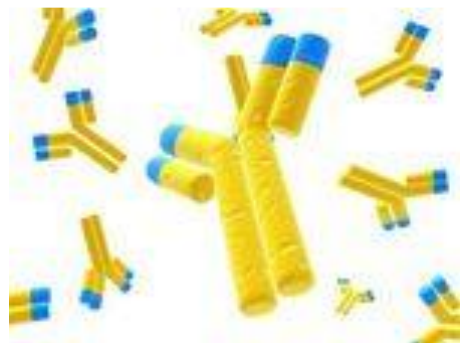
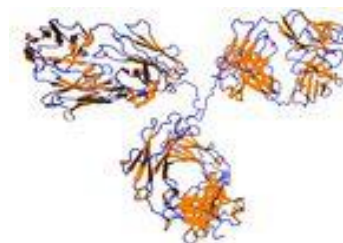


# Applications of HTRF and Tag-lite Assays for HTP Antibody Screening

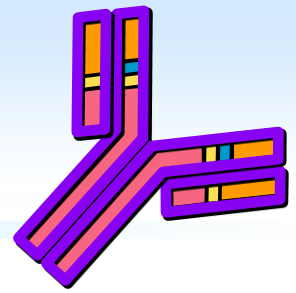


**Brigitte Devaux, PhD**  
Bristol Myers Squibb, Redwood City CA



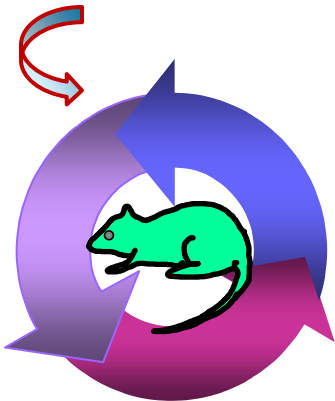
**HTRF Symposium**  
**April 25, 2013**

# Introduction



- Generate human therapeutic antibodies to large number of targets (soluble or membrane bound) for various therapeutic indications
- Also produce reagent antibodies for research/development/clinical applications
  - Hybridoma technology – Fusion of splenocytes with myeloma partner
- Need to develop HTP assays amenable to screening hybridoma supernatants from fusion plates
  - ◆ Create potential for screening on multiple antigens or cell types
    - Tagged antigen versus irrelevant Ag with same tag
    - human/cyno/mouse Ag target
    - Conjugated Ag versus un-conjugated Ag (with PEG for example)
    - Homologues from same family
    - Transfectants and cancer cells

Immunization



# Hybridoma Workflow

Human Ig Tg mice  
spleens

Balb/c spleens

**E-Fusion**  
Splenocytes + fusion partner

HuG/K Screening  
HTRF  
Automation-96

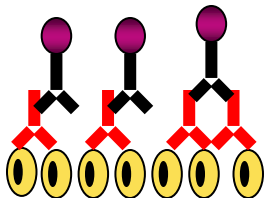
Antigen Screening

ELISA (soluble Ag)  
Hybridoma-96

HTRF (soluble Ag)  
AD-384

FACS (cells)  
Hybridoma-96

FMAT/Tag-lite (cells)  
AD-384



Ag specific Hit Picking  
Automation-24

Rescreen of  
positive lines

Subcloning – Expansion  
Purification

# HTP Antigen Screening Assays

---

## Requirements:

- ◆ Detect antibody binding to antigen (Soluble or Cell based)
- ◆ Robust & flexible assays
- ◆ No wash steps
- ◆ Work in conditioned media
- ◆ Sensitivity: Adequate for Ab concentration range from 10 ng-10  $\mu$ g /ml
- ◆ +/- response (no need for EC50 determination)
- ◆ Up to 10,000 samples in a run
- ◆ Time sensitive (24 hour turn around)
- ◆ Rapid assay development against a wide range of targets

# Presentation Outline

---

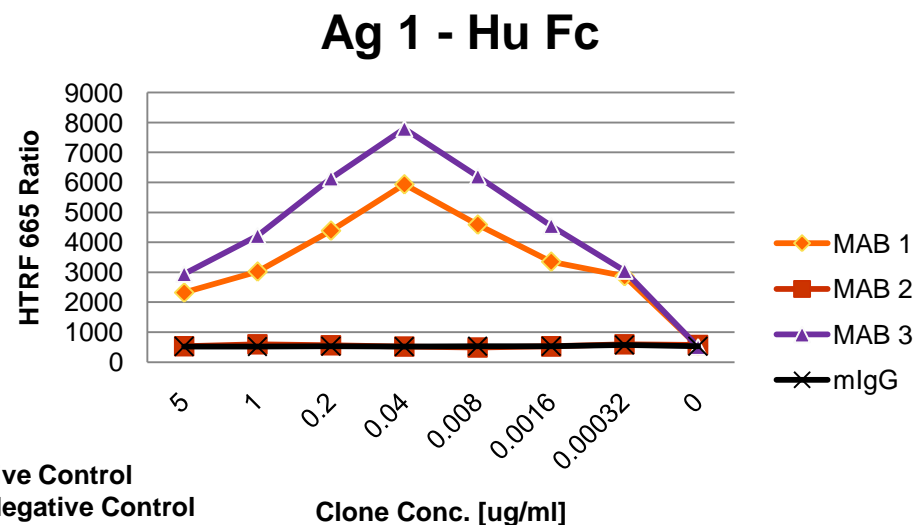
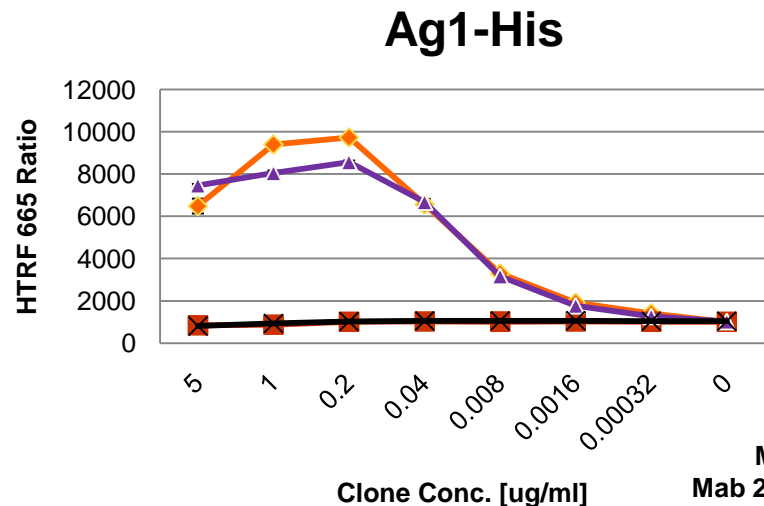
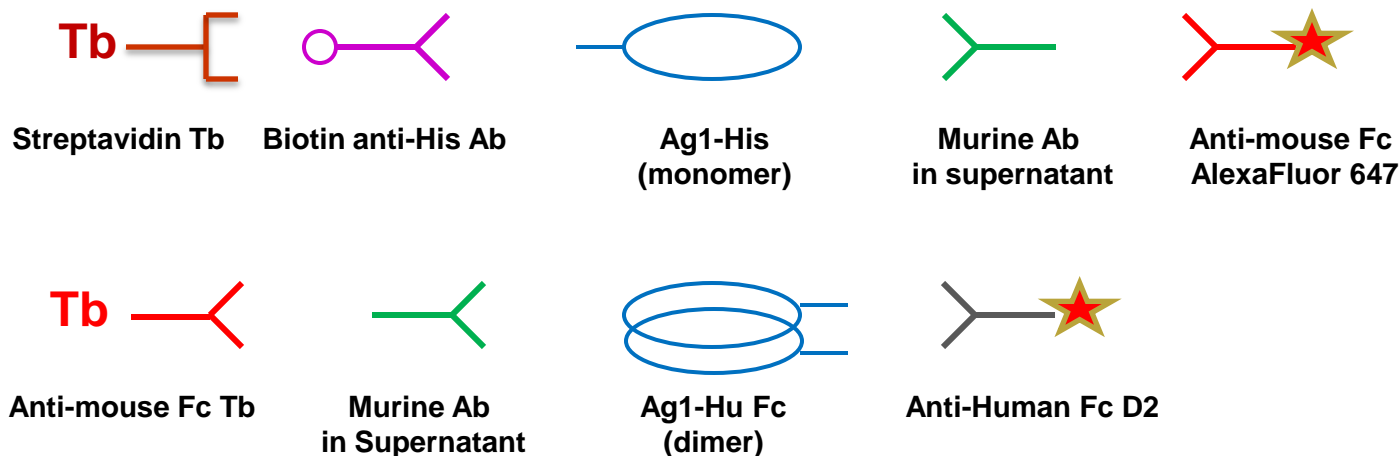
- **HTRF antigen screening assays (3 examples)**
  - HTRF with tagged antigens (e.g. His, Fc - Ag1 & 2)
  - HTRF with non-tagged antigen (Ag3) and labeled secondary reagents (non-blocking polyclonal)
- **HTRF with multiple labeled antigens in one well (duplexing, 2 examples)**
  - HTRF with Drug Ab IgG4 versus irrelevant IgG4
  - HTRF with PEG versus non PEG Ag
- **HTRF Blocking assay (1 example)**
- **HTRF Tag-Lite (2 examples)**
  - Tag-lite with GPCR1
  - Tag-lite with GPCR2

# Presentation Outline

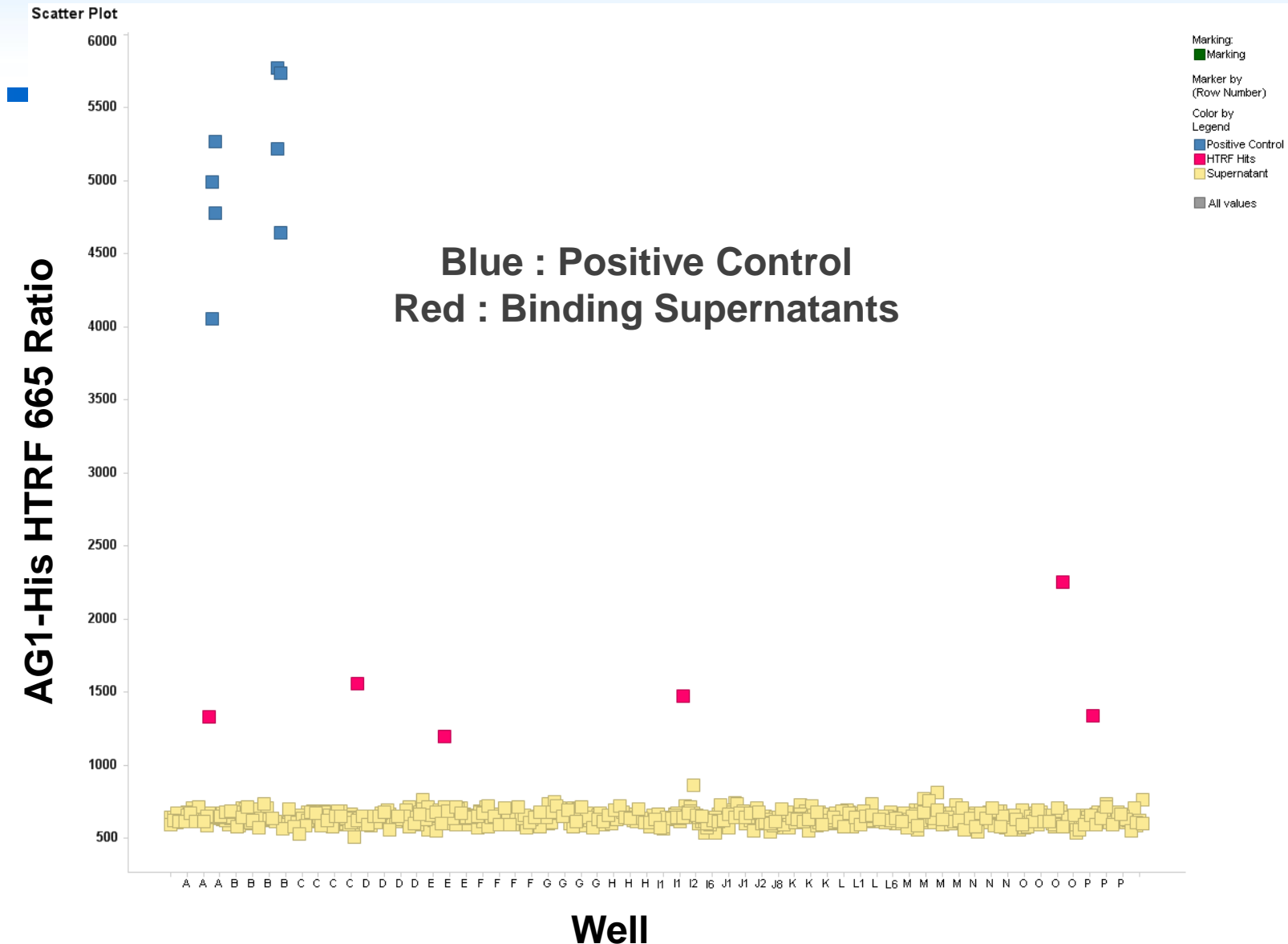
---

- **HTRF antigen screening assays (3 examples)**
  - HTRF with tagged antigens (Ag1 & 2)
  - HTRF with non-tagged antigen (Ag3) and labeled secondary reagents (non-blocking polyclonal)
- **HTRF with multiple labeled antigens in one well (duplexing, 2 examples)**
  - HTRF with Drug Ab IgG4 versus irrelevant IgG4
  - HTRF with PEG versus non PEG Ag
- **HTRF Blocking assay (1 example)**
- **HTRF Tag-Lite (2 examples)**
  - Tag-lite with GPCR1
  - Tag-lite with GPCR2

# HTRF with Tagged Antigens (His, Fc): Ag 1



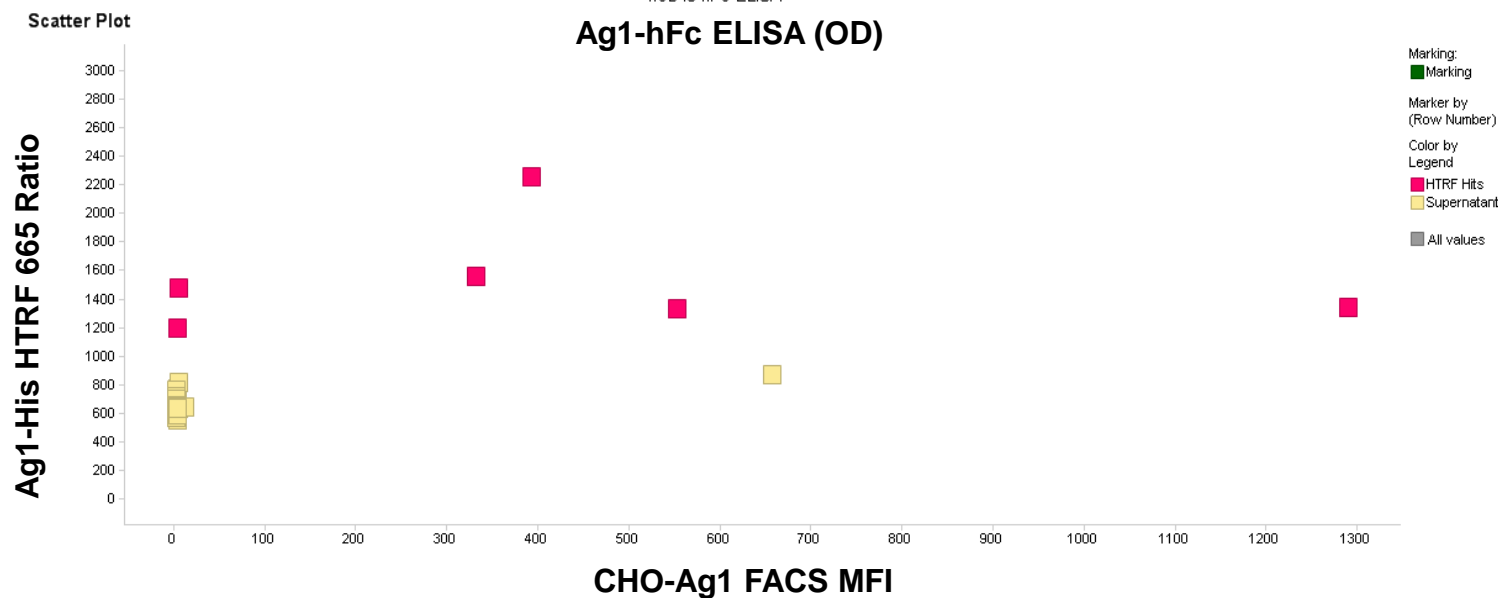
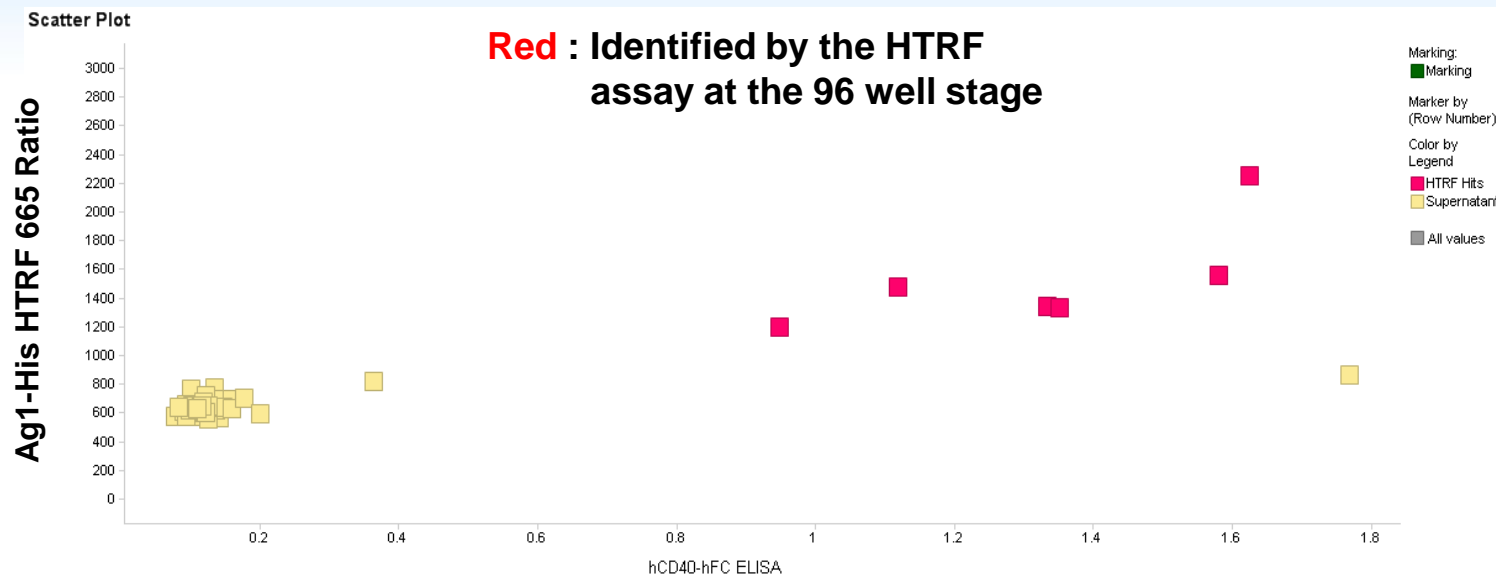
# HTRF Ag1-His Fusion Screening at 96 well stage



Identification of Six Ab Supernatants with Binding Activity



## Supernatants tested in Ag1 ELISA and FACS assays at 24 well stage



**Same hits in HTRF and ELISA assays**  
**Not all HTRF/ELISA hits bind by FACS**

# HTRF with Tagged Antigens (His): Ag 2

Antigen has a large number of leucine rich repeats on its extracellular surface

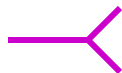
Primary screening: HTRF and ELISA in parallel

- HTRF binding in solution
- ELISA with Ag coated on plate

HTRF Format



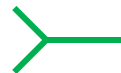
Anti-mouse Fc Tb



mouse anti-His



Ag2-His

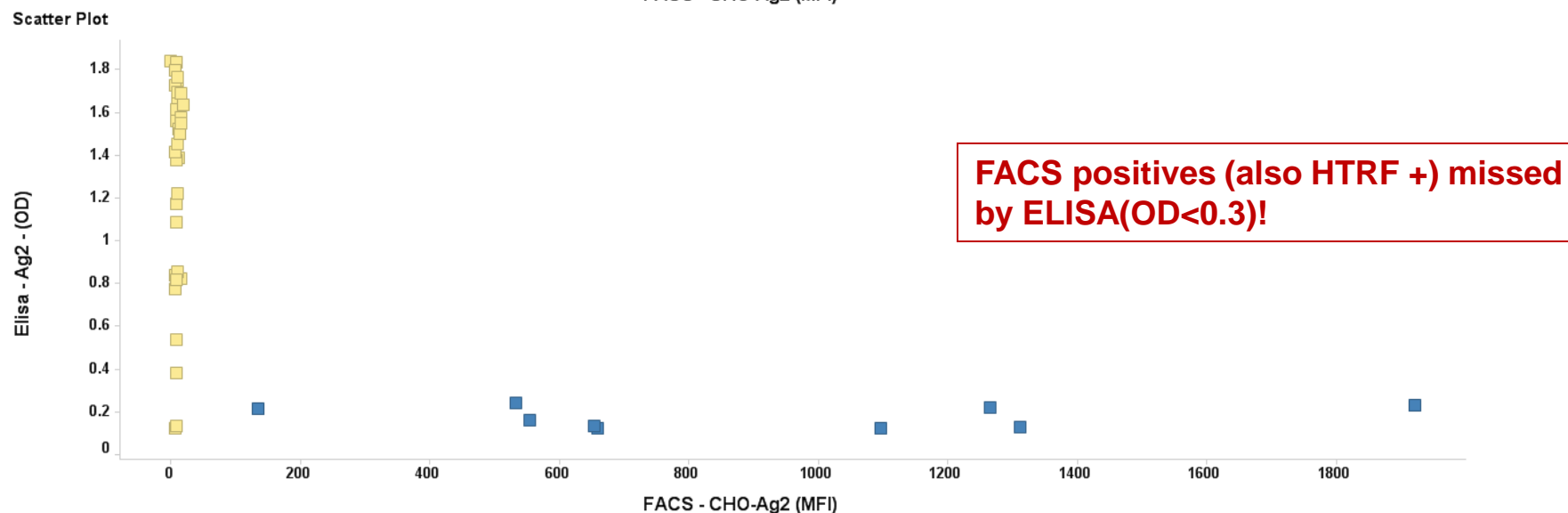
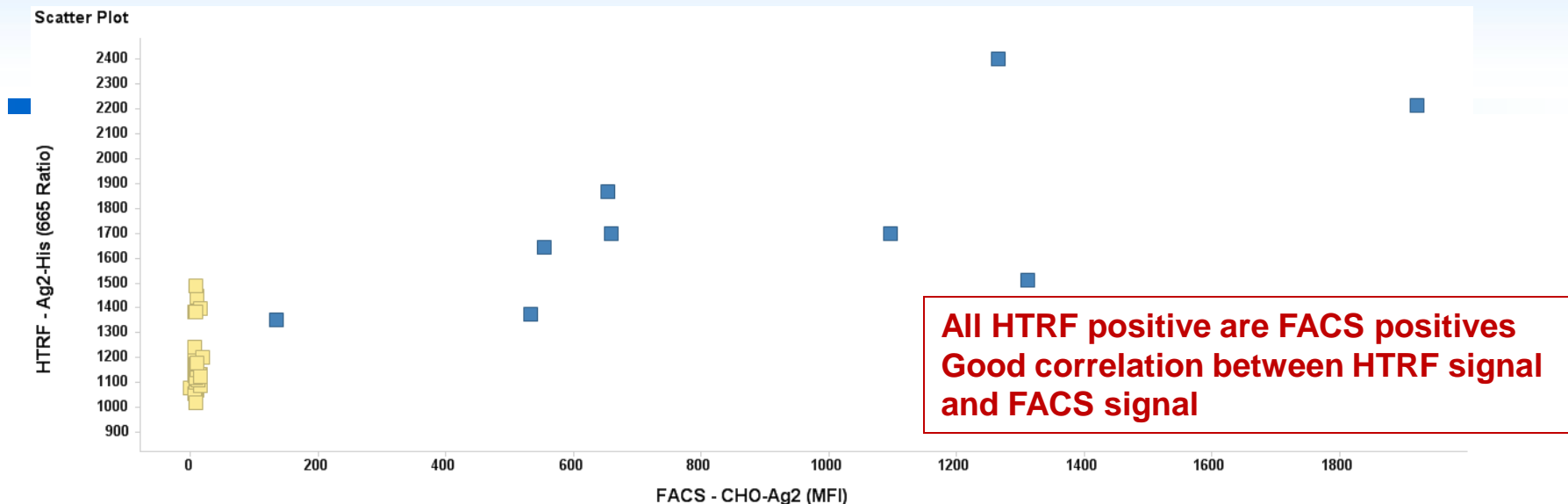


Human Clone  
in supernatant



Anti-Human Fc  
AlexaFluor 647

# Ag 2 - HTRF & ELISA vs. FACS

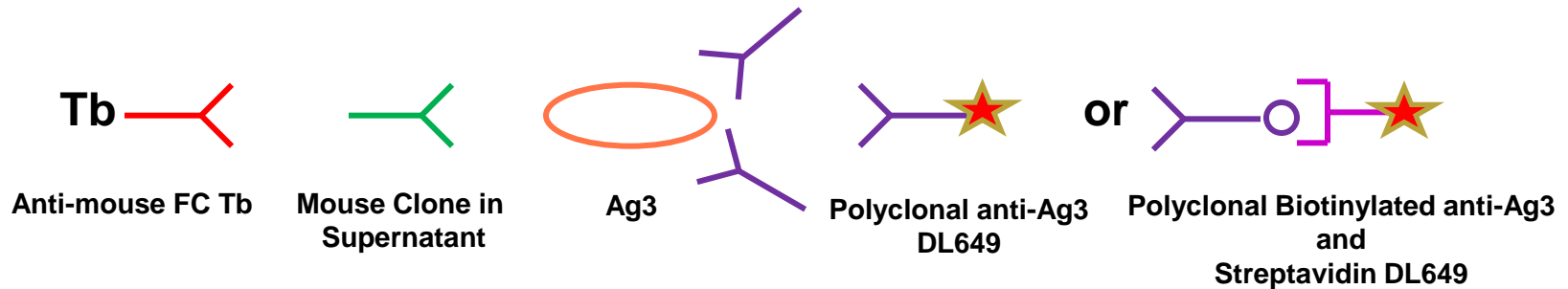


Blue are FACS positive - subcloned

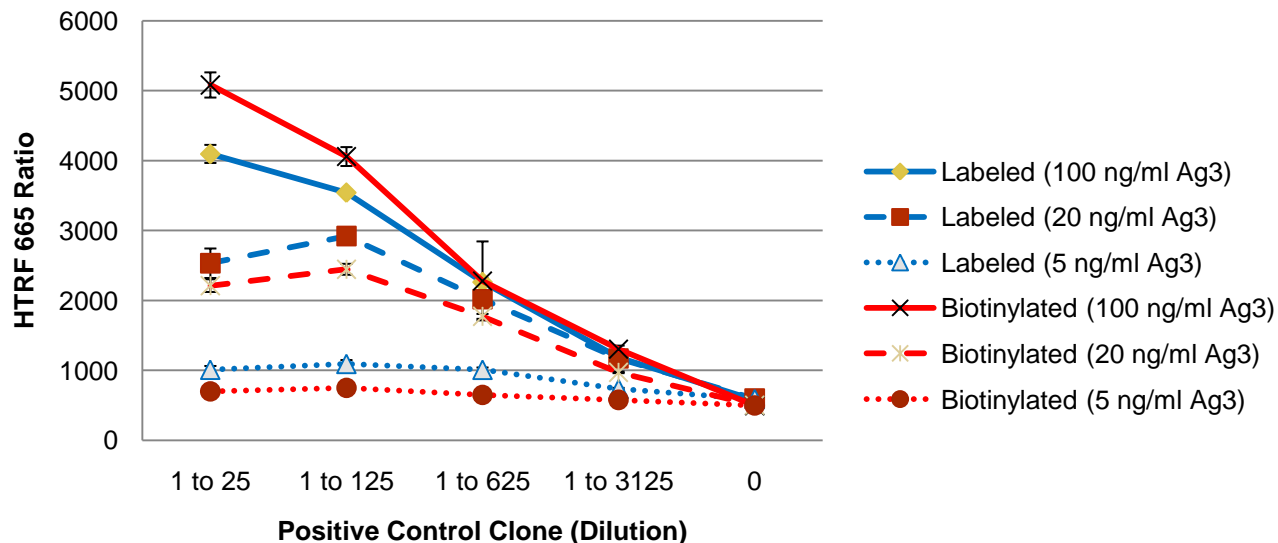


Bristol-Myers Squibb

# HTRF with Labeled Secondary Reagent Detection: Ag3



## HTRF - Ag3 - Labeled vs Biotinylated Secondary Antibody

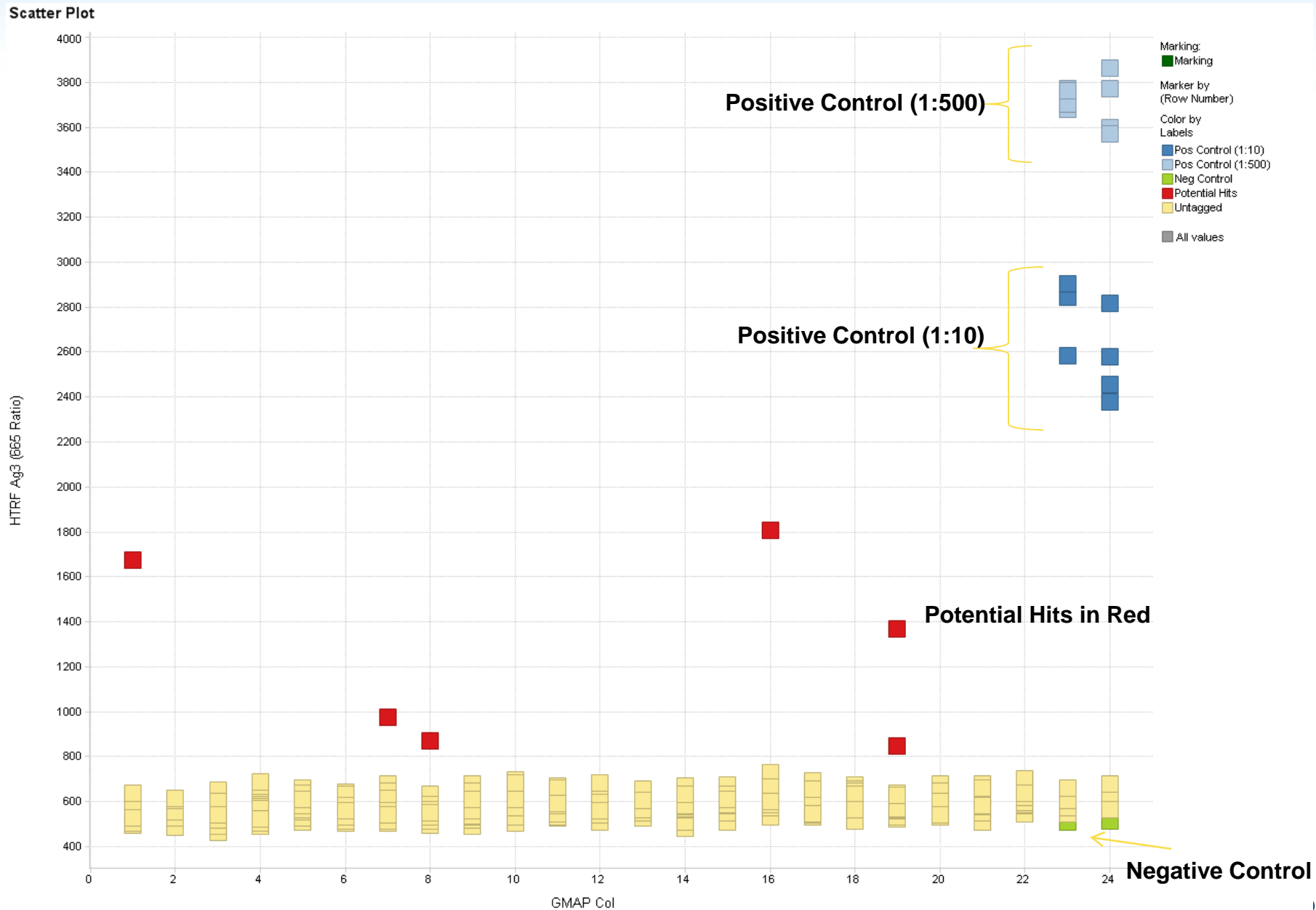


Direct and indirect labeling of polyclonal generated same results



Bristol-Myers Squibb

# Ag3 – HTRF 650 Ratio



# Presentation Outline

---

- **HTRF antigen screening assays (3 examples)**
  - HTRF with tagged antigens (Ag1 & 2)
  - HTRF with non-tagged antigen (Ag3) and labeled secondary reagents (non-blocking polyclonal)
- **HTRF with multiple labeled antigens in one well (duplexing, 2 examples)**
  - HTRF with Drug Ab IgG4 versus irrelevant IgG4
  - HTRF with PEG versus non PEG Ag
- **HTRF Blocking assay (1 example)**
- **HTRF Tag-Lite (2 examples)**
  - Tag-lite with GPCR1
  - Tag-lite with GPCR2

# HTRF with multiple Labeled Antigens in one well

## Duplexing: Drug Ab IgG4


Search for **anti-Id Abs**

The antigen is an Ab!

Do not want Abs binding to

Fc domain

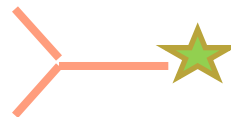
**“2 assays in 1”**

Tb   
Anti-mouse Fc Tb

  
Mouse Clone in  
Supernatant

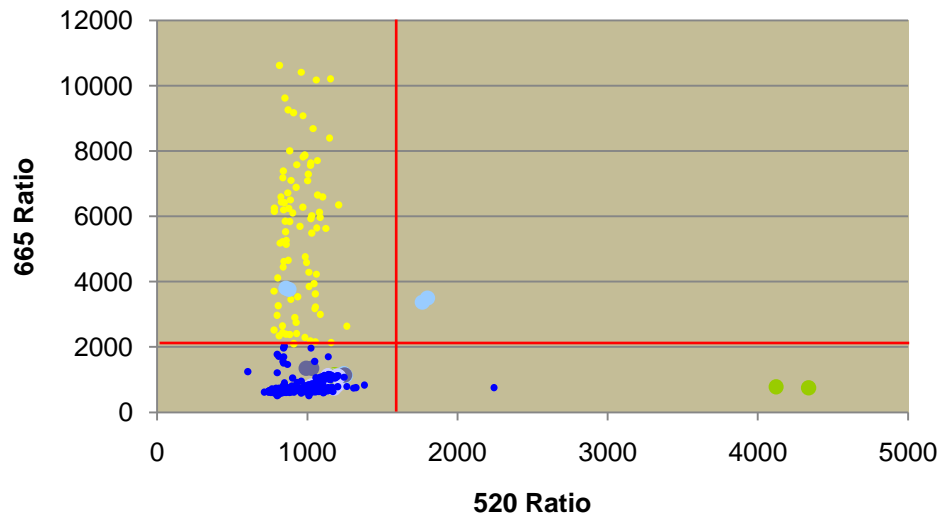
  
Drug Ab (IgG4) DL649

**AND**

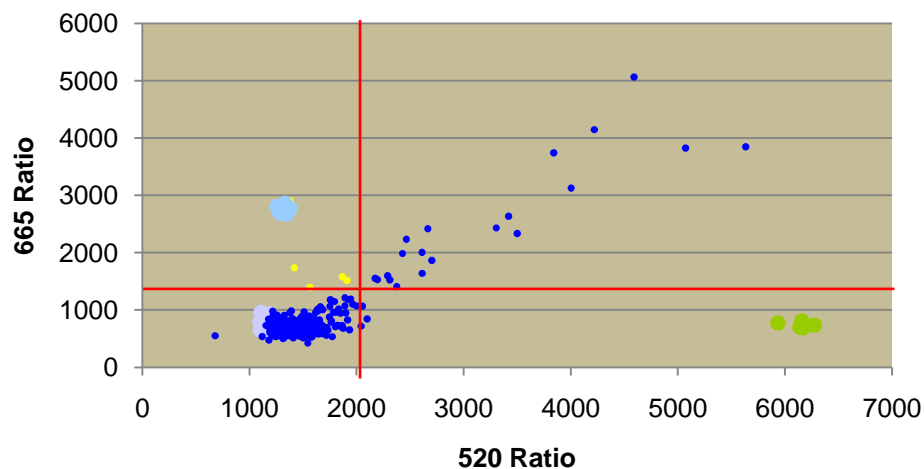
  
Irrelevant IgG4 DL488

Yellow : Selected hits  
Light Blue : Positive Control  
Dark Blue or Pink : Not selected

Drug Ab IgG4 DL 649



Immunogen :  
Drug Ab V domain only



Immunogen :  
Drug Ab IgG4

Irrelevant IgG4 DL 488

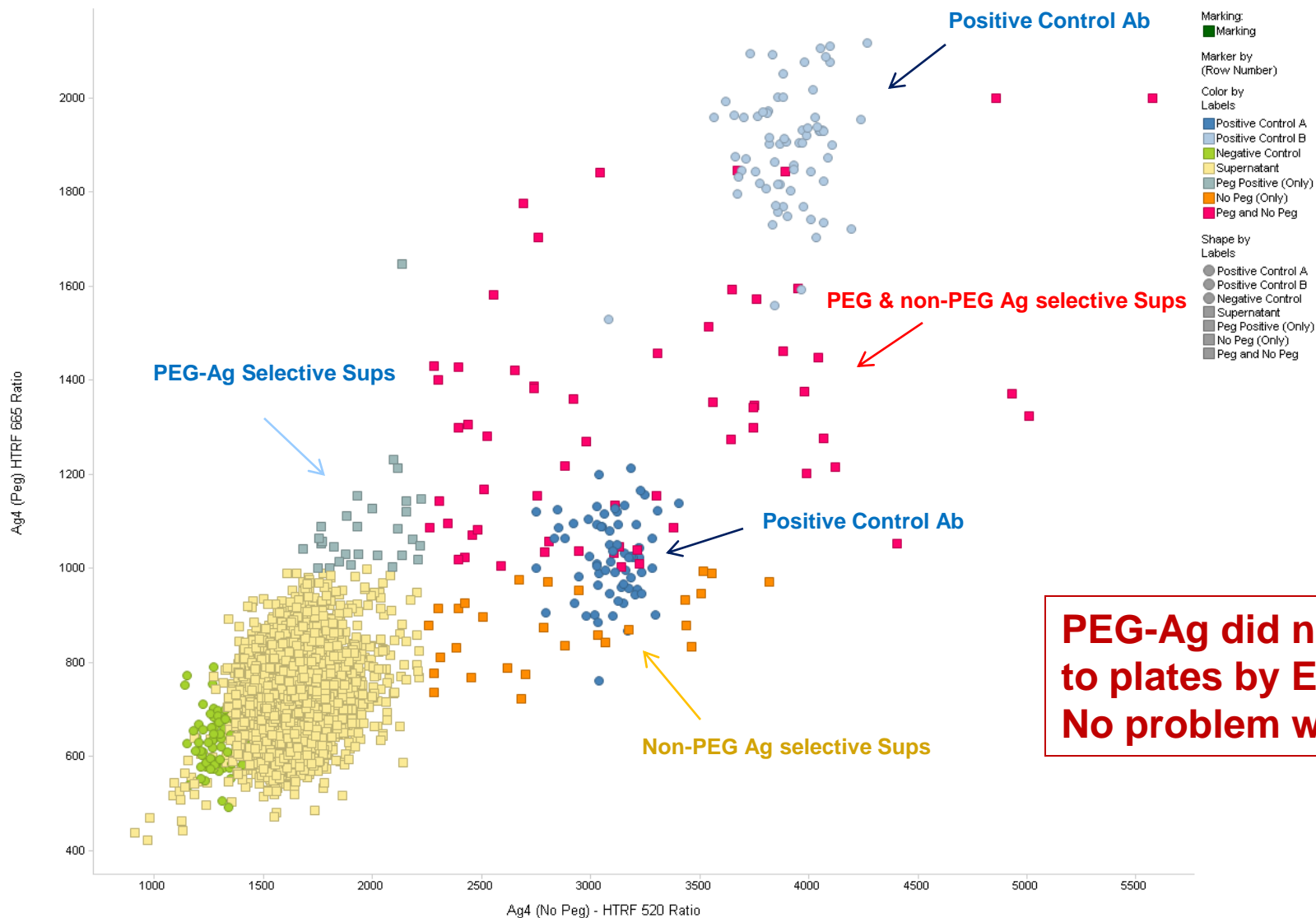
**More anti-Ids obtained when  
immunizing with V part of the Ab**



# HTRF with PEG-conjugated Antigen versus non-PEG Ag

## Duplexing: PEG-Ag versus non-PEG Ag

Scatter Plot



**PEG-Ag did not coat  
to plates by ELISA –  
No problem with HTRF**

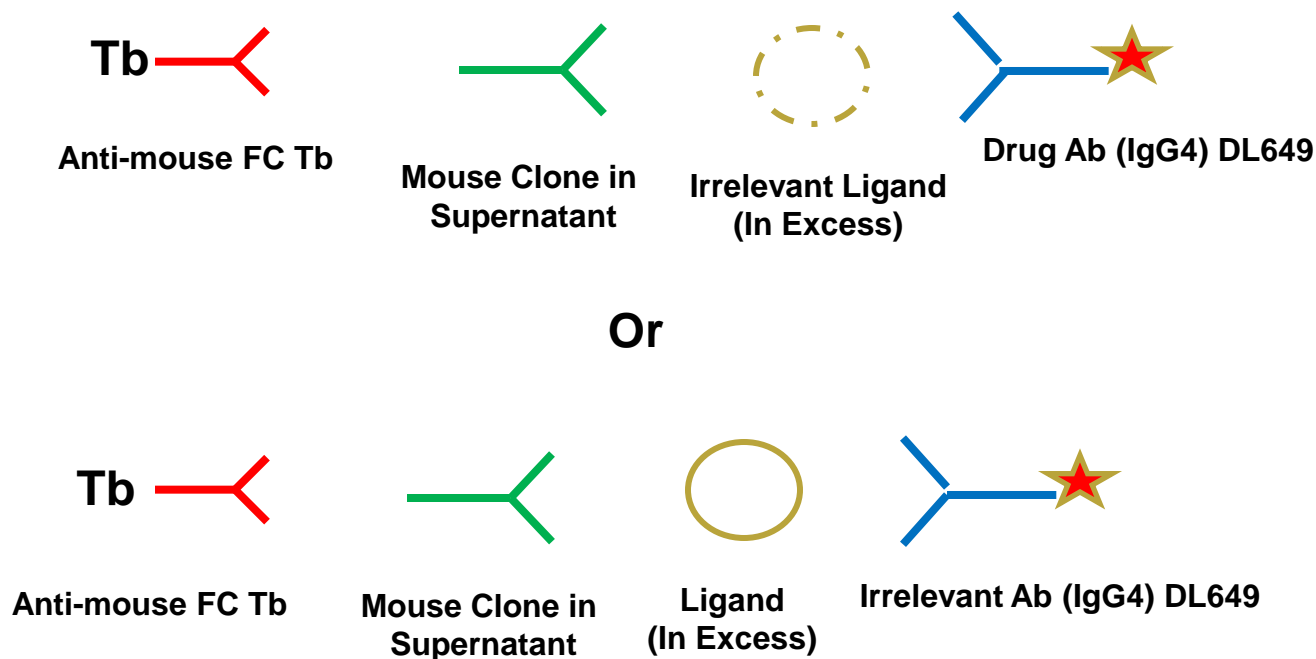
# Presentation Outline

---

- **HTRF antigen screening assays (3 examples)**
  - HTRF with tagged antigens (Ag1 & 2)
  - HTRF with non-tagged antigen (Ag3) and labeled secondary reagents (non-blocking polyclonal)
- **HTRF with multiple labeled antigens in one well (duplexing, 2 examples)**
  - HTRF with Drug Ab IgG4 versus irrelevant IgG4
  - HTRF with PEG versus non PEG Ag
- **HTRF Blocking assay (1 example)**
- **HTRF Tag-Lite (2 examples)**
  - Tag-lite with GPCR1
  - Tag-lite with GPCR2

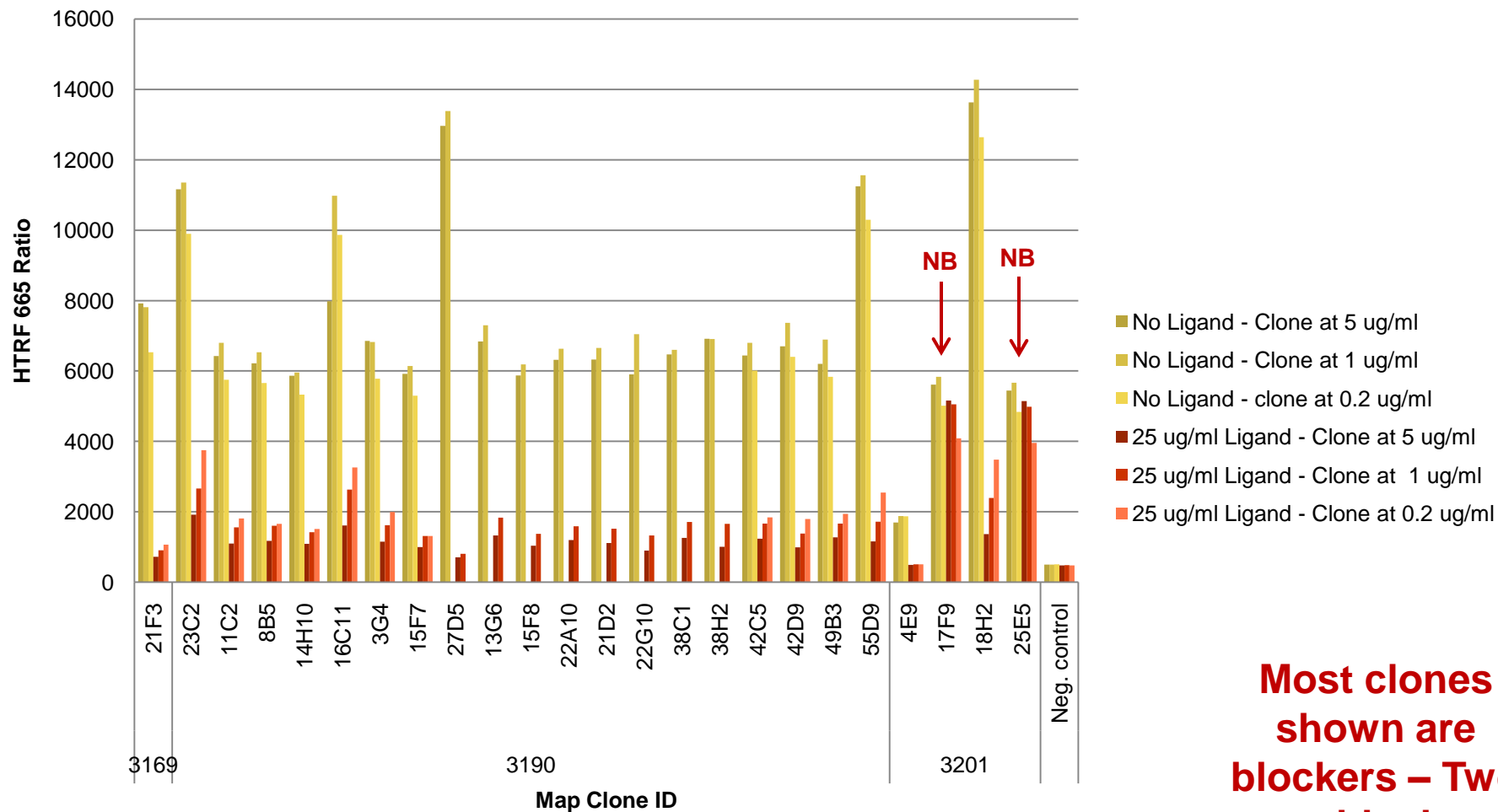
# HTRF Blocking assay

**Goal:** identify anti-Ids which bind to Ab when ligand is bound (non-blockers)  
or anti-Ids which interfere with ligand binding (blockers)



# HTRF Blocking Assay +/- Ligand

## HTRF Assay - Drug +/- Ligand - Anti-Id Ab Titration



**Most clones shown are blockers – Two non-blockers (anti-framework)**

# HTRF: Conclusions

## Advantages

- **Homogenous assay - No wash**
- **Sensitive** (5ng/ml or less)
- **Rapid** (1 hr incubation time)
- **Easily developed and automated**
  - ◆ Reagents stable
  - ◆ Low volume 384 well plates
  - ◆ Small reaction volumes
- **Binding in solution**
  - ◆ Binding antigen to plates for ELISA may block a subset of epitopes

## Disadvantages

- **Homogenous assay – Hook effect**
  - ◆ Can make distinguishing supernatants based on single point signals difficult
- **Requires antigen to be tagged or labeled**
  - ◆ May remove access to certain epitopes
- **Matrix (conditioned media) can interfere with HTRF ratio**
  - ◆ Ideally positive control in media



# Presentation Outline

---

- **HTRF antigen screening assays (3 examples)**
  - HTRF with tagged antigens (Ag1 & 2)
  - HTRF with non-tagged antigen (Ag3) and labeled secondary reagents (non-blocking polyclonal)
- **HTRF with multiple labeled antigens in one well (duplexing, 2 examples)**
  - HTRF with Drug Ab IgG4 versus irrelevant IgG4
  - HTRF with PEG versus non PEG Ag
- **HTRF Blocking assay (1 example)**
- **HTRF Tag-Lite (2 examples)**
  - Tag-lite with GPCR1
  - Tag-lite with GPCR2

# Tag-lite Technology: Applied to two GPCR Targets

Mix and Read ☺

Ligand or Ab binding to GPCR target

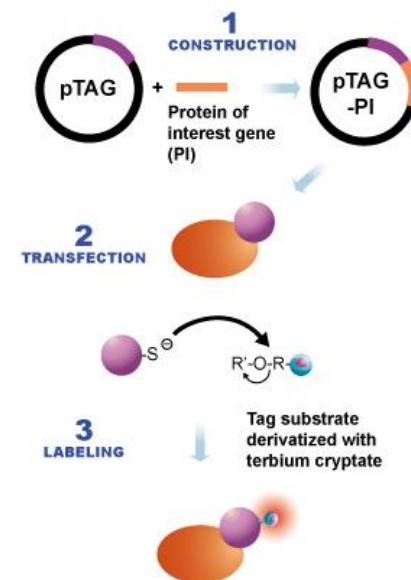
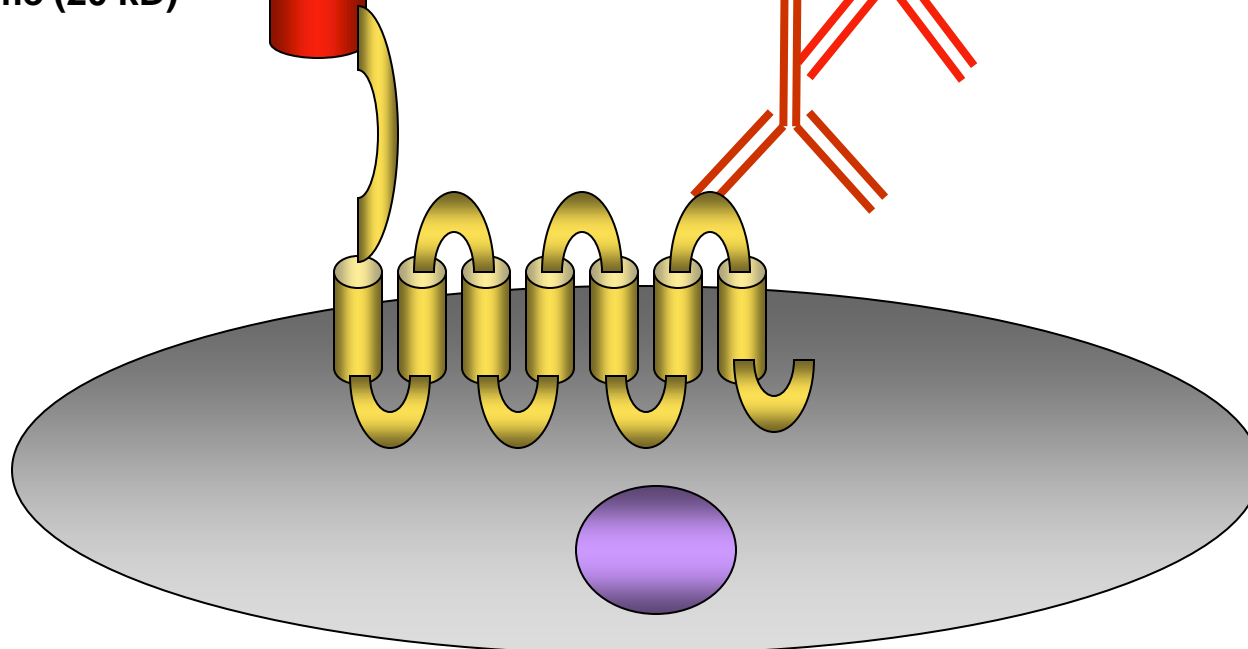
Alexafluor labeled Secondary Ab

Tb-Labeled Receptor



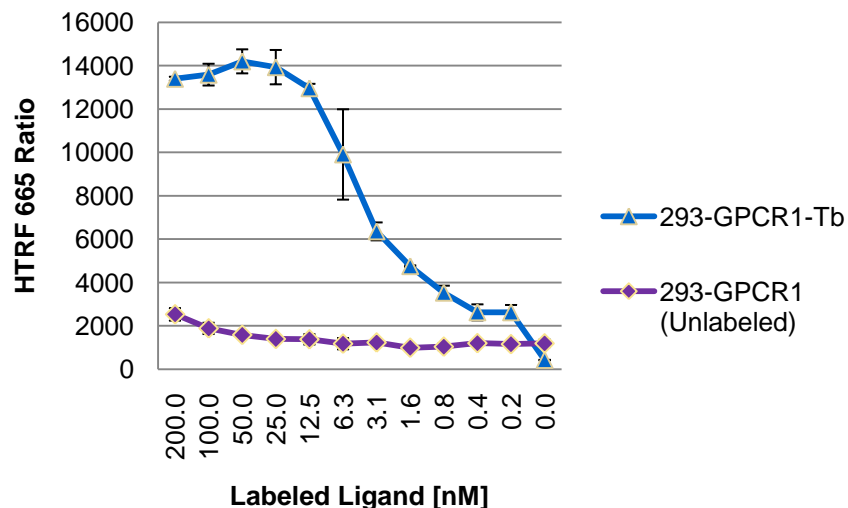
HTRF signal when binding occurs

Clip, Snap or Halo enzyme (20 kD)

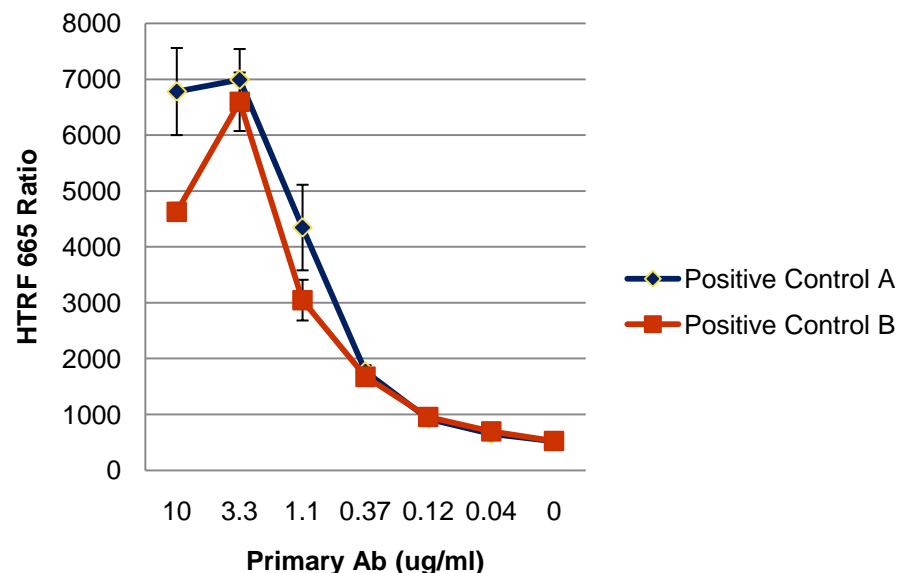


# GPCR 1 – HTRF Ligand & Ab Binding to Cells

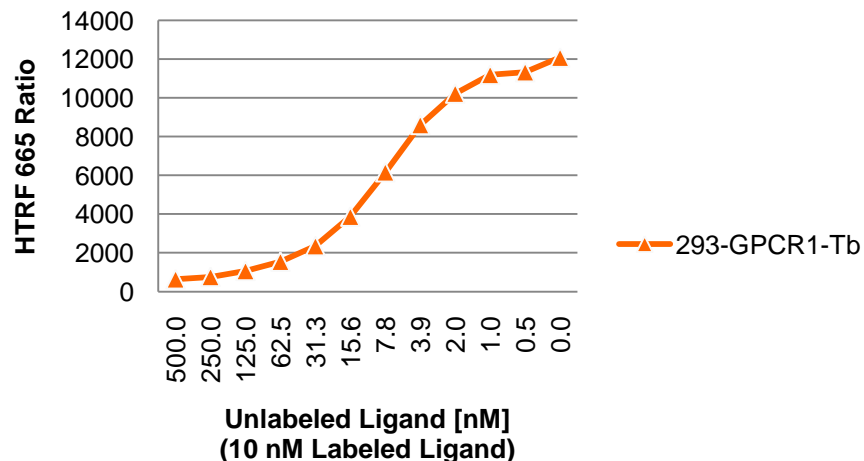
## Tag-lite - Cell Comparison



## Tag-lite - Primary Ab Titration



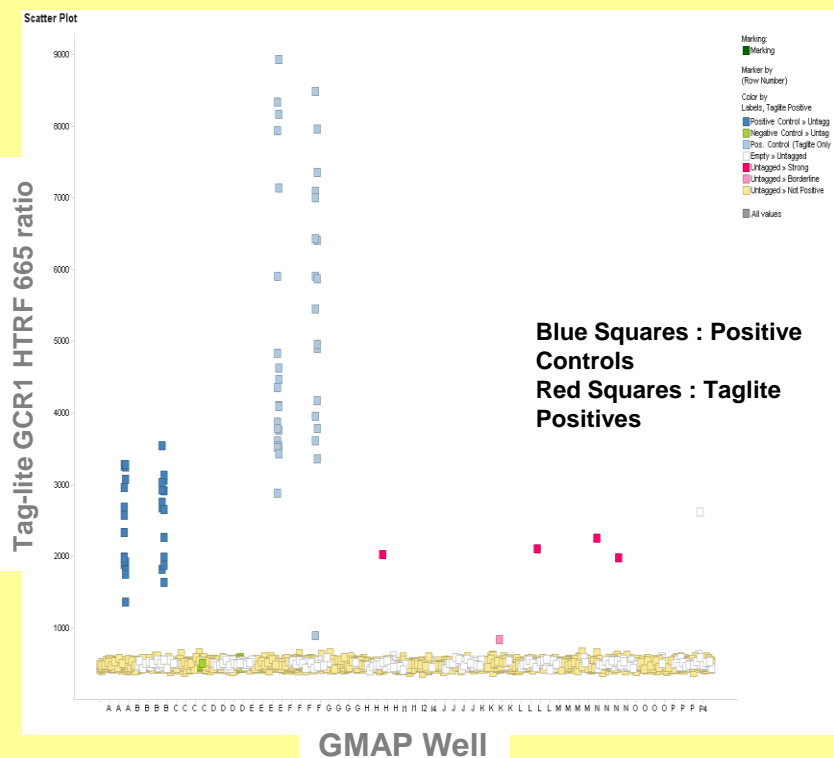
## Tag-lite - Unlabeled Ligand Titration



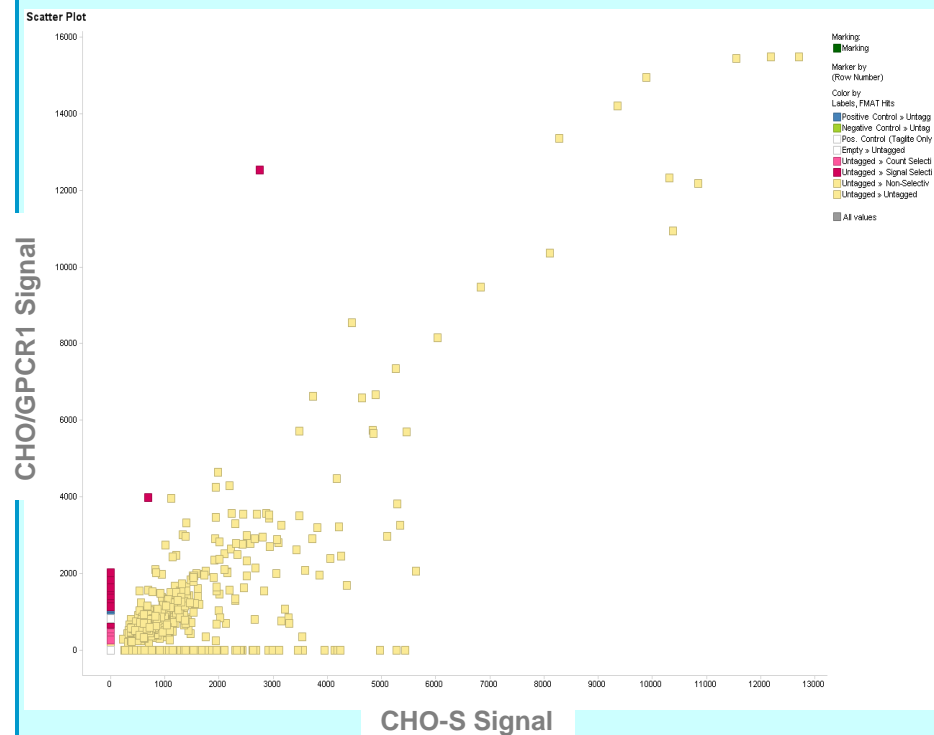


# Taglite – CHO/GPCR1 – Binding of Ab supernatants

## CHO/GPCR1 - Tag-lite GPCR1-Tb











## CHO-GPCR1 - FMAT



Many non-specific hits identified when FMAT was used  
Results much cleaner with Tag-lite



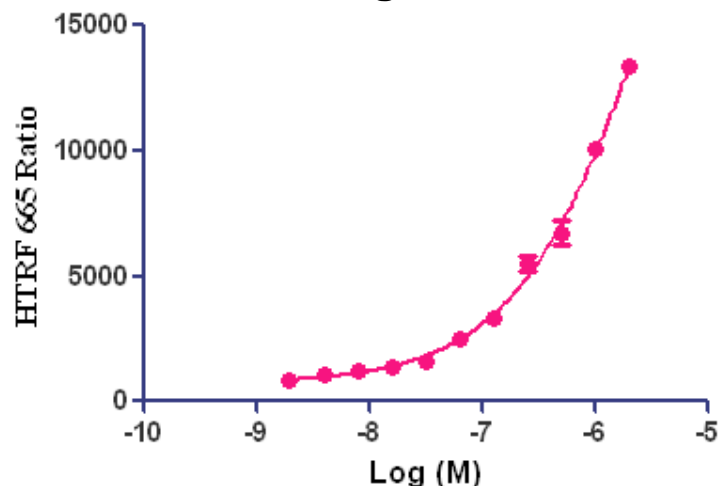
# Ab Binding on Various Cell Lines - FMAT

Clone	CHO-Ag	CHO	293-Ag	293	293Ag Snap	293Ag SnapTb	293-irrAg SnapTb
Pos.Control	+	-	+	-	+	+	-
MulG (-cont)	-	-	-	-	-	-	-
Ab Clone 1	+	+	-	-	-	-	-
Ab Clone 2		-		-			-
Ab Clone 3		-		-			-
Ab Clone 4	+	+	+	+	+	+	+

# GPCR 2 - HTRF Ligand Binding

No ligand binding cell assay available due to low affinity binding of ligand to GPCR receptor  
No possibility to assess blocking activity of antibodies or small molecule inhibitors

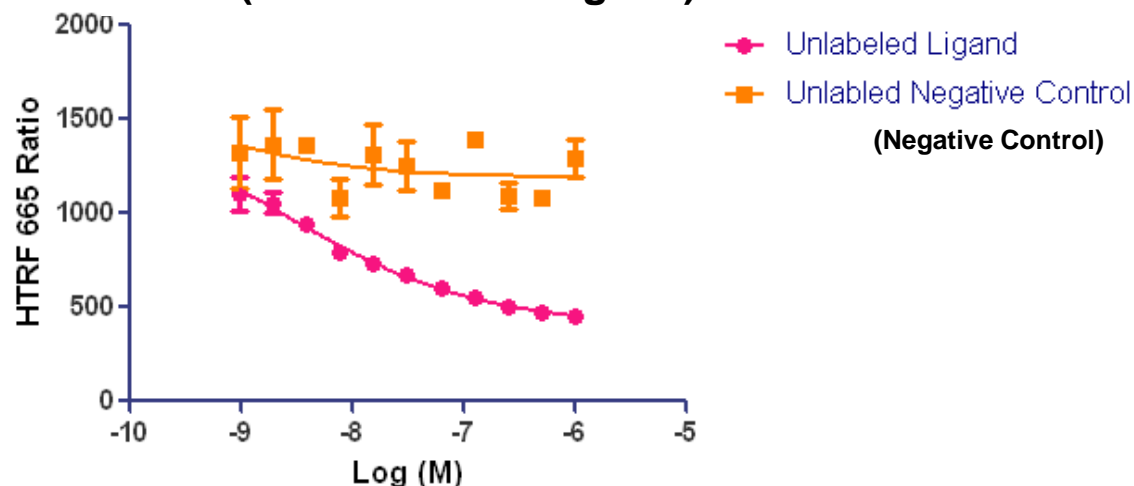
## Labeled Ligand Titration



50 nM ligand chosen for further inhibition experiments

(balancing the needs for maximizing the signal and minimizing ligand concentration)

## Unlabeled Ligand Competition (50 nM labeled Ligand)

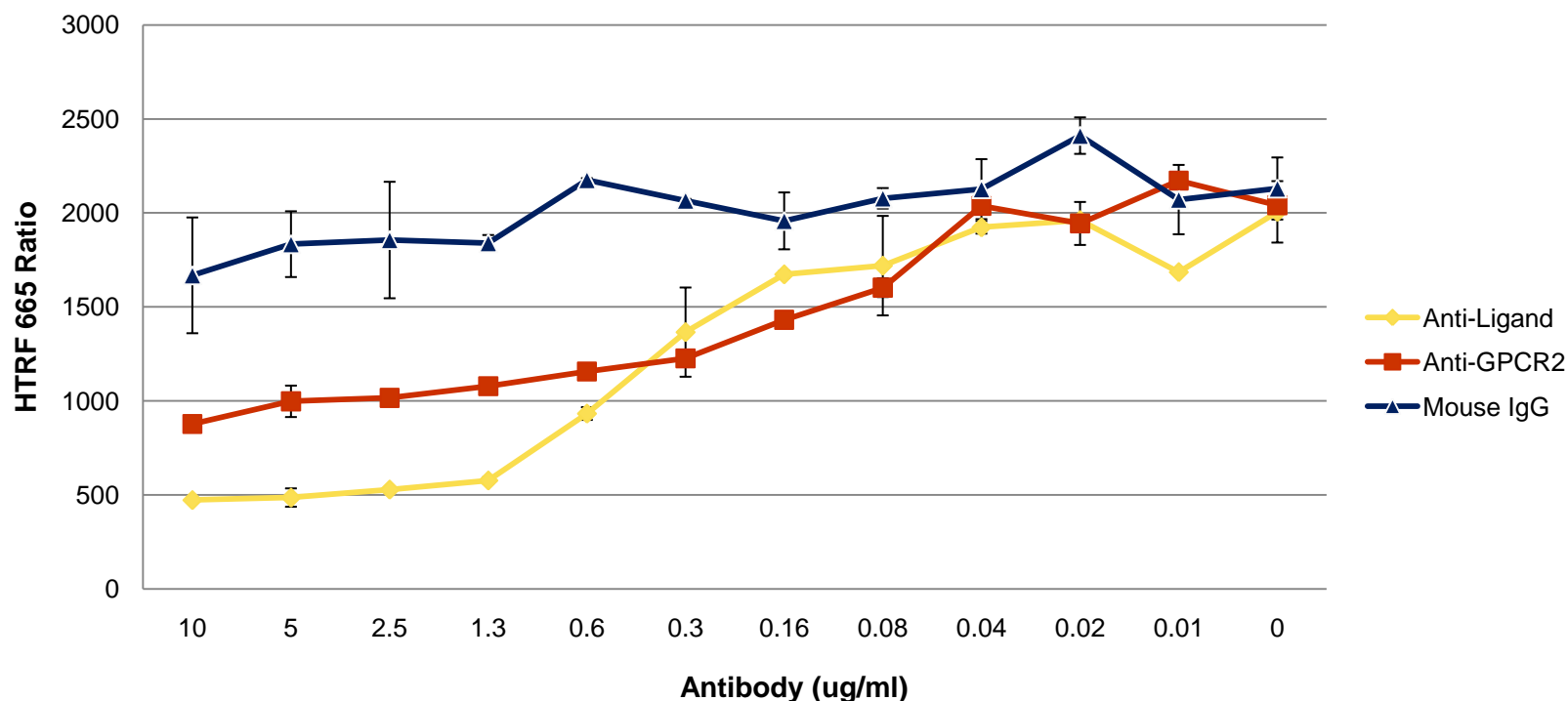


Cold ligand IC<sub>50</sub> = ~ 3nM

High concentrations of ligand are needed

# Inhibition by Blocking Antibodies

## Taglite - GPCR2 - Blocking Antibodies



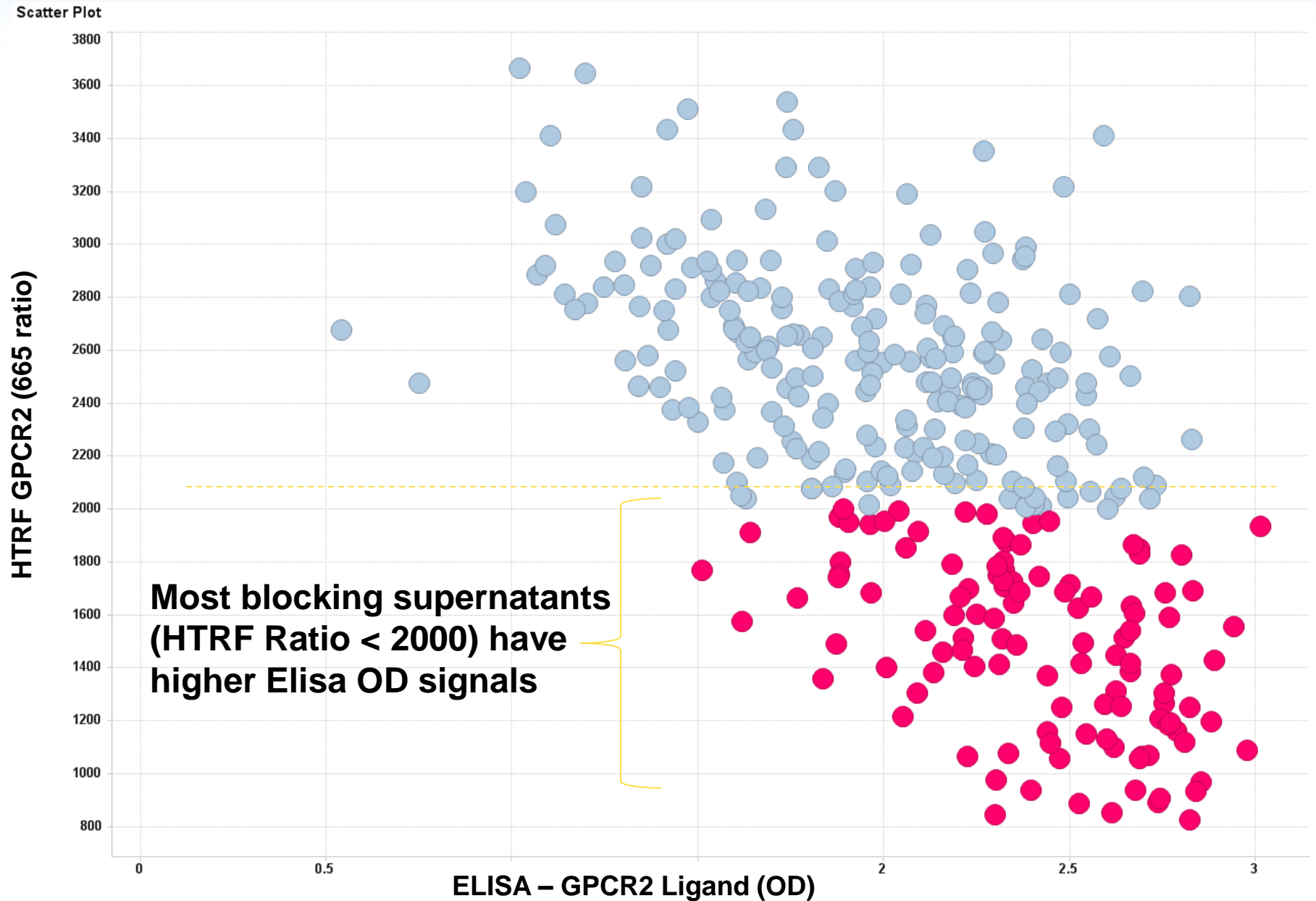
Anti-ligand Ab preincubated with ligand for 15 minutes  
Anti-GPCR2 and mIgG preincubated with cells for 15 minutes

**Both anti-ligand Ab and anti-GPCR2 Ab demonstrate significant inhibition of ligand binding, while irrelevant mouse IgG does not**

**Anti-GPCR2 does not appear to reach 100% inhibition, even when saturating (background is ~500)**

# Blocking with Anti-GPCR2 Ligand Supernatants

## ELISA (binding) versus Tag-lite (blocking)



# Taglite Conclusions

---

## Strengths:

- ◆ Easy, rapid, robust assay development
- ◆ Faster to execute than FMAT
- ◆ Less influenced by non-specific cell binding
- ◆ Can examine ligand/receptor interactions, even in cases where ligand binding assay could not be established using other means due to low affinity binding of ligand
- ◆ Can use same cells to look at binding and signaling

## Weaknesses

- ◆ Presence of Snap tag on the receptor can affect antibody binding
- ◆ Labeling ligand can alter its properties

# Thanks to

---

**Matthew Tomlinson**

**Scott Kurose**

**Edwin Haghazari &  
Na Tang From Cisbio**

**Paula So**

**Andrea Tatum**

**Anan Chuntharapai**

**Helen Cai**

**Erika Meaddough**

**Peter Brams – Keith Joho – Nils Lonberg**