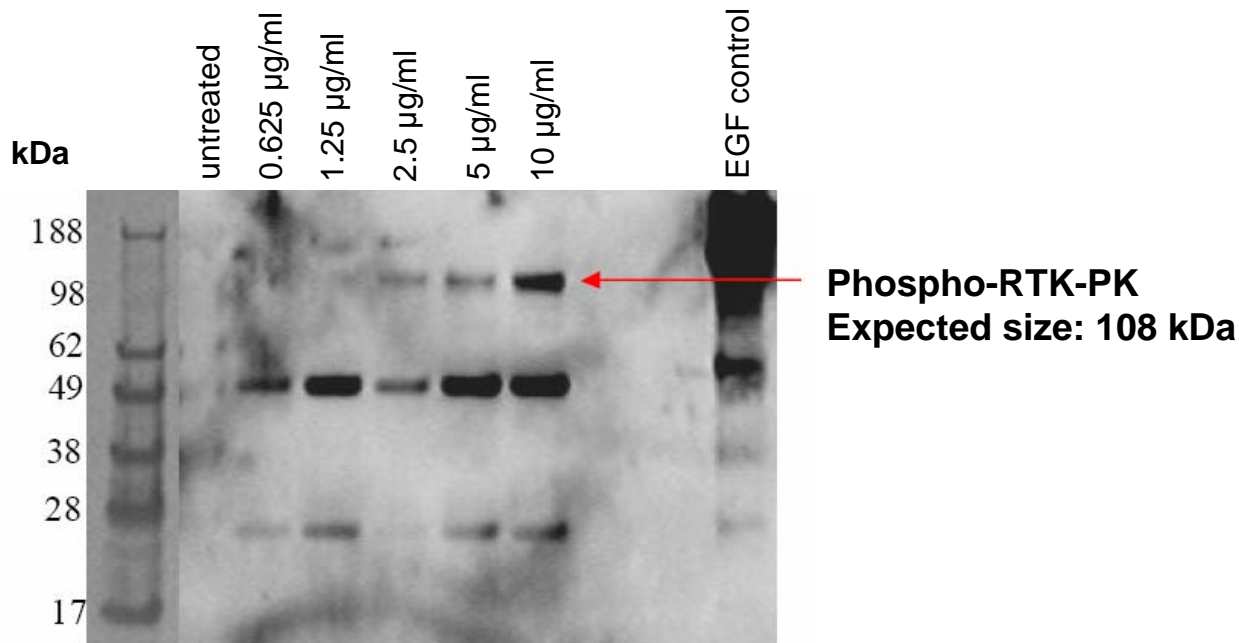
U2OS RTK/PTPN6
S/B = 5-6



Bayer HealthCare
Bayer Schering Pharma

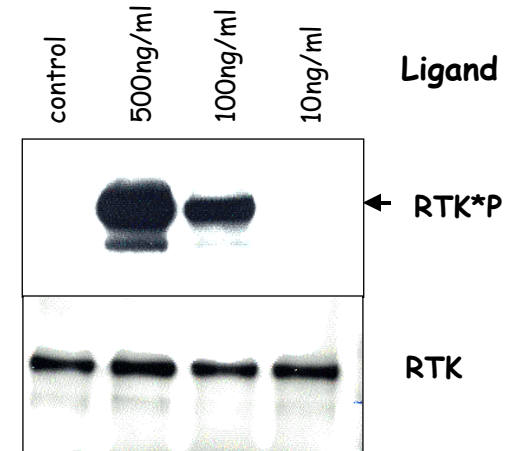
The data – Milestone 3

Aim: Show ligand-mediated auto-phosphorylation of RTK in Western Blot



IP: α -myc
IB: α -phospho-tyrosine

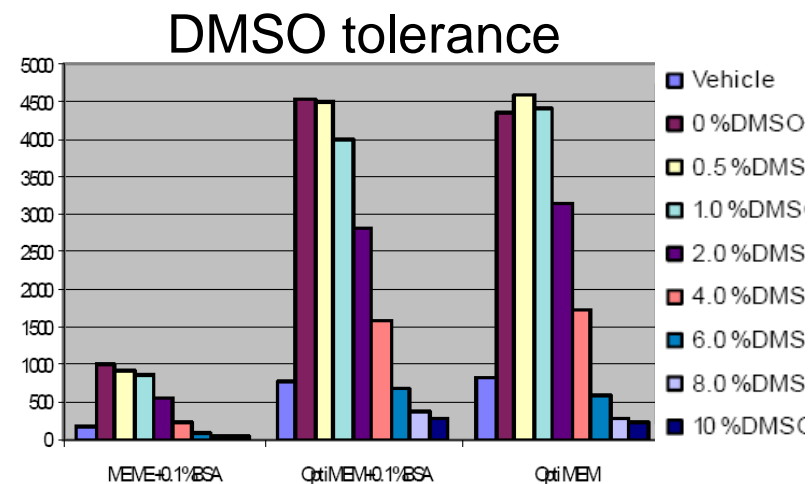
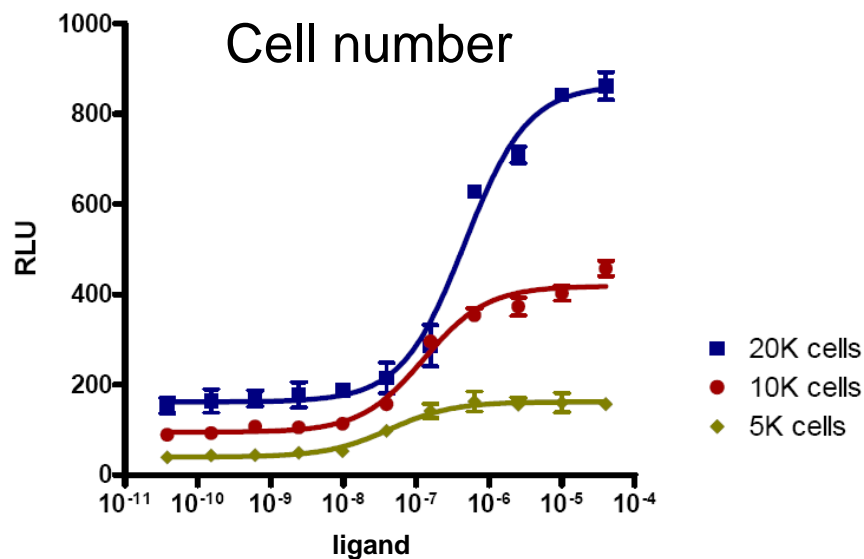
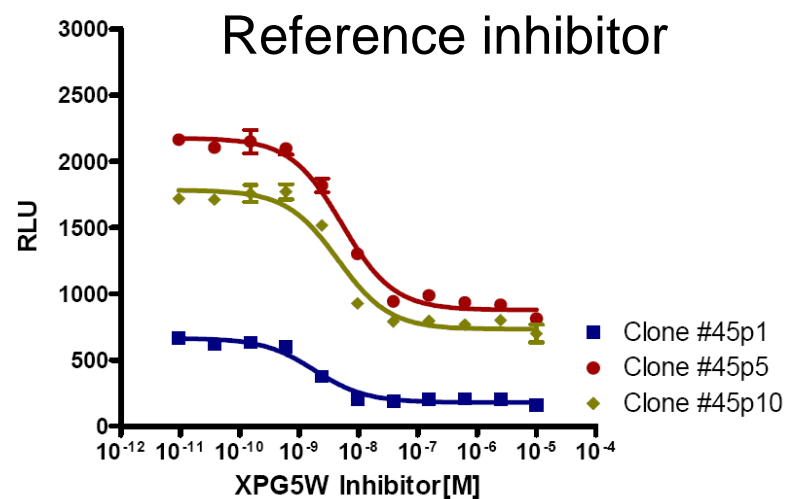
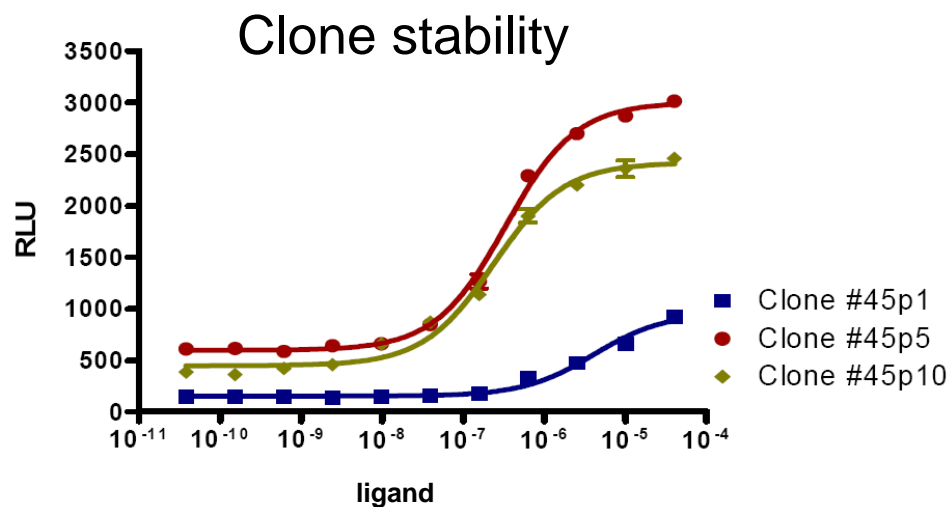
U2OS RTK/PTPN6



BSP data in CHO-cells

Data in line with expected outcome

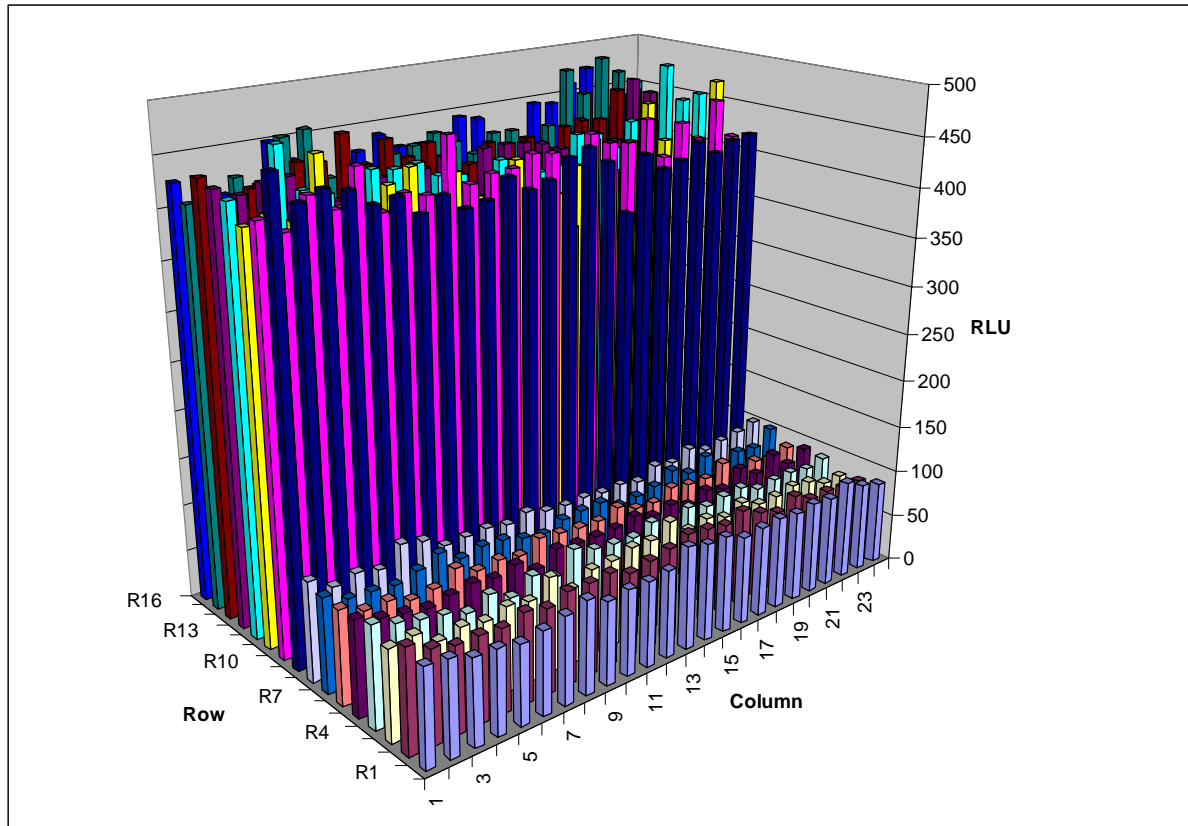
The data – Milestone 4



Bayer HealthCare
Bayer Schering Pharma

The data – Milestone 4 cont.

Aim: show Z'-factor > 0.55 for whole plate 384 well assay with 10.000 cells



	mean	SD	%CV
high	424.7	22.3	5%
low	89.7	6.5	7%

S/B	4.7
Z'F	0.74



Bayer HealthCare
Bayer Schering Pharma

Assay transfer, adaptation, optimization

1. Does the DiscoverRx protocol work?
2. Can we miniaturize ?
3. Can we use frozen cells ?



Bayer HealthCare
Bayer Schering Pharma

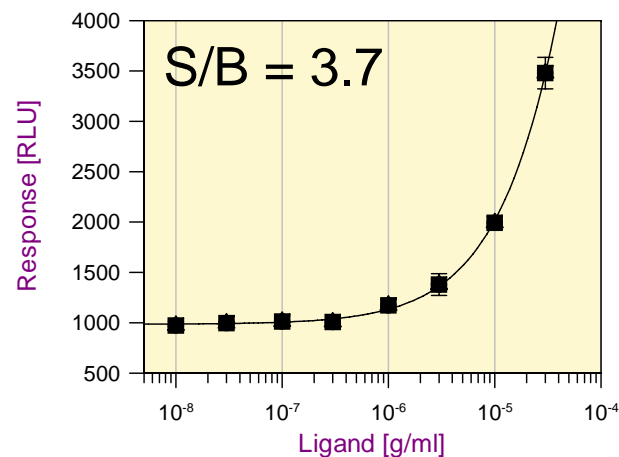
1. Does the DiscoverRx protocol work?

Steps	Volumes (384-well)
Step 1: Plate Cells & Incubate Overnight @ 37°C	Add 20 μ L of cells in each well at a preferred density of ~10,000 cells per well. Cells should be seeded in MEME +0.1% BSA.
Step 2: Treat Cells @ RT	Add 5 μ L of 5X concentrations of ligand made up in MEME (3-fold serial dilutions, highest final concentration = 40 μ g/mL). Treat for 1 hour at room temperature.
Step 3: Add CL Mix @ RT	Add 12 μ L of CL mix. Incubate at room temperature and in the dark for 60 minutes.
Step 4: Read Samples	Samples can be read on any standard luminescence plate reader.

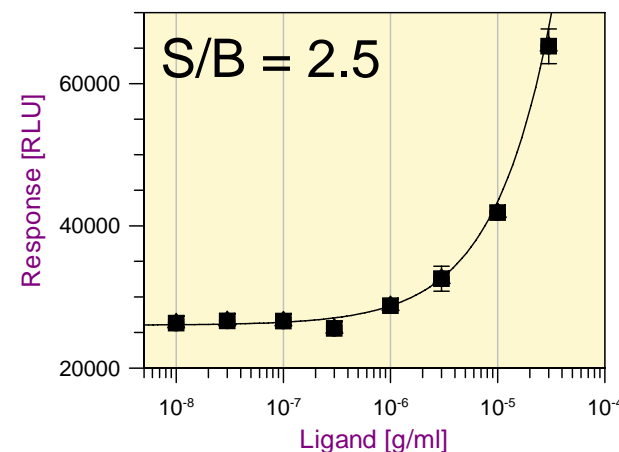
Yes, but...

...EC50 of ligand quite high, no saturation

PE ViewLux



BMG Pherastar

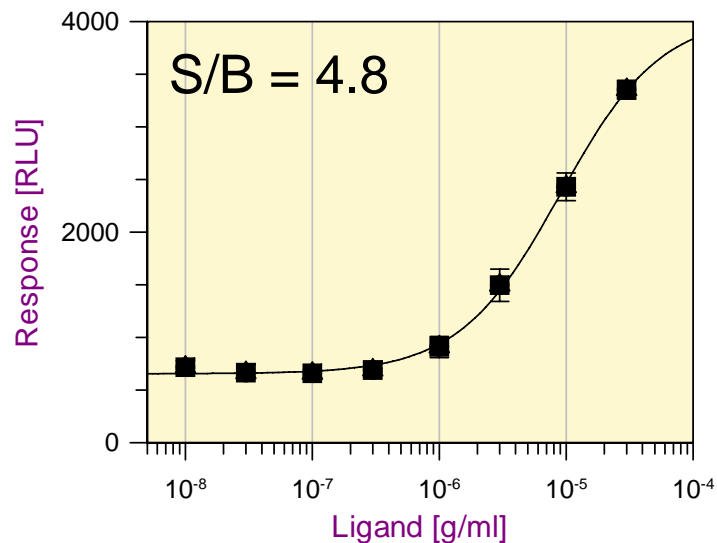


Bayer HealthCare
Bayer Schering Pharma

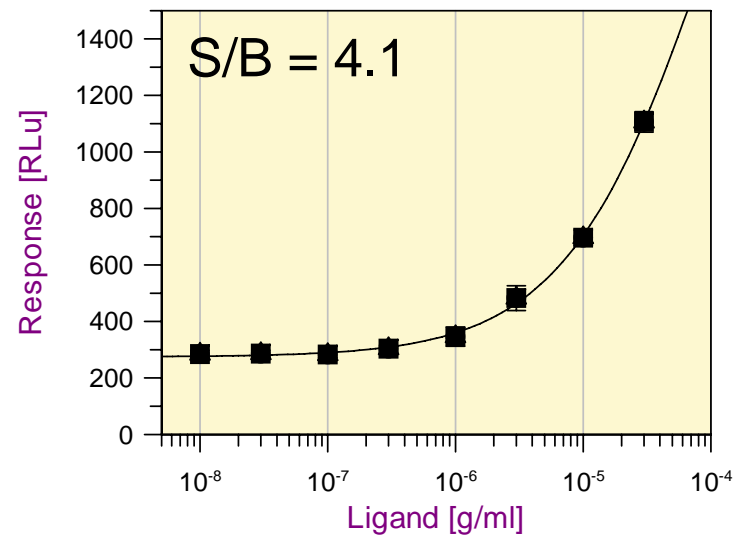
2. Can we miniaturize ?

1. 5 μ l cells (5000) into 384 SV tissue culture plates, ON 37°C
2. 2 μ l Ligand stimulation, 60 min RT
3. 3 μ l detection

Standard protocol, 10000 cells



Small volume (5000 cells)



Yes, we can...

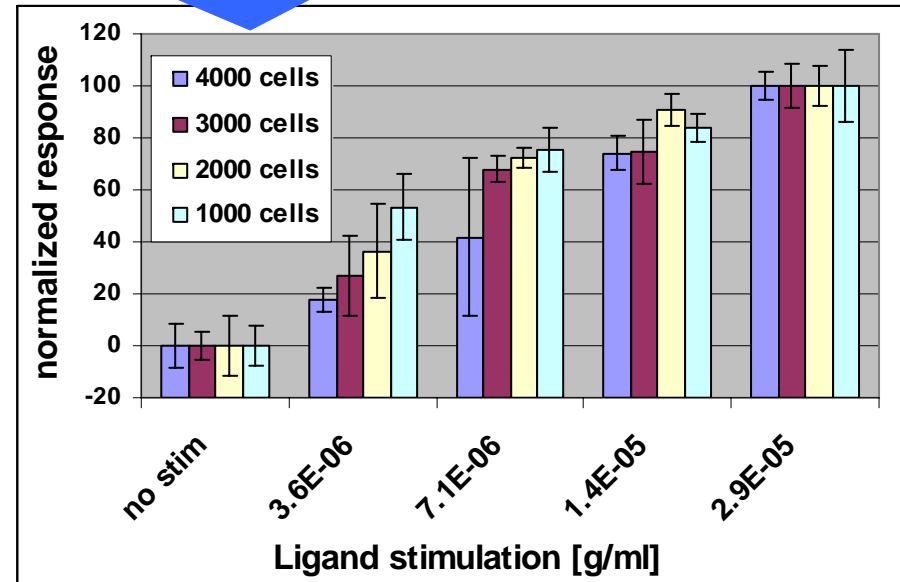
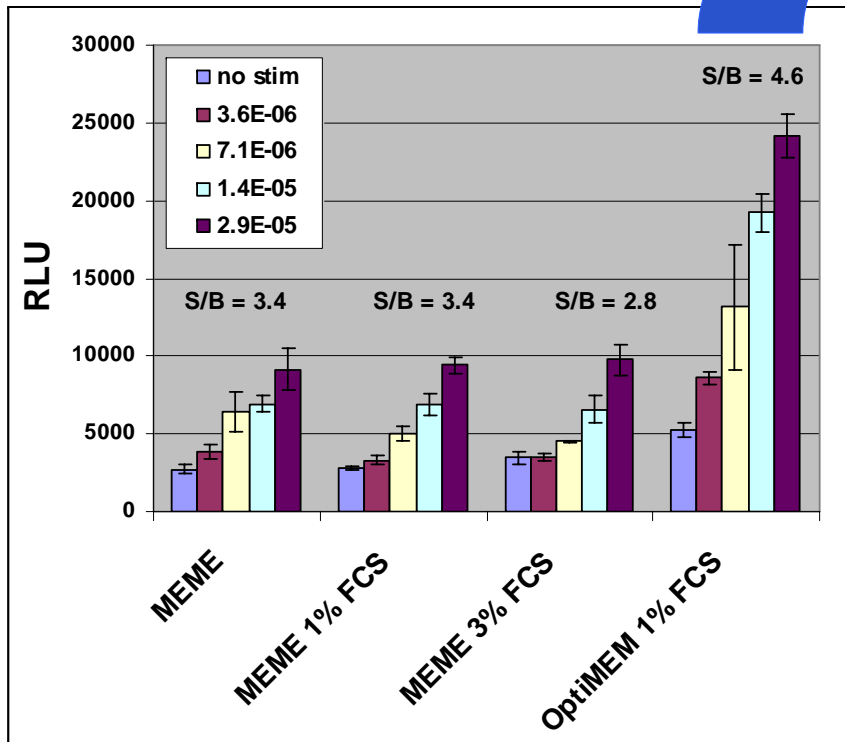
- less signal, but similar S/B
- again EC50 of ligand too high, no saturation



Bayer HealthCare
Bayer Schering Pharma

3. Can we use frozen cells ?

1. Thaw frozen cells and dilute in MEME + 0 / 1 / 3 % FCS or OptiMEM + 1% FCS
2. 5 µl cells (4000) into 384 SV tissue culture plates, ON 37°C
3. 2 µl Ligand stimulation, 60 min RT
4. 3 µl detection



cell #	4000	3000	2000	1000
S/B	4.6	4.5	3.9	3.8
Z'F	0.7	0.6	0.6	0.4
EC50 (µg/ml)	10	5.7	4.7	3.9

Yes, we can...

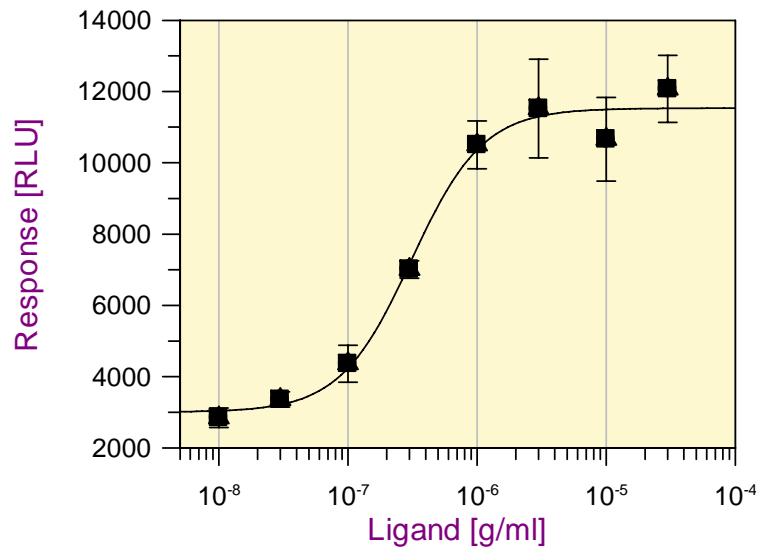
- FCS has no impact
- OptiMEM is enhancing signal and S/B

- Less cells in OptiMEM improves EC50

4. Optimized assay using OptiMEM

➔ OptiMEM without additives optimal for this assay

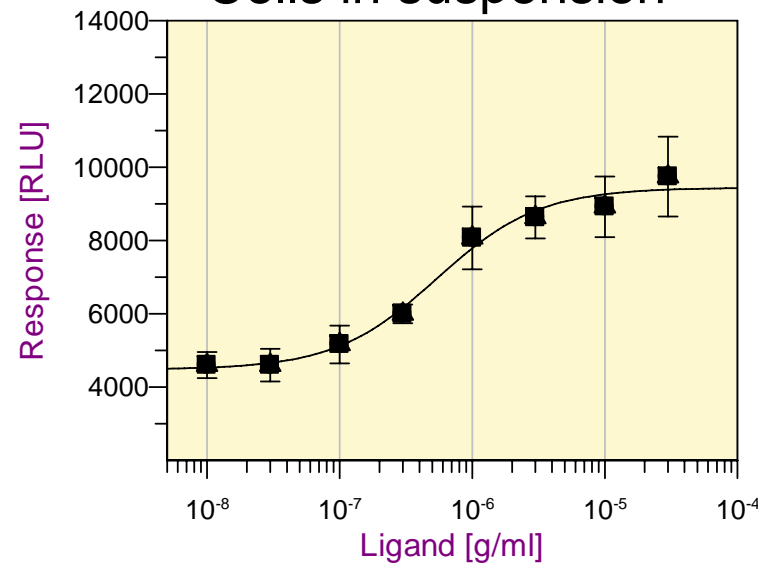
Cells adherent



S/B = 4.2

EC50 = 0.3 µg/ml

Cells in suspension



S/B = 2.2

EC50 = 0.5 µg/ml

- EC50 as expected
- Adherent cells with superior S/B

Summary and Outlook

- A cell-based, ligand inducible RTK-SH2 interaction assay was developed by DiscoverRx in an assay development project with Bayer Schering Pharma AG.
- The assay meets the requirements in terms of S/B, Z'-factor and IC50s of reference inhibitors
- The assay could be successfully miniaturized and applied to frozen cells
- Further work will aim at testing robustness in automation and 1536 well format

Acknowledgements

DiscoverX

Edwin Haghazari
Mimi Nguyen
Martine Ercken

Birgit Beyermann-Lopata
Sanj Kumar
Keith Olson

Bayer Schering Pharma

Dagmar Zeggert-Springer

Arndt Schmitz
Dieter Zopf
Verena Voehringer
Dirk Brohm
Heike Petrul
Sandra Bruder
Volkhart Li

Karsten Parczyk
Anke Müller-Fahrnow



Bayer HealthCare
Bayer Schering Pharma