

Exploiting HTRF for novel drug classes: Stabilizing 14-3-3 protein-protein interactions

The Lead Discovery Center (LDC)



- The Lead Discovery Center (LDC) was established in 2008 by the technology transfer organization of the Max-Planck Society, as a novel approach to capitalize on the potential of basic research for the discovery of new therapies for diseases with high medical need.
- As an independent company with an entrepreneurial outlook, the LDC closely collaborates with research institutions, universities and industry. Our aim is to transform promising early-stage projects into innovative pharmaceutical leads that reach initial proof-of-concept in animals.



MAX-PLANCK-GESELLSCHAFT

LDC's drug discovery network



MAX-PLANCK-GESELLSCHAFT

- 80 institutes
- 16.900 employees < 6.600 scientists
- >12.000 publications p.a.; 32 nobel laureates
- additional 7.700 young & guest scientists
- ~40 institutes with life science (biomedical) oriented research programs
- €1.73 Bio. annual research budget
- central tech transfer unit:

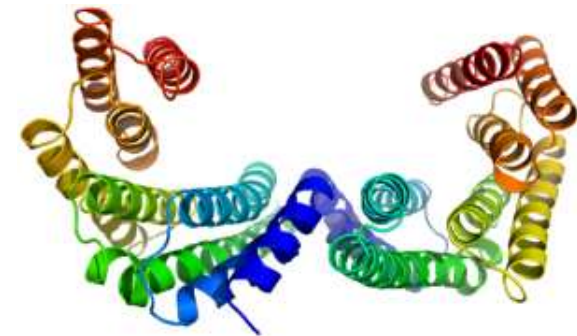


Max-Planck-Innovation

3.200 inventions
1.900 contracts
90 spin-offs

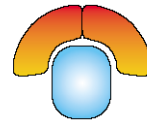
- ★ Max-Planck Institute
- ★ Academic Collaboration Partner
- ★ Pharma/Biotech Network

- ubiquitous in eucaryotes
- 7 isoforms in humans and higher plants
- highly conserved primary sequence
- physiological activity is mediated by direct protein-protein interactions
- key-regulators of signal transduction, cell cycle control, apoptosis, primary metabolism
- more than 500 interacting proteins described

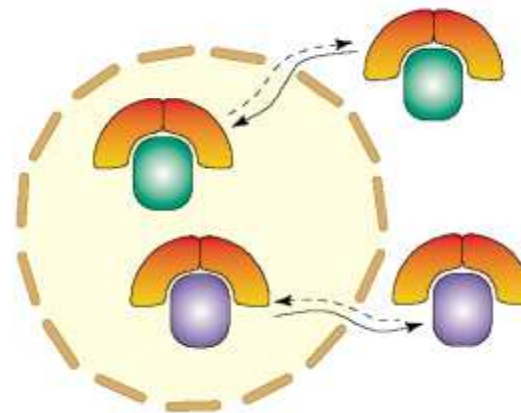


Modulation of the target activity by 14-3-3

1) enzymatic activity



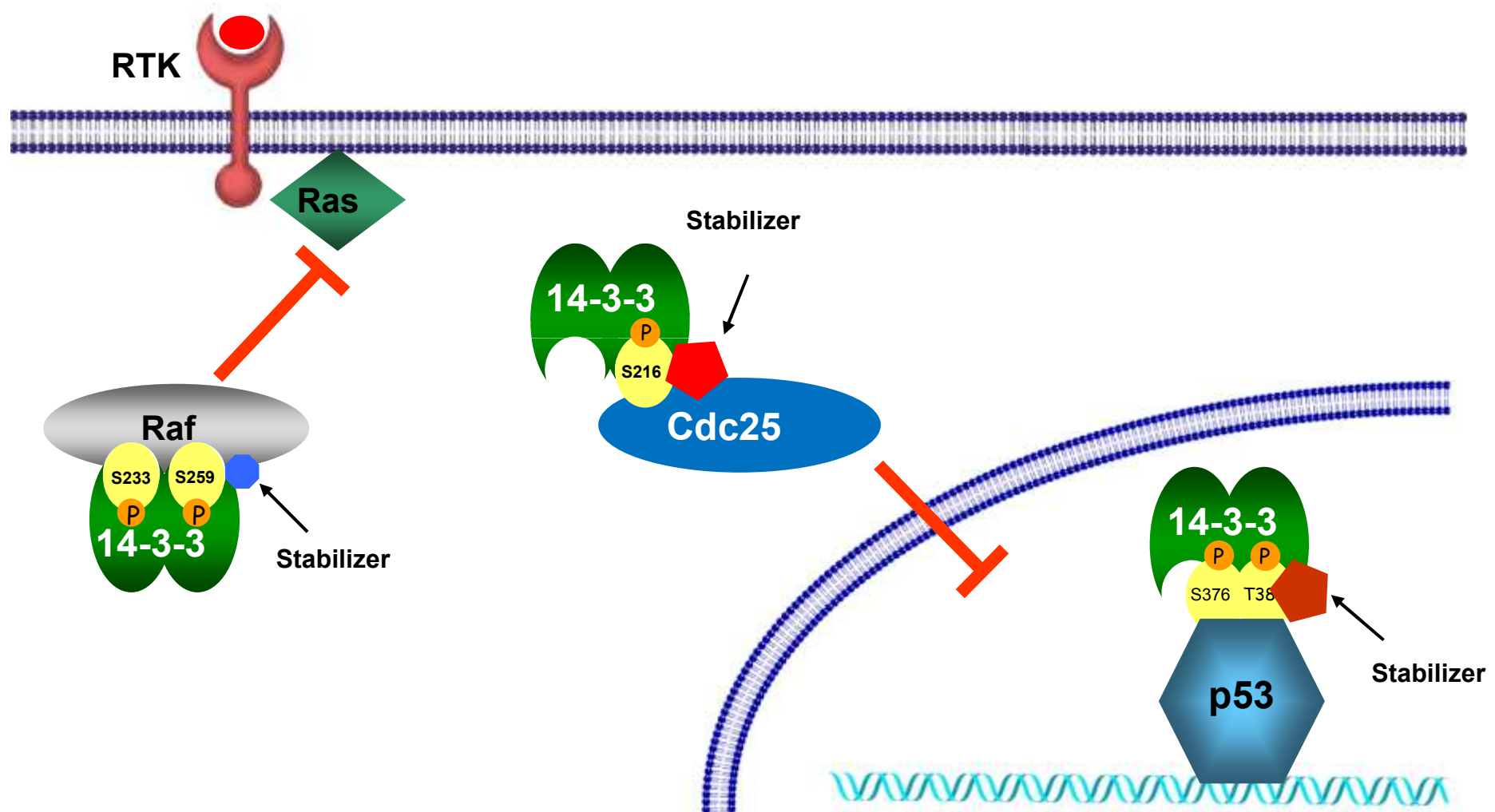
2) subcellular localization



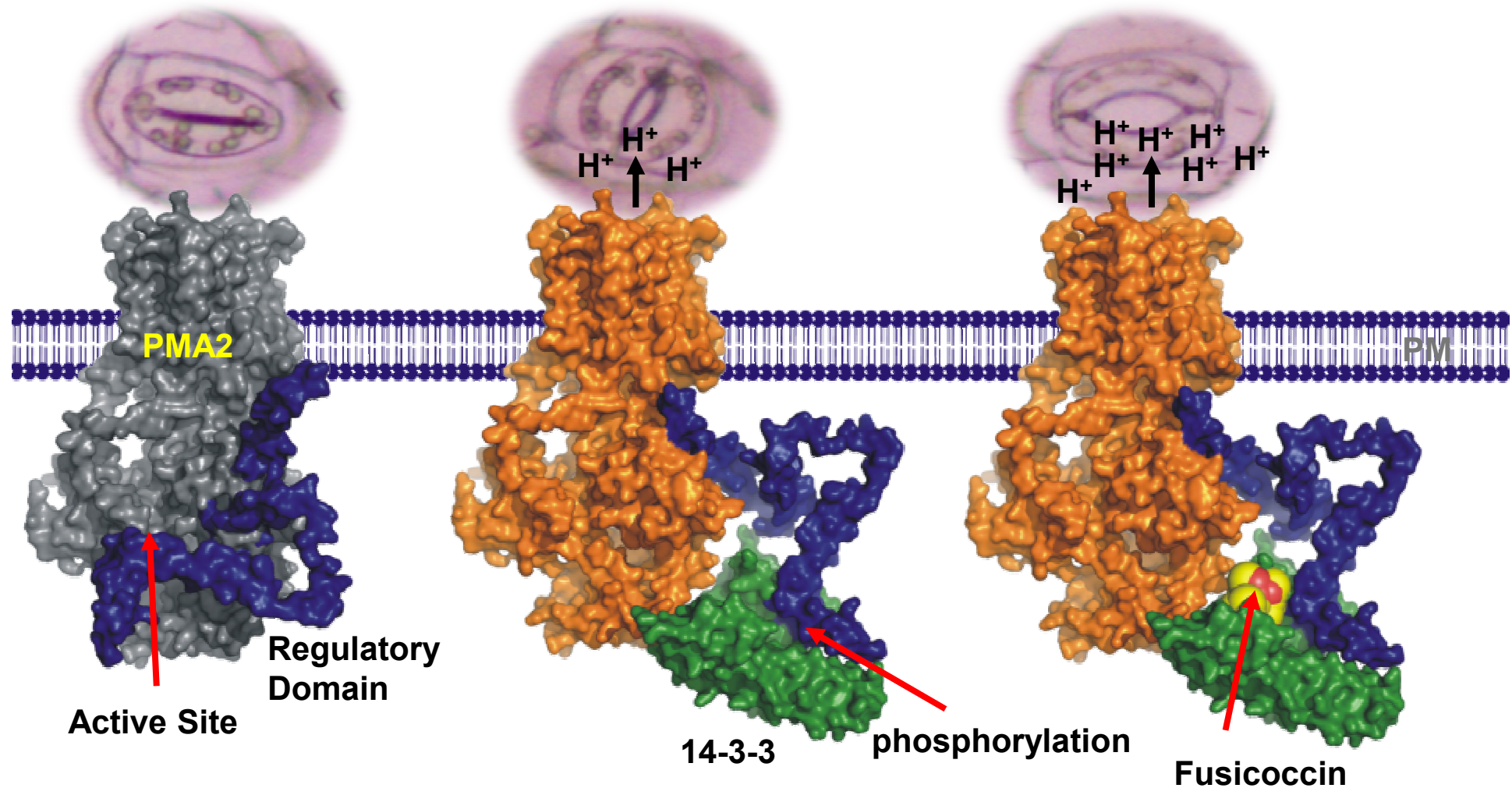
3) ability to interact with further protein partners



Why stabilize the 14-3-3 interaction



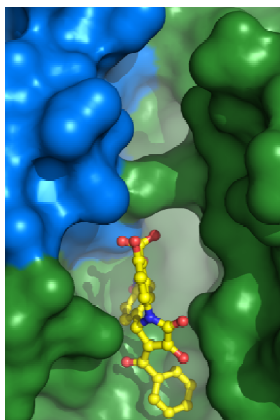
Nature's example: Activation of the H⁺-ATPase PMA2 by Fusicoccin



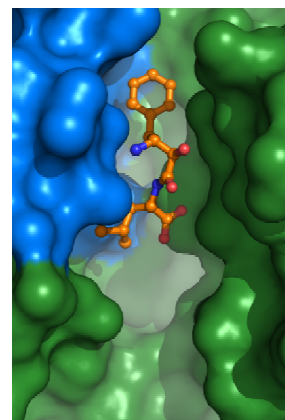
Structures courtesy by Christian Ottmann

A feasibility study

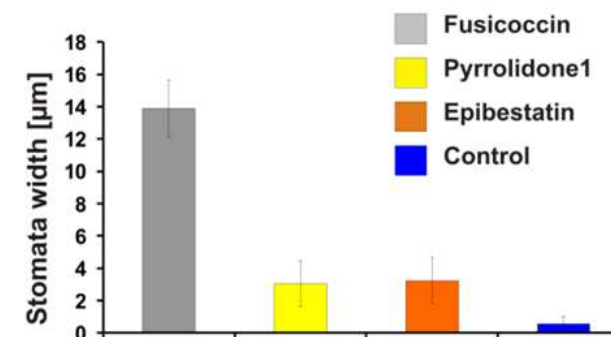
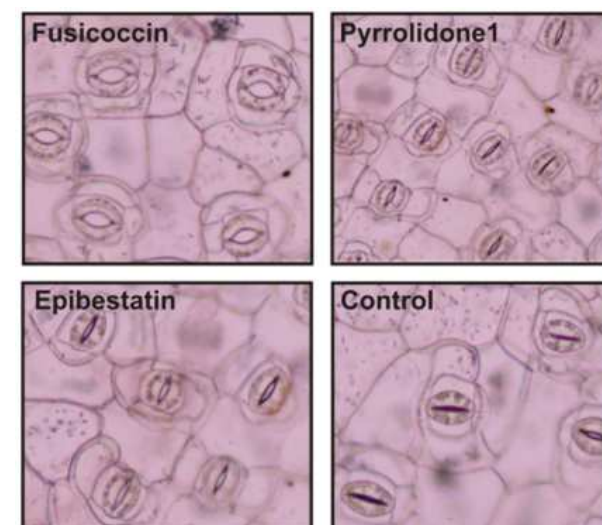
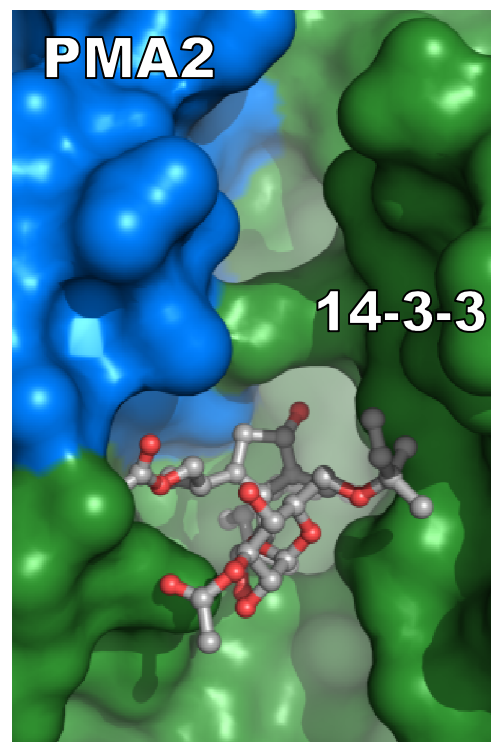
Pyrrolidone1



Epibestatin



Fusicoccin



Rose, R. et al: Angewandte 2010, 386, 913-919.

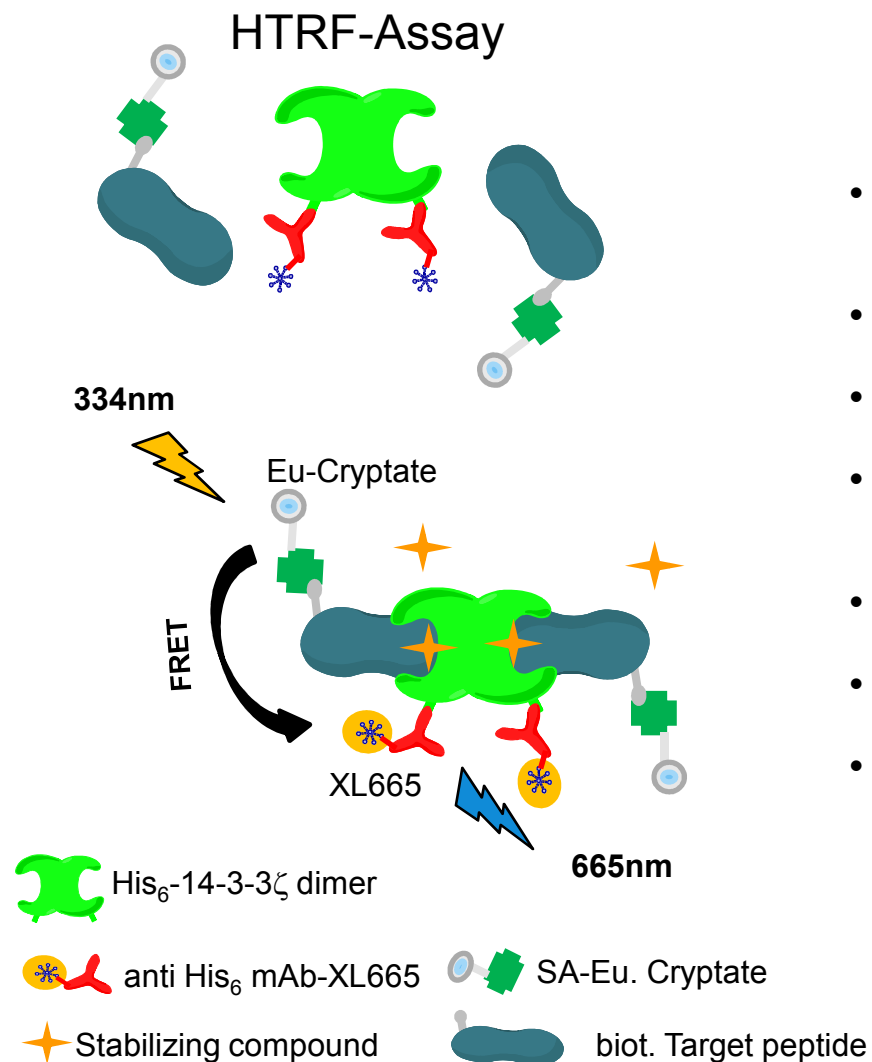
Prerequisites

- Homogeneous („mix and read“)
- Sensitivity
- Miniaturizable to 384well and 1536well format
- Compatible to Screening-Hardware

Common assay formats

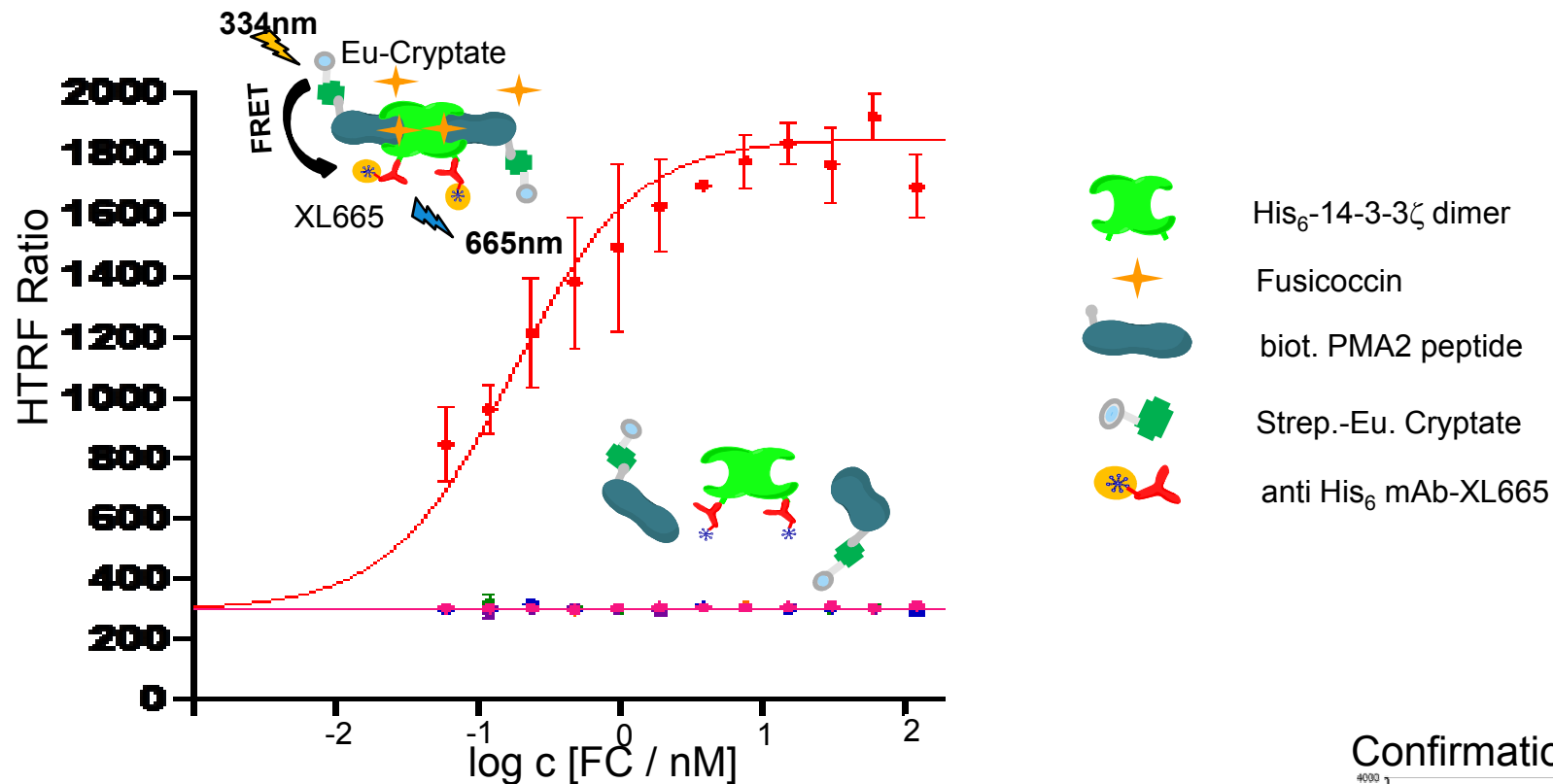
- **F**luorescence **P**olarization (**FP**)
- α -Screen
- **H**omogeneous **T**ime **R**esolved **F**luorescence (**HTRF**)

Assay development: Our choice HTRF



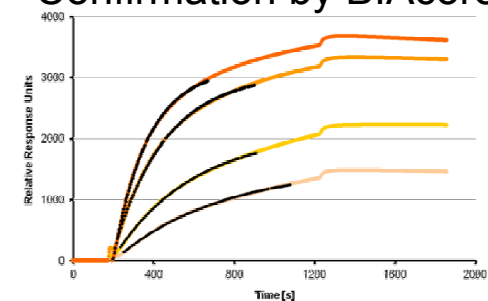
- HTRF toolbox reagents give a good flexibility for assay development
- Signal stability
- Sensitive readout in the far red region (665nm)
- Donor Em. at 620nm can be used as an internal reference
- Ratiometric readout (665/620nm)
- More than one distributor
- Good support (trouble shooting)

Assay development: HTRF assay for PMA2



- HTRF assay performed for: 14-3-3 / PMA2 \pm Fusicoccin
- K_d : < 3 nM
- Confirmation by BIAcore: K_d : 0.85 nM

Confirmation by BIAcore

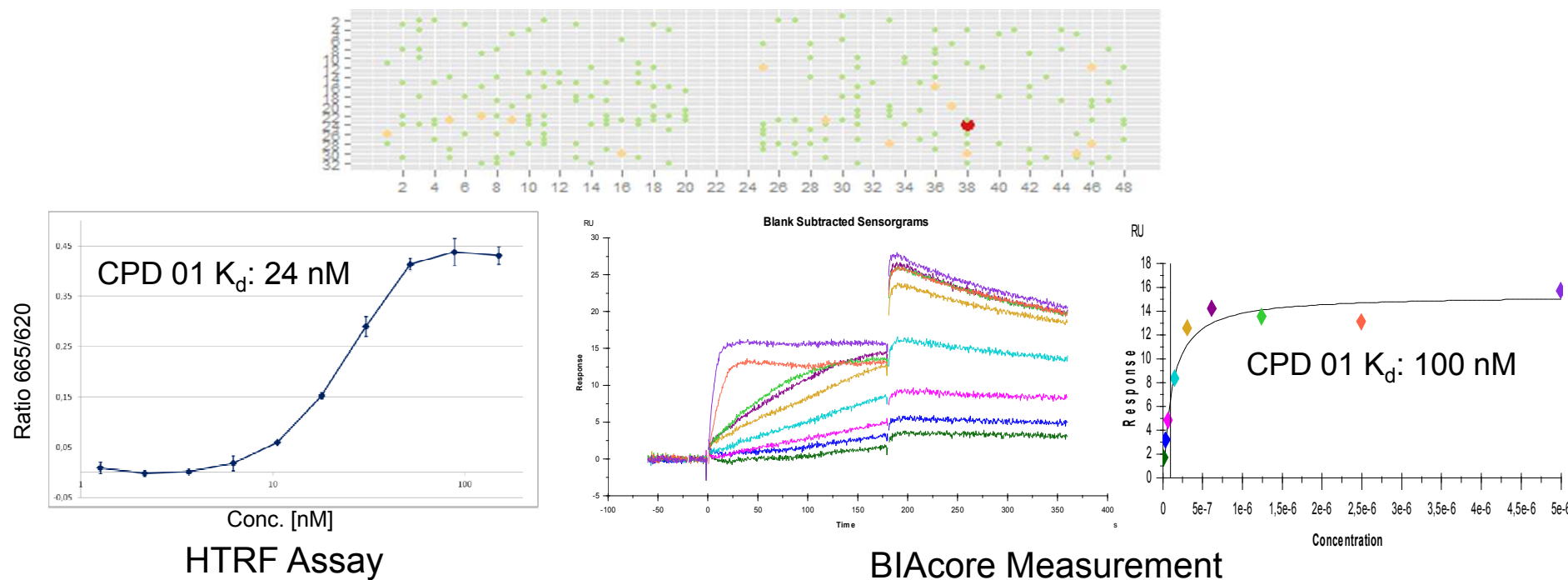


Assay development: Adaptation to screening platform



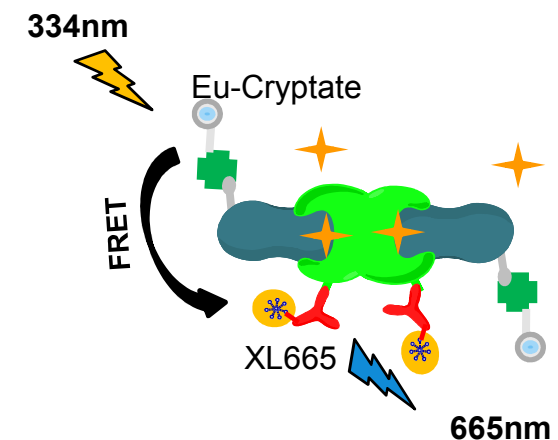
- **Scale down to 1536well format**
- **Total assay volume 8 μ l**
- **Signal / Background ratio: 3**
- **Signal / Noise ratio: 25**
- **Z'-factor: 0,76**

Identification of novel 14-3-3 PPI stabilizers



- ~150.000 compounds were screened
- 320 compounds were picked as primary hits
- 2 two scaffolds could be identified as new stabilizers for 14-3-3 PPI
- 1 scaffold was verified by BIAcore with a nM K_d

- **HTRF is a robust generic assay technology used in HTS environments**
- **The HTRF toolbox reagents proved to be a versatile instrument to set up PPIs assays**
- **Modulation of 14-3-3 PPIs by small molecules is possible**
- **Could lead to novel ways of pharmacological intervention by targeting 14-3-3 PPIs**
- **Goal: Development of a screening-toolbox to probe the extensive 14-3-3 interactome for novel PPI stabilizers**



Acknowledgement



Jan Eickhoff
Bert Klebl
Peter Nussbaumer
Axel Choidas
Malgorzata Skwarczynska
Carsten Degenhart
Uwe Koch
Matthias Baumann
Gerd Rüther
Corinna Lechleitner



MAX-PLANCK-GESELLSCHAFT



Bengt Hallberg



Luc Brunsveld

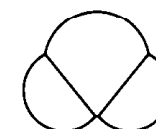


Isabelle Landrieu



CGC
Chemical Genomics Centre
of the Max Planck Society

Christian Ottmann
Lars Röglin
Philipp Thiel
Maria Bartel
Rolf Rose



mpi

Herbert Waldmann
Alfred Wittinghofer

14-3-3Stabs

IAPP Grant: 286418



Matthias Stein-Gerlach

Max-Planck-Innovation



General Assembly:

Jörn Erselius

Axel Ullrich

Herbert Waldmann



Ministerium für Innovation,
Wissenschaft, Forschung und Technologie
des Landes Nordrhein-Westfalen



MAX-PLANCK-GESELLSCHAFT



Thank you!

Lead Discovery Center GmbH

Otto-Hahn-Str. 15

44227 Dortmund

Telefon +49.231.97 42 70 00

www.lead-discovery.de

Industrial Advisory Board:



**Boehringer
Ingelheim**



Alexander Wolf
Assay Development and Screening
Lead Discovery Center GmbH
Otto-Hahn-Str. 15
44227 Dortmund
Germany

wolf@lead-discovery.de