

Science For A Better Life

Ultra HTS @ Bayer:

Use of IP-One and Tag-Lite assays in GPCR drug discovery

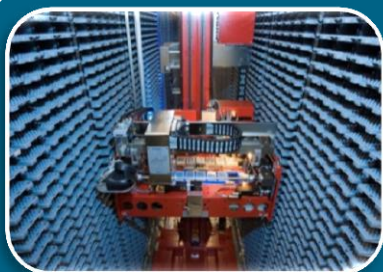
25th of April 2013

Katrin Nowak-Reppel, PhD, Lab Head

Agenda

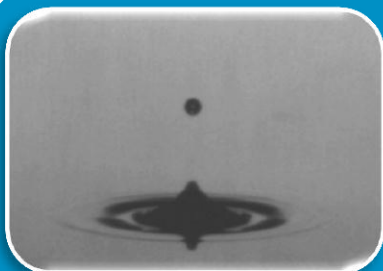
- Overview uHTS @ Bayer/Lead Discovery Berlin
- Highlights of cell-based screening
- 1536MTP format IP-One assay: development and screening
- Development of Tag-Lite binding assay in 1536MTP format
- Summary

Ultra HTS @ Bayer/Lead Discovery Berlin



Whole library screening

- ~3 million compounds for maximum of diversity
- Standard format is 1536MTP



www.labcyte.com

„Ready-to-assay“ plates

- Acoustic pre-dispersion of compounds in assay plates
- Flexible use for different assay days



Fully automated & Benchtop uHTS

- Addition-only homogenous assays
- Non-homogenous and kinetic assays (e.g. High Content Imaging assays)



Assay technology preferences for uHTS @ Bayer/Lead Discovery Berlin



Homogenous addition-only assay technologies

Endpoint assays for higher flexibility in time

Application for several target classes of different indications

Miniaturization potential to fit to 1536MTP format and reduce costs

Potential to use frozen cells and perform assay in suspension

➔ HTRF assay is a preferred assay technology



Cell-based uHTS @Bayer/LD Berlin

Use of frozen cells

- Reduce inter assay variance → Quality
- Thaw only these cells required for daily screening run → Flexibility
- Scale experiment corresponding to requirements → Efficiency

Evaluation of cell lines

- Physiological relevant cell lines (e.g. oncology – tumor type addressed)
- Endogenous expression of targets
- Avoid artificial expression of target

Phenotypical screening – High content analysis screening

- Whole pathway can be addressed
- Multiple parameter accessed at once
- High physiological relevance

GPCR screening project background

Background:

- Gq protein coupled receptor
- Member of β -subgroup of rhodopsin receptor subfamily
- Target in family with 2 relative receptors

Project goal:

Identification of specific antagonists interfering with agonist-induced GPCR activation



Development of an uHTS compatible IP-One Assay

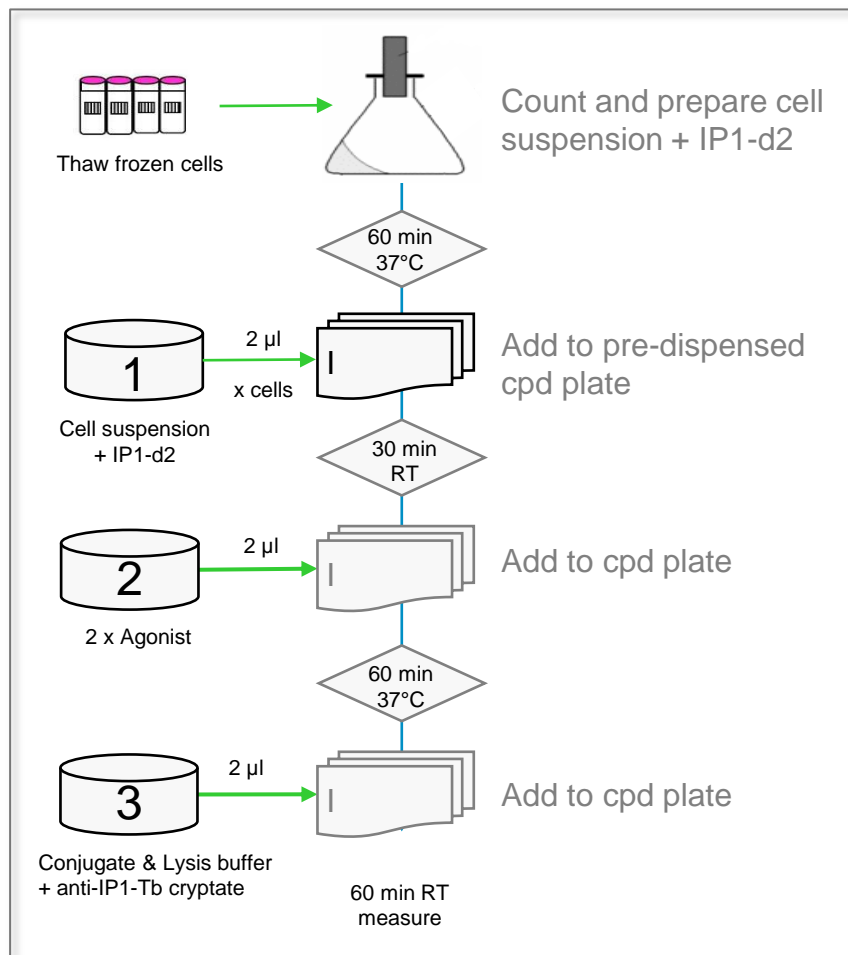


Start of assay development with standard IP-One HTRF Assay protocol.

Parameters tested during AD:

- Different cell lines and ligands
- Cell number/well
- Incubation time cells with compounds before ligand addition
- Incubation time and temperature of cells, compounds and agonist
- Volumina of addition steps
- Plate colour
- Dilution of IP1-d2 and anti-IP1-Tb cryptate
- ...

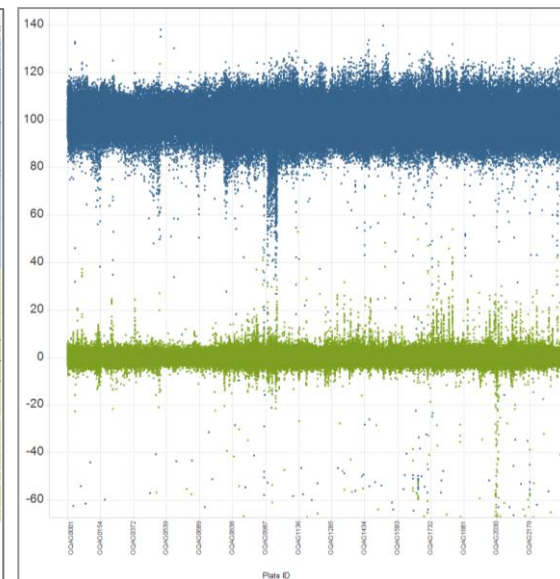
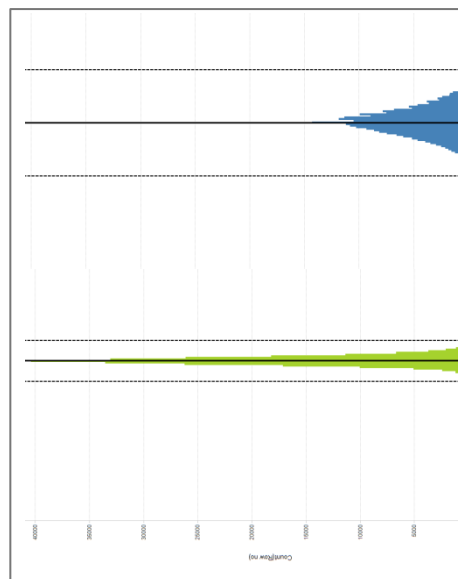
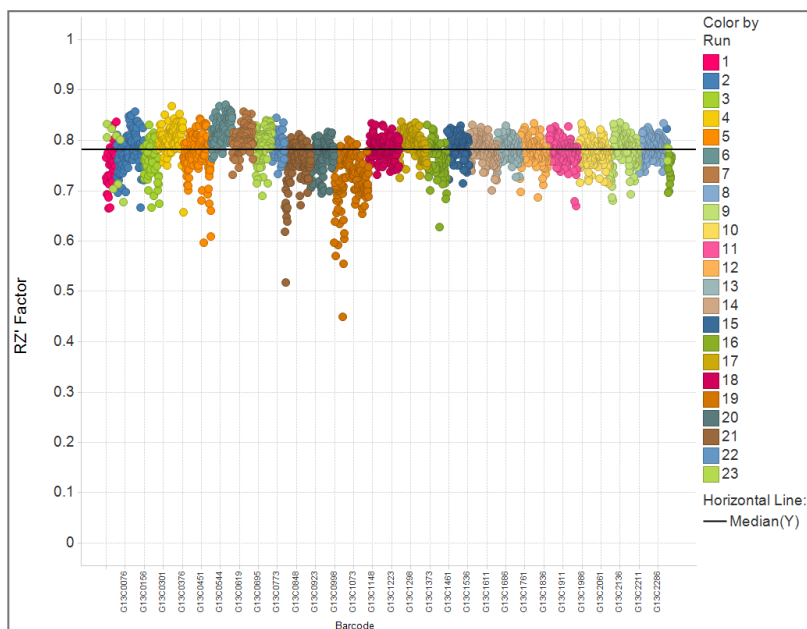
Setup of uHTS compatible IP-One assay



- Assay volume: 4µl
 - Detection volume: 6µl
 - Final compound concentration: 10µM
- Flexibility by use of frozen cells
 - Accuracy by use of ready-to-assay plates
 - Efficiency by reduction to 4µl volume and only three addition steps



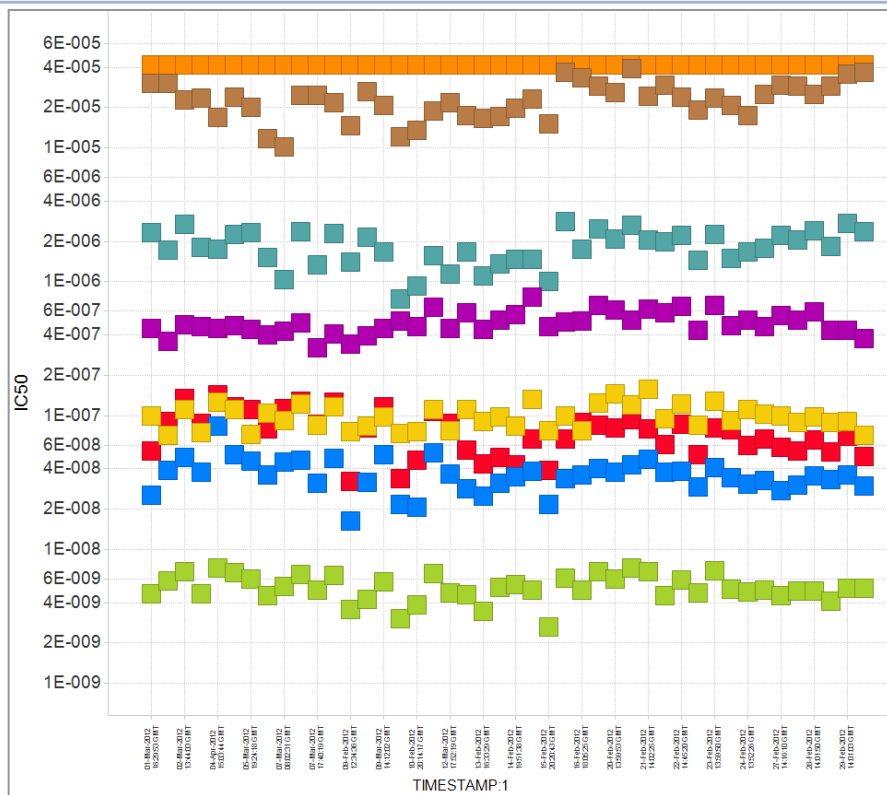
IP-One screen proves excellent assay quality



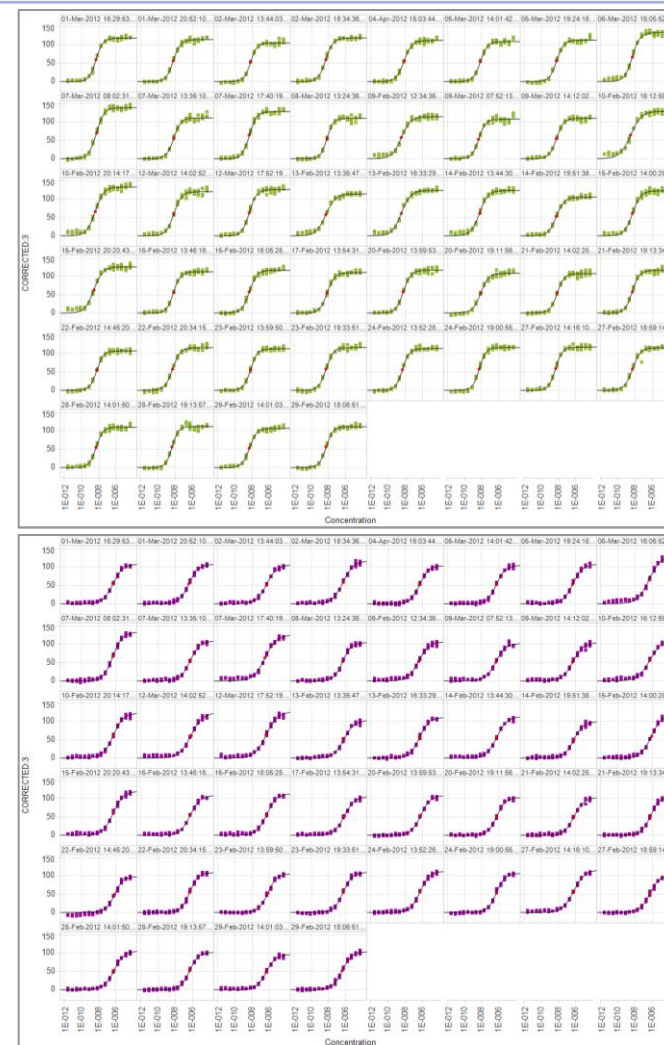
High robust Z' factors in every run

Broad separation of controls

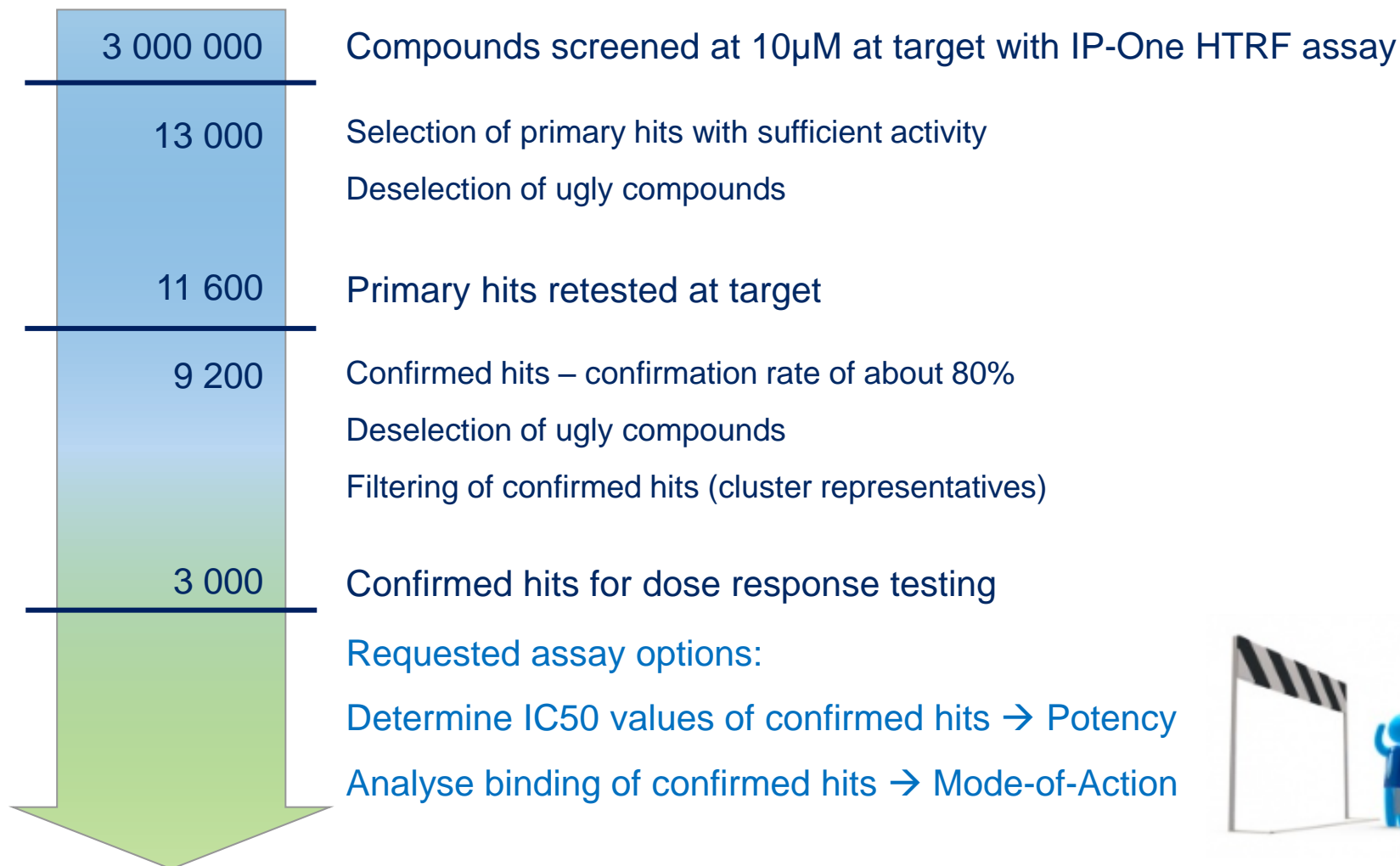
IP-One screen proves excellent assay quality



Constant IC50 values of reference compounds over whole screening



Screening cascade



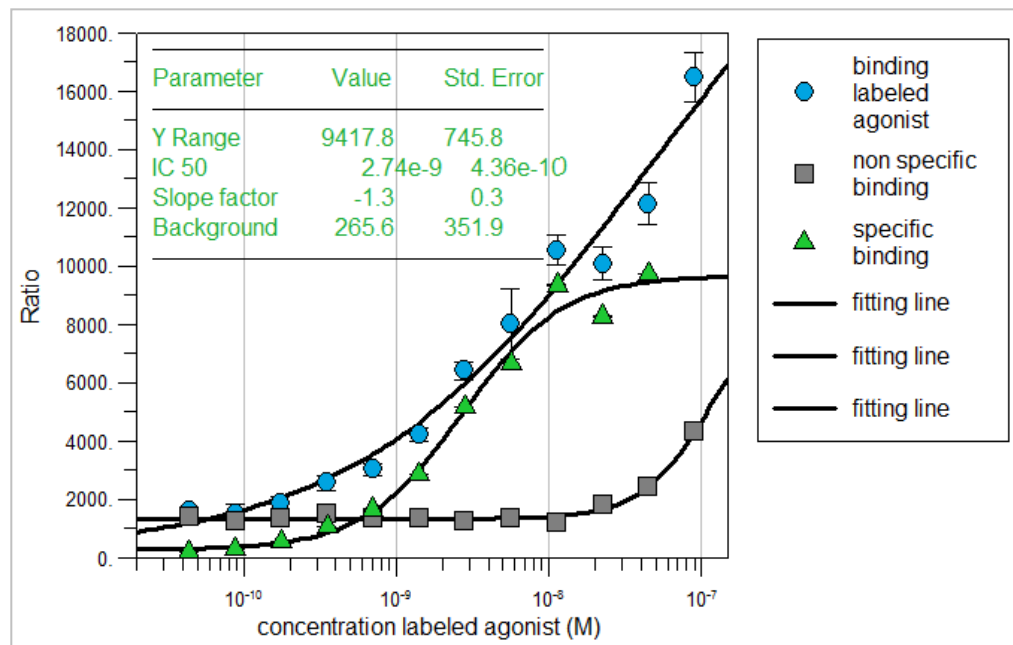
Transfer and establishment of available Tag-Lite binding assay



Starting point: Tag-Lite assay for target available as kit

(384MTP format, 20µl volume, scheduled for medium throughput)

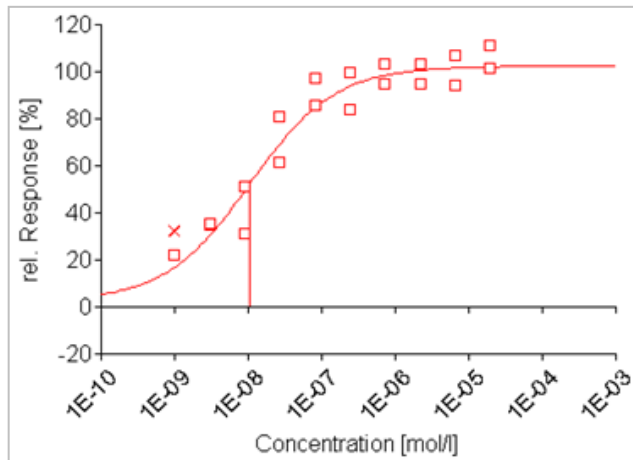
First inhouse experiment:



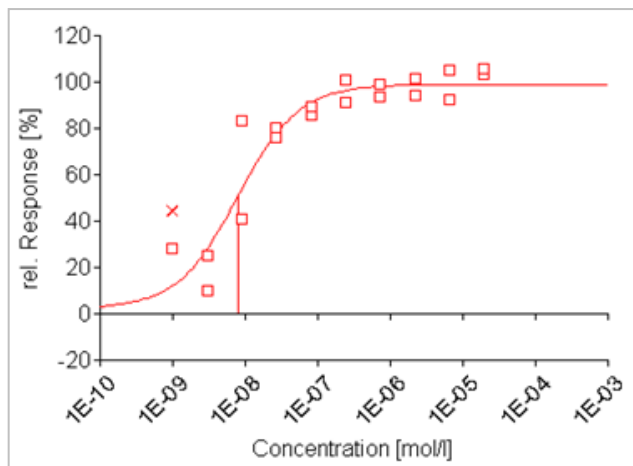
- 384MTP format
- Assay volume: 20µl
- Frozen cells
- Compound transfer

- Assay transferred successfully
- IC50 of labeled agonist confirmed

Tag-Lite competition experiment with pre-dispensed compounds

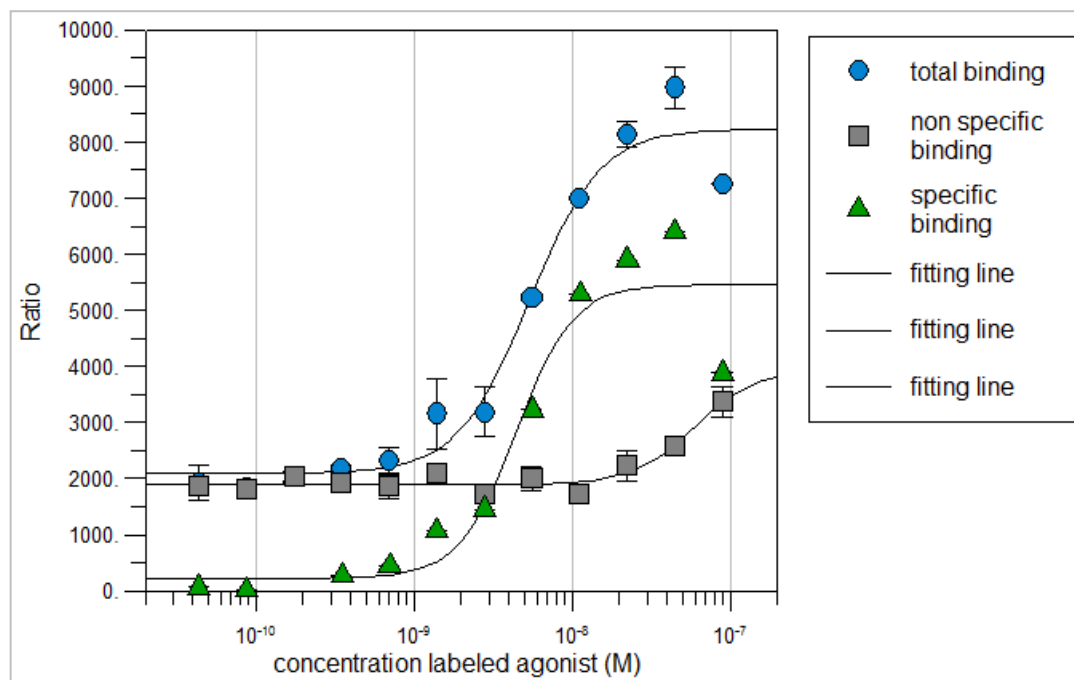


- 384MTP format
- Assay volume: 20µl
- Frozen cells
- 100nl compounds pre-dispensed
- Transfer of labeled agonist



- Excellent quality of competition experiment
- Use of pre-dispensed plates successful
- Data of orthosteric antagonists known from literature confirmed

Volume reduction of Tag-Lite assay



- 384MTP format
- Assay volume: 8 μ l
- Frozen cells

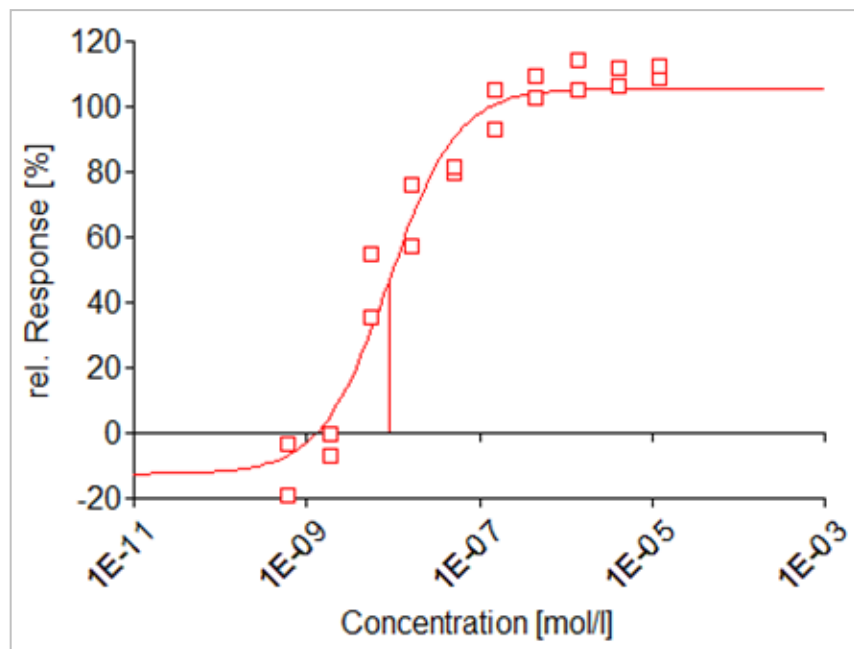
- Binding curve of labeled agonist reproduced in lower volume
- Reduction of assay volume successfully
- Assay performance in 1536MTP should be possible

Miniaturization of Tag-Lite assay

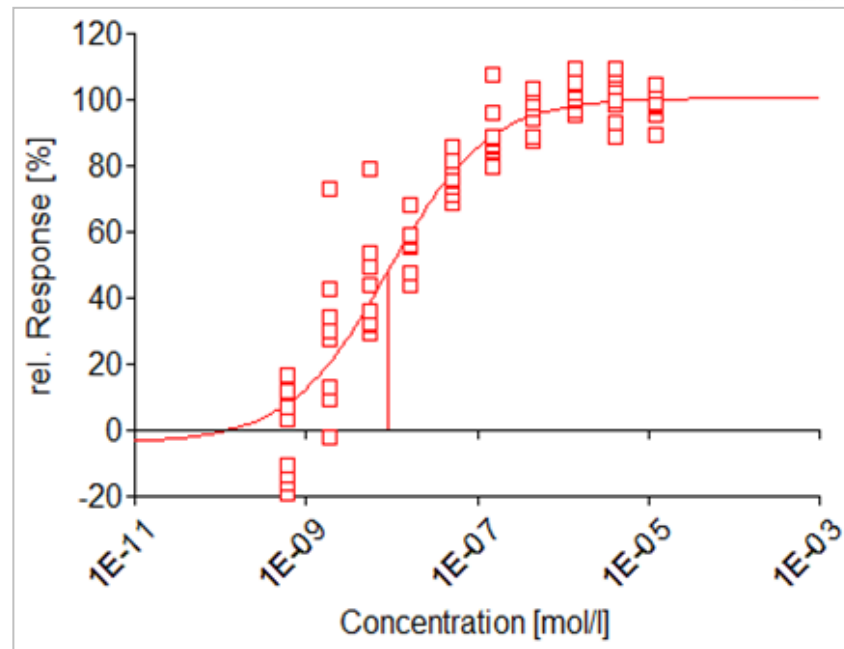
Competition dose response curve of known antagonist in



384MTP



1536MTP



➤ Transfer of miniaturized Tag-Lite assay to 1536MTP format successful

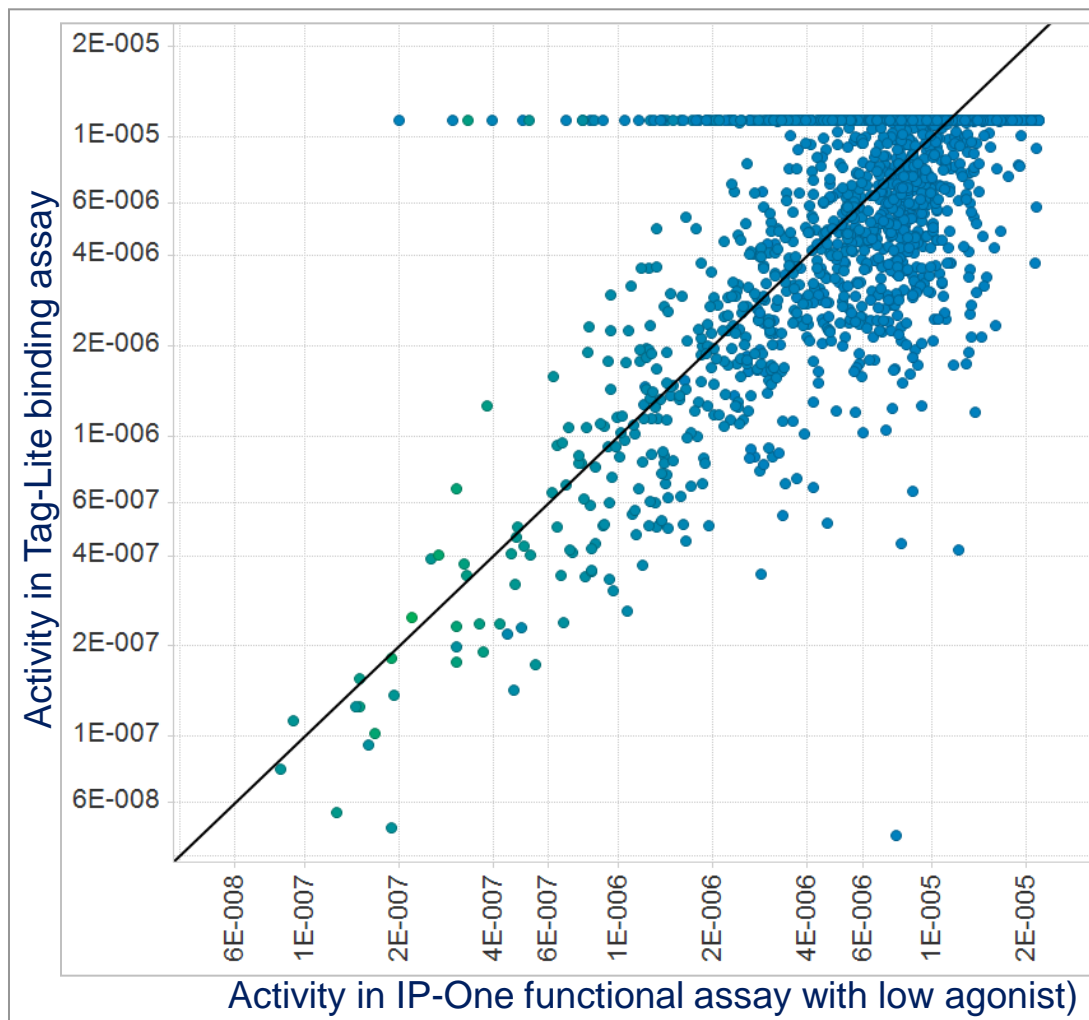
Automatization of miniaturized 1536MTP format Tag-Lite assay



Additional parameters tested and optimized:

- Automatized cell dispensation on compounds
- Dispensation of solution with labeled agonist
- Incubation before measurement
- Signal stability over time
- Screen batches of cells and labeled agonist

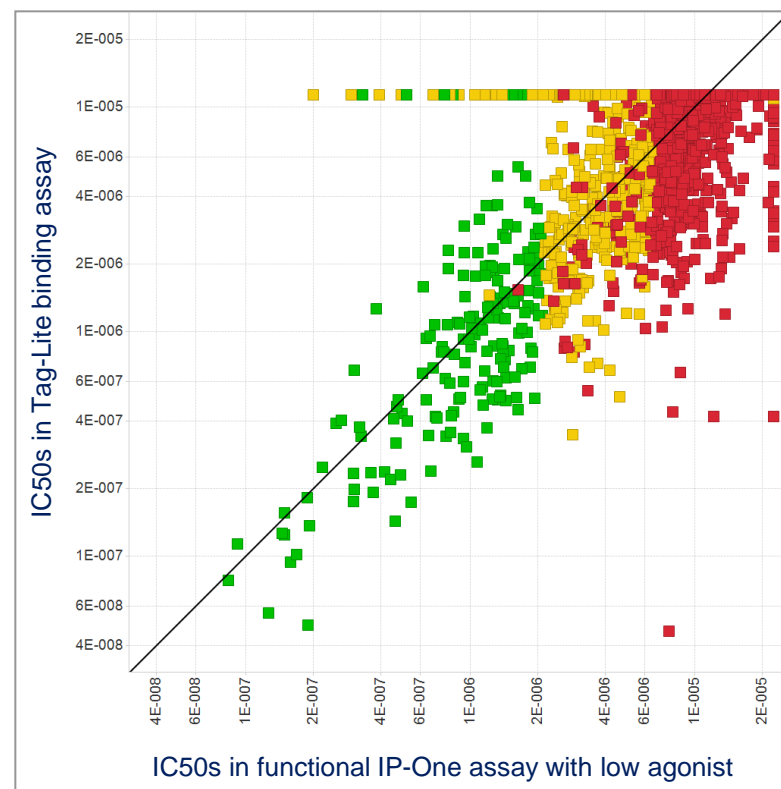
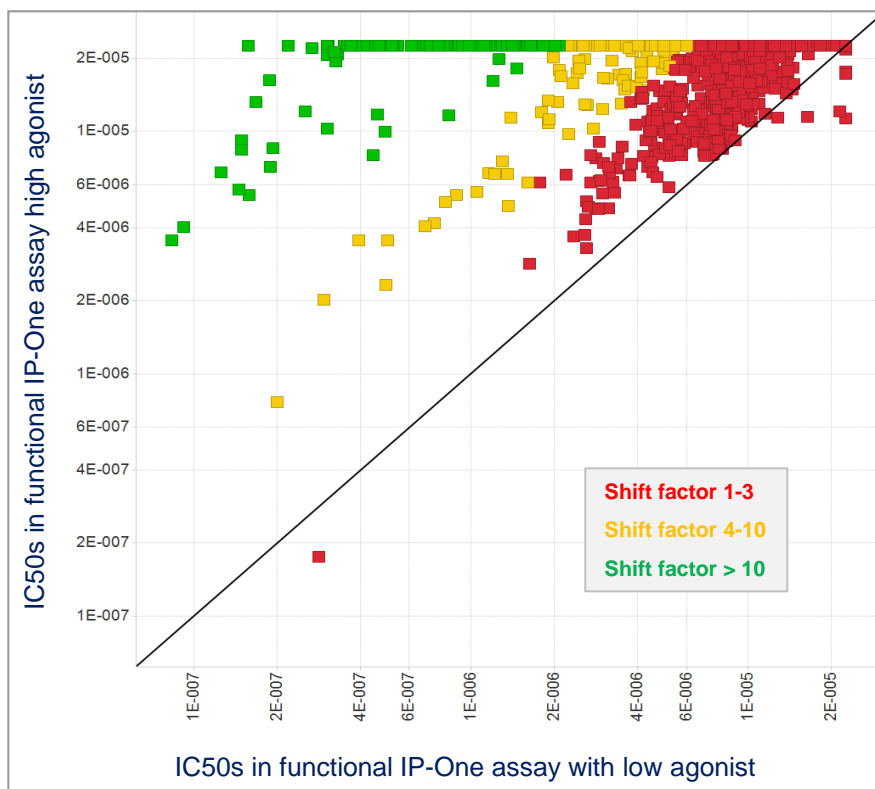
Correlation of data in binding and functional cell-based HTRF assays



- Dose response curves of ~3000 compounds
- IP-One functional assay
- Tag-Lite binding assay
- 1536MTP format
- Frozen cells

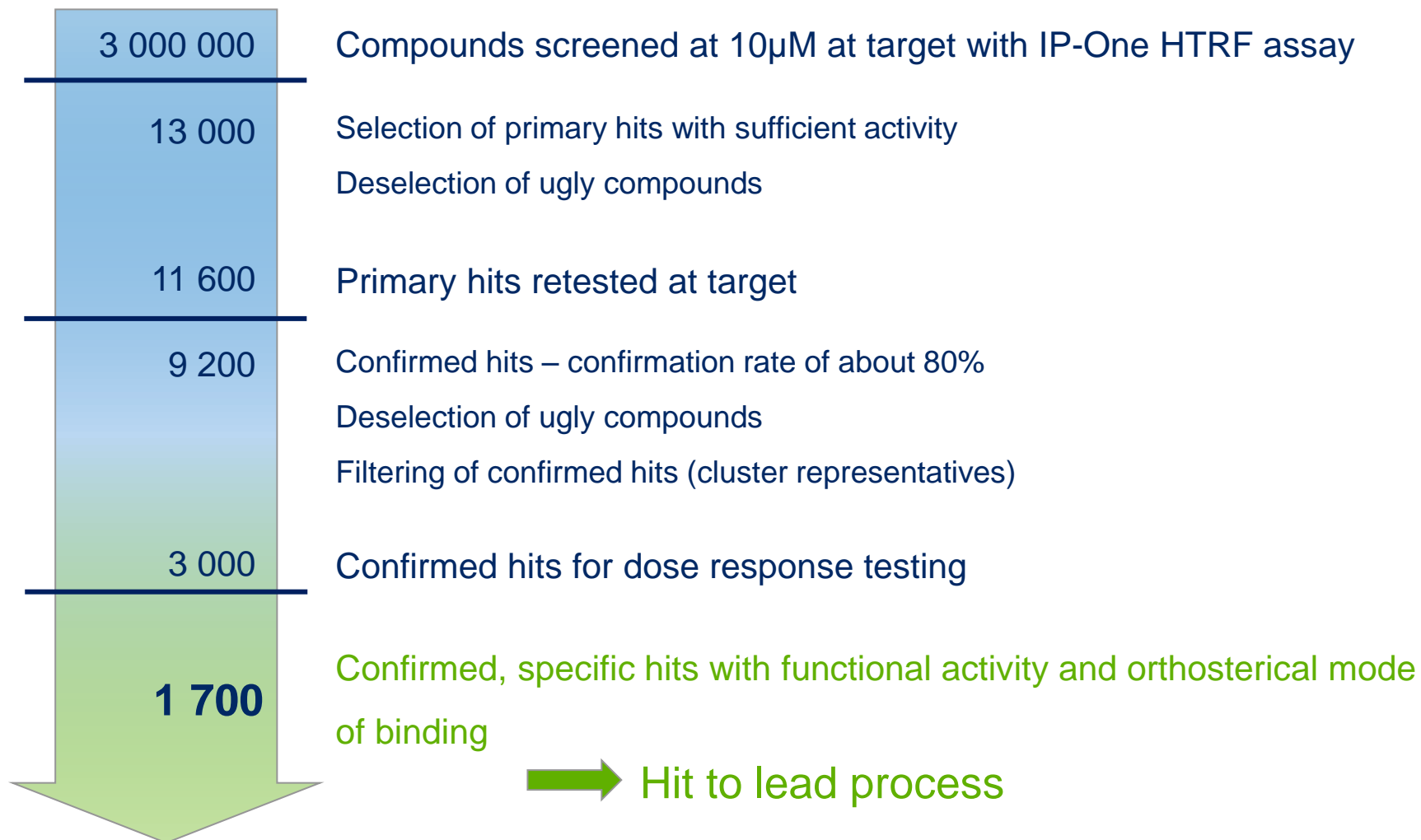
➤ Good correlation of functional and binding data

Comparison of results with different Mode-of-Action assay settings



- Compounds which could be shifted with increase of agonist display potent binding activity in Tag-Lite assay

Screening cascade





Summary

An IP-One HTRF assay in 1536MTP format and by using frozen cells was developed.

The uHTS campaign with this IP-One assay yielded in excellent quality and promising hits.

To address binding an available Tag-Lite assay was further optimized and miniaturized for performance in 1536MTP format.

Dose response curves of ~3000 compounds were performed in this Tag-Lite assay and displayed good correlation to functional data.

IP-One and Tag-Lite HTRF assays are suitable technologies for high-throughput cell-based assays and amenable for further automatization.

Acknowledgements

Jessica Schulze

Sabrina Reimann

Benjamin Bader

Karsten Parczyk



Marielle Mazille, Cisbio Bioassays

Reiner Hein, Cisbio Bioassays



Science For A Better Life

Thank you