

# Case Study: HTS of **GPR-54** with IP-One Assay

Marcie Glicksman, PhD

Senior Director of Leads Discovery

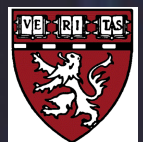
Laboratory for Drug Discovery in Neurodegeneration  
Brigham and Women's Hospital & Harvard Medical School

Assays and Cellular Targets

October 31, 2006



*Harvard Center for Neurodegeneration and Repair*



# Outline

---

- What is LDDN?
- GPR54 Background
- Why IP-One
- HTS protocol
- Summary



# Laboratory for Drug Discovery in Neurodegeneration



*Mission of the LDDN is to ...*

**Create a new model for drug discovery that integrates the best of industry and academics.**



**Discover chemical agents that can be used as lead structures in the development of drugs to treat neurodegenerative diseases.**



*Laboratory for Drug  
Discovery in  
Neurodegeneration*

# Laboratory for Drug Discovery in Neurodegeneration

## *Features of the LDDN model ...*

- Hypothesis-driven, screening-based approach.
- Managed by industry-seasoned professionals.
- Programs based on tight academic collaborations.
- Resourced for success – Commercialization of disease modifying therapeutics.
- Cover a broad range of targets in neurodegeneration.
- Forward-expansion – New collaborators and new disease areas.



---

# GPR54: Case Study for LDDN



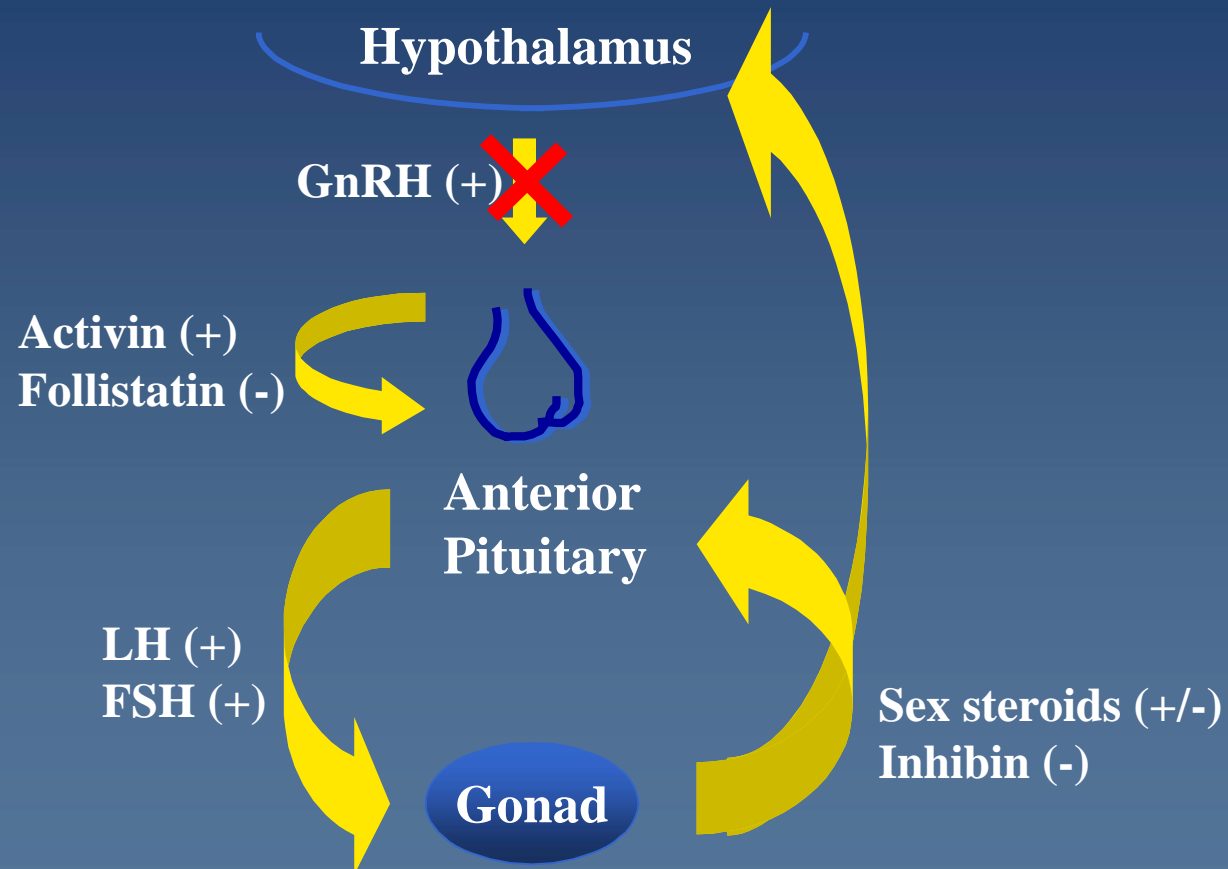
*Laboratory for Drug  
Discovery in  
Neurodegeneration*

# The Hypothalamic-Pituitary-Gonadal Axis

Aspartate, Leptin,

Dopamine, Glutamate, (+)  (-) GABA, Opioids

NE, NPY, **Kisspeptin**



# AIM

---

To identify GPR54 agonists, antagonists, and enhancers for therapeutic and research uses.

## **Agonists:**

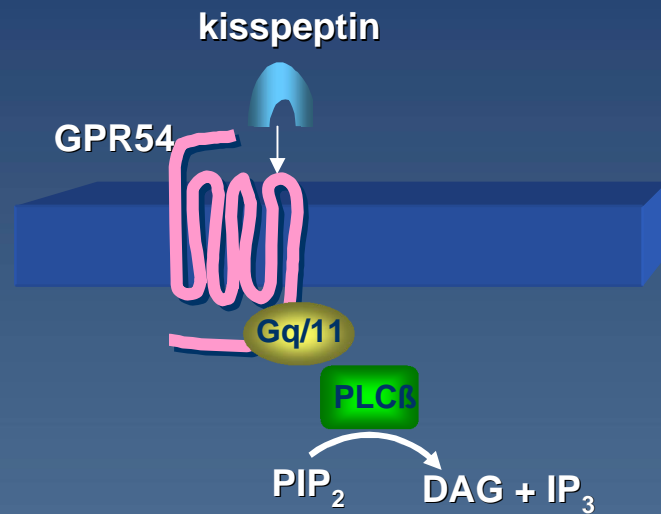
delayed puberty, infertility

## **Antagonists:**

hormone-dependent cancers, precocious puberty

# GPR54, A G Protein-Coupled Receptor

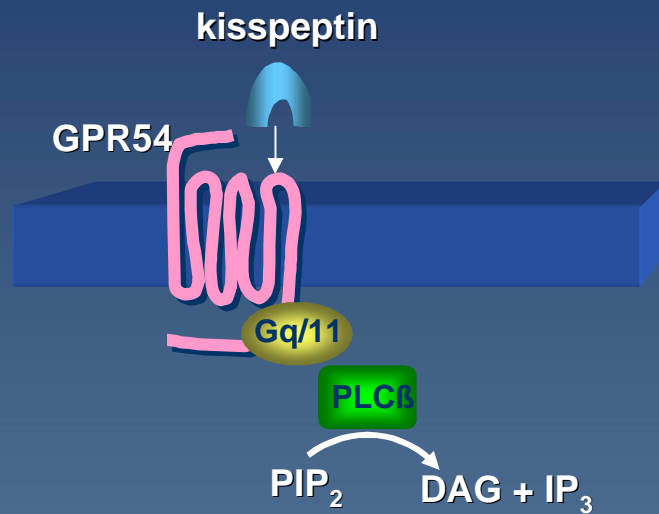
- GPR54 is expressed primarily in brain, pituitary, and placenta.



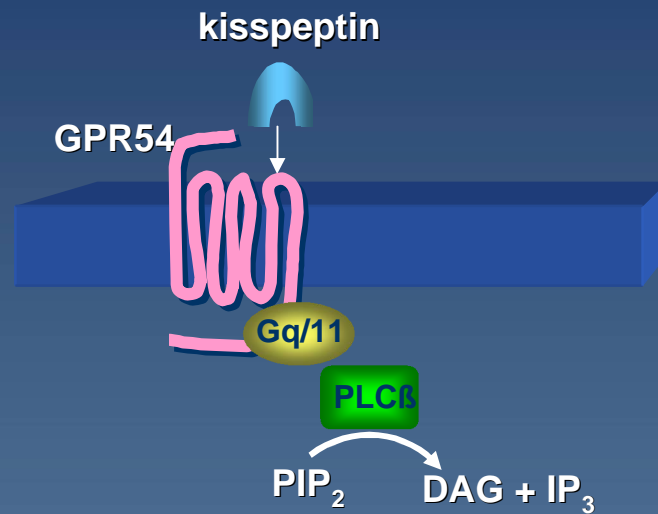


# GPR54, A G Protein-Coupled Receptor

- GPR54 is expressed primarily in brain, pituitary, and placenta.
- Natural ligand of GPR54 = kisspeptin, encoded by gene KiSS-1



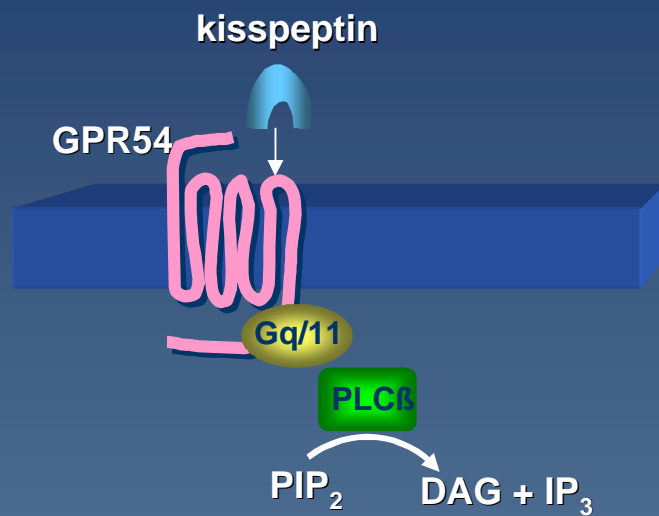
# GPR54, A G Protein-Coupled Receptor



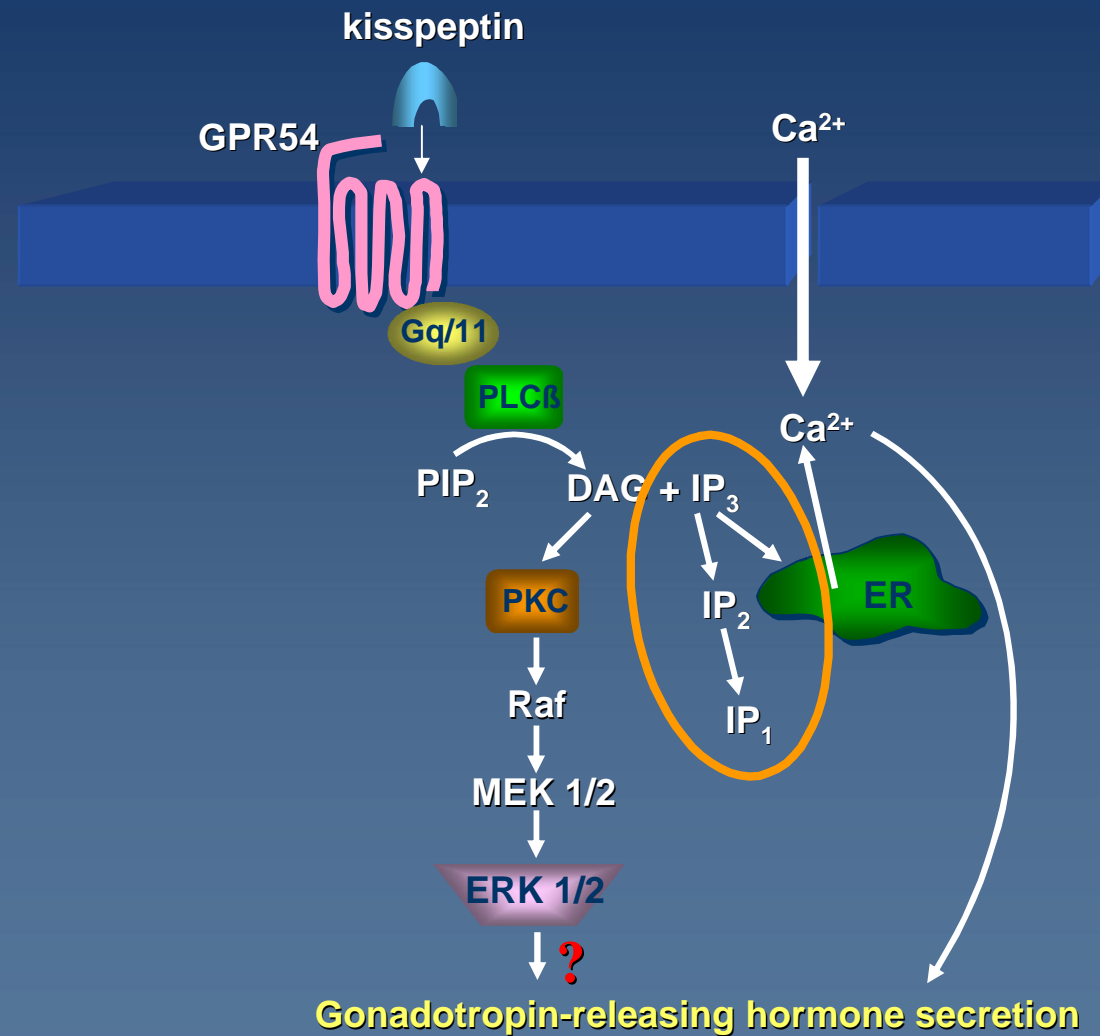
- GPR54 is expressed primarily in brain, pituitary, and placenta.
- Natural ligand of GPR54 = kisspeptin, encoded by gene KiSS-1
- KiSS-1 is expressed in hypothalamus and placenta.

# GPR54, A G Protein-Coupled Receptor

- GPR54 is expressed primarily in brain, pituitary, and placenta.
- Natural ligand of GPR54 = kisspeptin, encoded by gene KiSS-1
- KiSS-1 is expressed in hypothalamus and placenta.
- Kisspeptin-54 was named metastin for its ability to inhibit tumor metastasis.
- GPR54 knockouts do not undergo puberty and antibodies will block LH surge

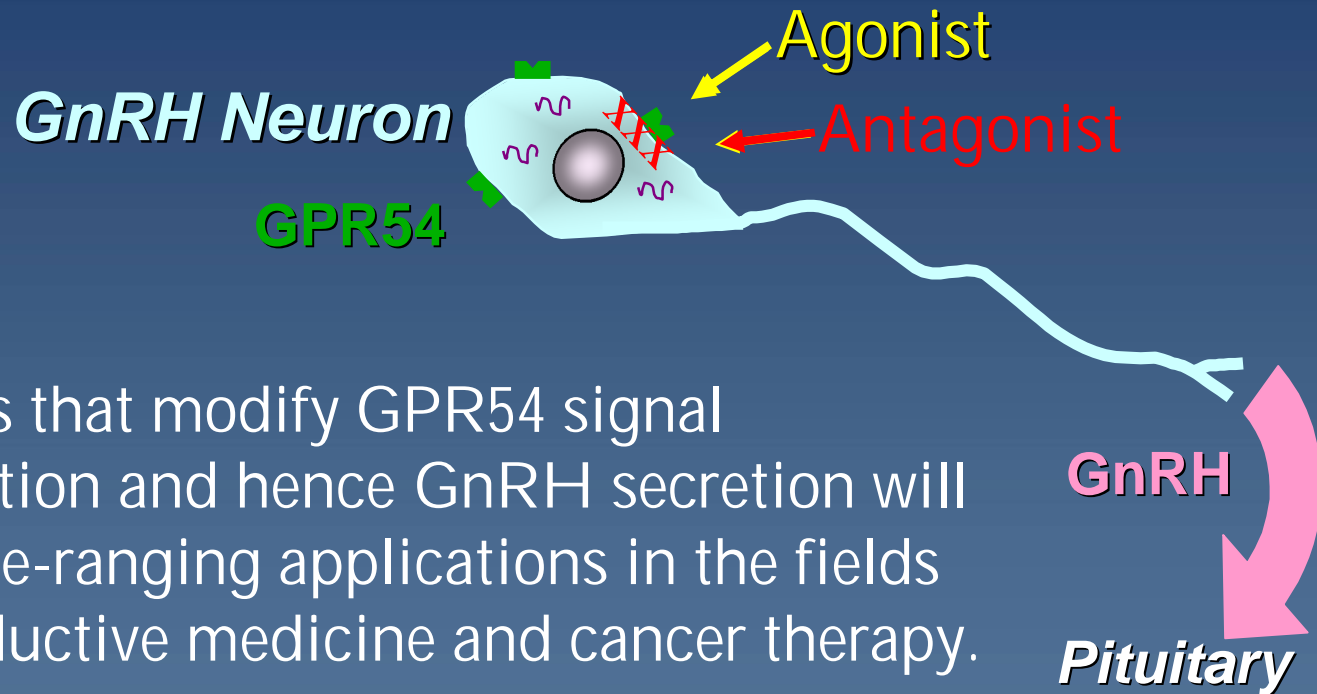


# GPR54-Coupled Signal Transduction Pathways



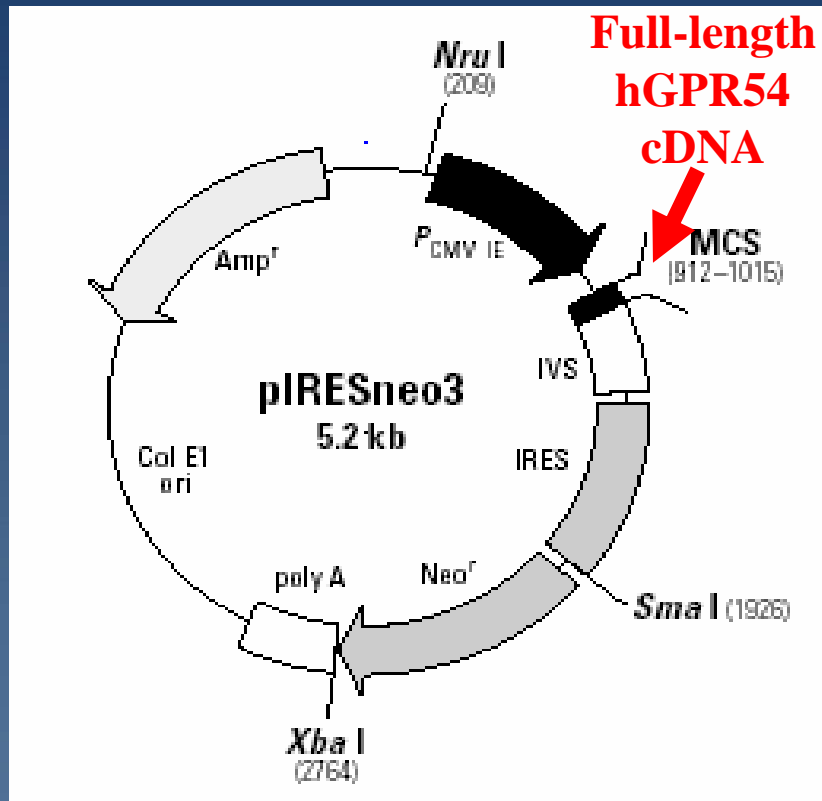
# HYPOTHESIS

Agonists and enhancers of GPR54 will promote GnRH secretion, while GPR54 antagonists will suppress GnRH secretion.



Reagents that modify GPR54 signal transduction and hence GnRH secretion will have wide-ranging applications in the fields of reproductive medicine and cancer therapy.

# Generation of Stably Transfected GPR54 Cell Lines



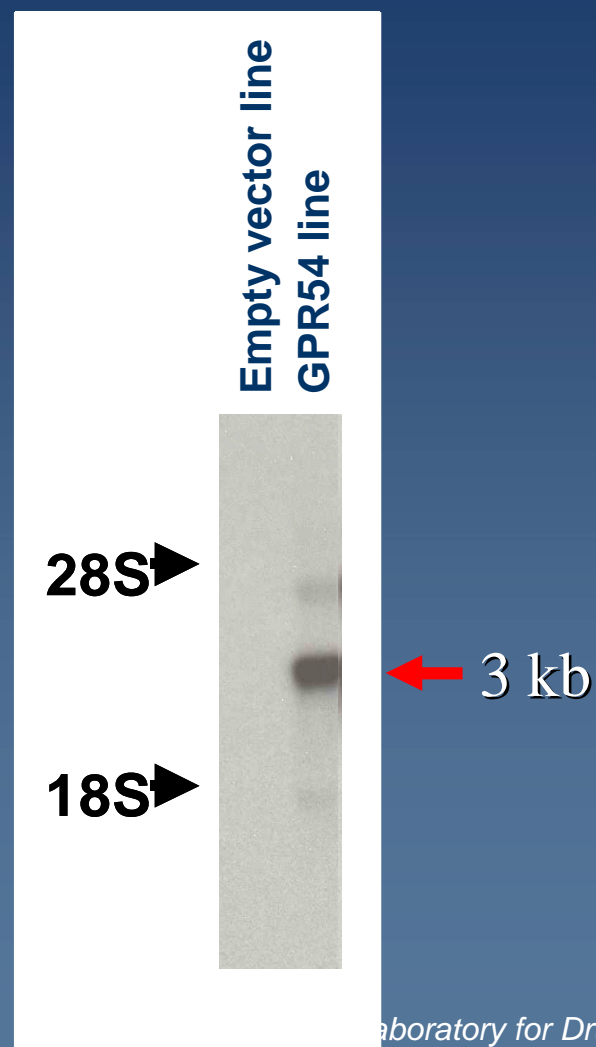
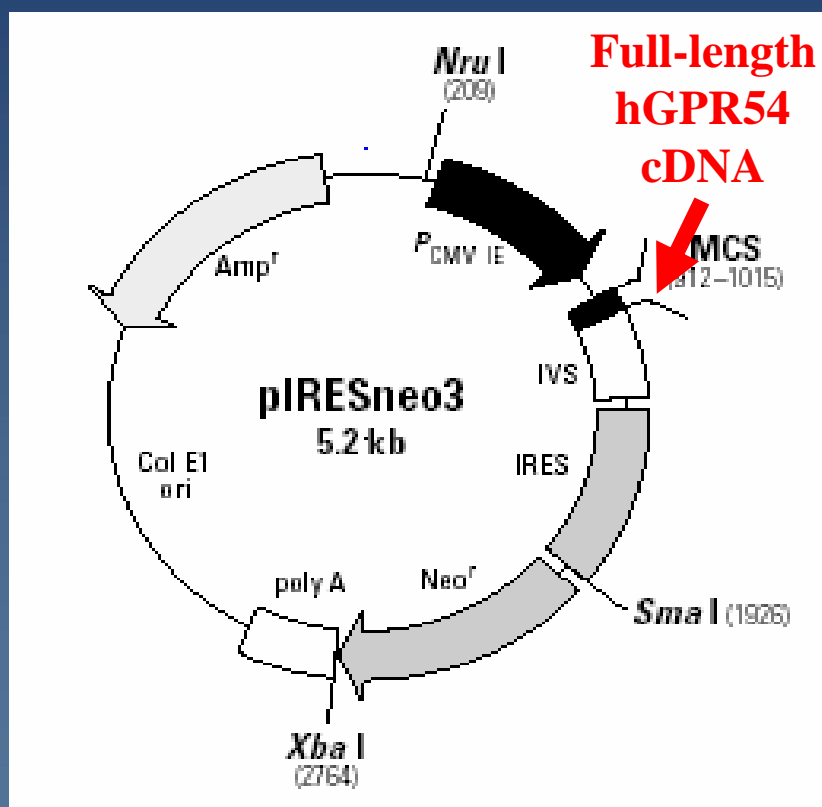
## ➤ Chinese hamster ovary (CHO) cell line

- Used in other published GPR54 stable cell lines
- Growth pattern simplifies clone selection

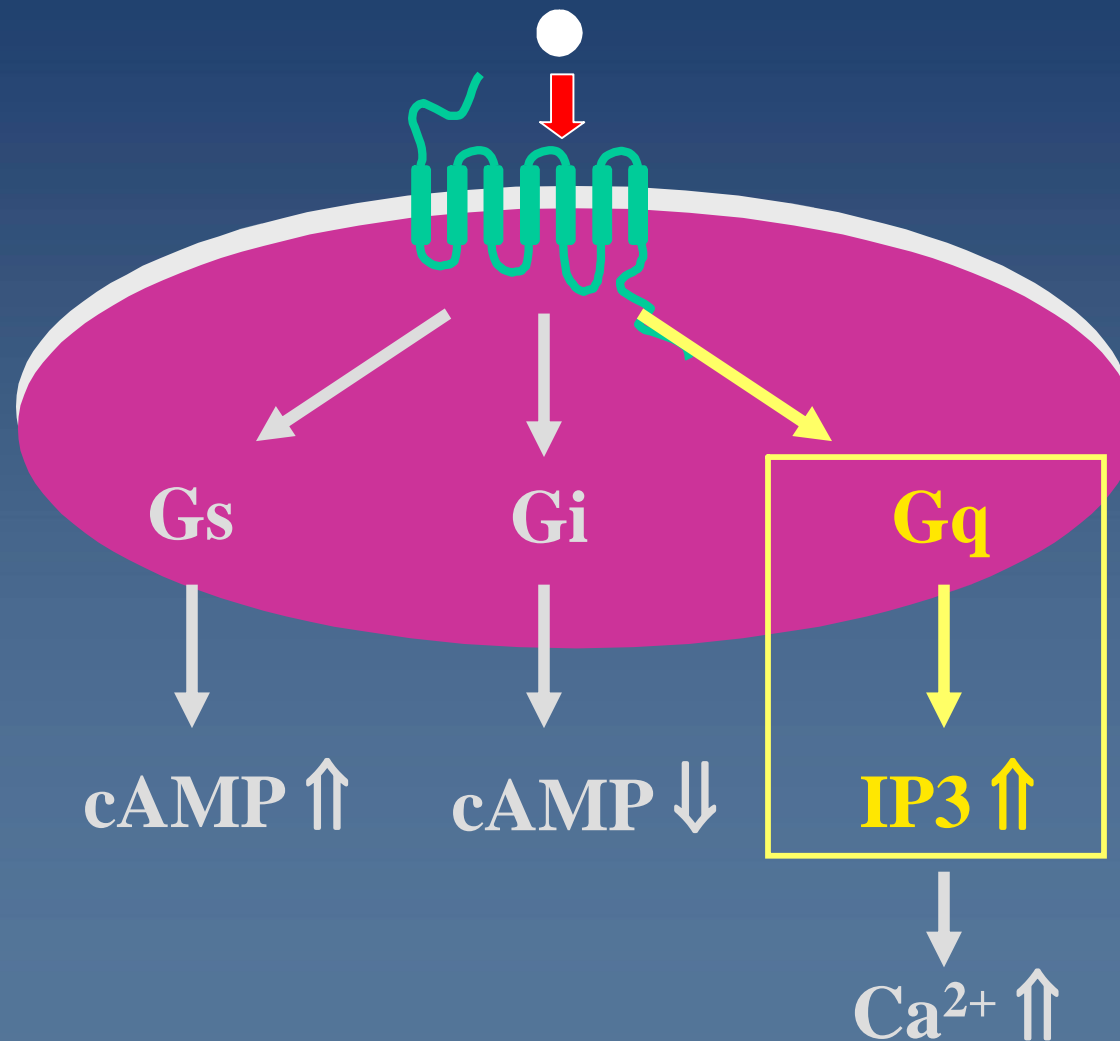
## ➤ Construct backbone – pIRESneo3 (Clontech)

- Bicistronic vector – gene of interest and selection marker are under control of same promoter, favoring selection of clones with higher expression of transgene

# Generation of Stably Transfected GPR54 Cell Lines

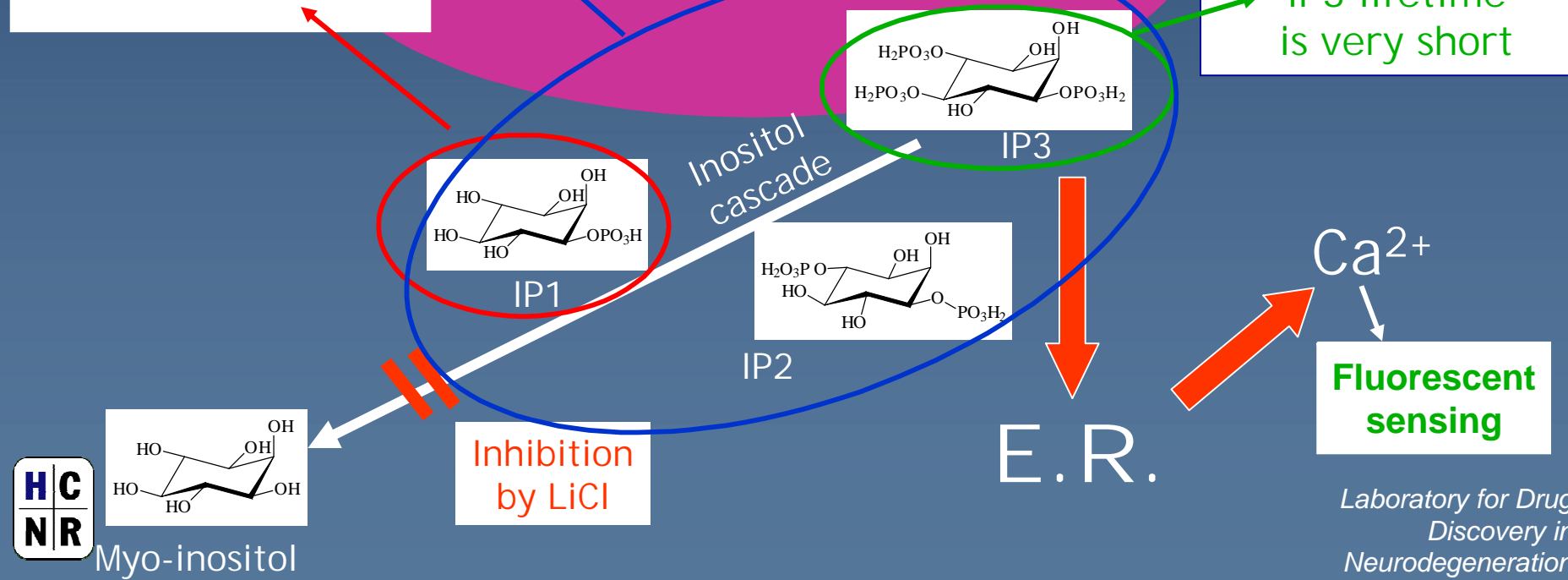
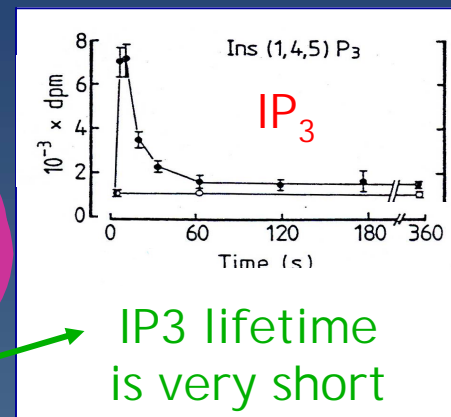
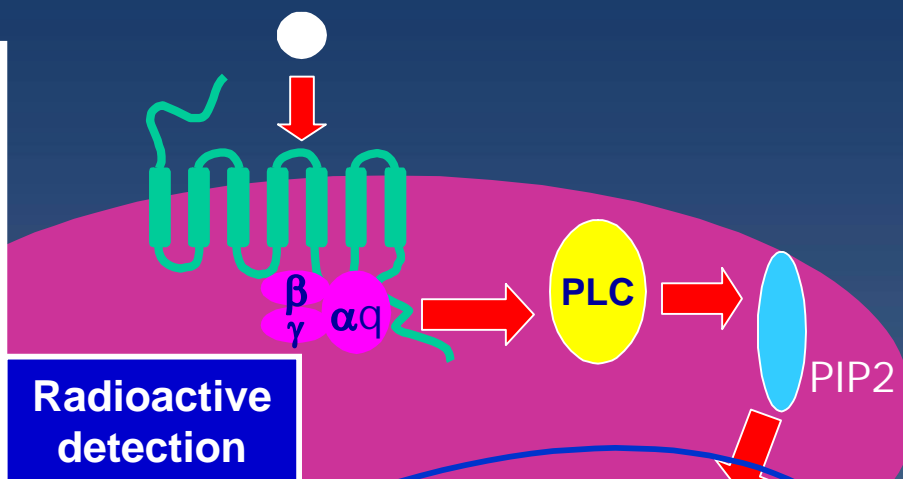
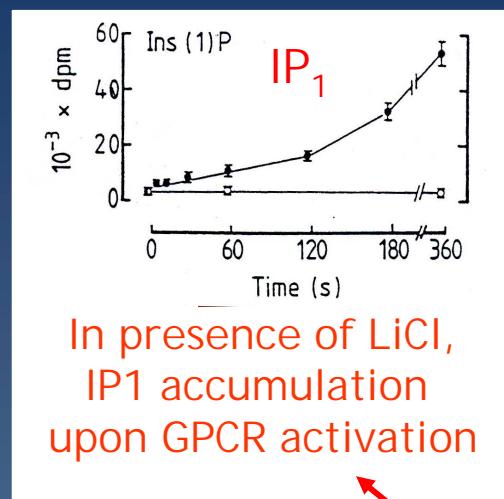


# GPR54 is a Gq coupled GPCR





# Gq coupled GPCR signalling pathway



# Why Choose IP-One for HTS

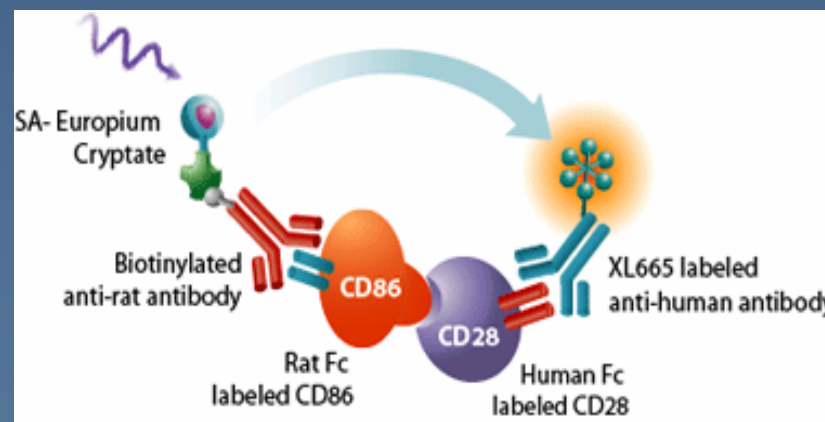
---

- Did not have access to a FLIPR for measuring  $\text{Ca}^{++}$  flux
- Miniaturization to 384-well format to reduce reagent consumption
- Minimal number of steps to facilitate automation and maximize speed and efficiency
- Non-radioactive detection for throughput, safety, and waste disposal considerations
- Must have satisfactory sensitivity, accuracy, and reproducibility (Z factor)



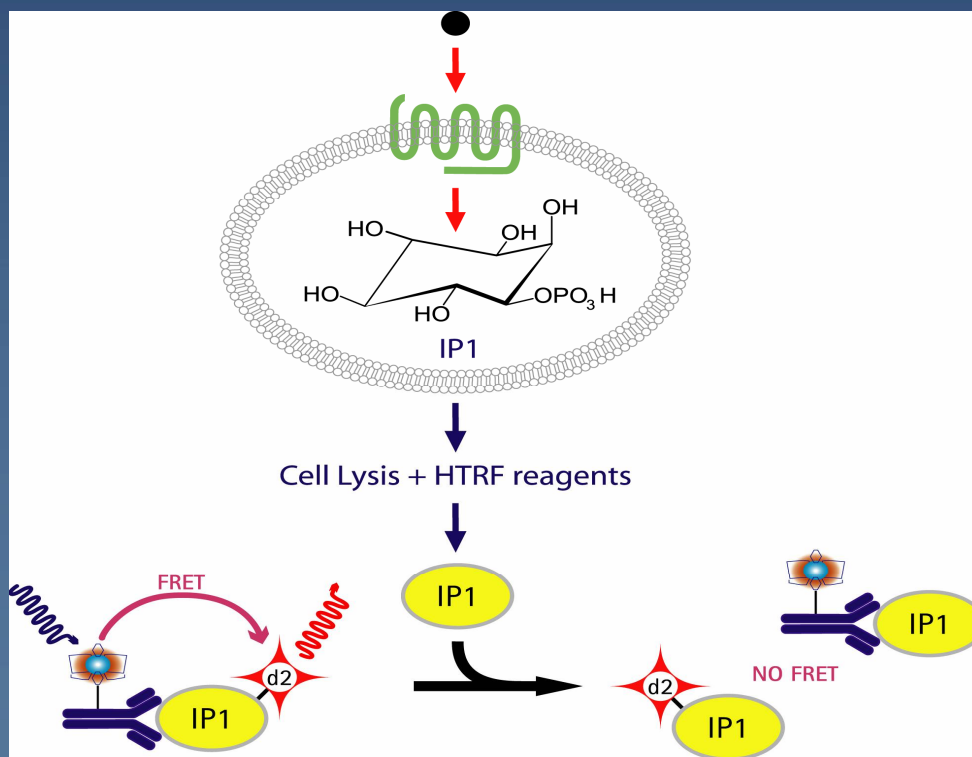
# Homogeneous Time Resolved Fluorescence

- HTRF® is a technology based on TR-FRET, a combination of FRET chemistry and the use of fluorophores with long emission half-lives.
- HTRF uses lanthanide with an extremely long half-life (Europium), conjugation of  $\text{Eu}^{3+}$  to cryptate, an entity which confers increased assay stability
- Use of a ratiometric measurement that allows correction for quenching and sample interferences.
  - Simplified assay miniaturization
  - Tolerant of additives e.g. DMSO & EDTA
  - Cell-based functional assay



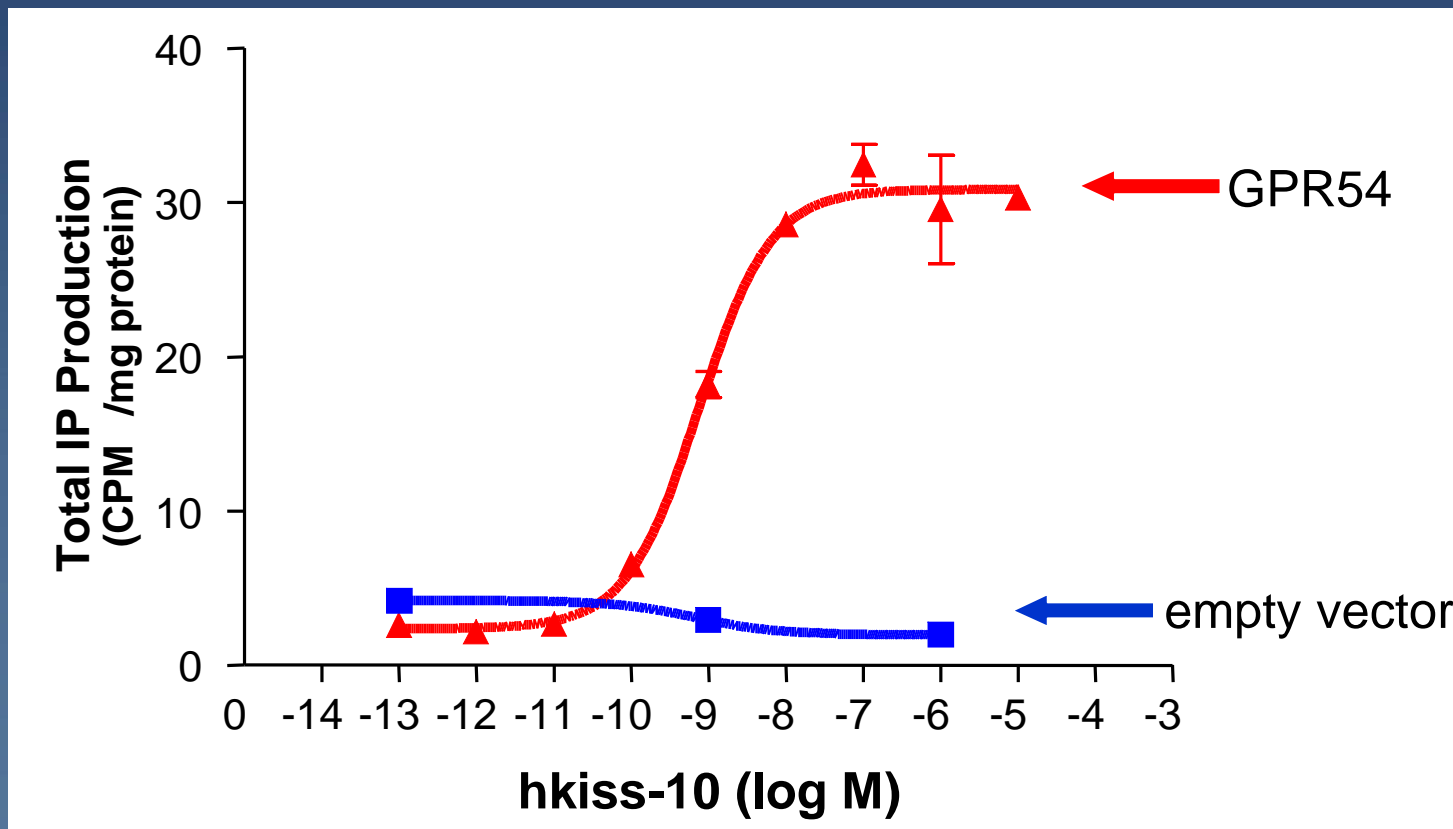
# The IP-One assay

- Competitive immunoassay with cryptate-labeled anti-IP1 monoclonal antibody and d2-labeled IP1.
- LiCl is used causing the accumulation of IP1 upon receptor activation.
- The assay can be run in a single microplate and requires only a single 1 hour incubation following cell stimulation.
- No cross-reactivity with 50 $\mu$ M of (phospho) inositides phosphates



# Kisspeptin Stimulation

- Kisspeptin Stimulates IP Accumulation In a Dose-Dependent Manner in GPR54 Stably Transfected Cells

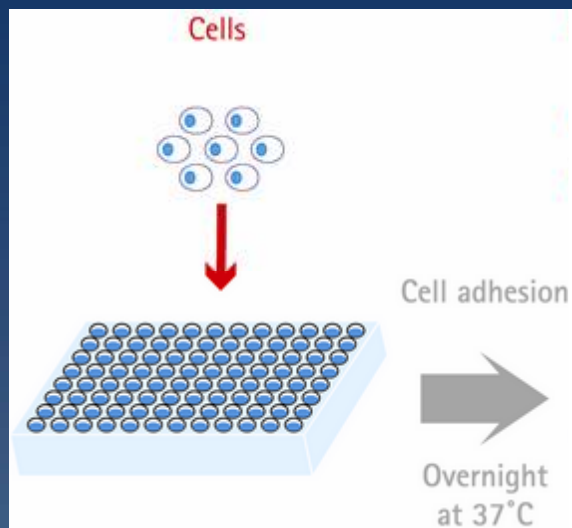


# IP-One assay miniaturization

---

Plate type	96	96 <sup>1/2</sup>	384	384 sv
Cell number	80,000	40,000	15,000 to 30,000	8,000
Total volume (μl)	100	50	20 to 40	10

# IP-One assay protocol

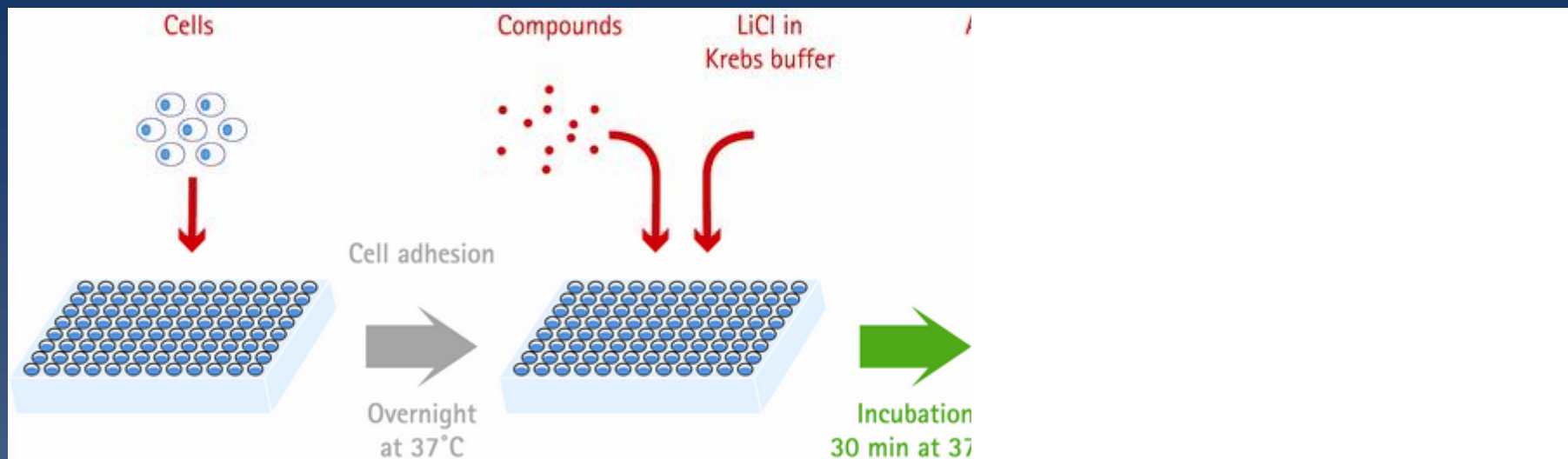


30,000 cells/40ul/well in  
white Nunc 384-plates



Seal with aeroSeals

# IP-One assay protocol



30,000 cells/40ul/well in  
white Nunc 384-plates



Seal with aeroSeals

Remove media



Add compound together  
with stimulation buffer



(add Kisspeptin for  
antagonist screen)

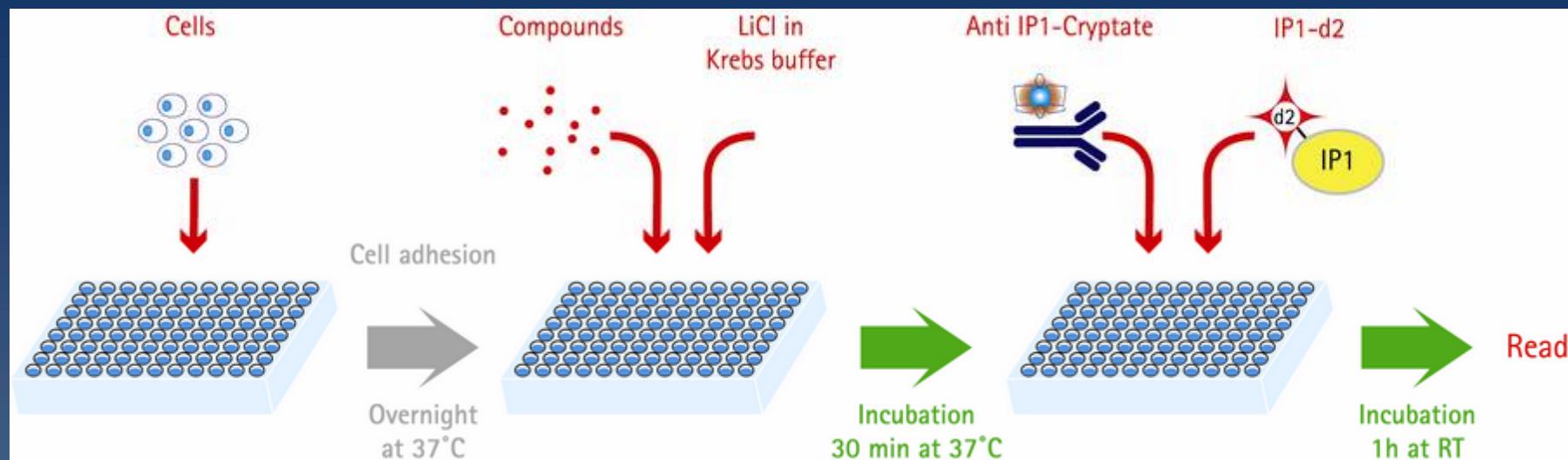


Seal with aeroSeals





# IP-One assay protocol



30,000 cells/40ul/well in  
white Nunc 384-plates



Seal with aeroSeals

Remove media



Add compound together  
with stimulation buffer



(add Kisspeptin for  
antagonist screen)



Seal with aeroSeals

Add IP-One kit D2



Add IP-One Cryptate  
reagent



Seal with aluminum seals

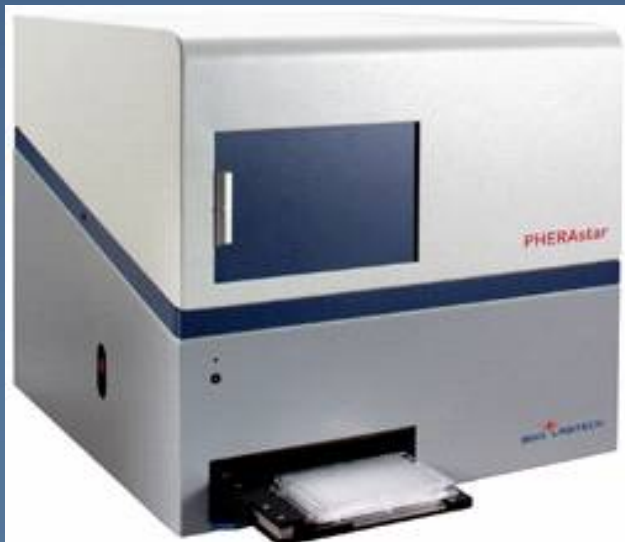


1 h-24 h



Laboratory for Drug  
Discovery in  
Neurodegeneration

# Robotics Used for the Screen



- Beckman FX core system
- BMG PHERAstar or PerkinElmer Envision
- High energy xenon flashlamp excitation
- Simultaneous dual wavelength measurement for HTRF
- Integrated into robotic system



# Compound Library (110,000 compounds)

➤ Computational filters applied to select compounds with an increased probability of oral bio-availability and blood brain barrier penetration

3000 Peakdale

7000 Bionet

5000 CEREP

16,000 Maybridge

50,000 ChemDiv

6000 Enamine

6000 IF Lab

- All small molecules adhere to Lipinski's rules.
- Low proportion of toxicophores
- Low proportion of unwanted functionalities.
- Maximization of molecular diversity.

1670 FDA-approved drugs from Prestwick

480 purified natural products

4,100 N-Ac-tetrapeptide amides

1570 compounds from academic organic chemists

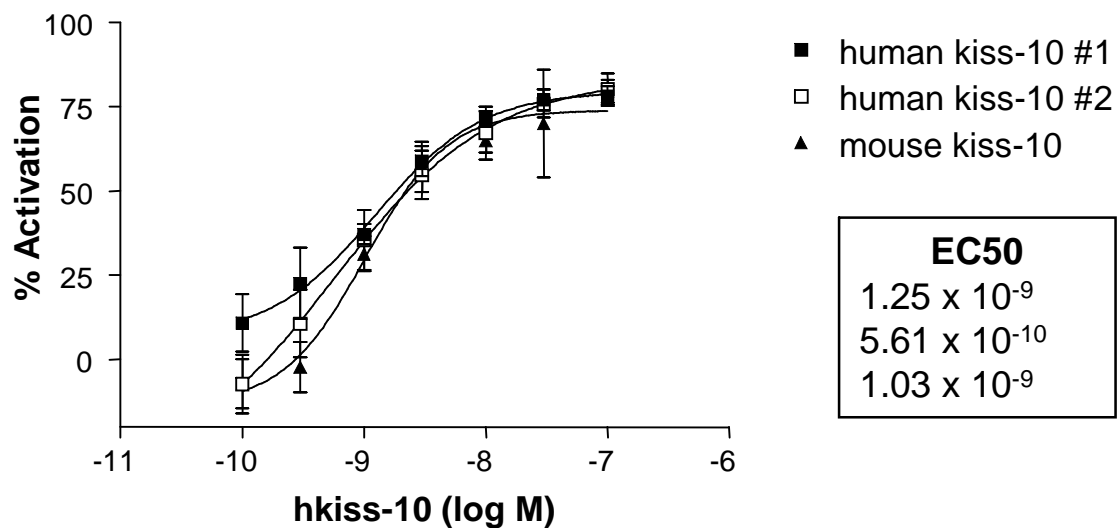
800 proprietary compounds synthesized by LDDN

- *Plus . . .* Program to collect additional compounds from academia.
- Additional compounds being purchased.



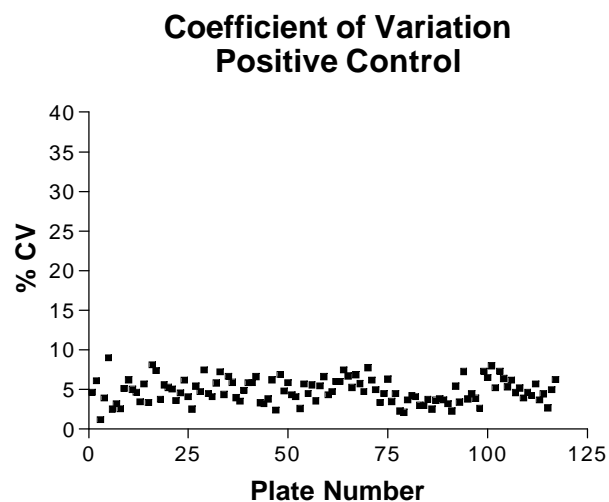
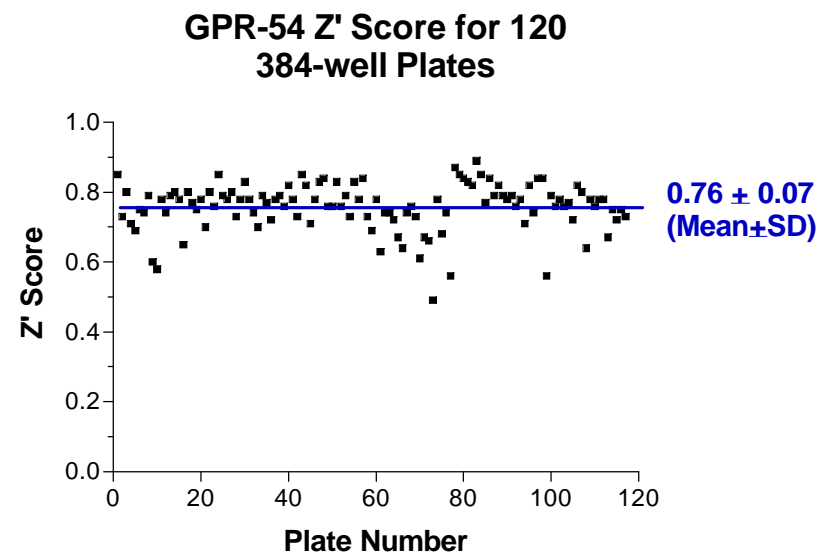
# Control for Automation

## ➤ % Activation with Kisspeptin Dose-Response

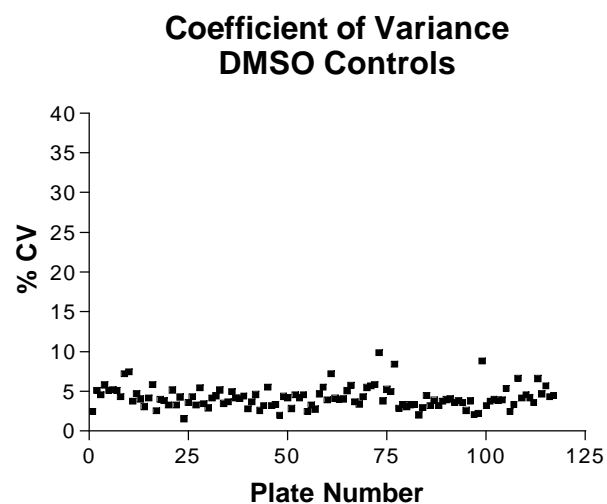


# Screening Statistics

$Z' = 0.7$   
S:B = 4-5



Maximal kisspeptin



DMSO

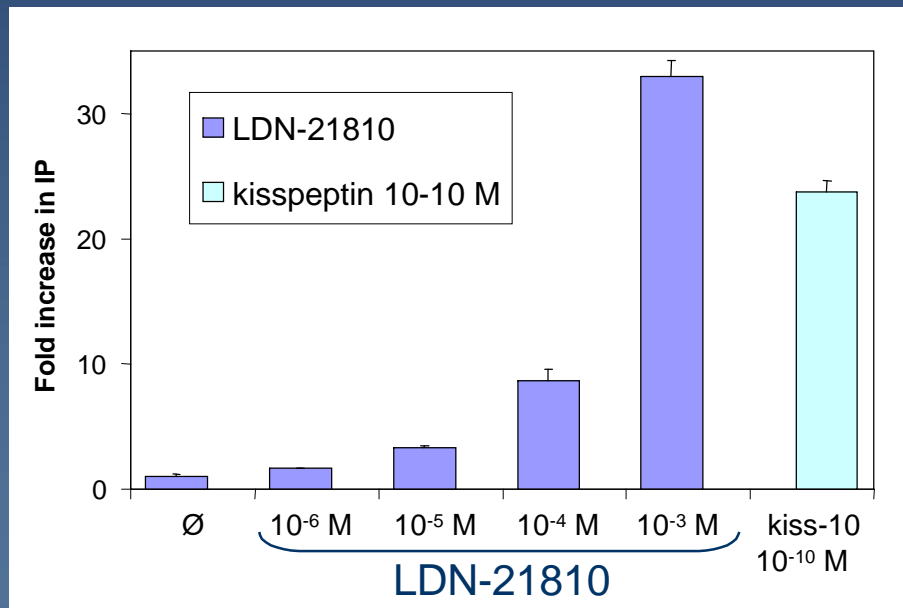
Laboratory for Drug  
Discovery in  
Neurodegeneration



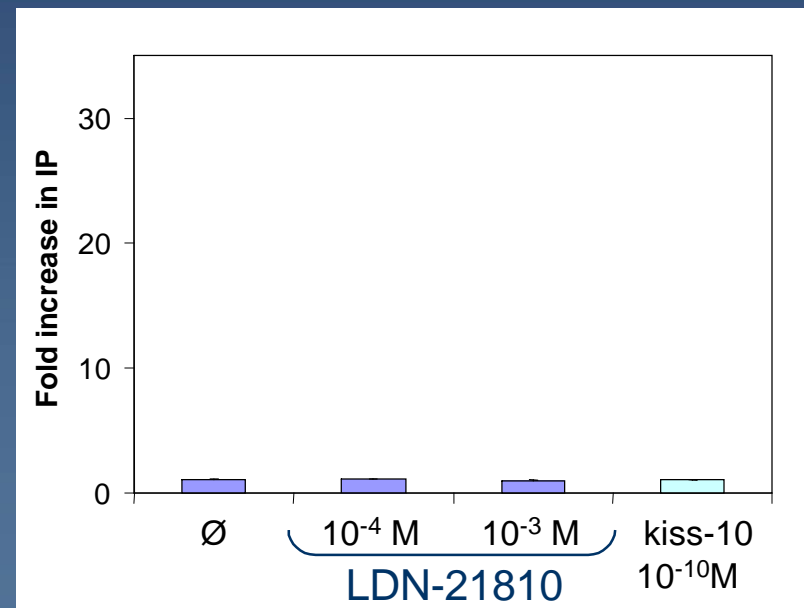
# LDN-21810 In Follow-up Assays

- **LDN-21810 Stimulates IP Production in a GPR54-Specific and Dose-Dependent Manner**

## GPR54 Line



## Empty Vector Line



# Agonist Screening Results

---

- The LDDN small molecule library of 110,000 compounds has been screened for GPR54 agonists.
- Currently characterizing one of the hits
  - LDN-21810.
- Unfortunately, HTS did not identify many hits. This is not surprising with peptide-ligand receptors.

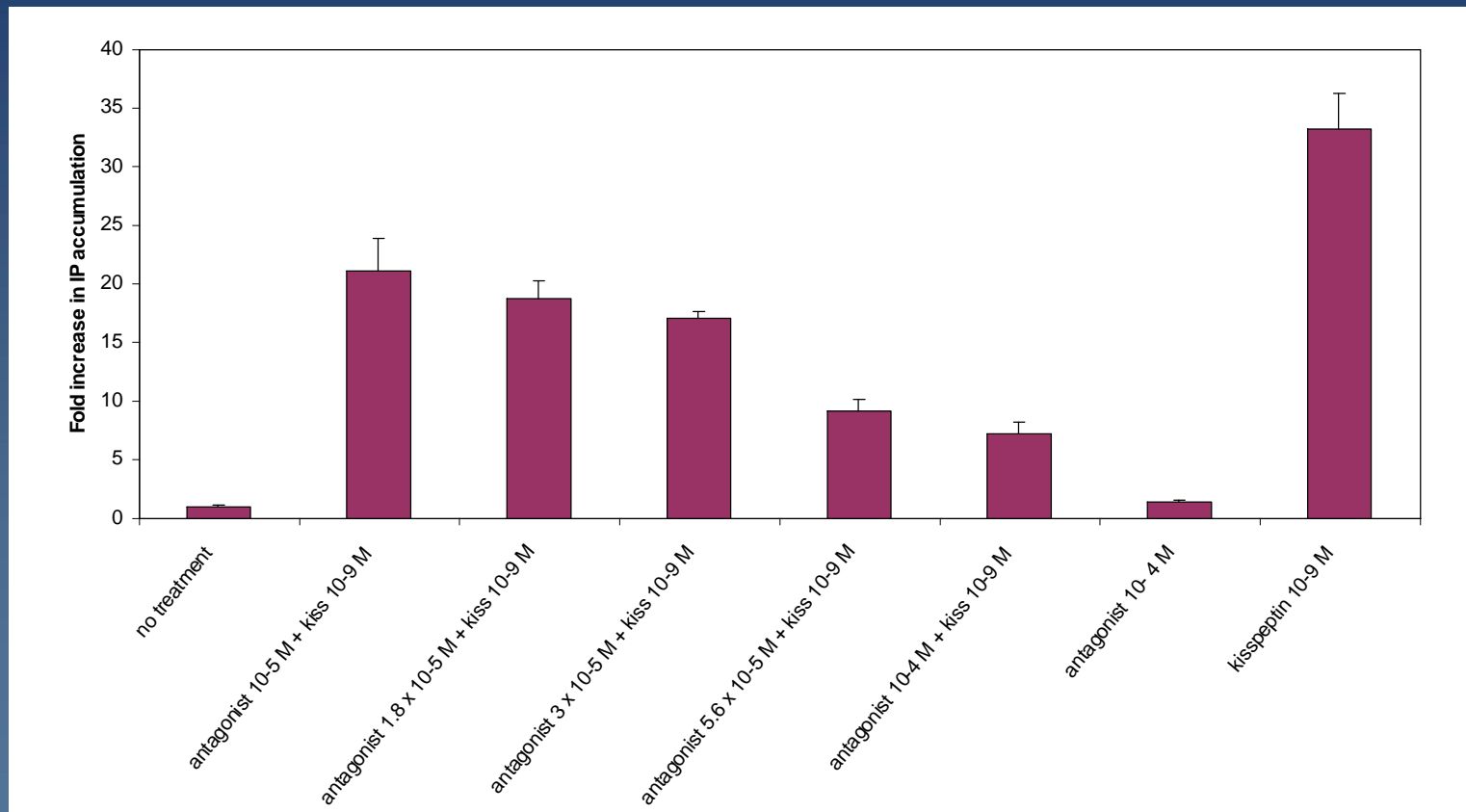
# Strategy for Antagonist Screening

---

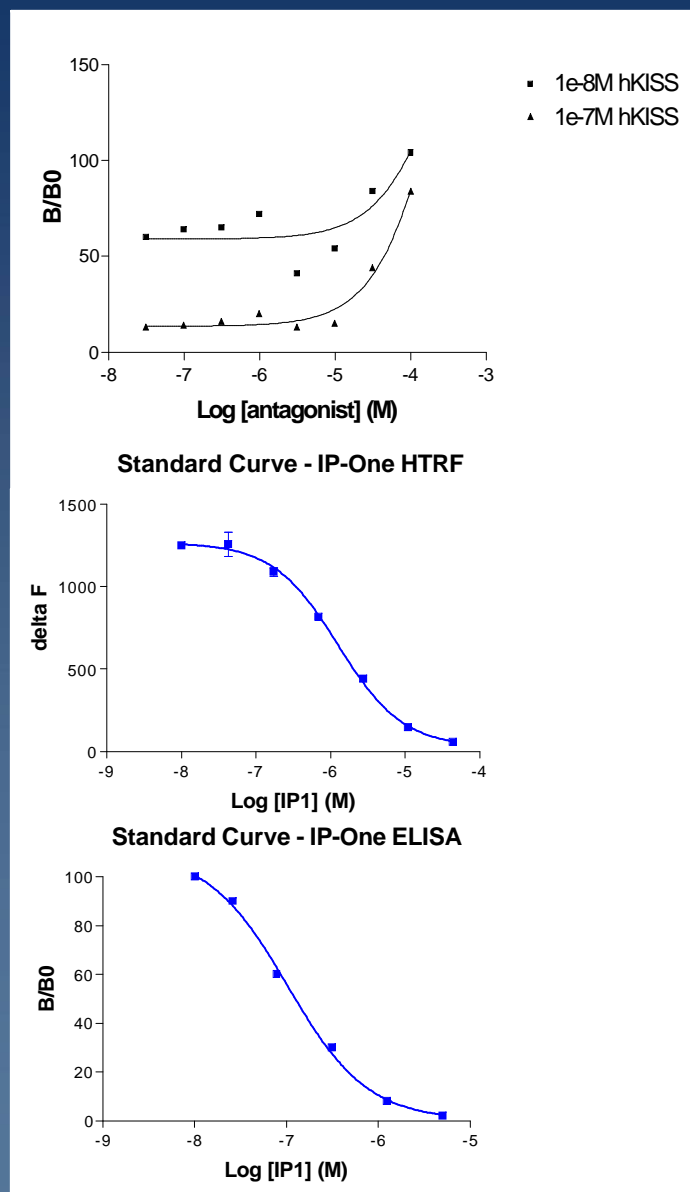
- The PLC inhibitor U-73122 used as a positive control.
- Negative controls: kisspeptin without U-73122 and no D<sub>2</sub> wells
- Currently undergoing optimization



# Radioactive IP assay with kisspeptin and U-73122 (antagonist)



# IP-One ELISA assay of GPR54 with kisspeptin



► ELISA is more sensitive than HTRF

► Data very similar

► Allows labs without fancy readers the ability to use IP-One

## SUMMARY

---

- A GPR54 stably transfected cell line was generated for screening and was found to respond appropriately in functional assays.
  - IP accumulation
  - ERK phosphorylation

## SUMMARY

---

- A GPR54 stably transfected cell line was generated for screening and was found to respond appropriately in functional assays.
  - IP accumulation
  - ERK phosphorylation
- An assay based on measurement of IP1 accumulation was optimized for high-throughput screening.

## SUMMARY

---

- A GPR54 stably transfected cell line was generated for screening and was found to respond appropriately in functional assays.
  - IP accumulation
  - ERK phosphorylation
- An assay based on measurement of IP1 accumulation was optimized for high-throughput screening.
- We completed an initial screen of an entire small molecule library for GPR54 agonists.

## SUMMARY

---

- A GPR54 stably transfected cell line was generated for screening and was found to respond appropriately in functional assays.
  - IP accumulation
  - ERK phosphorylation
- An assay based on measurement of IP1 accumulation was optimized for high-throughput screening.
- We completed an initial screen of an entire small molecule library for GPR54 agonists.
  - Positive hits from ligand screening were selected and verified using the screening assay (IP-One) and secondary functional assays.
  - The compound verified on secondary assays is now being tested in an animal model.

## SUMMARY

---

- A GPR54 stably transfected cell line was generated for screening and was found to respond appropriately in functional assays.
  - IP accumulation
  - ERK phosphorylation
- An assay based on measurement of IP1 accumulation was optimized for high-throughput screening.
- We completed an initial screen of an entire small molecule library for GPR54 agonists.
  - Positive hits from ligand screening were selected and verified using the screening assay (IP-One) and secondary functional assays.
  - The compound verified on secondary assays will then be tested in an animal model.
- Repeat screening of the library for GPR54 antagonists and enhancers is underway.



# Laboratory for Drug Discovery in Neurodegeneration

Ross Stein, PhD, *Director* LDDN

Marcie Glicksman, *Sr. Director* Leads Discovery

Greg Cuny, *Sr. Director* Chemistry

- Jake Ni
- April Case
- Mickey Huang

*In collaboration with:*

**Brigham and Women's Hospital**

- Wendy Kuohung
- Ursula Kaiser
- Deepa Mukhtyar

**Massachusetts General Hospital**

- William Crowley, Jr.
- Stephanie Seminara
- Samuel Posner



*Harvard Center for Neurodegeneration and Repair*