

Miniaturized Whole Cell-based GPCR cAMP Assay Using a Novel Detection System

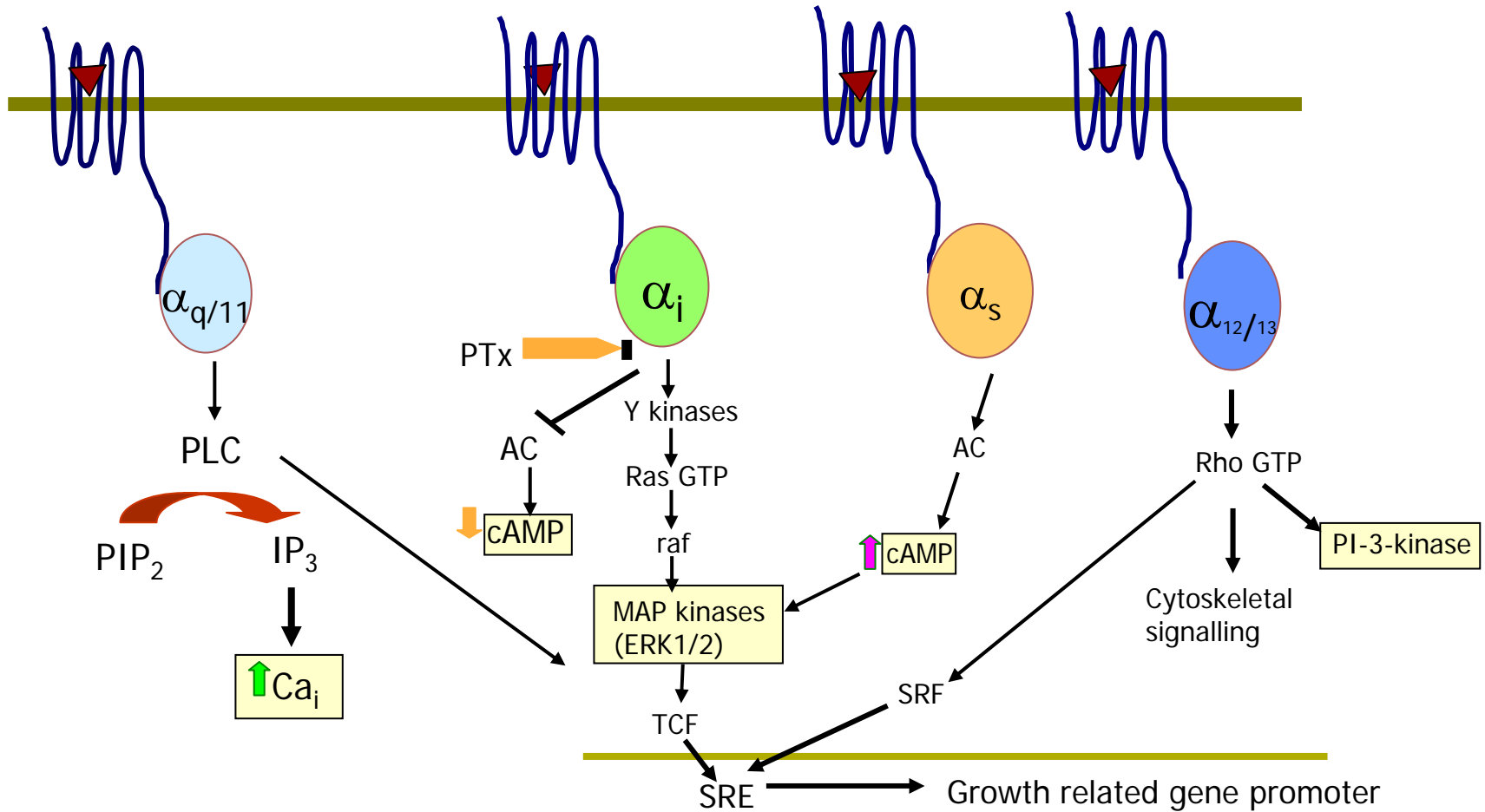
Geetha Shankar, Ph.D.
Associate Director
New Lead Discovery
Exelixis, Inc
South San Francisco, CA



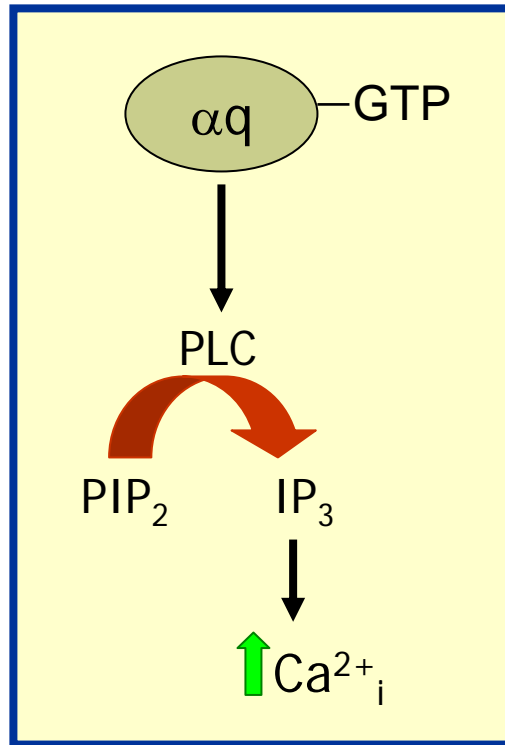
Overview

- GPCR HTS assays
- ACT:One cAMP biosensor
- Exelixis screening platform
- Case study: CB1 HTS campaign
- Summary

GPCR Signaling Pathways

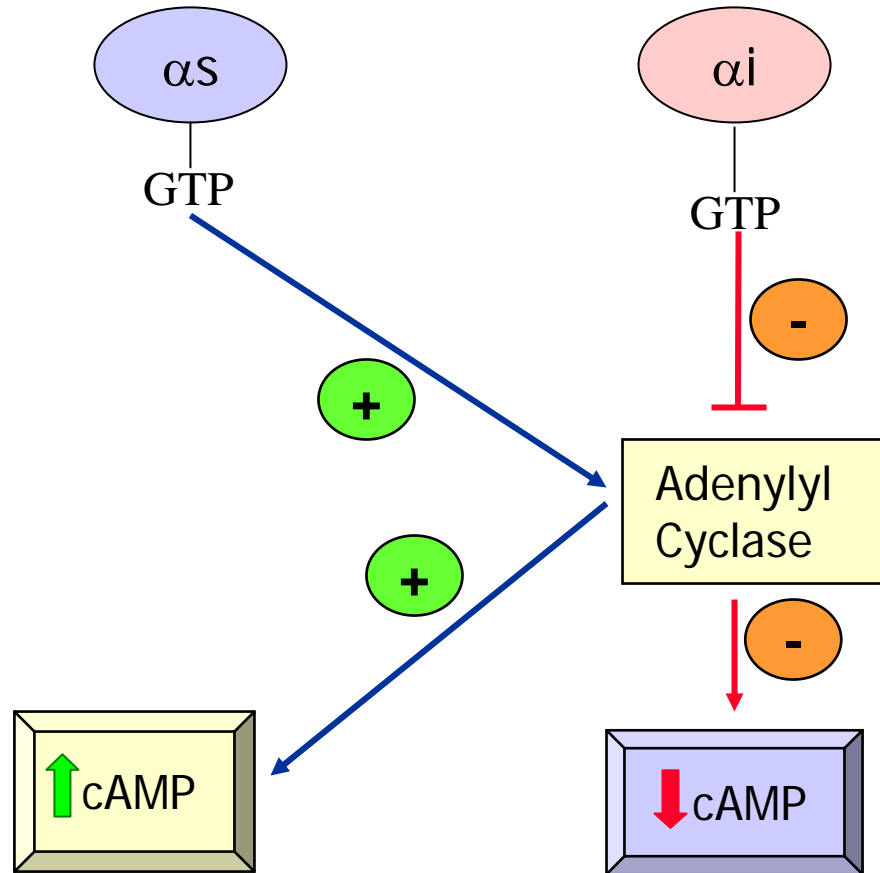


Gq and Calcium Signaling



- Fluorescence Imaging Plate Reader
- Load cells with calcium sensitive dye
- Changes in fluorescence correlate directly with changes in intracellular calcium
- 384-well format

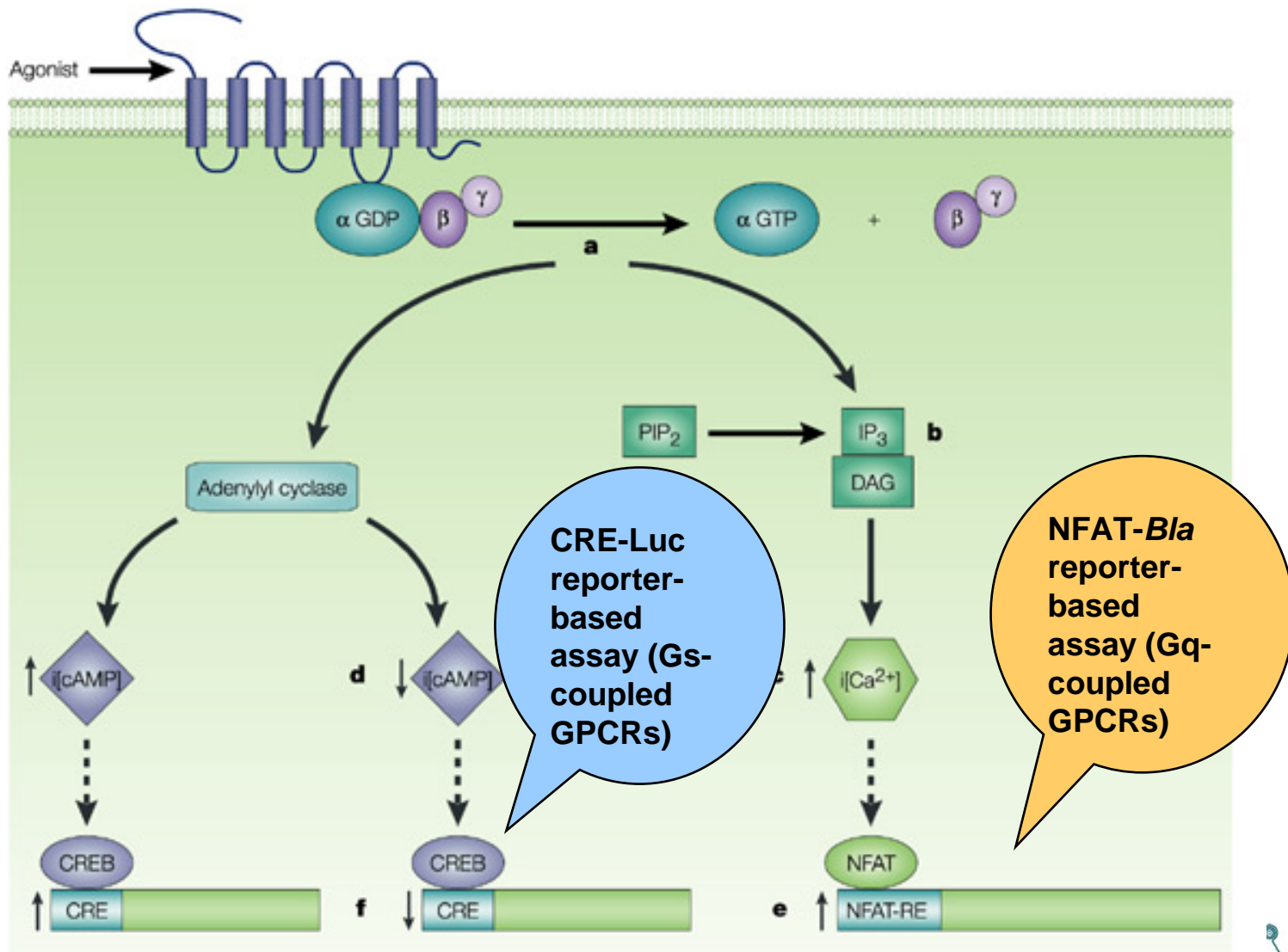
Gi and Gs-coupled receptors



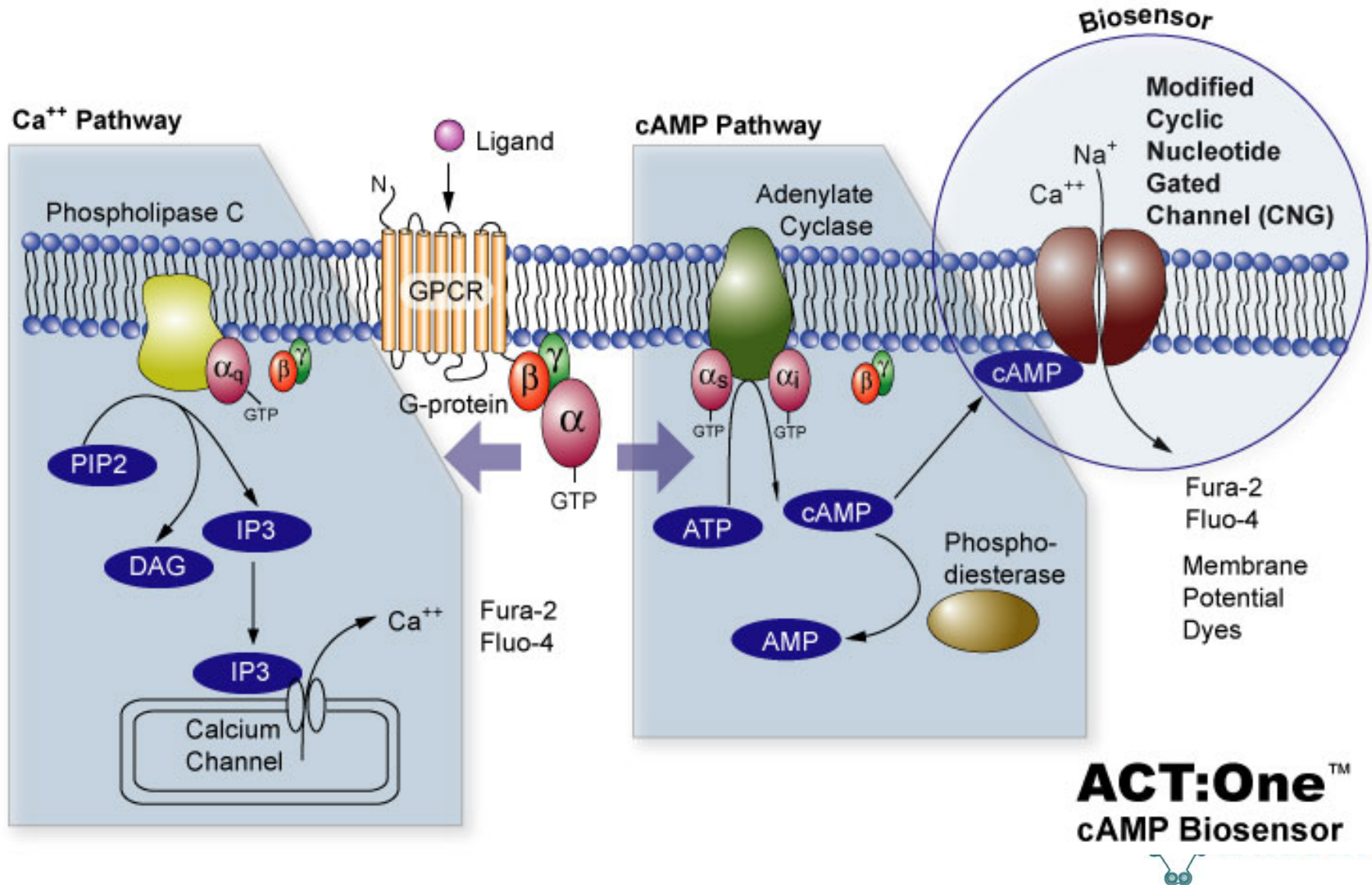
HTS strategies for non-Gq coupled receptors

- Chimeric $G\alpha$ proteins: Gqi5, Gqs, Gqx
 - ◆ Couple to $G\alpha$ as normal, but signal like Gq (via calcium)
 - ◆ FLIPR adaptable
- Promiscuous $G\alpha$ proteins
 - ◆ $G\alpha$ 15/16: Overexpress to drive signaling through calcium pathway
- cAMP assays for G_i and G_s coupled
 - ◆ Radioactive methods
 - ◆ FP
 - ◆ AlphaScreen
 - ◆ Enzyme complementation assays

Reporter-based assays for GPCRs



ACT:One Biosensor



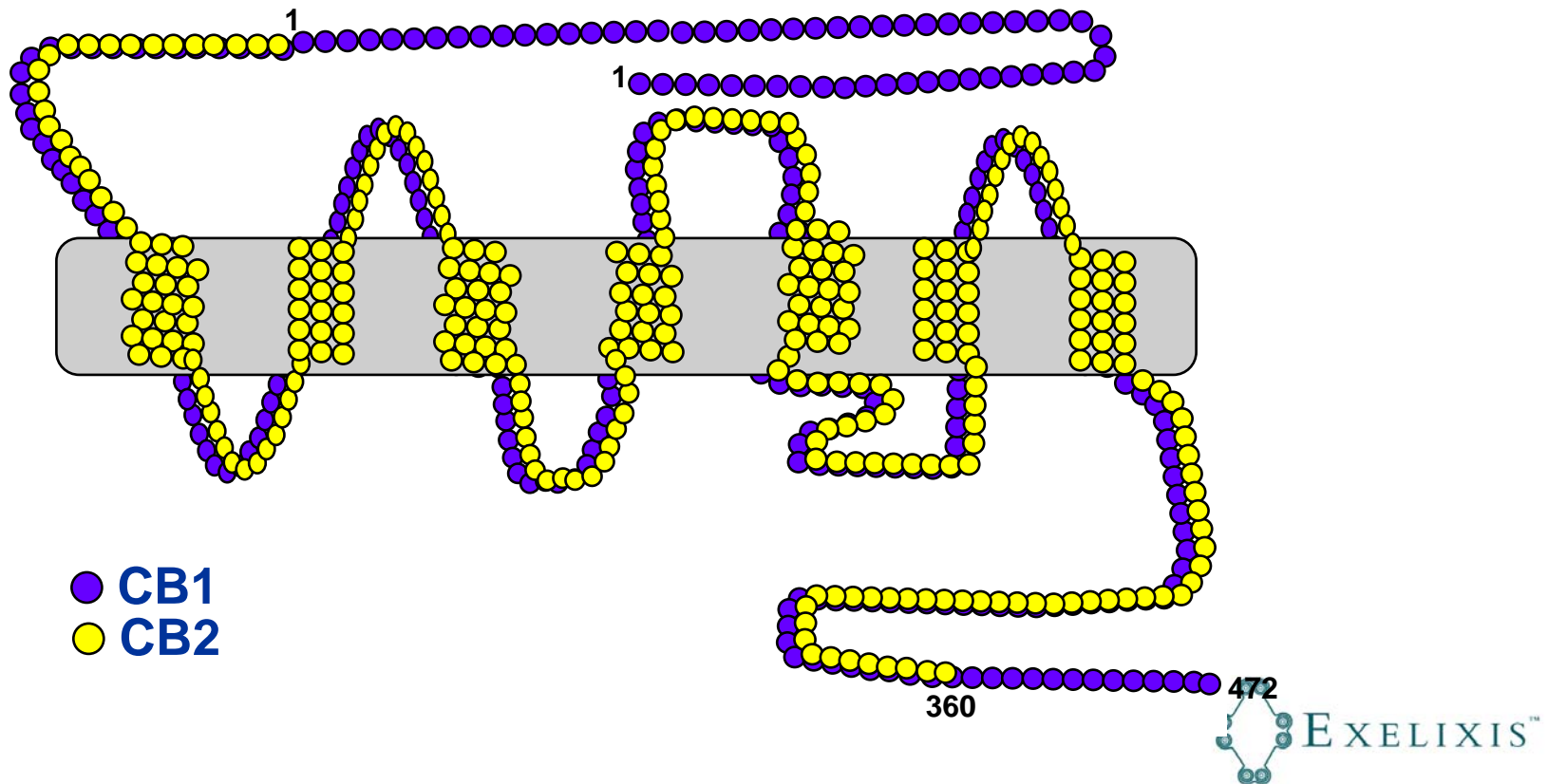
Biosensor based cAMP assay

- No modification of GPCRs required
- Endogenous and overexpressed receptors
- Gi and Gs coupled GPCRS
- Agonists and antagonists
- No radioactivity; fluorescence read-out (MP dye)
- End-point or kinetic measurements
 - ◆ Stable signal
 - ◆ Automation friendly
 - ◆ No special equipment/reader

Example: CB1 HTS campaign using CNG biosensor technology

Cannabinoid Receptor 1 Project

- Goal: Discovery of a novel, orally-bioavailable CB1 antagonist for the treatment of obesity
- Clinically-validated target – Rimonabant™
 - ◆ Phase 3 (Sanofi-Aventis)



The EC System Is a General Stress-Recovery System and Is Overall “Silent”; It Becomes Transiently Activated To:

Relax

reduction of pain and anxiety; modulation of body temperature, hormone production, smooth muscle tone, and blood pressure

Rest

inhibition of motor behavior and sedation

Forget

extinction of aversive memories

Protect

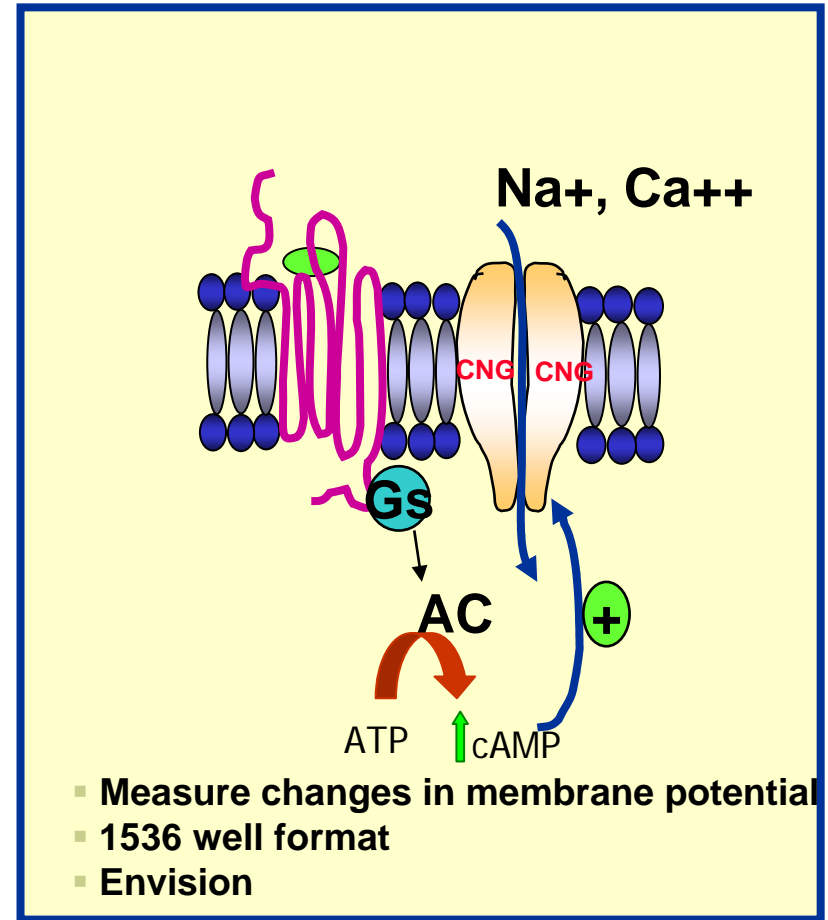
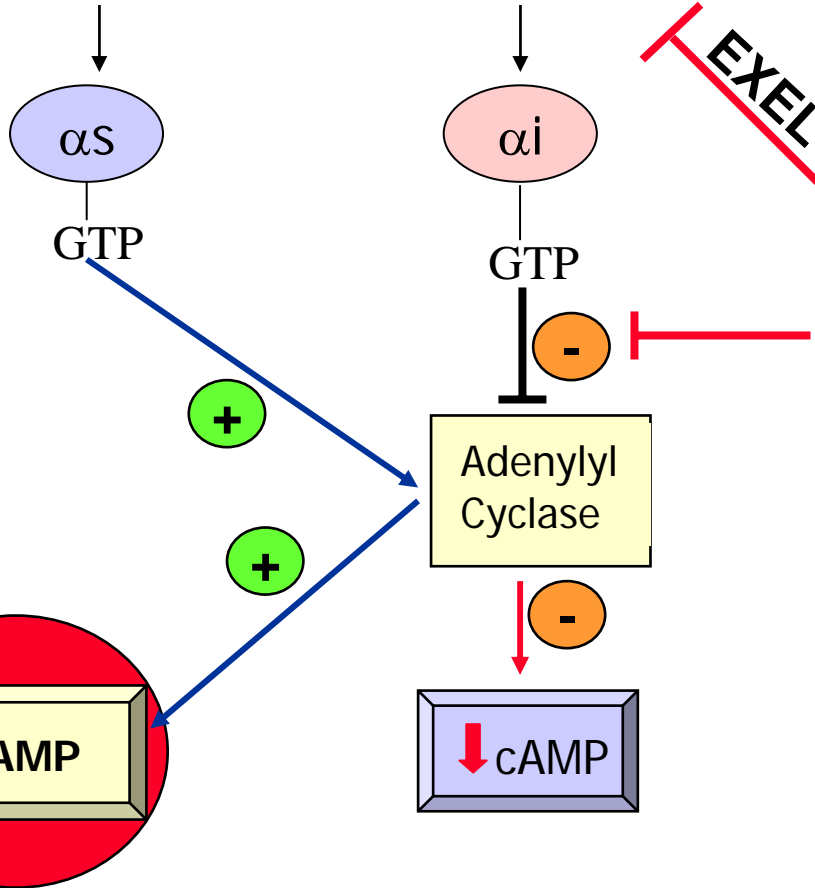
at both the cellular and emotional levels

Eat

appetite-inducing and reward-reinforcing effects

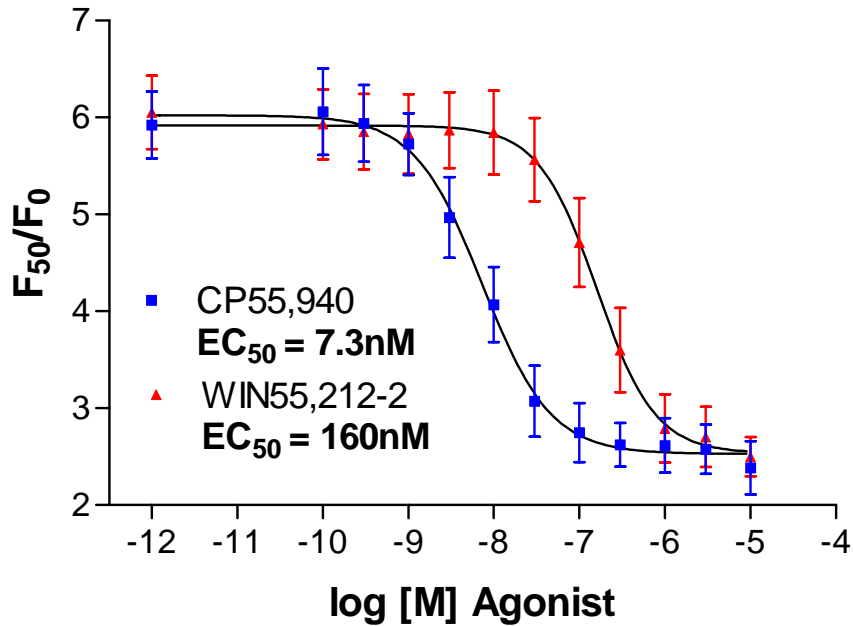
CB1 – Gi-coupled receptor

Isoproterenol CB1 (CP 55,940)

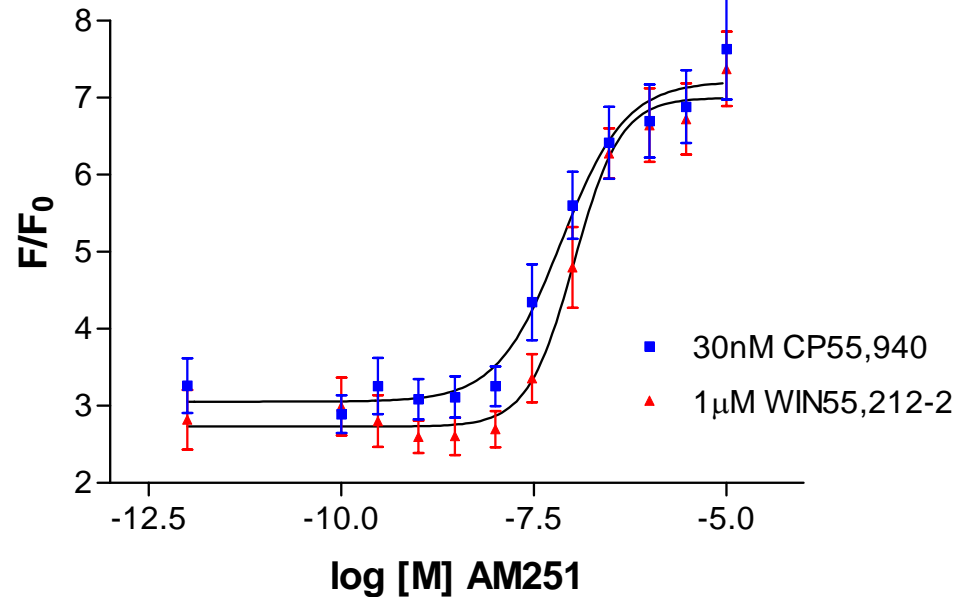


CB1 Pharmacology: Act:One cAMP Assay

1536-well format



Agonist stimulation



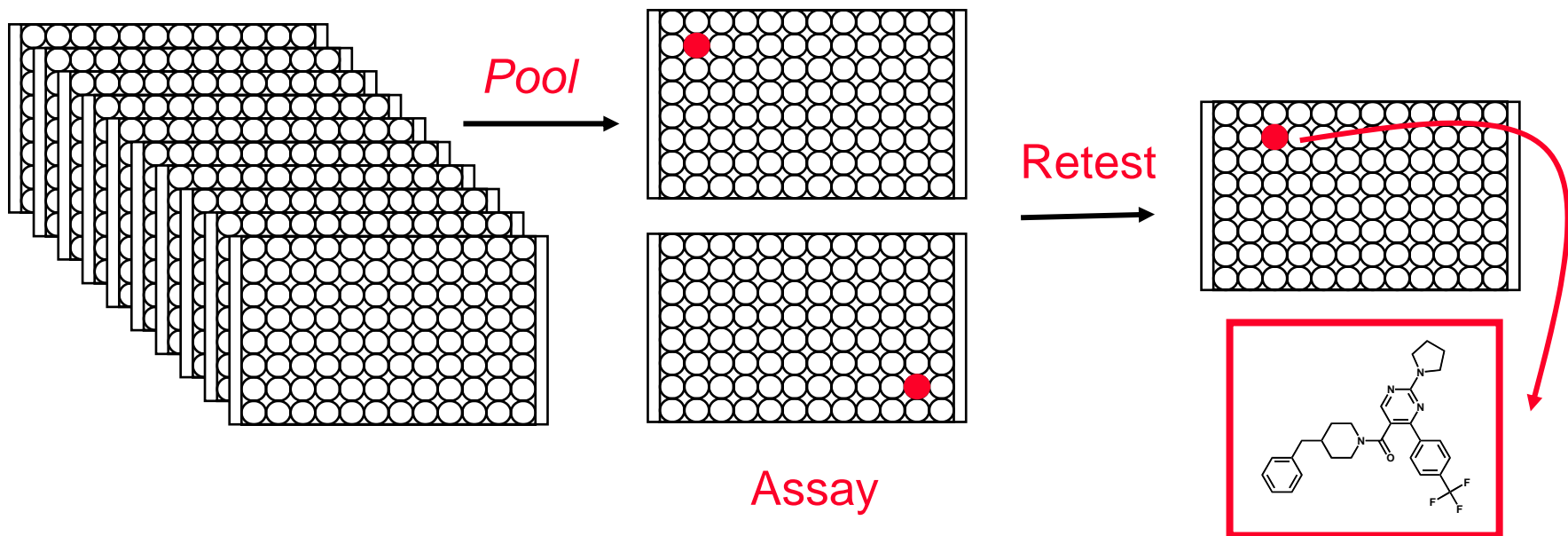
Inhibition by AM251

High Throughput Screening Platform

- HTS is not capacity limiting
 - Routine screening of 4+ MM cmpds
- Dynamic technology
 - 384/1536 microtiter plate formats
- Continued increase in library size
 - Complete HTS in 1-2 weeks



Orthogonal Compound Pooling



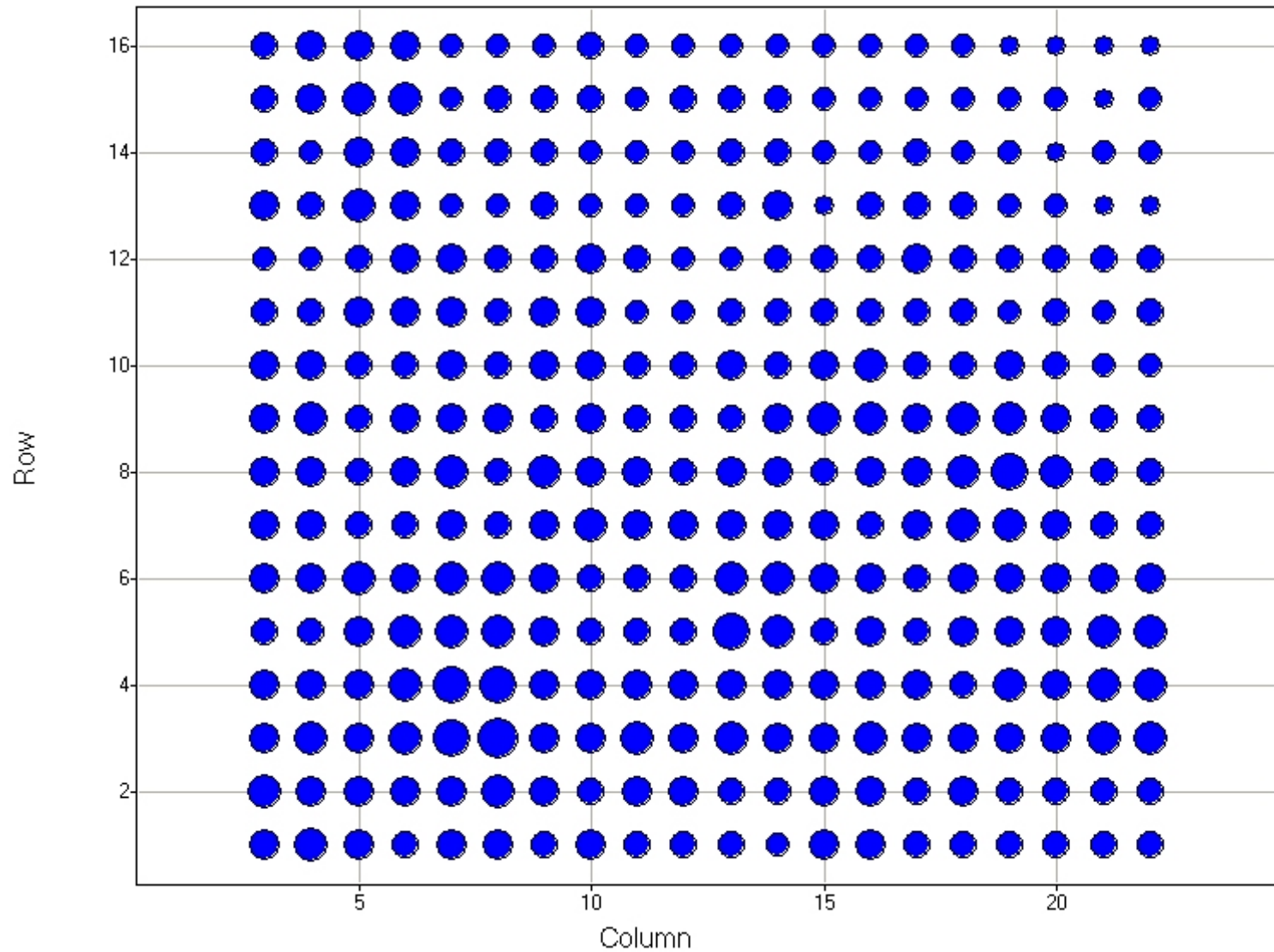
- 10 Compounds/well in two sets of mixtures
- Deconvolute replicates and confirm as single compounds
- Low false positive and false negative hit rate

Five-fold increase in HTS efficiency

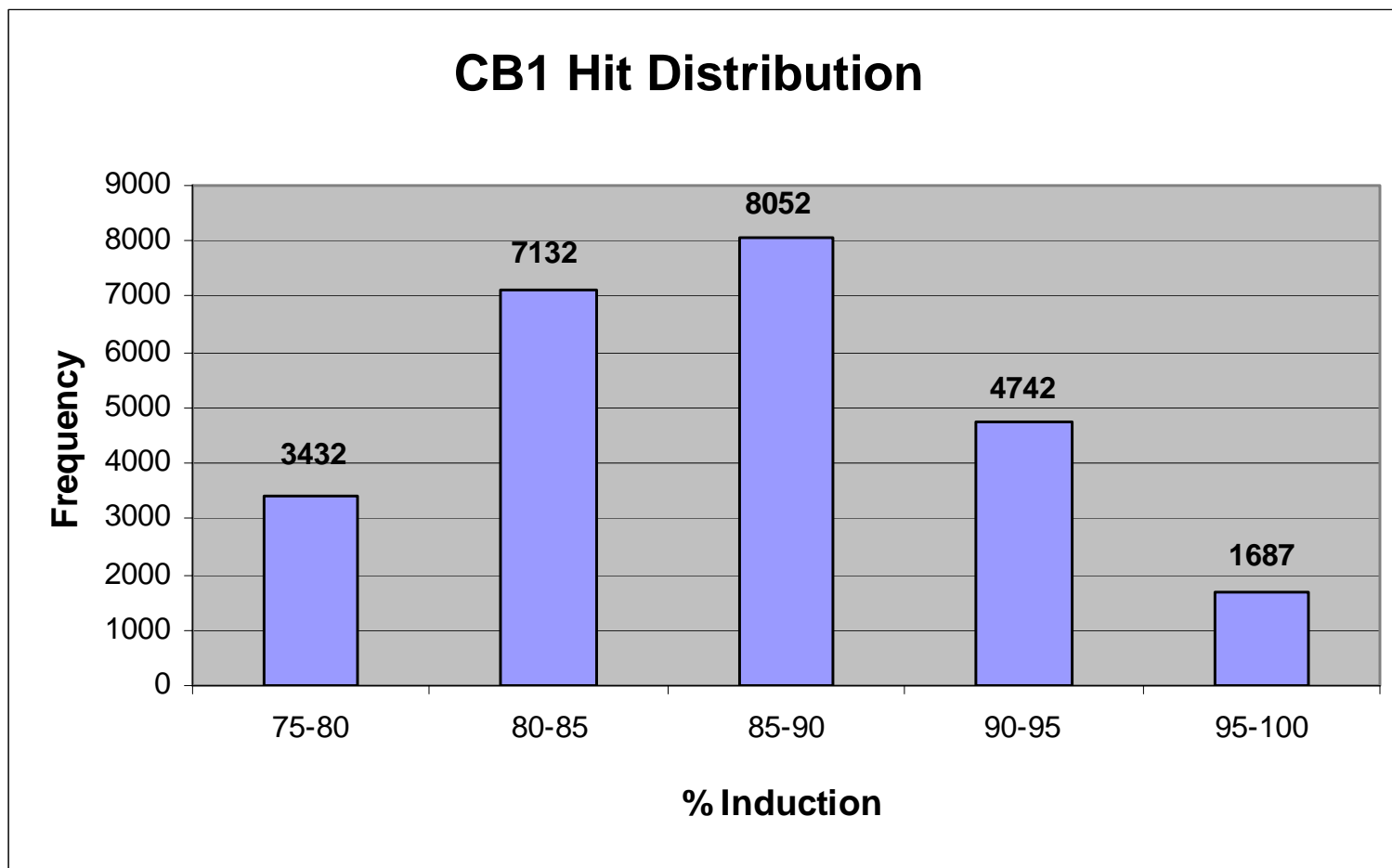
HTS Assay Performance

- 3.6 M Compounds screened at 1.7 μ M
- 1536-well format
- 4000 cells/well; 50 minute incubation; end-point fluorescence read
- Novel, robust assay format for G_i -coupled receptors
 - ◆ $Z' = 0.4 - 0.6$
 - ◆ Mean S/B = 2.2
 - ◆ # hits = >25,000 (75-100%)
- Identified multiple potent antagonists

Hit distribution



Hit distribution



Confirmation Summary

- 667 Compounds confirmed upon single point retesting
- 208 Compounds evaluated in dose-response studies
 - ◆ 4 < 10 nM, 73 <100 nM, 123 <500 nM IC₅₀ (cAMP)
- Eleven scaffolds confirmed via resynthesis
 - ◆ IC₅₀ values 1-200 nM
 - ◆ >60% derived from internal combi-chem synthesis
- PD studies used to prioritize lead series

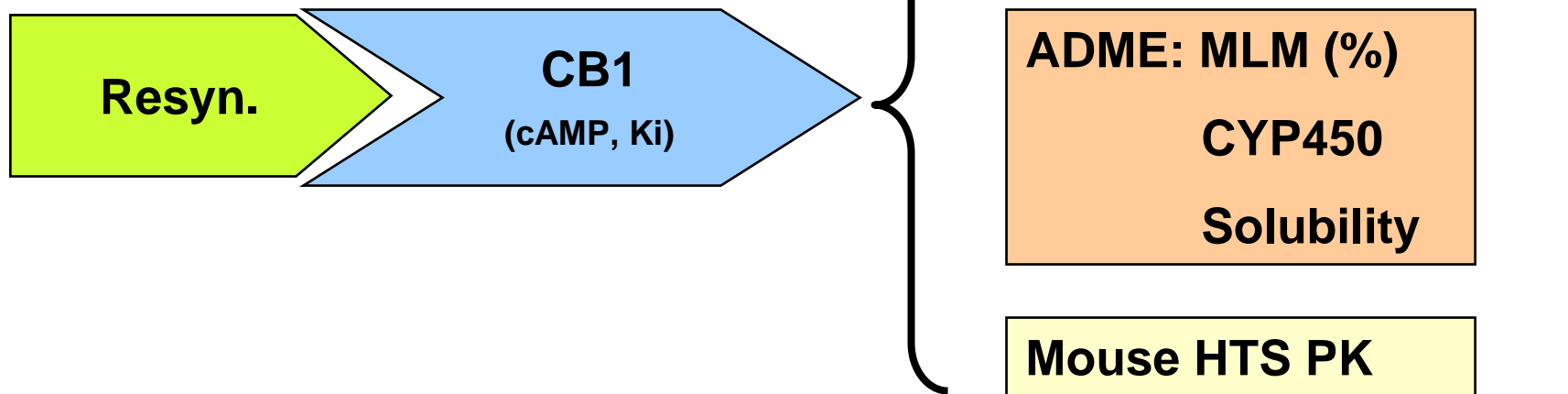
Summary of Confirmed Actives

Compound (# examples)	CB1 – cAMP (nM)	CB1 – Ki (nM)	PD (100 mpk) % reversal
Rimonabant	4	3	80 (30 mpk)
EXEL-3627 (85)	1.5		79 (300 mpk)
EXEL-9628 (17)	3.2	>100	8 (300 mpk)
EXEL-5412 (20)	3.8	-	-
EXEL-4933 (62)	4	50	36
EXEL-3124 (45)	8	-	-
EXEL-2355 (52)	13		31 (300 mpk)
EXEL-7594 (1)	23	76	3
EXEL-5450 (13)	77		-40 (300 mpk)
EXEL-5649 (8)	100	-	-
EXEL-4197 (10)	107	-	-
EXEL-6921 (6)	183	-	

Lead Validation

■ Compound Progression

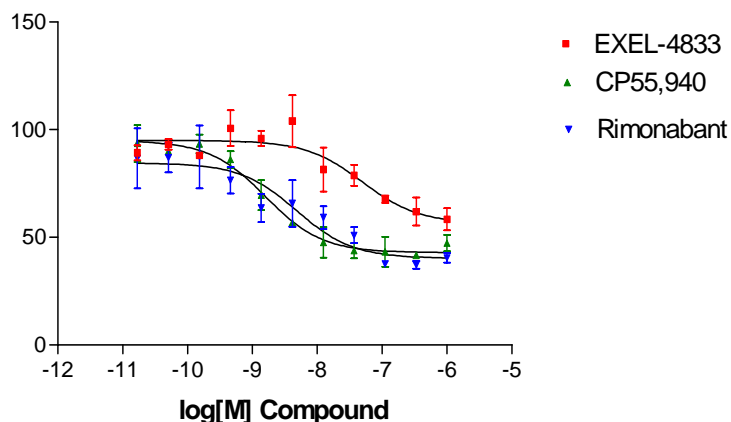
- ◆ Active resynthesized HTS compounds
 - cAMP, Displacement binding (<10 nM)
- ◆ Pharmacodynamic assay
 - Reversal of CB1-agonist induced hypothermia (p.o.)



In vitro displacement binding

SAR-dependent binding observed

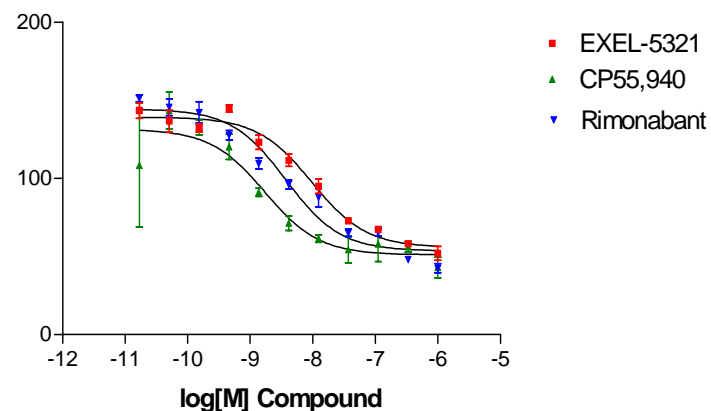
EXEL-4933 with 3H CP55940



EC₅₀: 8 nM

Ki: 50 nM

EXEL-5321 with 3H CP55940



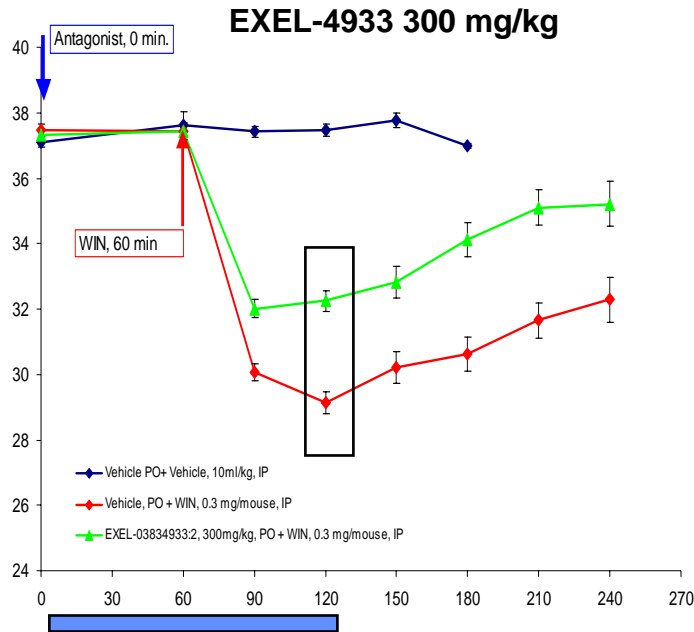
EC₅₀: 5 nM

Ki: 10 nM

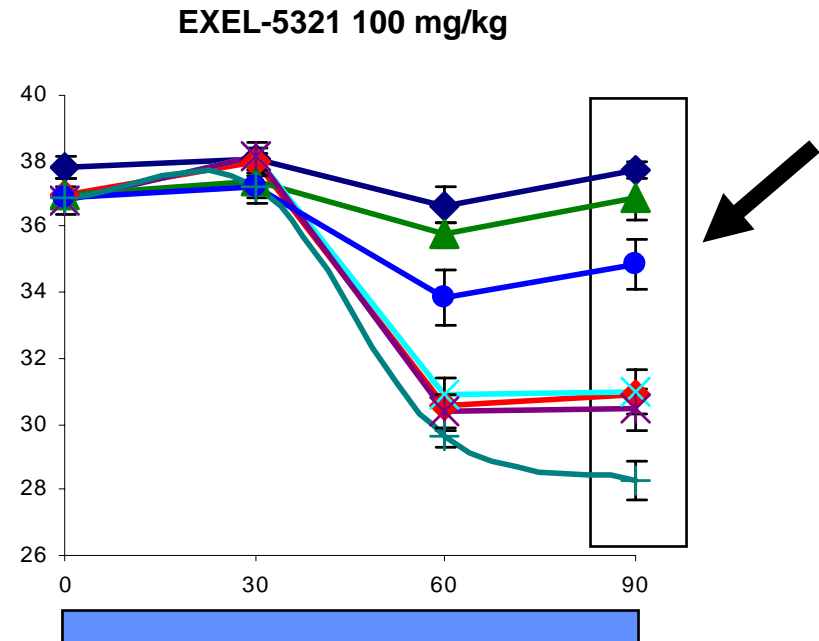
Correlation with in vivo efficacy

EXEL-4933 mediated prevention of hypothermia

- Validated model mediated by the cannabinoid system



37% inhibition of hypothermia



68% inhibition of hypothermia

Summary

- CNG channel technology offers a novel detection method for cAMP measurements
- Technology has wide applicability to Gi and Gs coupled GPCRs
- Significant cost-savings with miniaturization
- Stable, robust signal allows for detection of potent hits

Acknowledgments

- GPCR

- ◆ Scott Ogus
- ◆ John Lesnick
- ◆ Nicole Moore

- HTS

- ◆ Shaun Nguyen

- Kirk McMillan, VP New Lead Discovery