



Novartis Institutes for BioMedical Research  
Rochdi Bouhelal  
SBS Sept. 2005

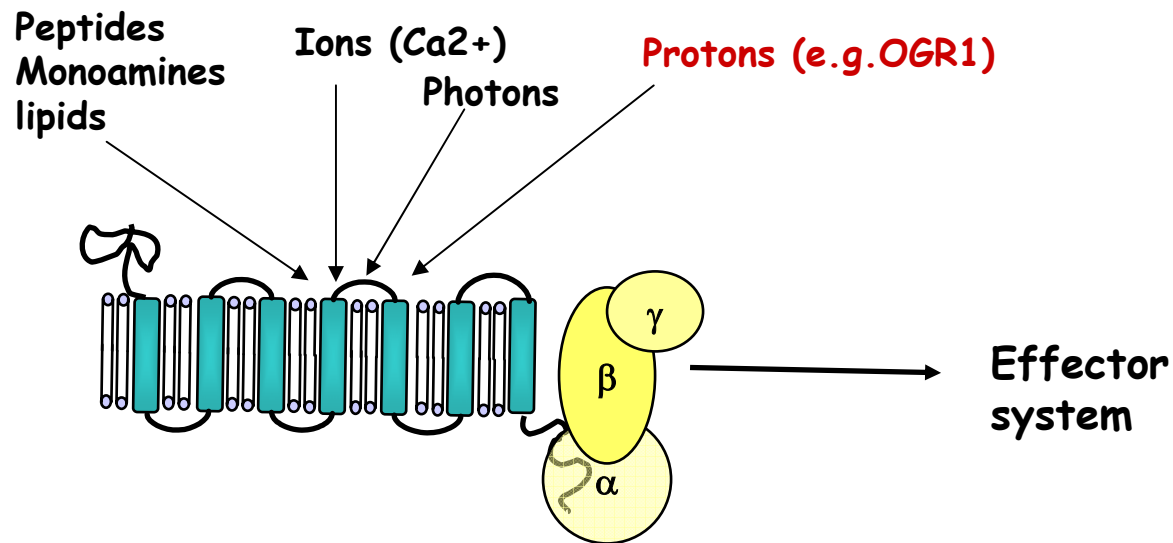


**A new HTRF inositol phosphate assay to monitor G<sub>q</sub> coupled GPCRs responses**

– Comparative study calcium mobilisation / IP1

Ina Hammerl, Stéphane Martinez, Monique Amoravain and Rochdi Bouhelal

## GPCRs: A universal communication system



### Signal Transduction mechanisms

- cGMP PDE,
- Adenylate cyclase (cAMP) & Protein kinase A
- Phospholipase C , IP3 , calcium & Protein kinase C,
- MAP kinase pathways,
- Ion channel channels conductance

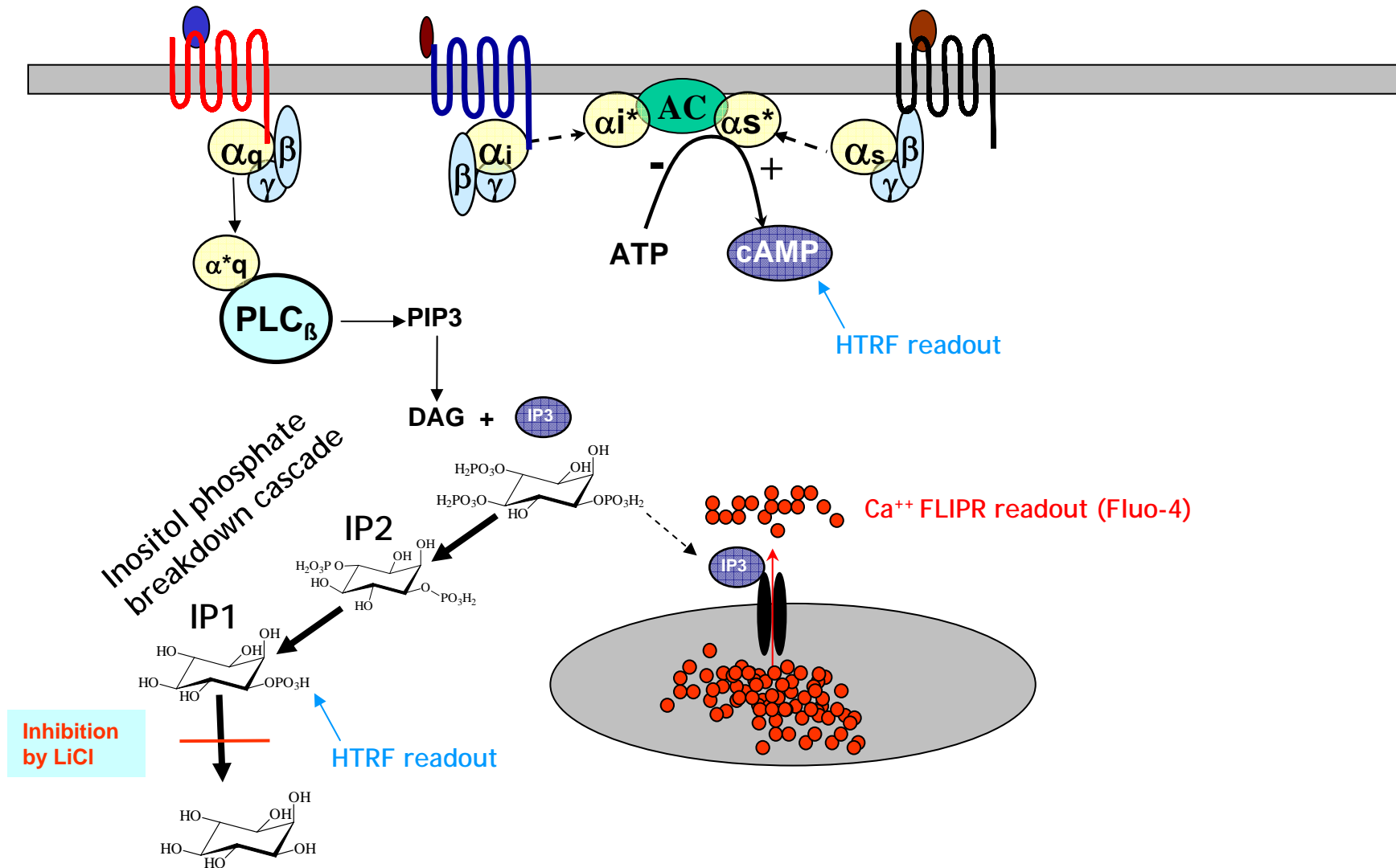
## Need for a HTS assay format for IP formation

- Several assay formats exist to monitor all events in the GPCR activation cascade with a high throughput.
- While such assays to monitor **cAMP** and **calcium** levels were developed the past 10 years and are now widely used in the pharmaceutical industry, current IP technologies are limited by their low throughput and safety issues.
- Need of assays amenable to HTS or MTS in the lead discovery process.
- Recently, a homogeneous HTRF assay was developed by Cisbio which measures IP1 the last component of the PIP2 degradation pathways.
- A study was initiated at Novartis with the objective to validate this novel assay format and to evaluate its usefulness in our discovery processes.

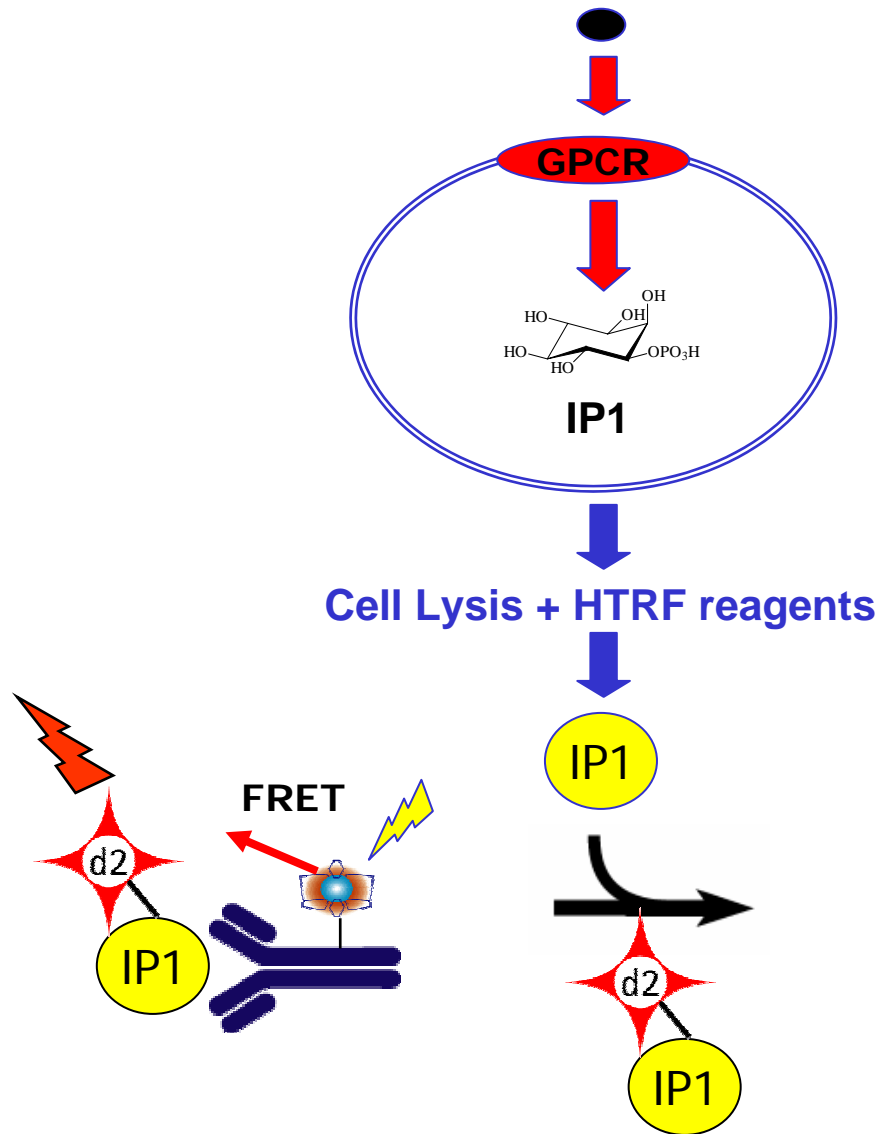
## Study description

- The assay was assessed in various cell systems
  - HEK293 and CHOK1 cell expressing native GPCRs
  - CCL39 cell expressing the recombinant human parathyroid calcium sensing receptor (HupCaR).
- Pharmacological characterization of GPCRs with agonists and antagonists.
- In addition, Novartis libraries are tested in the HupCaR calcium mobilization assay using the FLIPR technology and the HTRF IP1 accumulation assay using an imaging reader (Viewlux)
- Aim: Comparison of hit rates obtained in the two assays and the suitability of the IP1 assay as a primary or secondary assay for GPCR screening.

# Gq signaling pathways

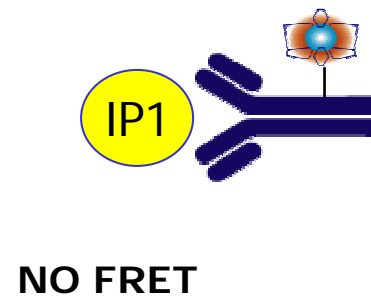


## IP-one assay principle



based on unique and proprietary reagents

- A fluorescent analog of IP1
- A monoclonal antibody against IP1



## Methods: HTRF & Calcium assays

### HTRF / IP1

#### Cell seeding

- 10 000 cells/well/50 ul media in a white tissue culture treated 384 MTP
- 24 h incubation at 37 °C and 5 % CO<sub>2</sub>

#### Cell stimulation

- Cell media removal
- Cells stimulation with agonists / compounds
- diluted in HBS Buffer with (LiCL 50 mM) for 30 min
- addition of HTRF reagents and incubation for 1 h

#### HTRF Reading:

- read plate in Viewlux at 665 and 620 nm

### FLIPR Fluo4 calcium

#### Cell seeding

- 10 000 cells/well/50 ul media in a black tissue
- culture treated 384 MTP with clear bottom.
- 24 h incubation at 37 °C and 5 % CO<sub>2</sub>

#### Cell loading:

- After removing cell media, load cells with Fluo4
- incubate 1 h at 37°C and 5 % CO<sub>2</sub>

#### Cell washing:

- Remove remaining Fluo4 by washing plates

#### Cell stimulation & Calcium Reading

- Read plate in FLIPR at 525 nm during compound injection





## Endogenous P2YR and M3R in HEK293 cells

# Endogenous Muscarinic M3R and P2YR in HEK293 cells

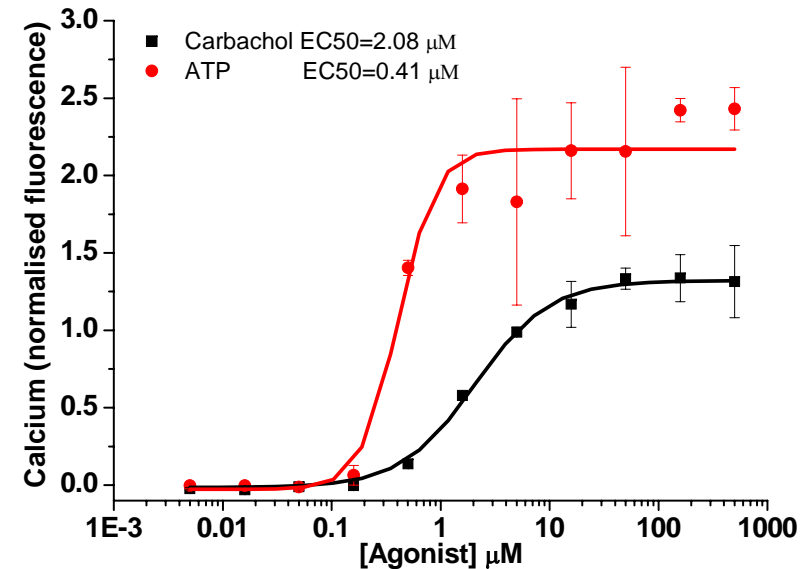
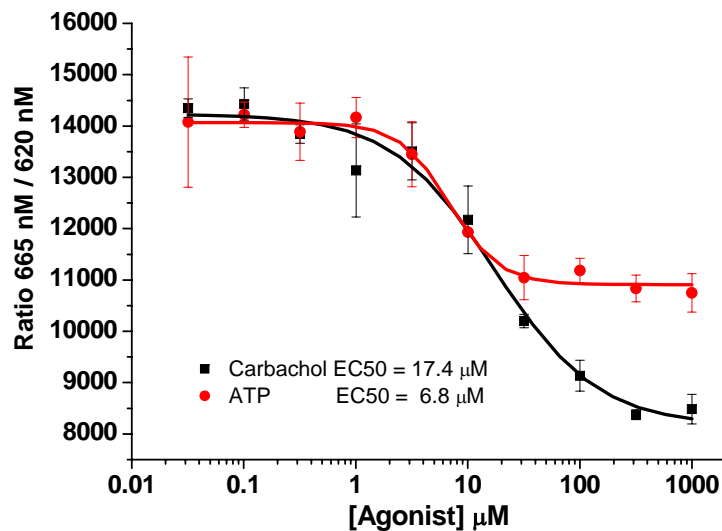
HTRF / IP1

60,000 cells, 24 h

**Agonist effects**

FLIPR Fluo4 calcium

10,000 cells, 24h

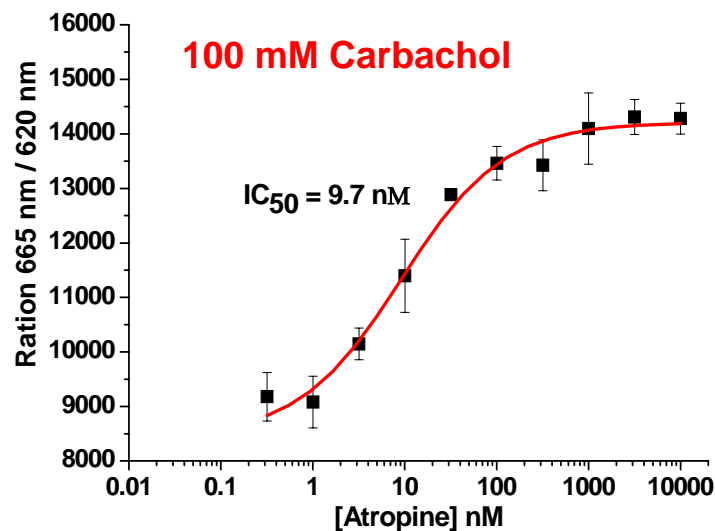


- Carbachol and ATP are more potent in the calcium assay
- Curve shift factors 8 and 16 for carbachol and ATP
- Differences between efficacies in the two assay systems are also reflected.
- Calcium mobilisation assay is more sensitive (amplification step) in particular with low receptor expression (endogenous receptors)

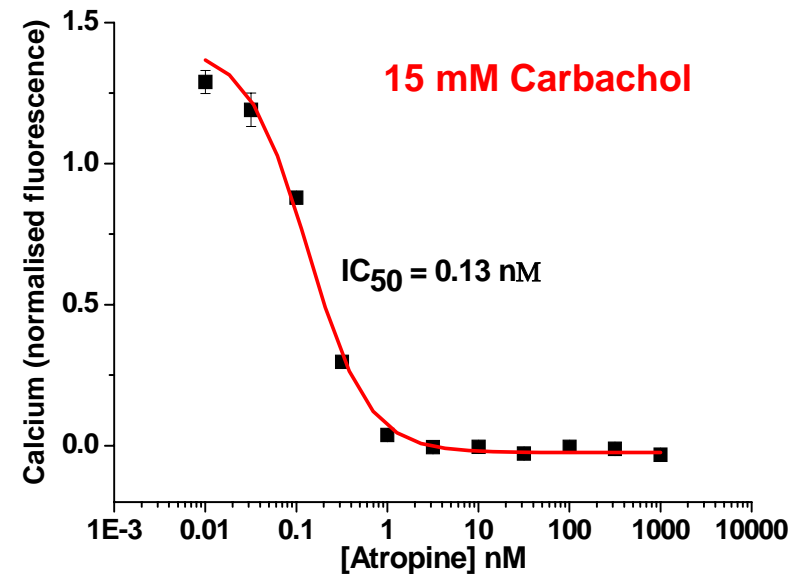
# Endogenous Muscarinic M3R in HEK293 cells

## Antagonist effects

### HTRF / IP1



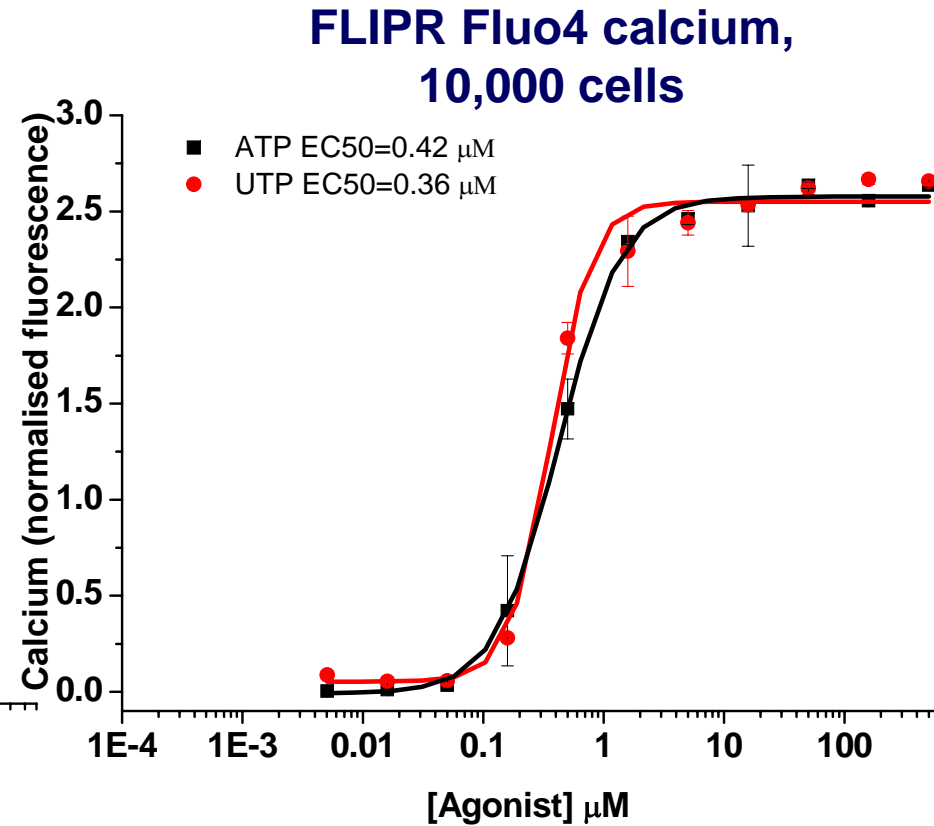
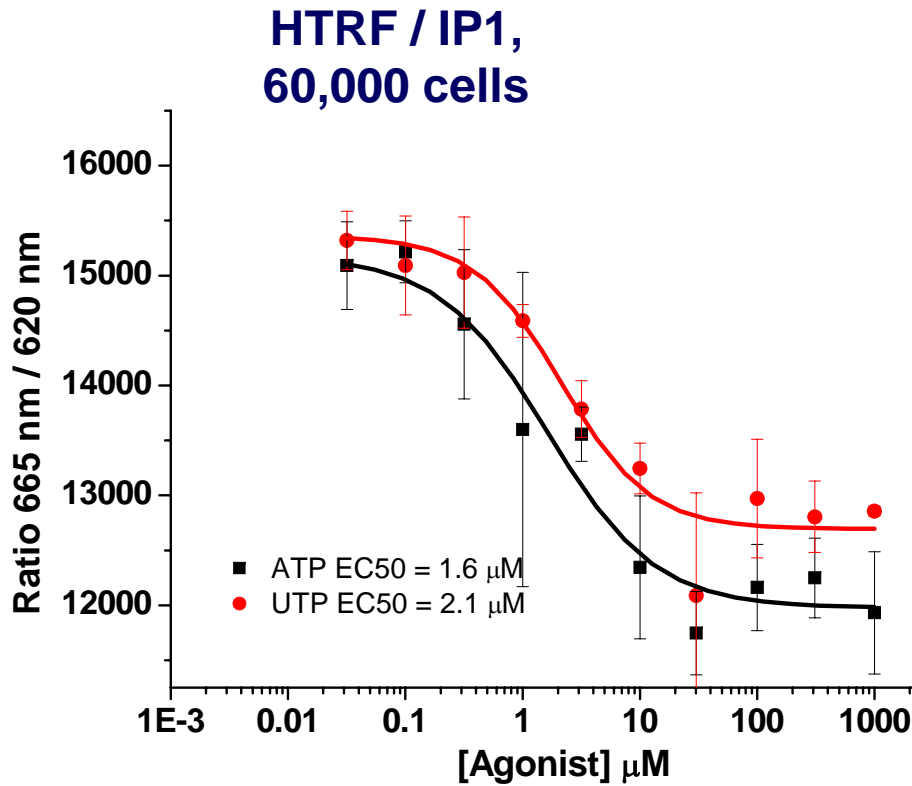
### FLIPR Fluo4 calcium



- Atropine potency higher in calcium assay (75 fold)
- Reflects differences of agonist concentrations and receptor number used  
 Carbachol 100  $\mu$ M in IP1 / 15  $\mu$ M in Calcium  
 60,000 cells in IP1 / 10,000 in calcium

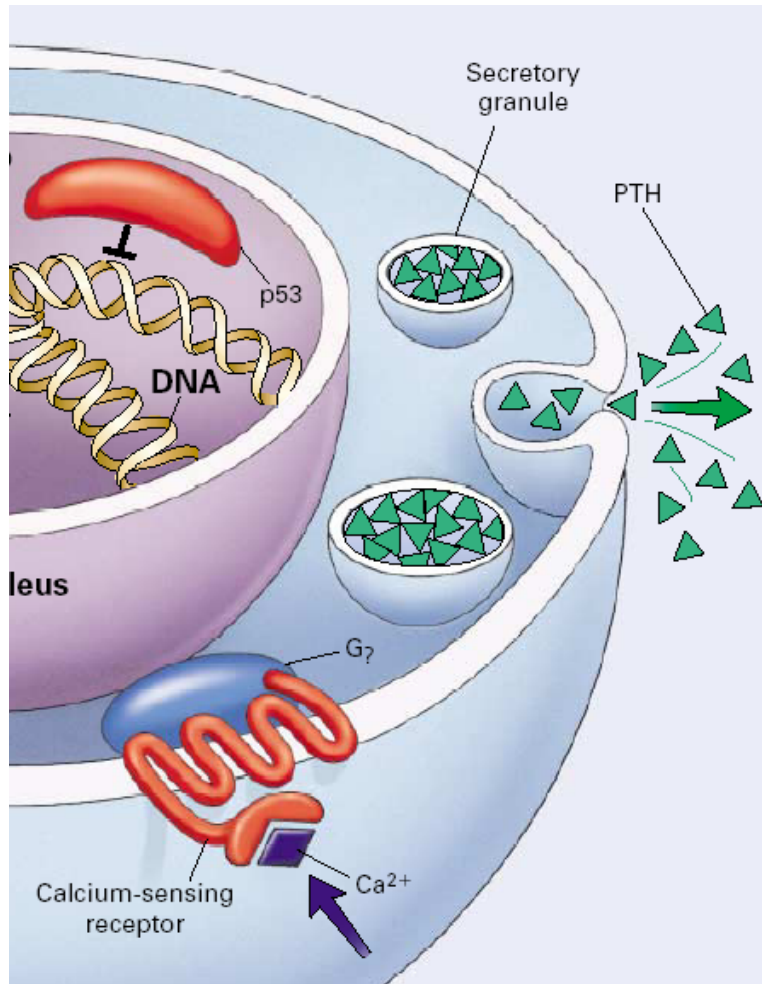
**Conclusion:** No major difference in antagonist potency expected

## Endogenous P2Y receptors in CHOK1 cells: Agonist effects



**Recombinant human calcium sensing receptor  
(HupCaR) in CCL39 cells**

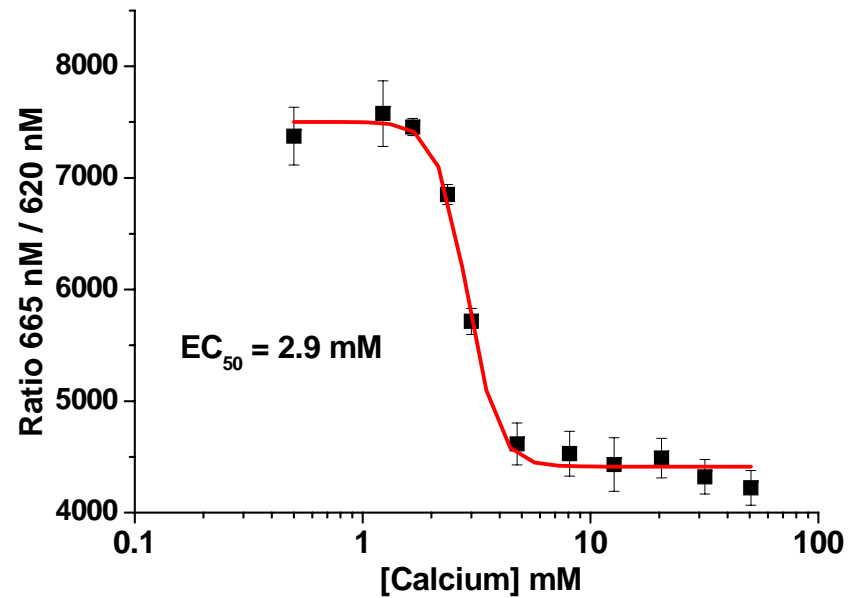
## The parathyroid calcium sensing receptor



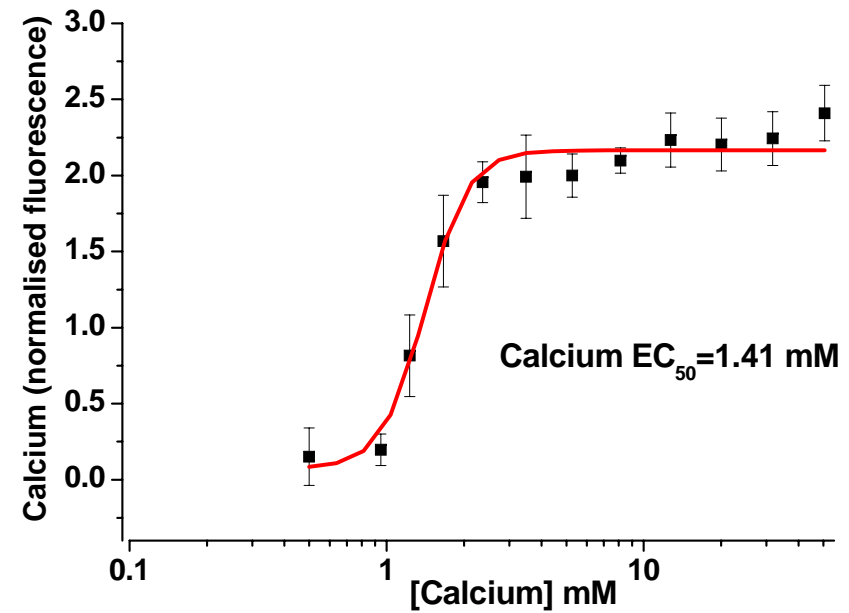
- A Gq coupled receptor highly expressed In the parathyroid gland
- Uses circulating calcium as an “agonist”
- Can be blocked by allosteric negative modulators
- Controls PTH release
- Role in bone formation

## Effect of calcium on IP1 & Calcium mobilisation

### HTRF / IP1



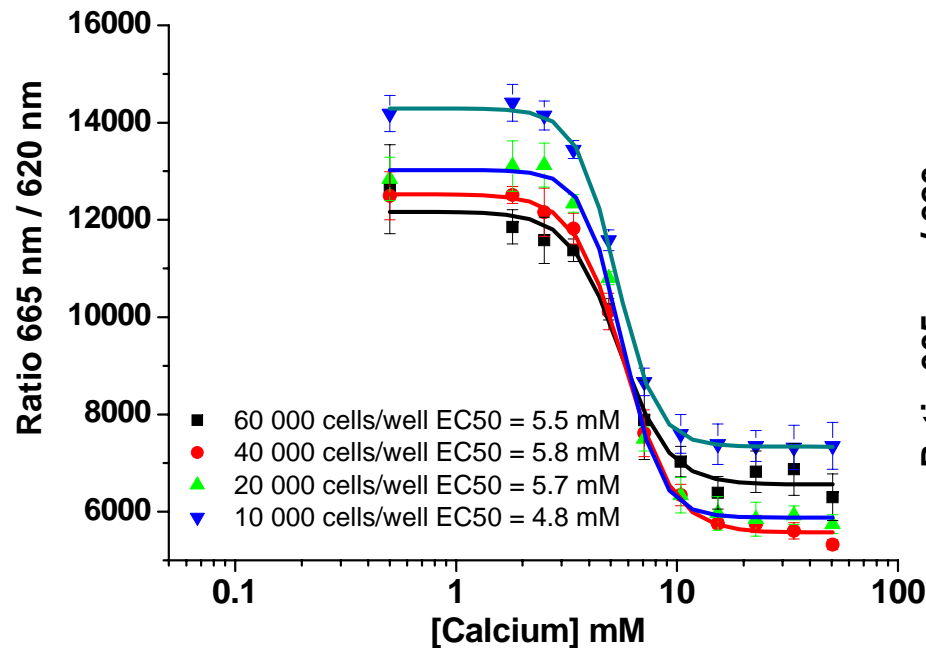
### FLIPR Fluo4 calcium



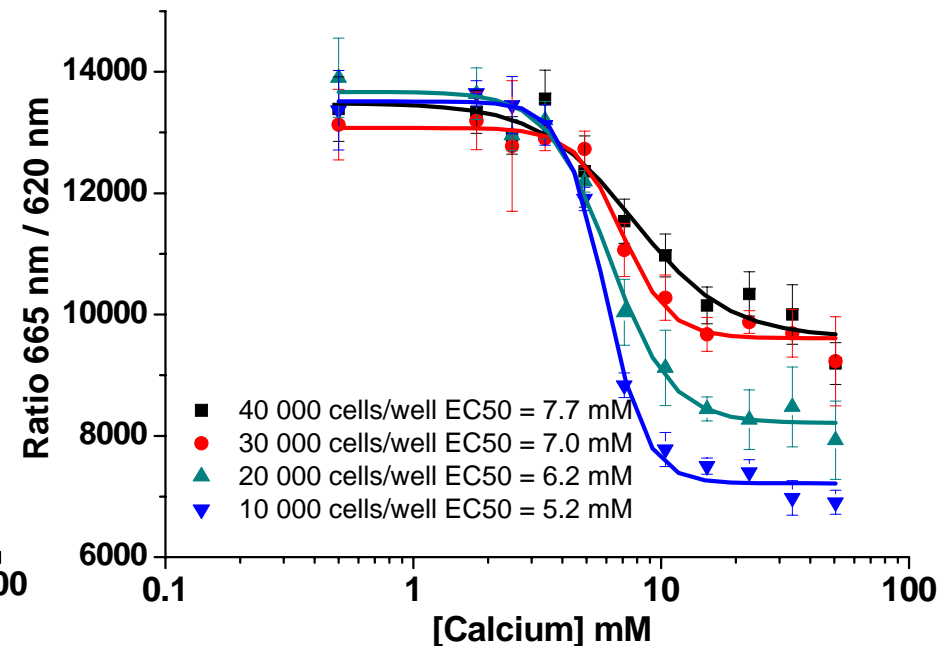
10000 cells/well, 08.07.05

## HupCaR: Evaluation of HTRF IP1 assay parameters: cell density

### Cell growth in plate over 24 h



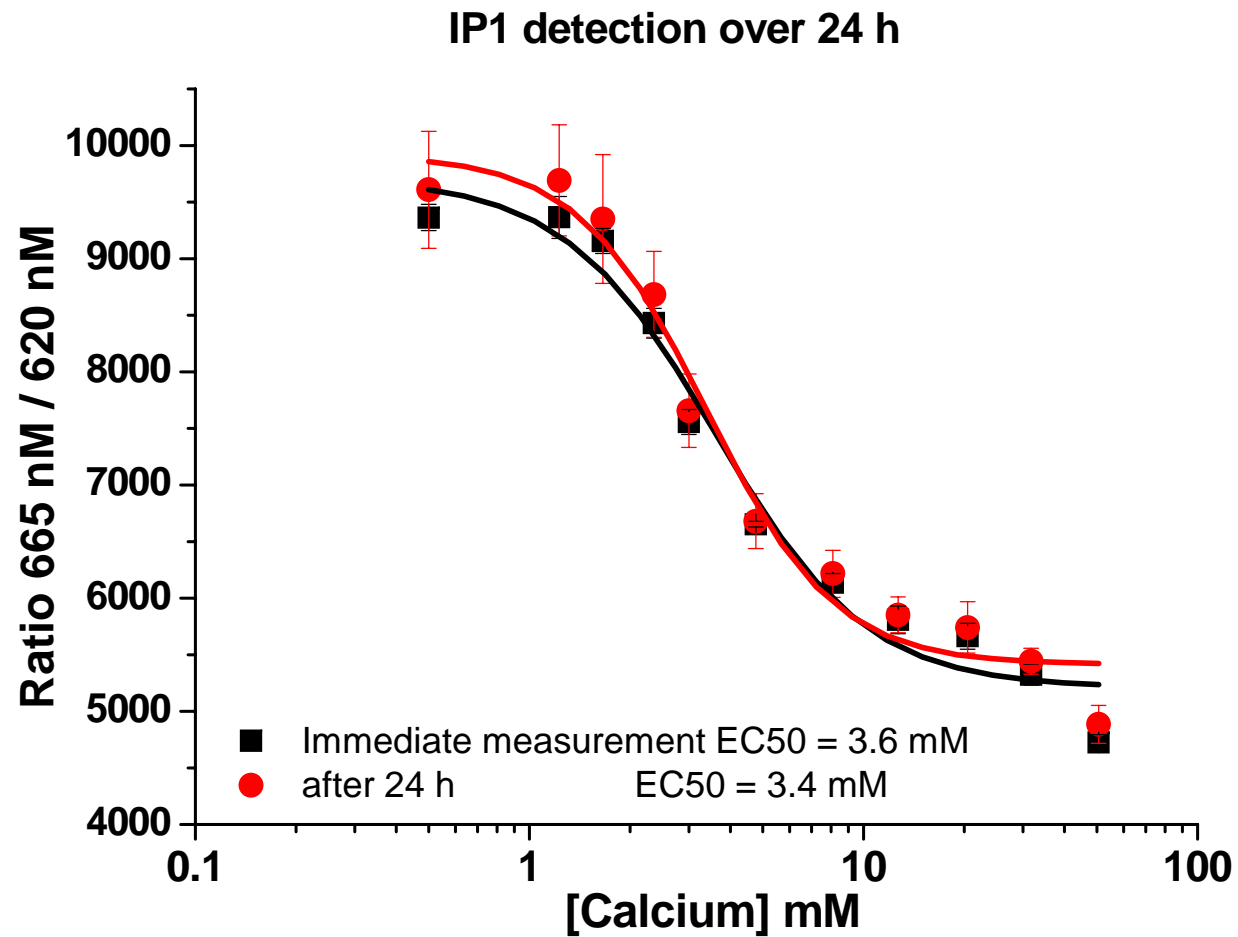
### Cell growth in plate over 48 h



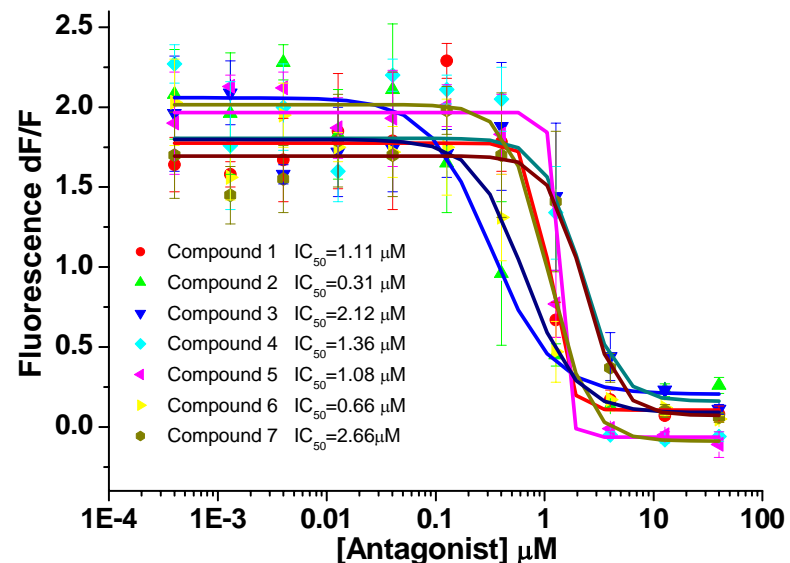
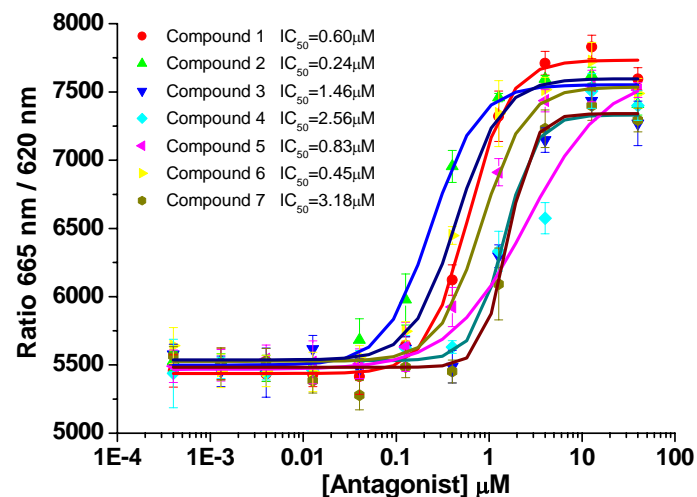
- Potencies unchanged
- Cell density can be reduced to 5000-10000 cells /well over 24 or 48 h



## HTRF signal stability over time

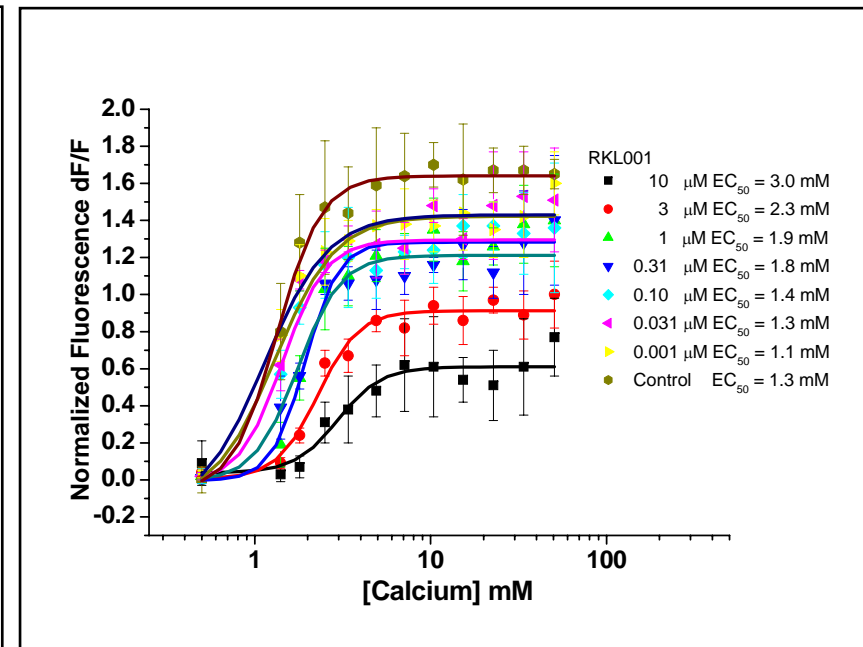
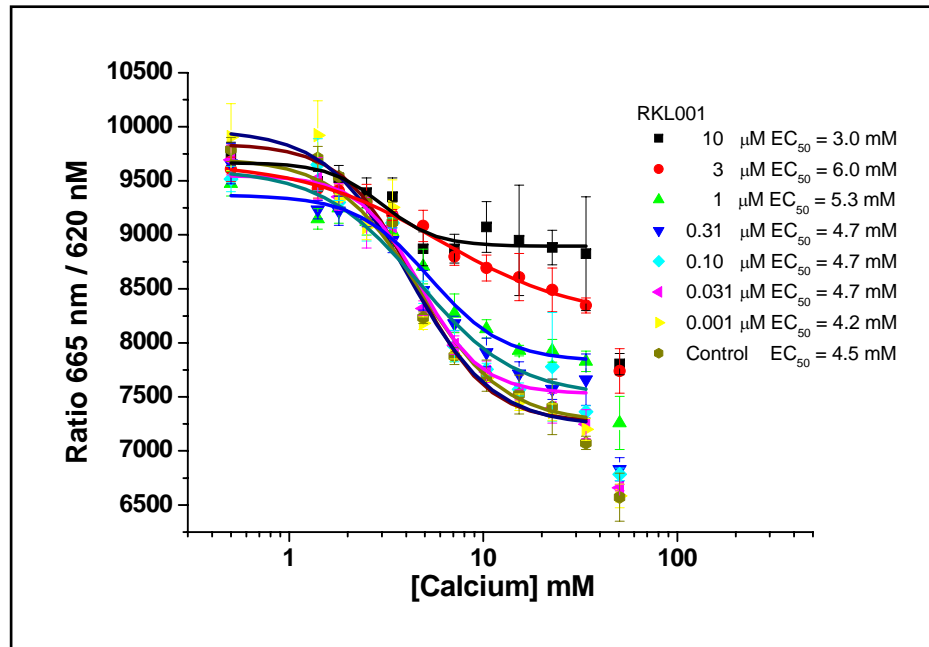


# Antagonist effects in the HTRF & Calcium assays



Compound ID	IP1 (HTRF)	Calcium (FLIPR)
	IC50 ( $\mu M$ ) n = 3	
1	0.58 $\pm$ 0.02	0.74 $\pm$ 0.32
2	1.00 $\pm$ 0.67	0.66 $\pm$ 0.51
3	1.23 $\pm$ 0.38	1.75 $\pm$ 0.84
4	2.91 $\pm$ 0.81	1.98 $\pm$ 0.54
5	0.61 $\pm$ 0.21	1.32 $\pm$ 1.04
6	2.51 $\pm$ 1.87	1.32 $\pm$ 1.43
7	1.88 $\pm$ 1.13	3.06 $\pm$ 1.14

## Antagonist effects: RKL001



- Varying blocker concentration depresses maximal activity without major changes in calcium  $\text{EC}_{50}$

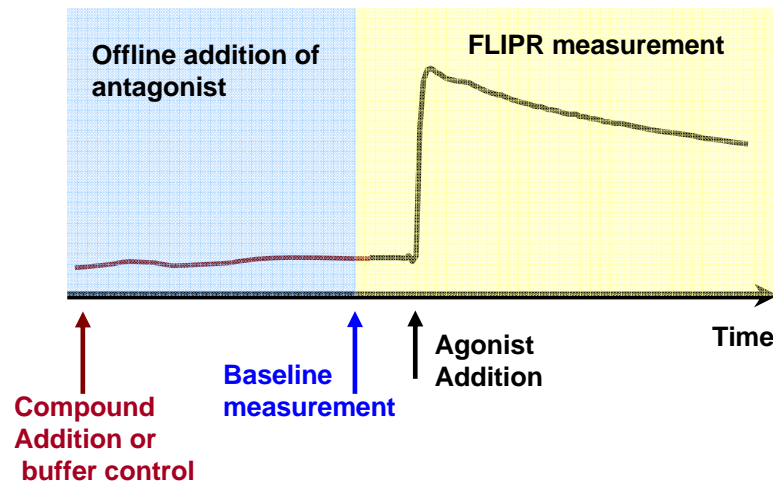
## IP1 in HTS

### Study aim

- Use the IP one assay in a productive screening campaign
- Comparison to calcium mobilization screen
- 7744 natural product tested

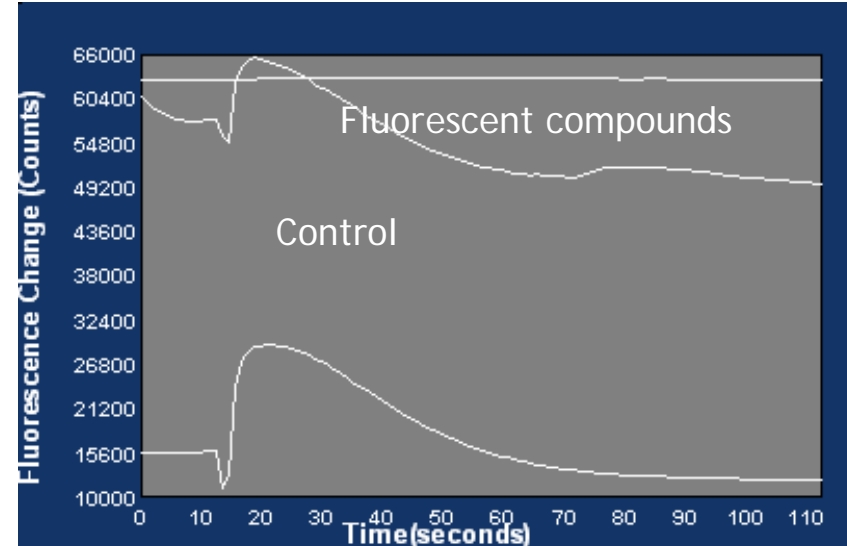
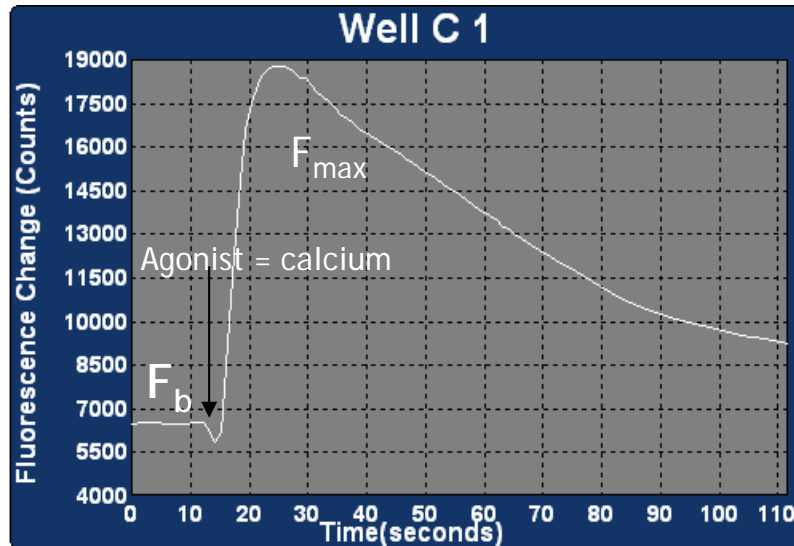
## FLIPR assay types: Data Analysis

### Antagonist FLIPR assay set-up



- 
- $F_{\text{basal}}$  or  $F_b$  = Fluorescence before agonist injection
  - $F_{\text{max}}$  = Fluorescence maximum or peak
  - $dF/F$  = Normalised Fluorescence =  $(F_{\text{max}} - F_b) / F_b$
  - $F_{b \text{ sample}} / \text{average } F_{b \text{ controls}}$  = high ratio indicates autofluorescence/ toxicity of compound

## FLIPR data handling for antagonist assays



Two values of fluorescence calcium responses are exported

$F_b$  corresponding to the value prior to agonist injection,  $F_{max}$ , the fluorescence at the signal peak.

From these the values two parameters were then calculated and exported

Calculate  $dF/F = F_m - F_b / F_b$  for High controls H, Low controls (L, Buffer) and Samples (S).

Activity (A) expressed as a percent of the maximal stimulation induced by agonist .

—  $A (\%) = (S-H) / (H-L) * 100$

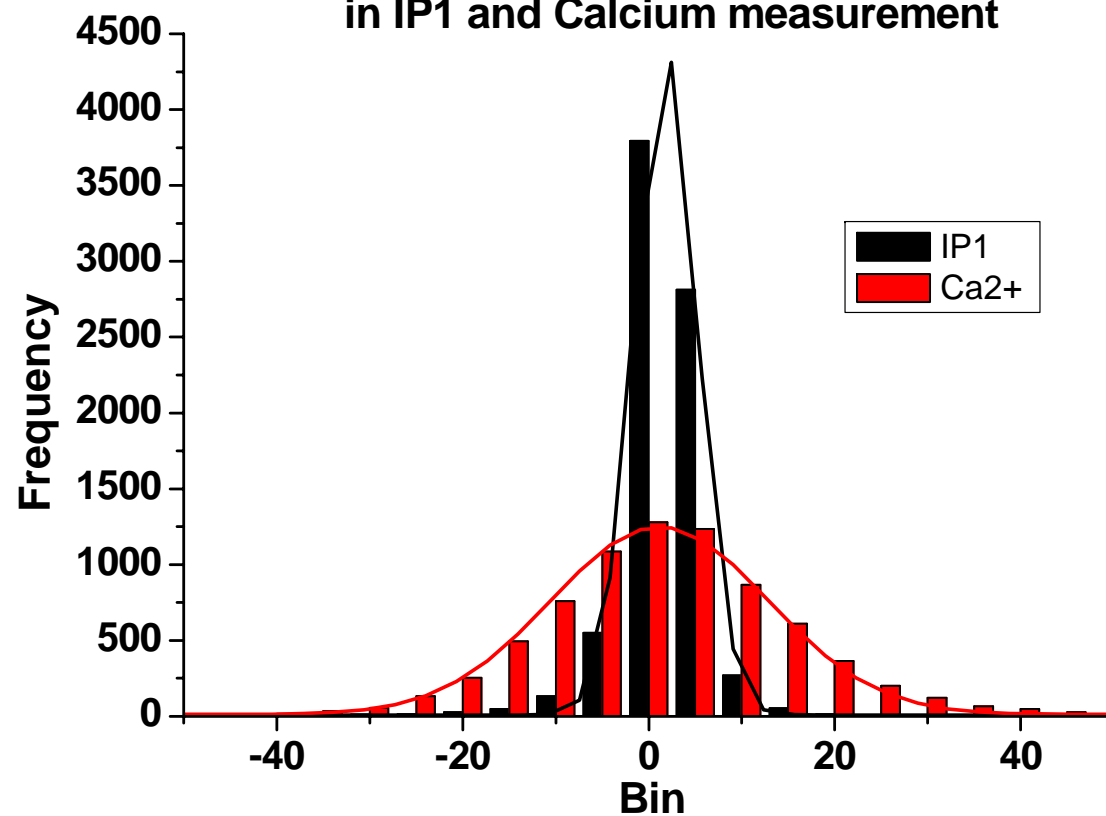
— A quality parameter ( $R_b$ ) which indicates whether the compounds acted through a physiologically relevant mechanism)

$R_b = F_{b,s} / F_{b,H}$

— This parameter is used to reject compounds showing a sustained activity (toxicity / autofluorescence)

## Calcium receptor HupCaR miniscreen - HTRF & Calcium assay

Analysis of 7744 compounds from Napr collection  
in IP1 and Calcium measurement



Threshold		-30	-40	-50	-60	-70
Hit number	IP1	17	4	2	2	1
	Ca <sup>++</sup>	165	75	50	37	28

## Miniscreen: Data quality

	IP1	Ca <sup>2+</sup>
EC <sub>50,Ca</sub> [mM]	4.94 ± 3.39 n=9	1.97 ± 0.53 n=7
IC <sub>50, Anta</sub> [μM]	0.48	1.3

	IP1	Ca <sup>2+</sup>
mean % change	-0.45	-0.57
sd	4.74	15.02
3*sd	14.2	45.0
Z'	0.85 ± 0.05	0.56 ± 0.10
Z	0.86 ± 0.05	0.66 ± 0.06



## IP1 and Calcium: Compound interference

Compound	IP1	Calcium	
	% change	% change	Ratio Rb
1	6	-59	1.34
2	-4	-64	2.03
3	-3	-85	2.13
4	-3	-102	3.65
5	0	-102	4.07
6	-3	-100	4.09
7	-3	-102	3.96
8	0	-102	4.57
9	-2	-102	4.04
10	1	-93	3.8
11	5	-102	4.38
12	3	-100	3.02
13	-1	-93	2.33
14	-4	-72	1.45
15	-1	-72	1.58
16	-3	-97	1.77
17	-12	-102	2.59
18	-3	-91	2.42

– Frequent hitters are compounds scoring positive in all FLIPR screening campaigns due to their toxicity, their fluorescence at 488 nM or their interaction with common pathways.

– These compounds are not detected in the HTRF IP one assay

## IP1 and Calcium: Common hits

Compound	IP1	Calcium	
	% change	% change	Ratio Rb
1	<b>-37</b>	<b>-78</b>	0.75
2	<b>-31</b>	<b>-93</b>	0.82
3	<b>-22</b>	<b>-50</b>	0.70
4	<b>-35</b>	<b>-78</b>	0.73
5	<b>-34</b>	<b>-68</b>	0.73
6	<b>-30</b>	<b>-74</b>	0.74
7	<b>-38</b>	<b>-65</b>	0.77

— Hits interfering with the receptor of common mechanism

## IP1 and Calcium: PS hits in calcium

Compound	IP1	Calcium	
	% change	% change	Ratio Rb
1	-9	<b>-54</b>	0.91
2	1	<b>-88</b>	1.00
3	13	<b>-67</b>	0.86
4	3	<b>-55</b>	0.71
5	7	<b>-63</b>	0.93
6	2	<b>-58</b>	1.24
7	1	<b>-55</b>	1.02
8	-1	<b>-64</b>	0.71
9	-2	<b>-72</b>	0.70
10	3	<b>-62</b>	0.63
11	4	<b>-86</b>	0.66
12	11	<b>-67</b>	0.70
13	3	<b>-54</b>	0.78
14	-2	<b>-55</b>	0.85
15	-1	<b>-88</b>	0.81
16	-3	<b>-56</b>	0.84
17	-2	<b>-70</b>	0.62
18	-1	<b>-61</b>	1.05
19	2	<b>-71</b>	0.98
20	-5	<b>-72</b>	0.54
21	13	<b>-53</b>	1.05
22	12	<b>-52</b>	1.21
23	6	<b>-73</b>	0.71

– A number of compounds are found only in the calcium assay without modification of the calcium baseline

Common FLIPR hits, i.e. found in several FLIPR assay (

False positives in calcium assay; Known for FLIPR assays

Other mechanisms ?

## Conclusion / outlook

- IP1-one is a useful HTS assay using IP1 as a reporter for IP3
  - Homogenous assay. Some parameters were optimised (Cell density & culture condition)
  - Similar data obtained compared to calcium FLIPR with a slightly higher sensitivity for calcium (amplification mechanisms)
  - Hit rates lower
    - Can be used for secondary screening to exclude false positives and FLIPR specific hits
  - More robust. Better assay quality in productive screening set-up
- Further studies
  - More studies with controlled conditions with endogenous GPCRs.
  - More data to be produced with other recombinant system (SMW agonist GPCR assay with competitive antagonists. Potential calcium interference with IP1 assay ?)

## Acknowledgements

Novartis

K. Seuwen

D. Gabriel

Cisbio France

E. Trinquet

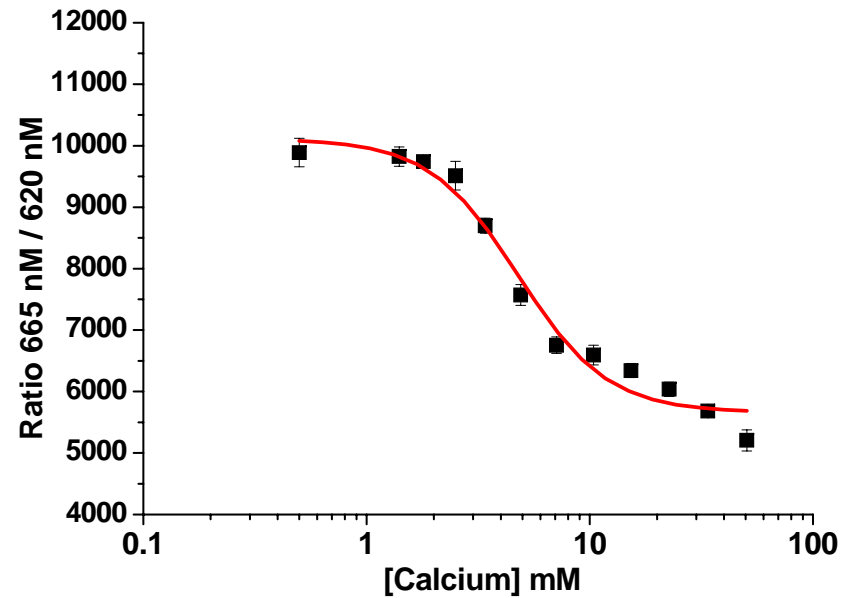
M. Fink

P. Seguin

J-L Tardieu

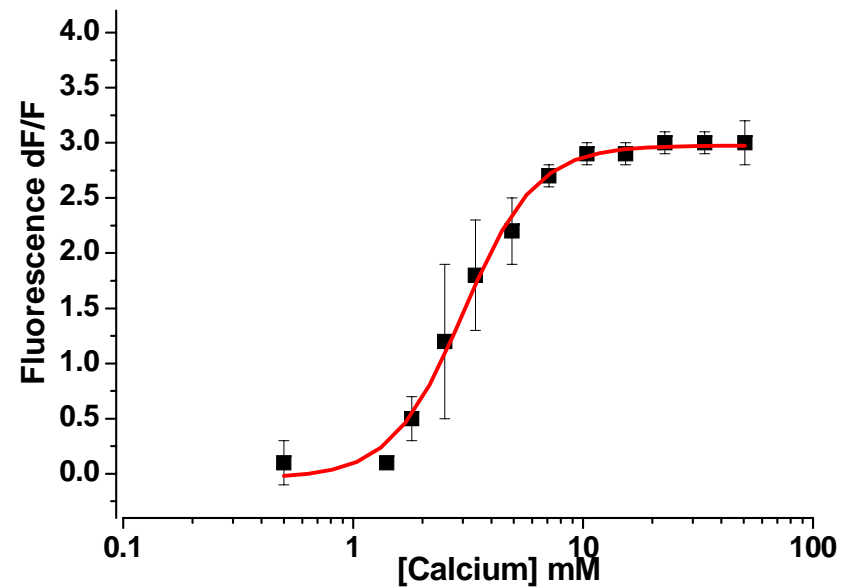
# The calcium sensing receptor HupCaR

## HTRF / IP1



$EC_{50} = 4.79 \text{ mM}$

## FLIPR Fluo4 calcium



$EC_{50} = 3.0 \text{ mM}$