



Novartis Institutes for BioMedical Research

Novel Functional Assay Approaches for GPCR Ligand Discovery

R. Bouhelal,

Centre for Proteomics Chemistry

Lead Discovery Platform

Basel, Switzerland

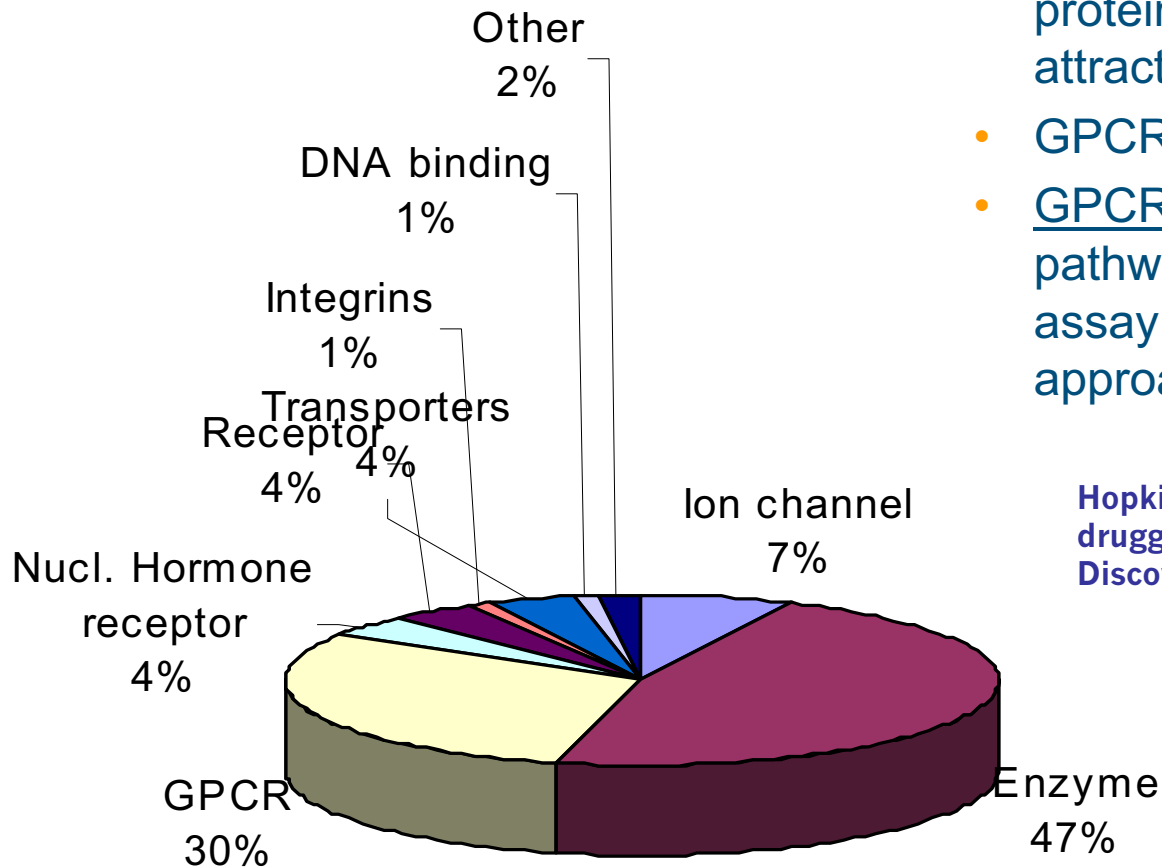
Screening Europe Barcelona
February 20-21, 2007

These slides were taken, courtesy of Rochdi Bouhelal, from a presentation given at Screening Europe in February 2007.



Drug targets for existing medicines

Novartis Institutes for BioMedical Research



- GPCR the most studied membrane protein and amongst the most attractive drug targets
- GPCR's déjà vu ?
- GPCR's renaissance: Novel pathways, molecular properties and assay technologies and HTS approaches are increasingly studied

Hopkins et al NRDD 1, 727-730 (2002). The druggable genome": Nature Reviews, Drug Discovery, Vol.1, September 2002

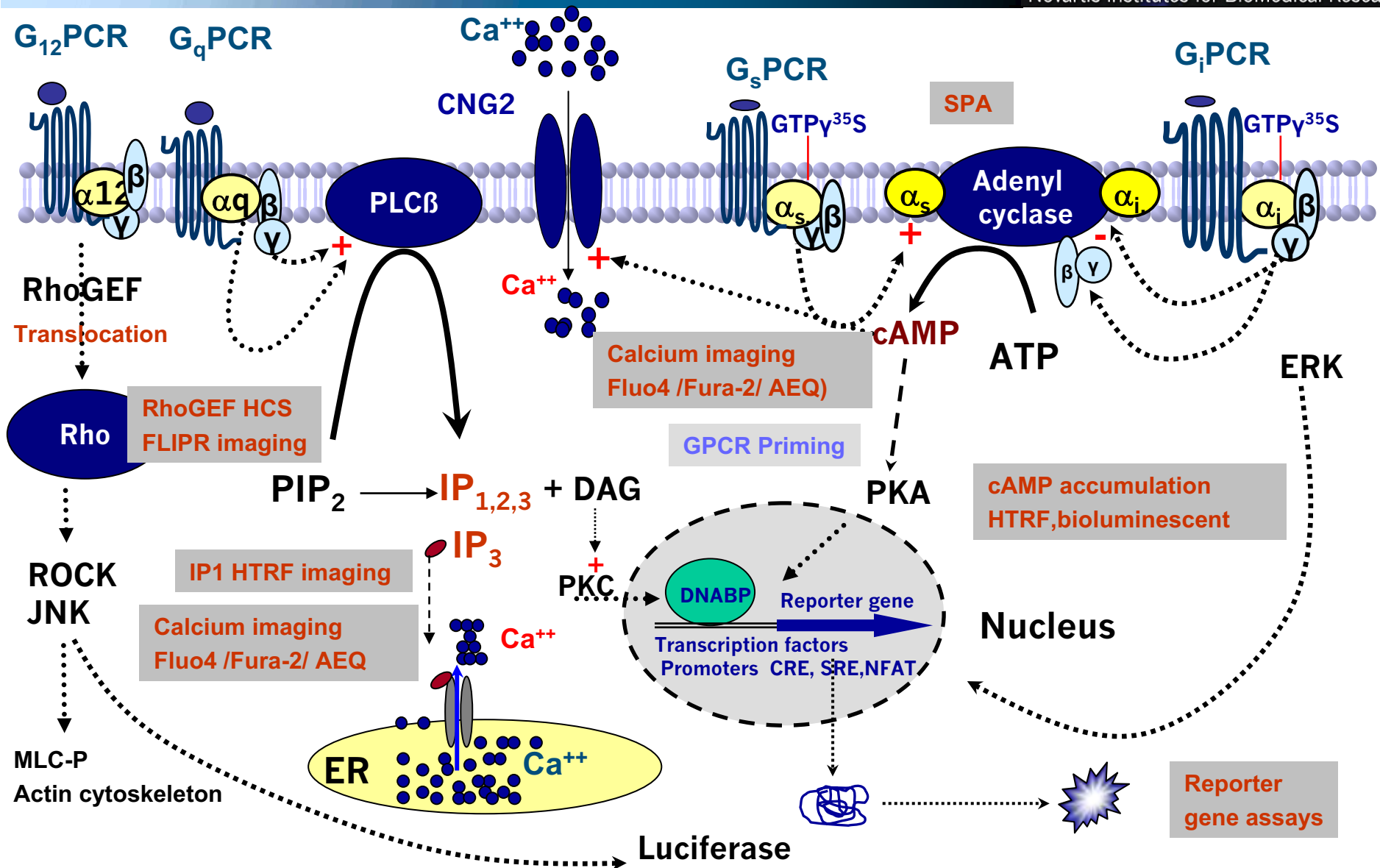
Monitoring GPCR activation: some facts & basic principles

Novartis Institutes for BioMedical Research

- **Functional & ligand binding GPCR assays**
 - The best approach to ligand discovery ?
 - **Ligand binding assay**
 - ⇒ in general straight forward but poor information content
 - ⇒ In general 1st choice for medicinal chemists
 - **Functional assays**
 - ⇒ Less simple to configure, high hit rates, false positives
 - ⇒ Access to the GPCR biology , Uncovers novel compound mechanisms
 - ⇒ May offer rich information output
- **Try to remain proximal to the receptor**
 - However reporter gene assays can be used in some circumstances
- **Imaging is used to enhance throughput**
 - One single screening platform for orphan receptors and for HTS
- **Use non-invasive assays**
 - Try to bring all receptors to signal through calcium
 - ⇒ Assay multiplexing (selectivity, several pathways)
 - ⇒ Allows to align more easily HTS data
 - Rich pharmacological data output including kinetic data analysis
 - ⇒ Helps excluding non-specific compounds

GPCR signaling pathways and applied technologies is lead discovery

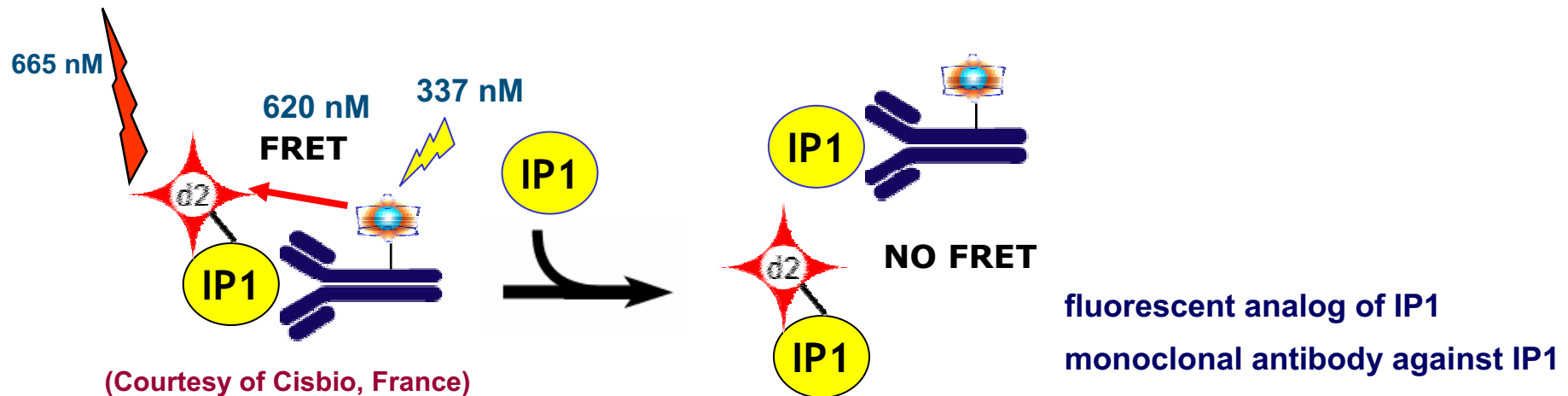
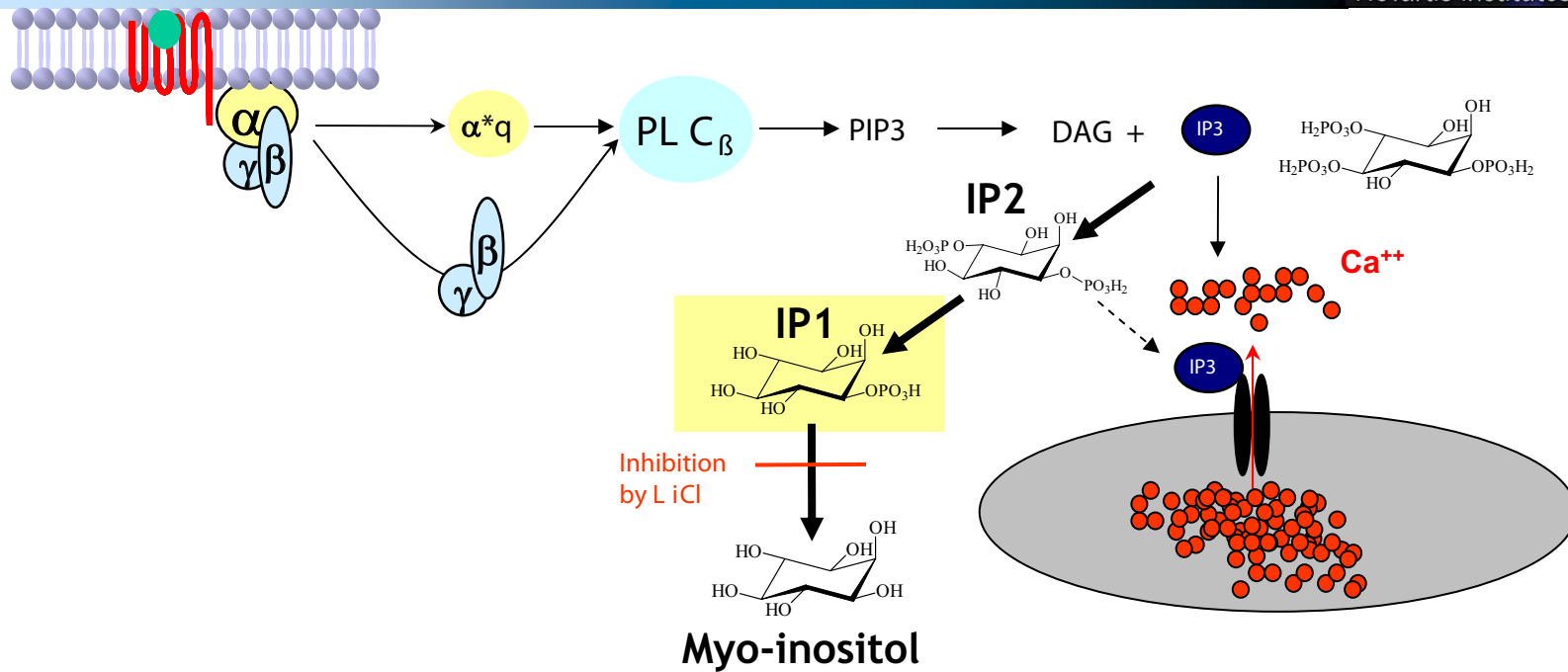
Novartis Institutes for BioMedical Research



Inositol phosphate 1 for activation Gq coupled receptors

IP1 accumulation HTRF assay

Novartis Institutes for BioMedical Research

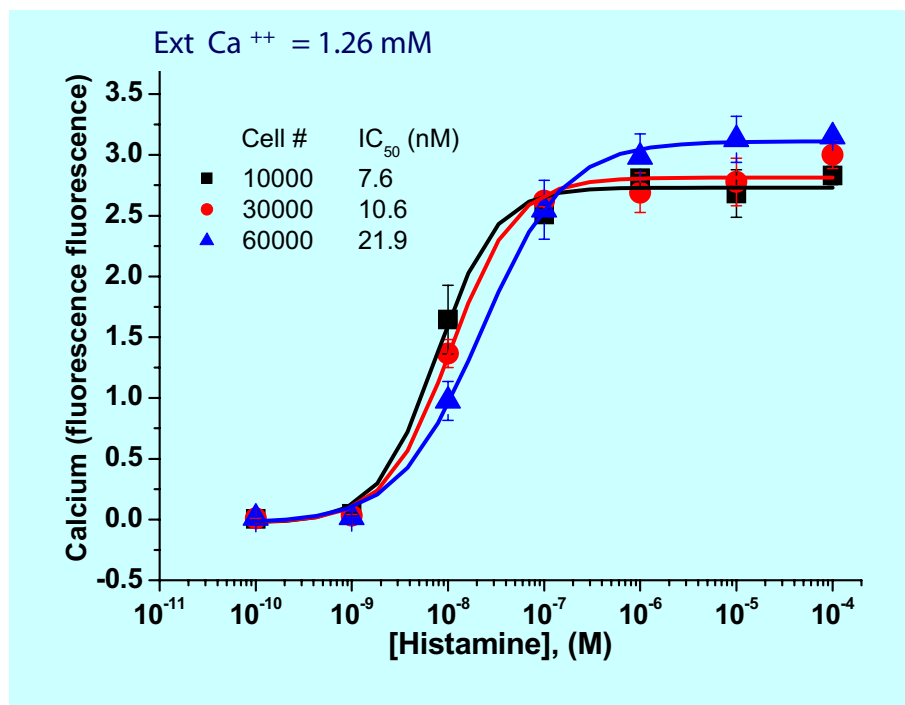


(Courtesy of Cisbio, France)

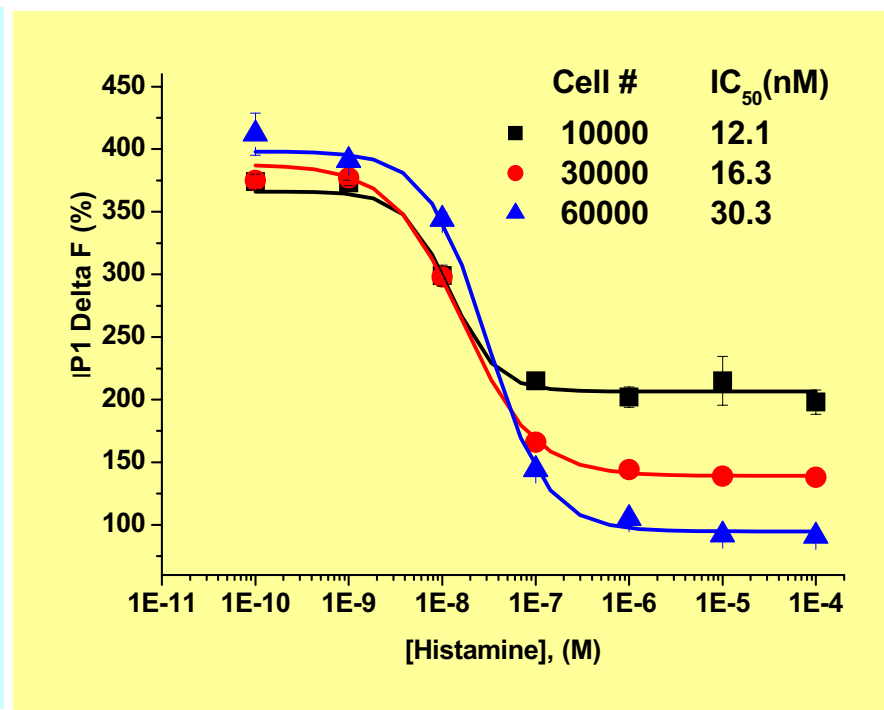
IP1 accumulation: Histamine responses in CHOK1-H1R

Novartis Institutes for BioMedical Research

Calcium FLIPR



IP1 HTRF

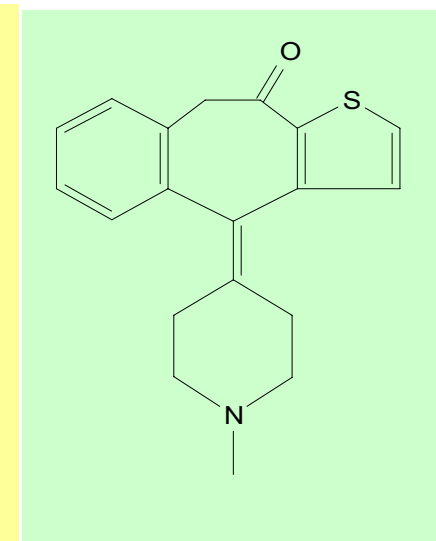
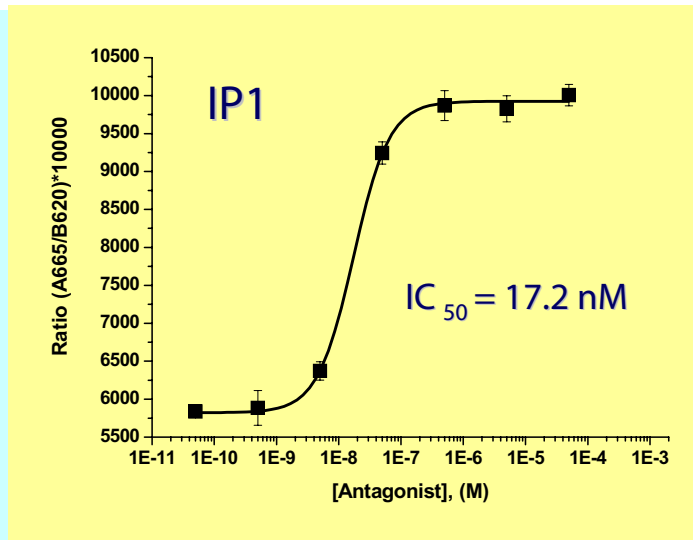
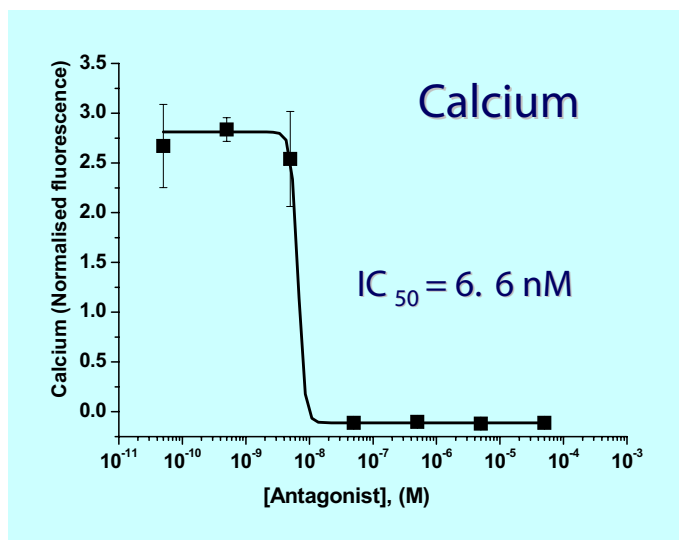


IP1 accumulation: Histamine H1 receptor in CHOK1 cells: antagonist effects

Novartis Institutes for BioMedical Research

Compound	IC ₅₀ (nM) Ca ⁺⁺ (FLIPR)			IC ₅₀ (nM) IP1 (HTRF)			IC _{50, Ca++} / IC _{50 IP1}
	exp 1	exp 2	mean	exp 1)	exp 2	mean	
Cetirizine	925	1450	1188	86	167	127	9.38
Ketotifen	6.4	6.6	6.5	5.2	17.2	11	0.58
Astemizole	460	458	459	375	865	620	0.74
Loratadine	3910	2350	3130	635	939	787	3.98
Clemastine	48.5	24.4	36	15.7	46.3	31	1.18
Doxepin	13.0	33.6	23	7.5	13.4	10	2.23
Mirtazapine	12.4	7.9	10	7.0	12.4	10	1.05

Antagonism by Ketotifen



IP1 accumulation for Secondary Screening & HTS

Novartis Institutes for BioMedical Research

- Alternative format for G_q-coupled receptors evaluated recently (see review June 06)
- Homogeneous format (HTRF)
- Applied for secondary screening of GPR40

← FLIP R / Ca⁺⁺ → ← HTRF / IP1 →

	CHOK1-C4-GPR40 Ca ⁺⁺		CHOK1-C4-GPR40 Ca ⁺⁺		CHOK1-C4-V1a Ca ⁺⁺		CHOK1-C4-GPR40 IP1	
Agonist	C18:2		synthetic agonist		Arg-vasopressine		NVP-BJX437	
Compound	IC ₅₀ (μM)	Max. inh. (% of control)	IC ₅₀ (μM)	Max. inh. (% of control)	IC ₅₀ (μM)	Max. inh. (% of control)	IC ₅₀ (μM)	Max. inh. (% of control)
RBAS1	>30	20	>30	20	>30	0	>30	31
RBAS2	>30	0	>30	0	>30	0	>30	0
RBAS3	3.6	100	3.0	100	>30	0	5.6	116
RBAS4	6.8	100	2.6	100	>30	0	7.2	109
RBAS5	2.1	60	4.8	62	6	32	>30	0

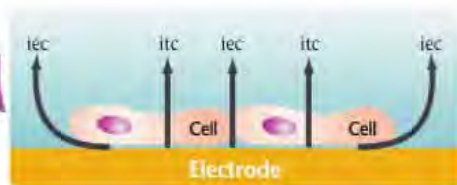
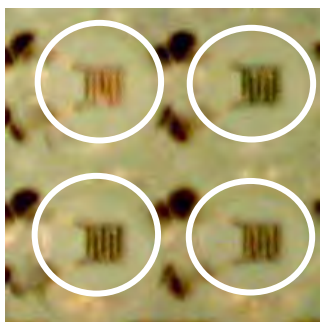
- Good correlation obtained between Ca⁺⁺ and IP1 data
- “Frequent hitters” from FLIPR screens readout can be excluded

CellKey™ System: Assay Principle

Novartis Institutes for BioMedical Research



- **Impedance based measurements**
- **Real-time, kinetic measurements across all 96 wells**
- **Automated fluid handling**
- **Simultaneous compound addition and read of all wells**
- **Monitors cell shape changes following receptor activation**
- **May constitute a phenotype assay reflecting cell migration**



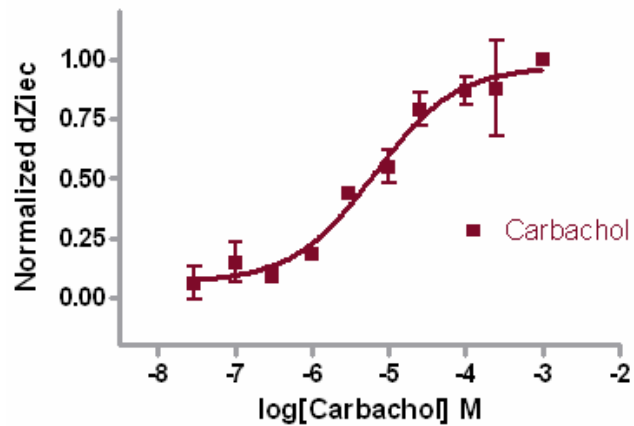
$$Z \approx V/I$$

→ Transcellular (Z_{itc})
→ Extracellular (Z_{iec})

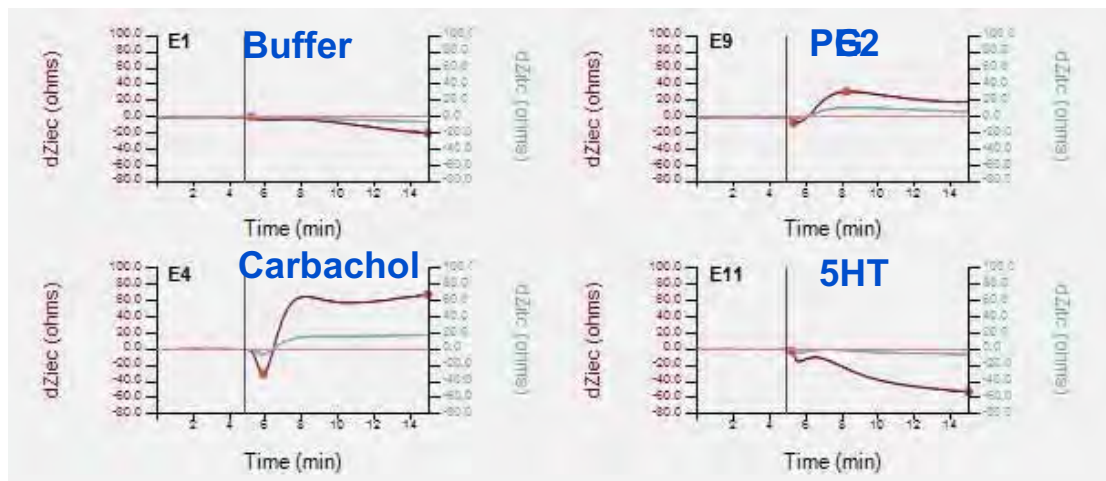
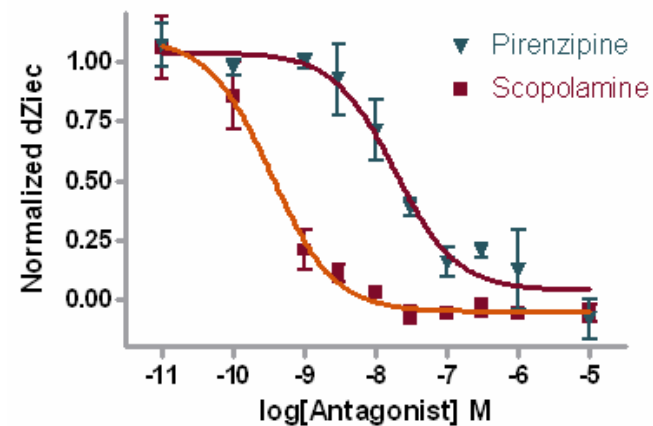
CellKey™ System: CHOm1 Muscarinic Agonist & Antagonist Effects

Novartis Institutes for BioMedical Research

060329 CHOm1 Agonist CRC



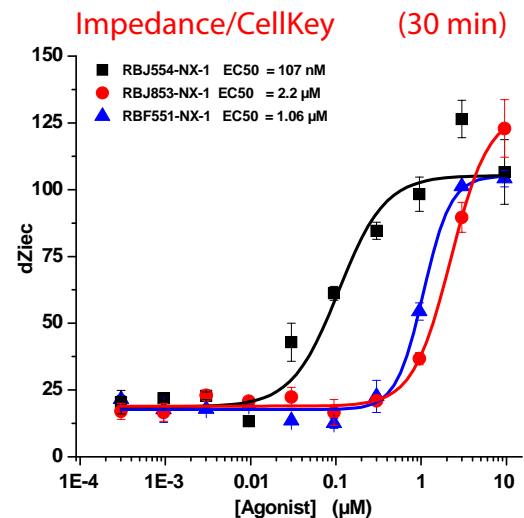
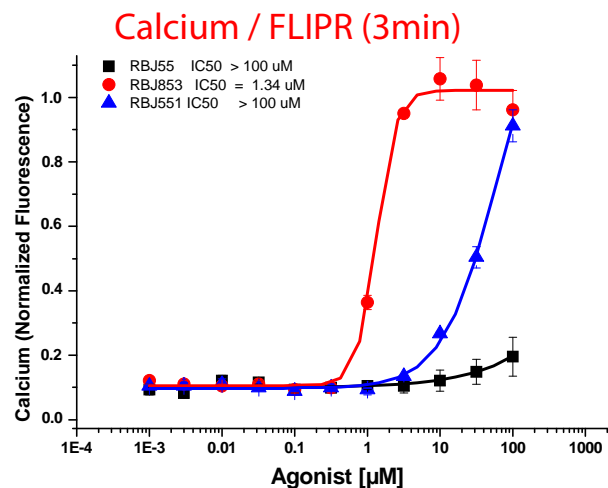
060329 CHOm1 Antagonists



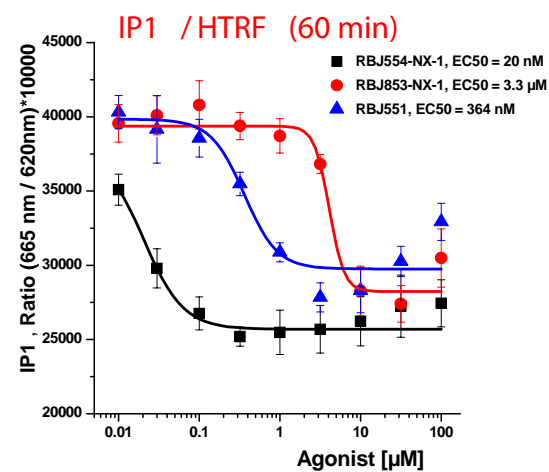
- Monitors activity and kinetic shape
- Shape determines GPCR coupling mode
- Good correlation with FLIPR data

Different potencies with three technologies

Slow acting agonists in three GPCR assay formats



Compound	Calcium (FLIPR)	IP1 (HTRF)	Impedance (CDS)
RBJ551	>100	0.36	1.1
RBJ554	>100	0.020	0.11
RBJ853	1.3	3.3	2.2



Conclusions / summary

Coupling	Ligand binding	G-protein activation	Signaling	2 _{nd} mess.	technology / Instrument	Comment
FUNCTIONAL ASSAYS						
G_q	SPA	-	PLC-?	Ca ⁺⁺	Fluo4 /FLIPR	multiplexing possible
			PLC-?	IP1	HTRF / Viewlux imaging	Temporal multiplexing
G_s	SPA	SPA GTP?S	AC activation	cAMP	HTRF / Viewlux imaging	Temporal multiplexing
G_s			AC activation	cAMP / Ca ⁺⁺ CNG2	Fluo4 /FLIPR	cAMP duplex mode possible, G _s /G _q duplex possible
G_s			PLC-?	Ca ⁺⁺	Fluo4 /FLIPR	Ca ⁺⁺ obtained via GPCR priming
G₇ 16 & chimeric				Ca ⁺⁺	Fluo4 /FLIPR	valid for ca 70 % Gi or G _s coupled receptors
G_i	SPA	SPA GTP?S	AC inhibition	cAMP	HTRF / Viewlux imaging	
G_i			PLC-?	Ca ⁺⁺	Fluo4 /FLIPR	Ca ⁺⁺ obtained via GPCR priming
G_{12/13}			RHO			
G_s, G_q, G_i, & G_{12/13}			PKC / PKA ?	role unknow	Cell key	Non-invasive free-label technology

- Novel generic assays being developed:
 - DiscoverX arrestin assays & sensigen assays