

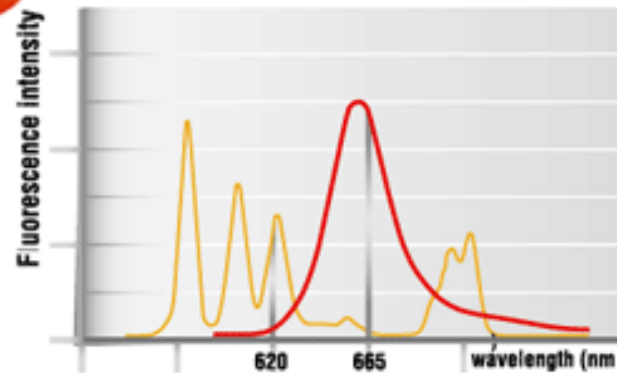
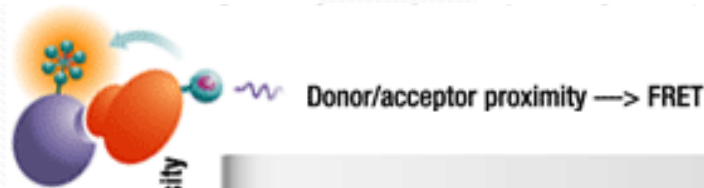
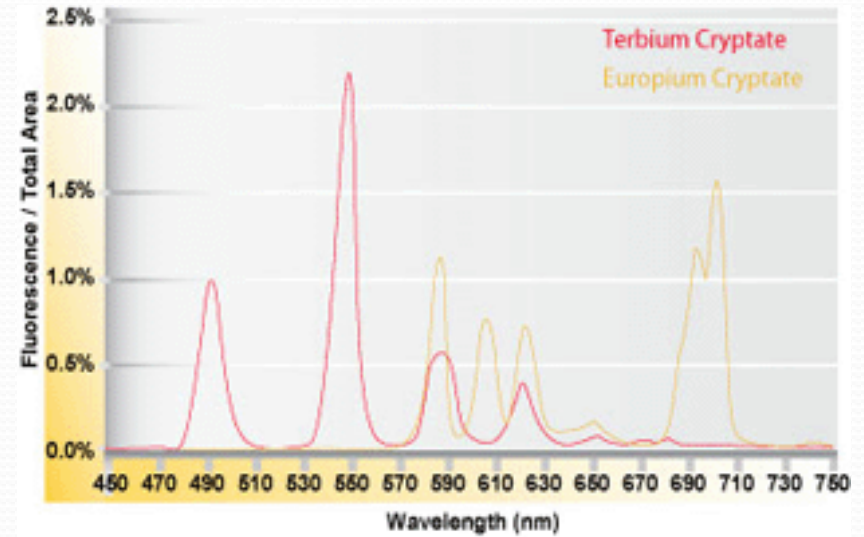
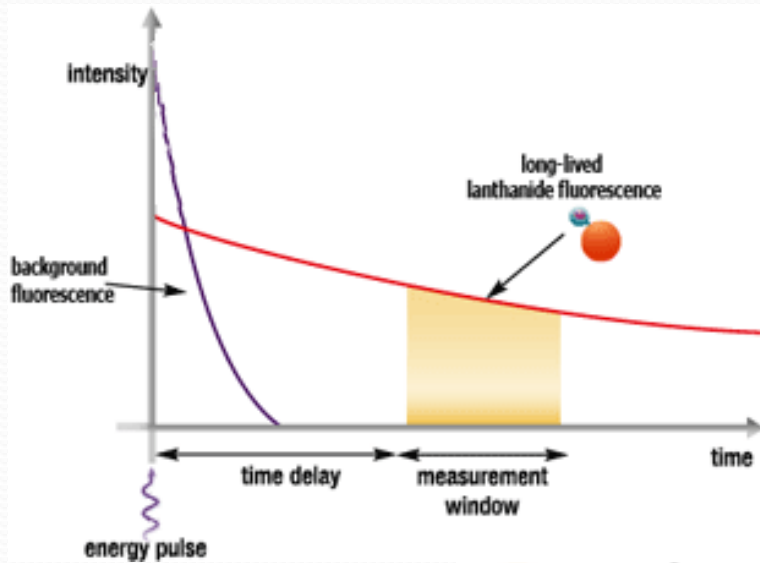
# HTRF: One technology, many uses in a pre-clinical laboratory setting.

SLAS 2012

# Outline of Discussion

1. Brief Review of HTRF Technology
2. Our Capabilities/Desires for the use of HTRF in drug development research.
3. Uses of HTRF Technology:
  - Binding Assay
    - Receptor Tyrosine Kinase ECD Dimer Competition Assay
    - Assay development discussion
  - Quantitation Assays
    - Cell line characterization
    - Serum samples
  - Cellular Activity Assay
    - pAkt and pERK evaluation
- Summary and Discussion of Additional Planned Uses of HTRF

# Homogenous Time Resolved Fluorescence:



# Benefits and Considerations of HTRF:

- Advantages:
  - Homogeneous: No plate binding, washing or detection development
  - Quick assay development.
  - Low matrix effects
  - Practical cost/well
  - Slow bleach rates of lanthanides allow multiple reads (kinetic runs)
- Disadvantages:
  - Requires a good set of reagents
  - Potential for “hook” effect

# General Methods:

- Biotek Synergy 2 Instrument

- Equipped with:

- 330 ± 40 Excitation Filter
    - 620 ± 5 Emission Filter
    - 665 ± 4 Emission Filter
    - Xenon flash lamp as excitation source

- Data Acquisition:

- Gen5 software
    - Sequential Mode: 665nm emission then 620nm emission
    - 100µs delay; 300µs integration
    - 10 – 20 reads/well
    - Gain set to highest expect signal source to maximize dynamic range

- Corning #3673 half-volume, 384-well white plates

- Reaction volumes 20µl

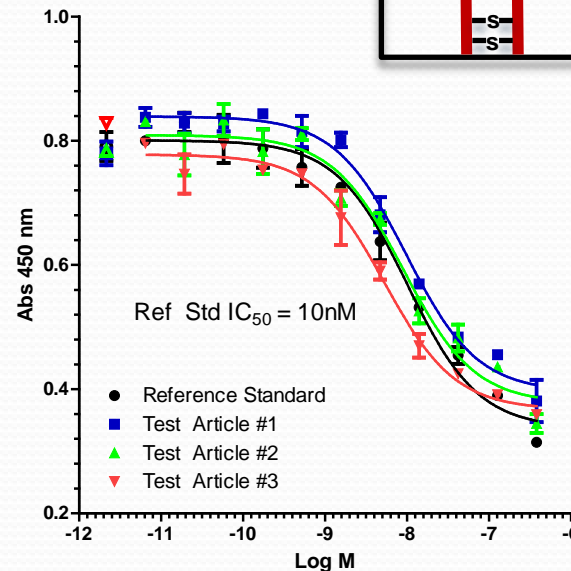
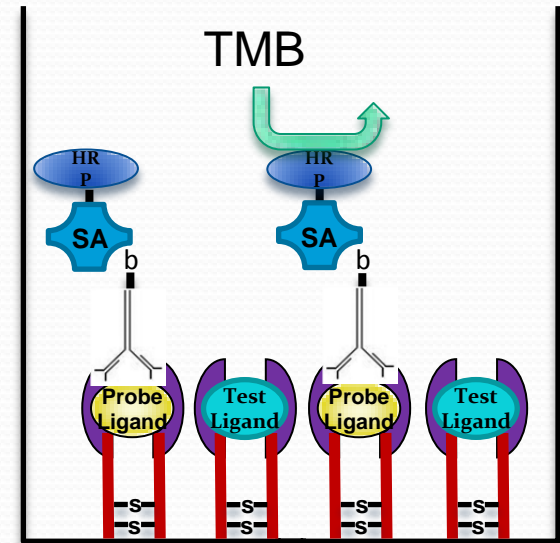
- Data analysis:

- $$\text{HTRF Ratio} = \frac{665_{em}}{620_{em}} \times 10,000 \text{ or } \Delta F\% = \frac{(HTRF_{Test} - HTRF_{Control})}{HTRF_{Control}}$$

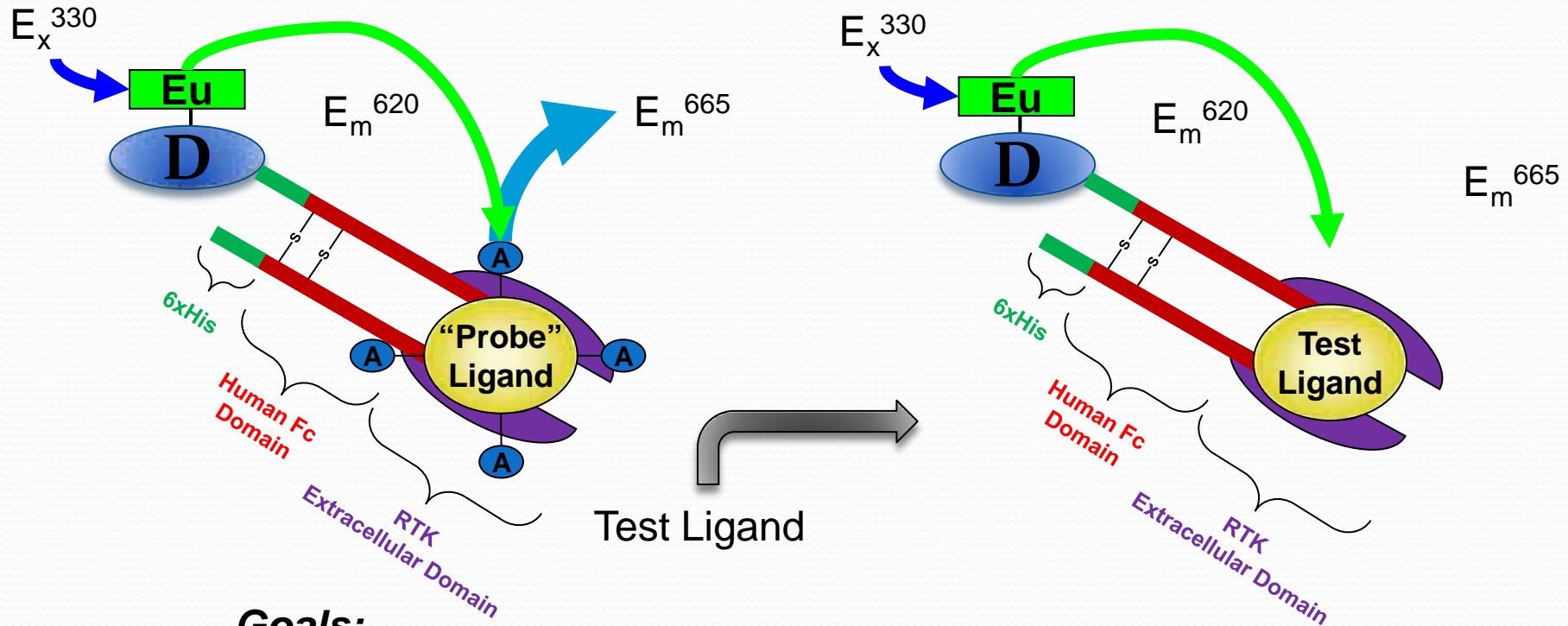


# Plate Bound Competition Assay:

1. Coat o/n with RTK-ECD (18h)
2. Wash
3. Block (1h)
4. Perform binding with probe/test ligands (1h)
5. Wash
6. Label with biotinylated detection antibody (1h)
7. Wash
8. Label with streptavidin-HRP (0.5h)
9. Wash
10. Incubate with TMB (0.5h)
11. Read in plate reader



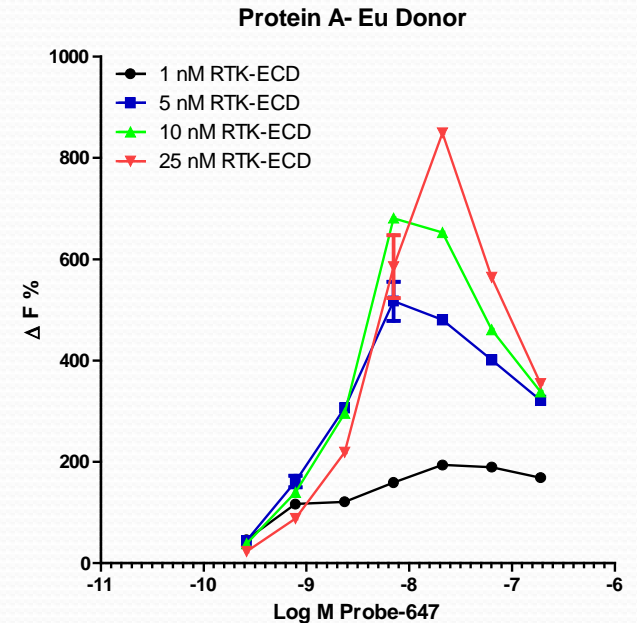
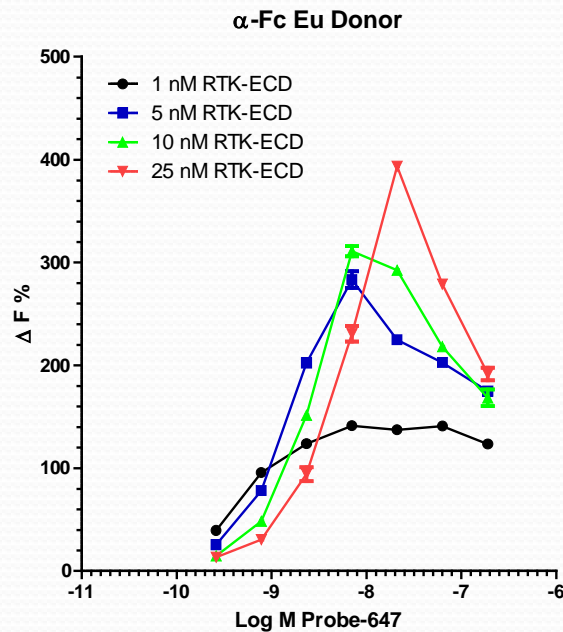
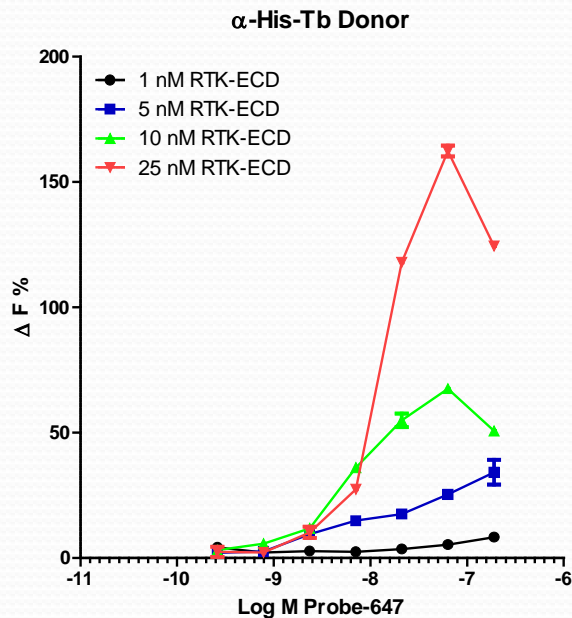
# HTRF Binding Assays: Competition Format



## Goals:

1. True estimate of  $IC_{50}$  of test ligands
2. Non-radioactive
3. Low (<10%) ligand depletion conditions
4. Works within the pM – nM range
5. Highly reproducible
6. Measure dozens/hundreds of variants at a time

# Identification of Donor Molecule:

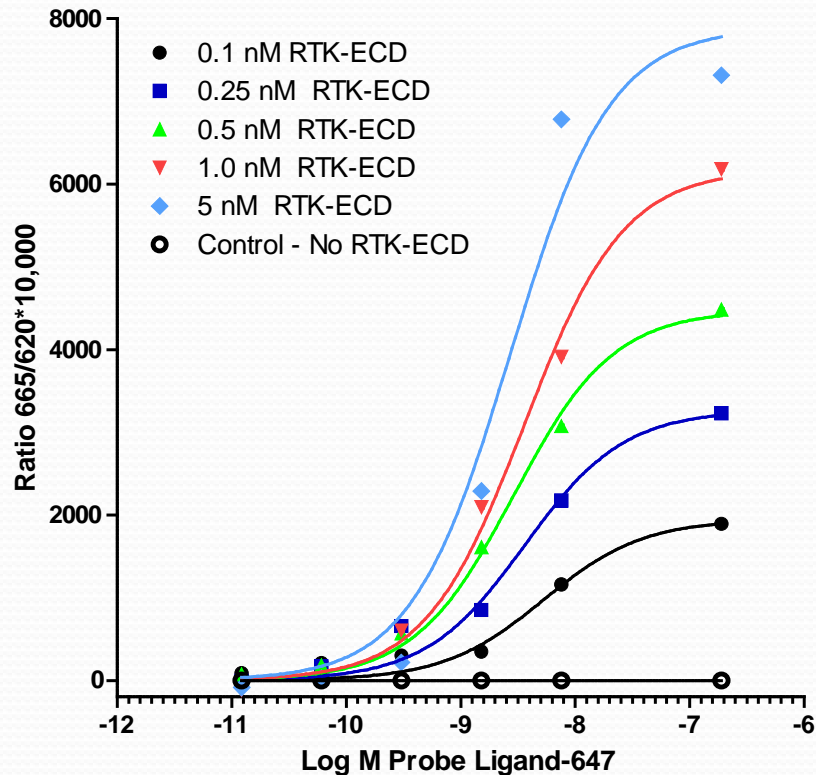


- Alexa-647 labeled Probe Ligand
- Donors at 1:400
- 2.0 h Incubation with all components

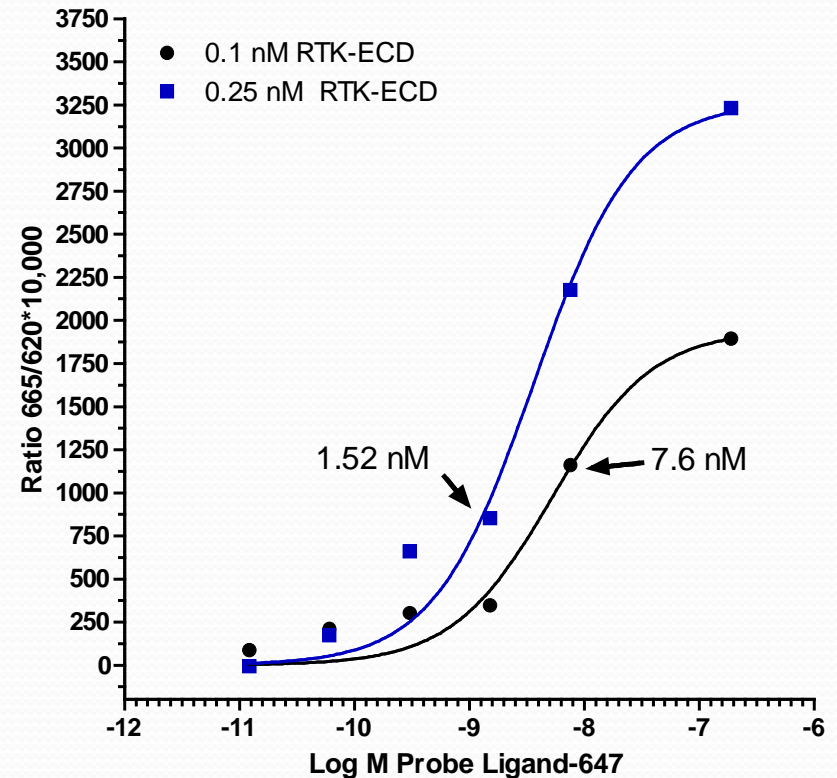


# Determining Practical Concentrations of RTK-ECD and Probe Ligand:

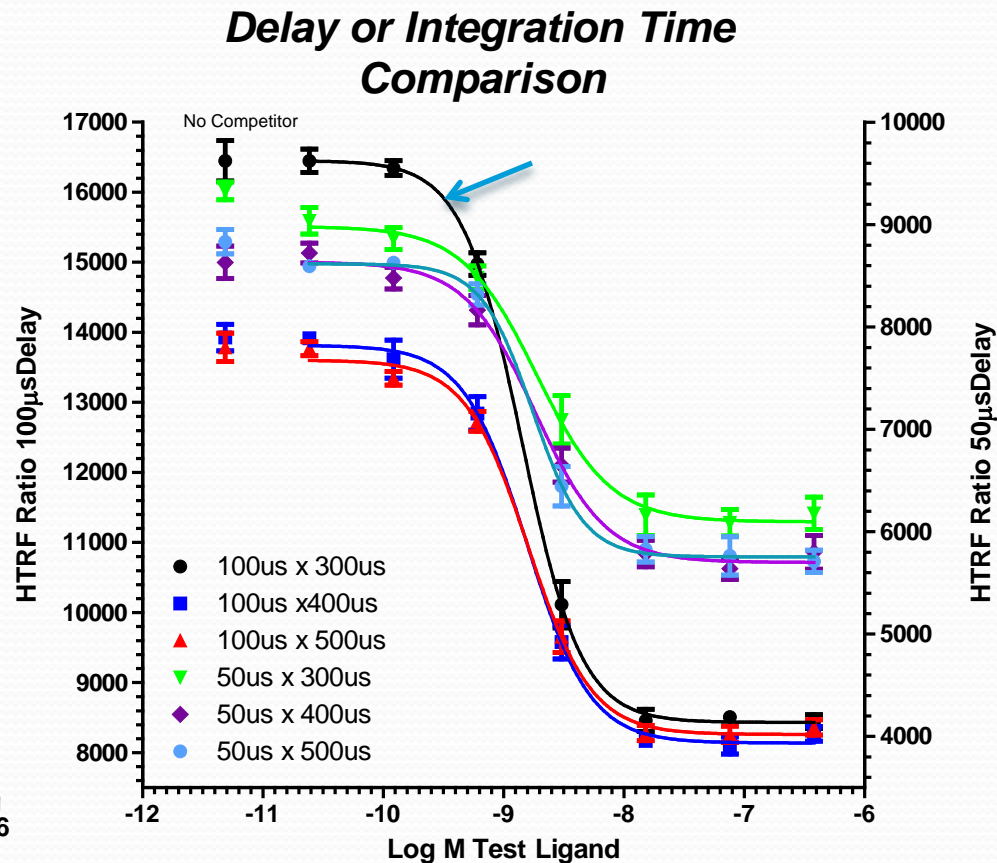
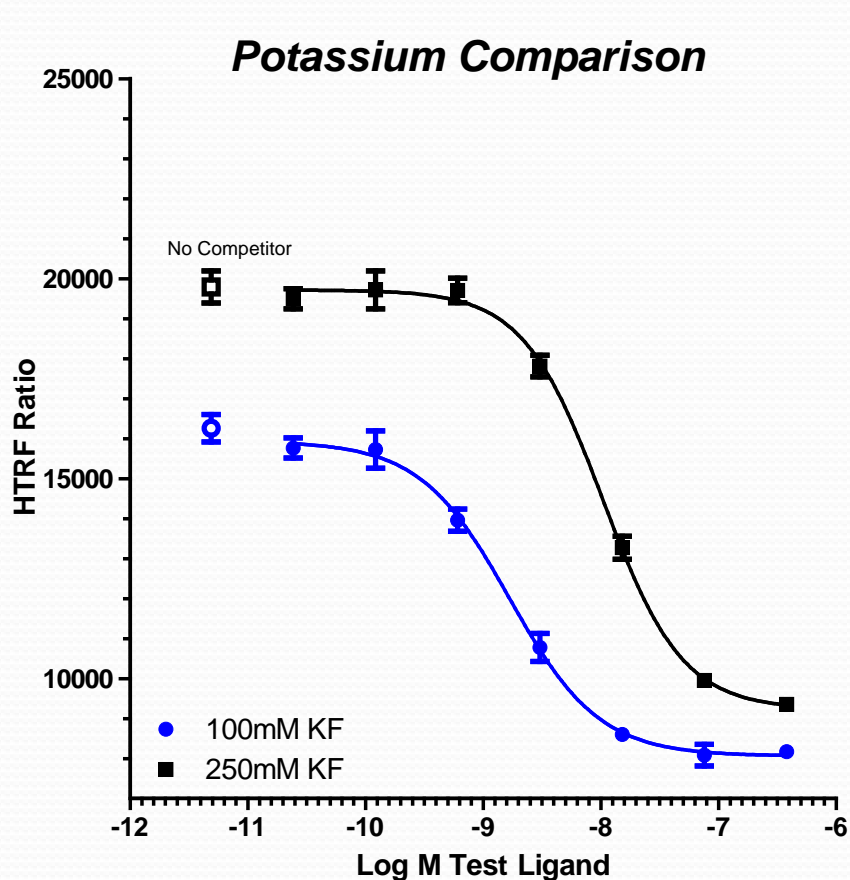
RTK Binding Assay  
Detection anti-Fc-Eu + Ligand-647



RTK Binding Assay  
Detection anti-Fc-Eu + Ligand-647



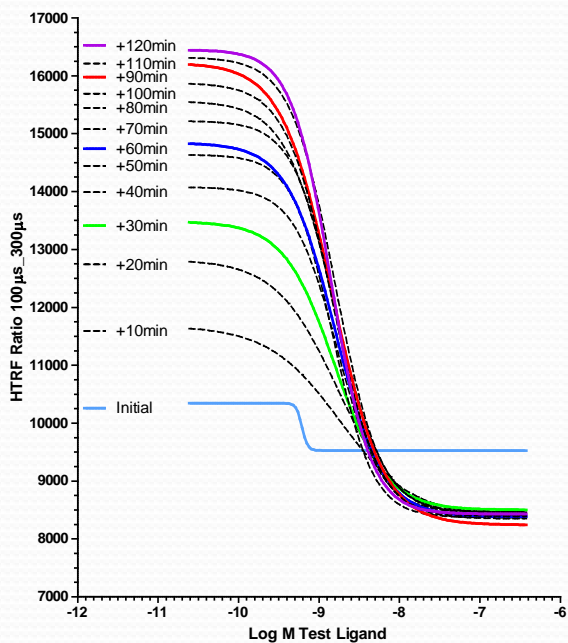
# Effects of KF and Read Times on Assay Performance:



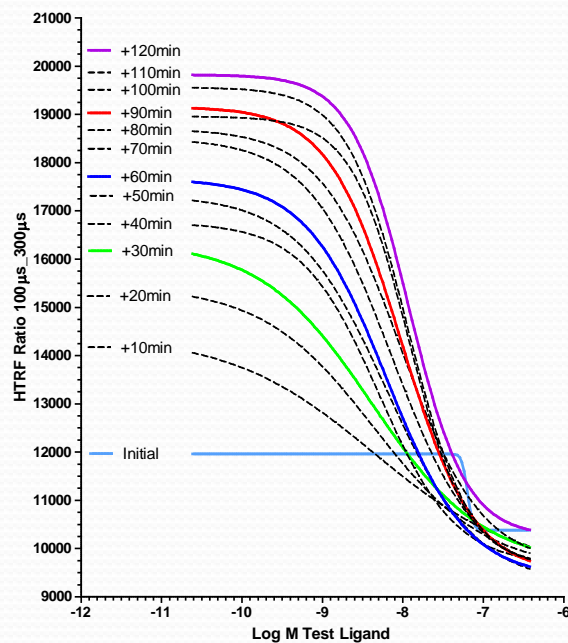
- biotinylated Probe Ligand (7.6nM) + 0.1nM RTK Dimer
- Strept-D2 + anti-Fc-Eu (1:400)
- 2.0 h Incubation with all components

# Time Course of Binding Equilibrium

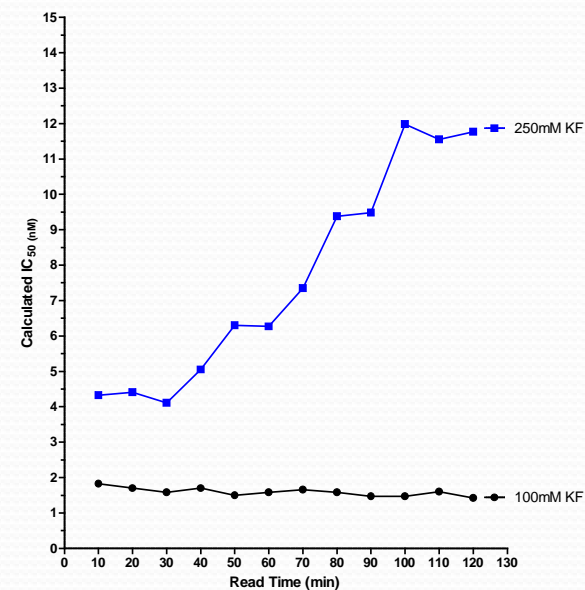
## Time Course 100mM KF



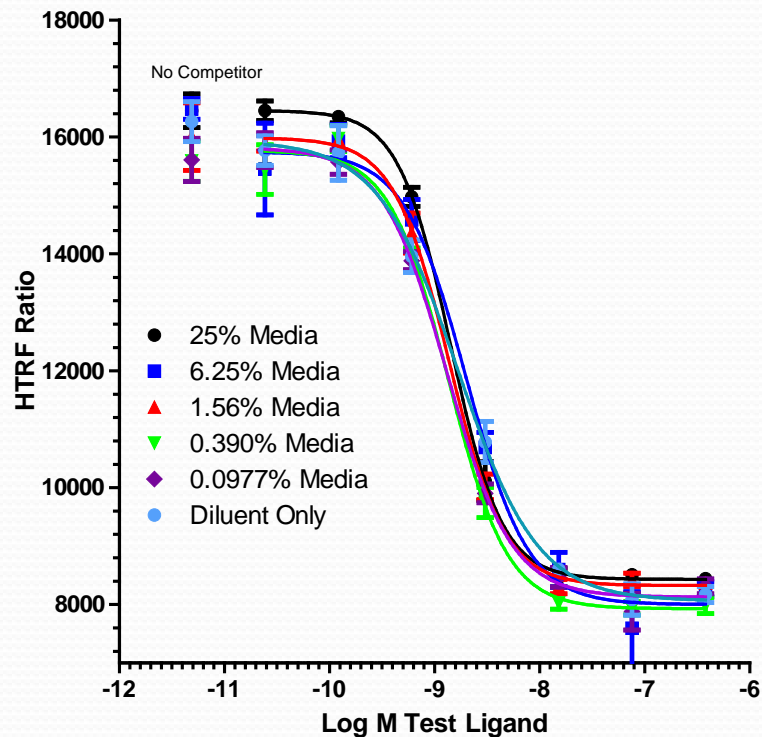
## Time Course 250mM KF



## Calculated $IC_{50}$ s



# Effect of Matrix Concentrations on Assay Performance:

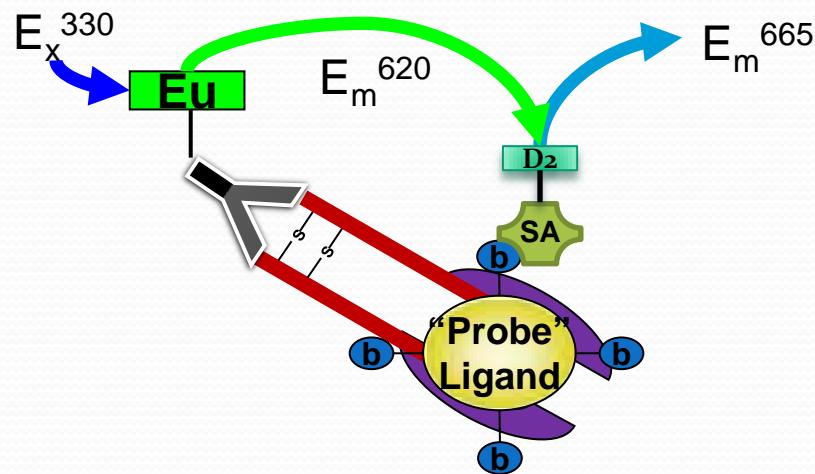


Column1	25% Media	6.25% Media	1.56% Media	0.390% Media	0.0977% Media	Diluent Only	Mean	Standard Dev
Bottom	8431	8000	8324	7928	8128	8063	8145.7	194.4
Top	16450	15745	15984	15770	15816	15934	15949.8	262.2
Log IC50	-8.847	-8.697	-8.861	-8.867	-8.882	-8.775	-8.8	0.07
HillSlope	-1.762	-1.464	-1.639	-1.546	-1.415	-1.141	-1.5	0.21
IC50	1.4E-09	2.008E-09	1.4E-09	1.357E-09	1.31E-09	1.7E-09	1.53E-09	2.70E-10
Span	8019	7745	7660	7842	7688	7871	7804.2	134.0

# Summary of HTRF RTK-ECD Competition Assay

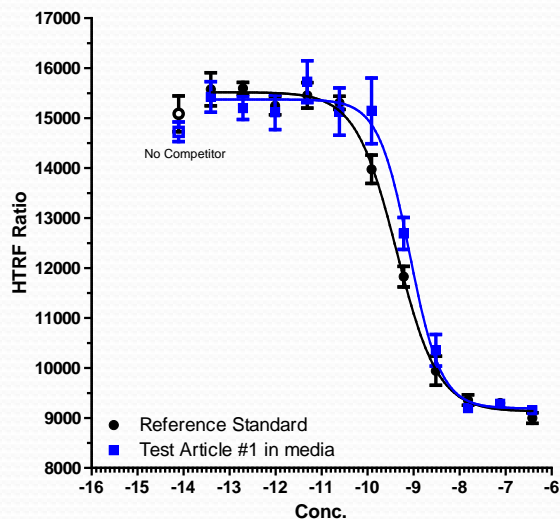
## Procedure:

1. RTK-ECD at 4x (5 $\mu$ l/well)
2. Reference Standard/Test Article in diluent + Probe Ligand at 2x (10 $\mu$ l/well)
3. Detection Reagents in diluent + KF at 4x (5 $\mu$ l/well)
  - StreptAv-D2 and anti-Fc-Eu
4. Read with HTRF settings

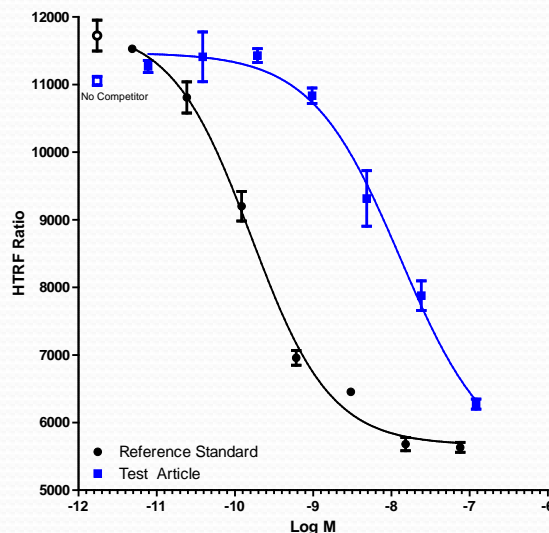


# HTRF Competition Assay in Practice:

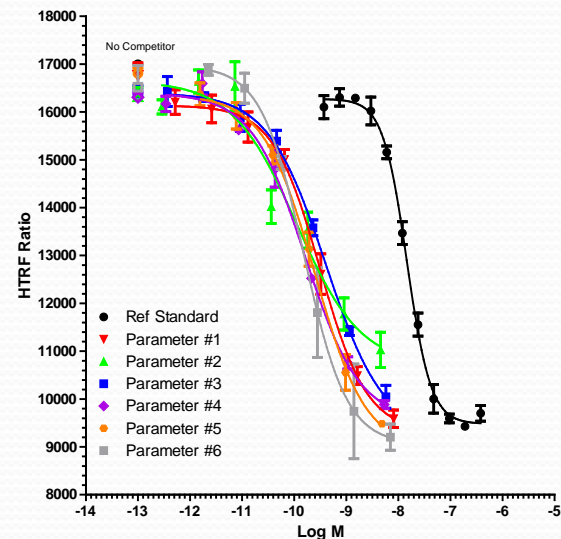
## Two Related Molecules



## Two Related Molecules



## Same molecule, different production parameters



**Plate Bound Assay**

Reagent	Cost	Amt	cost/well
RTK-ECD	\$200	50 ug	\$0.20
anti-Probe Abs	\$380	50 ug	\$0.02
Probe Ligand	\$295	10 ug	\$0.06
Strept-Av-HRP	\$695	0.5 ml	\$0.01
High Bind Plates	\$242	100 each	\$0.03
TMB	\$135	600 ml	\$0.02
<b>Total</b>			<b>\$0.34</b>

**HTRF Assay**

Reagent	Cost	Amt	cost/well
RTK-ECD	\$200	50 ug	\$0.02
Strept-D2	\$793	1 ml	\$0.04
anti-Fc-EU	\$790	1 ml	\$0.04
White Plates	\$900	100 each	\$0.09
<b>Total</b>			<b>\$0.19</b>

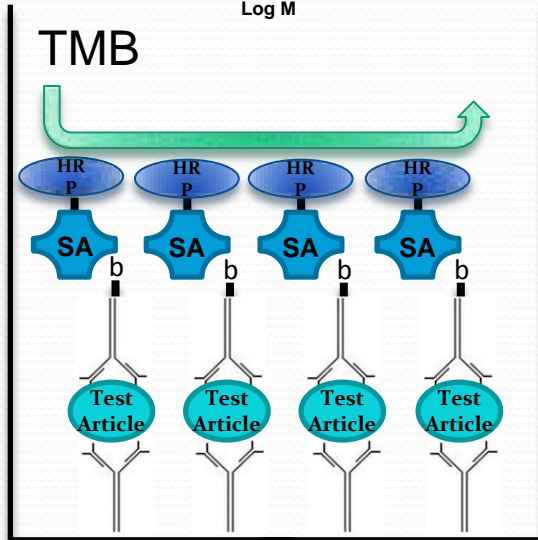
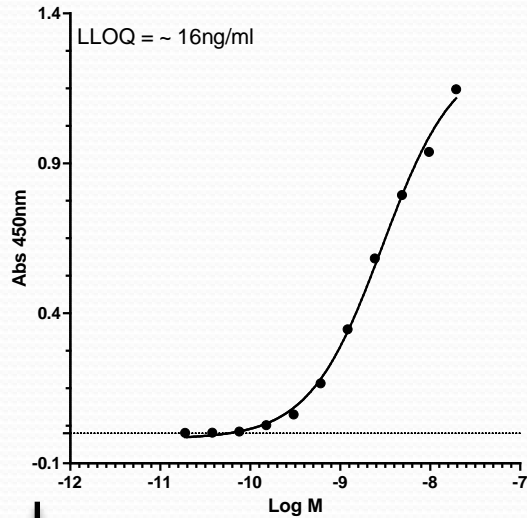
# HTRF Quantitation Assays:

- Applications:
  - Evaluating production amounts in conditioned media.
    - Both transient and clonal cell lines
  - Identifying best producing cell clone in cells under selection
  - Evaluating drug product present in biological test samples.

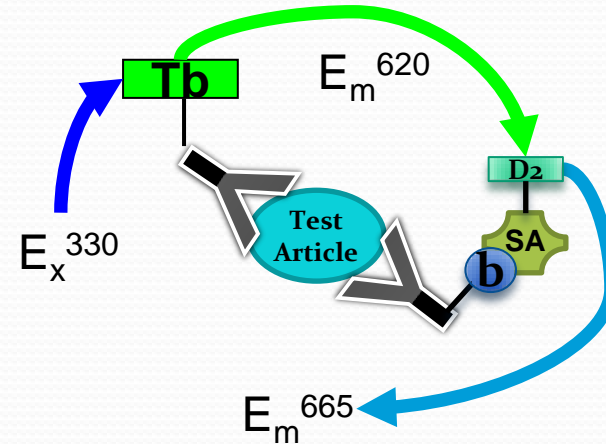
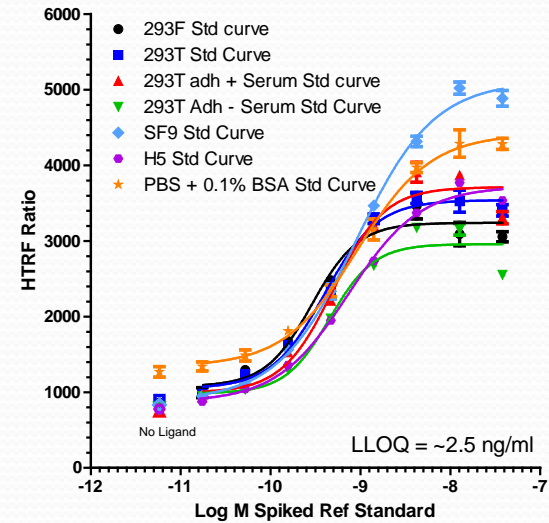
# Examples of Quantitation Assays:

## Sandwich Format

### Sandwich ELISA

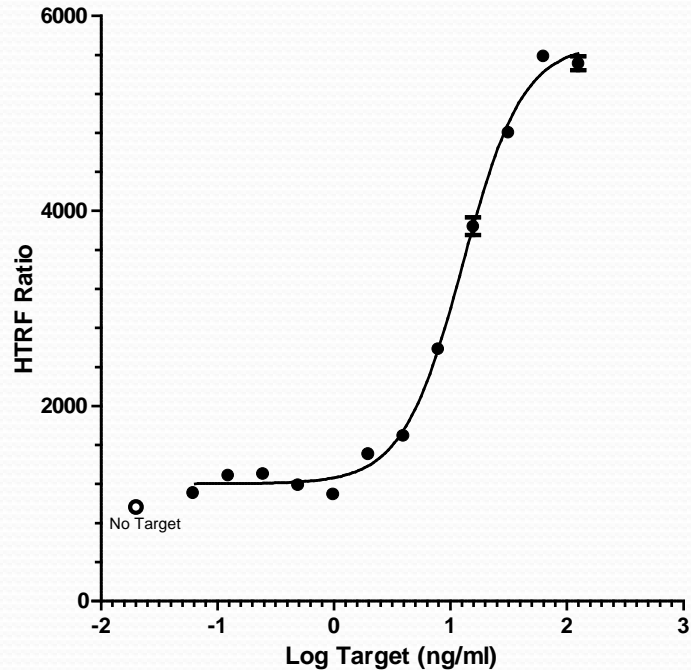
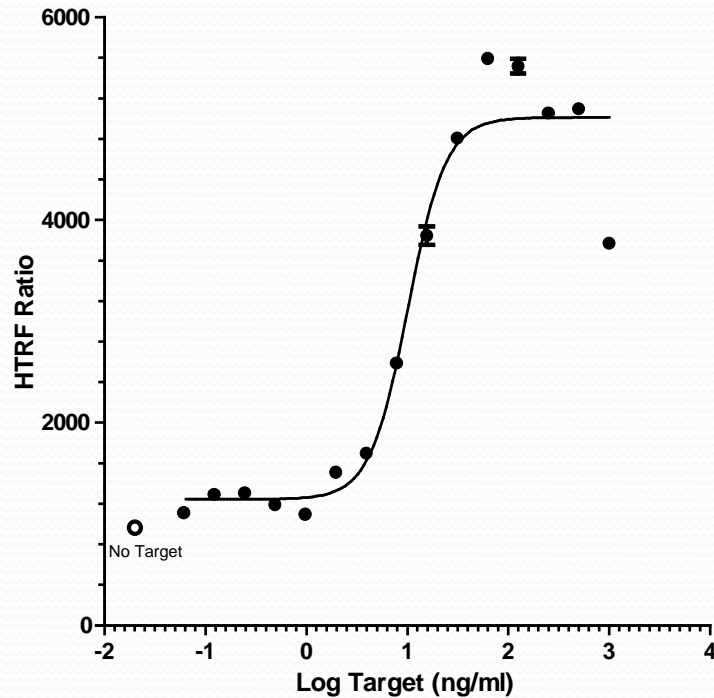


### Sandwich HTRF





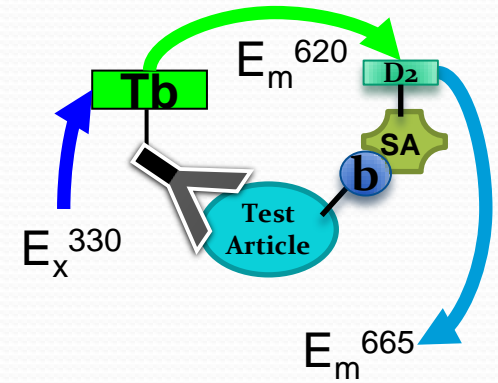
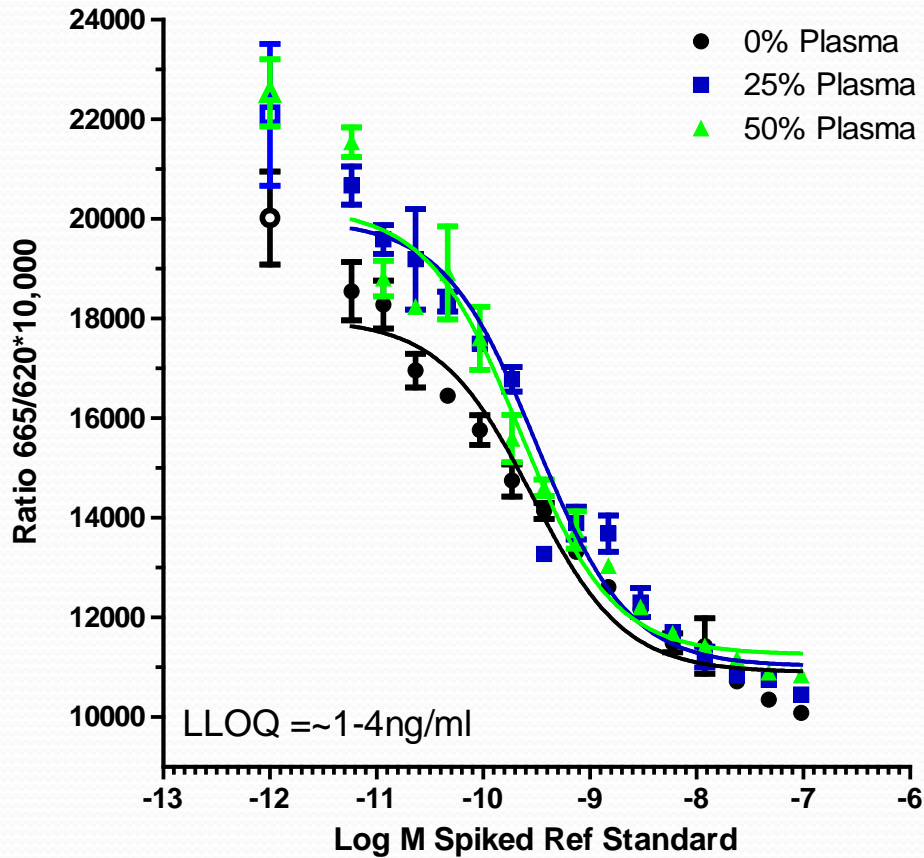
# Considerations with Quantitation:



Multiple dilutions of a test sample must be tested.

# Examples of Quantitation Assays:

## Competition Format



# Cellular Activity Assay:

## tAkt/pAkt-Ser<sup>473</sup> Cell Based Detection Kit

## pAkt-Ser<sup>473</sup> or Cellul'ERK-Cisbio

Activate 5-30 min

Fix 20 min

Wash (3 x 5min)

Quench 20 min

Wash (3 x 5min)

Block 1h

Wash (3 x 5min)

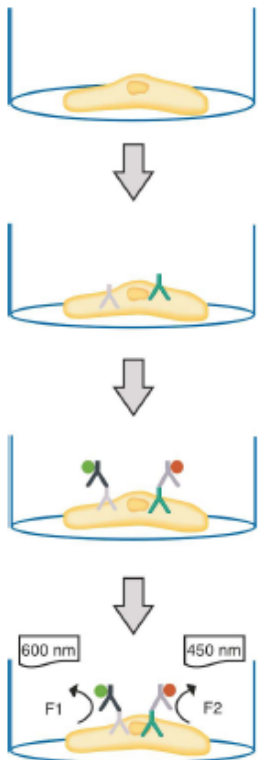
Label 1<sup>o</sup> 16h

Wash (3 x 5min)

Label 2<sup>o</sup> 2h

Wash (4 x 5min)

Develop ~30min



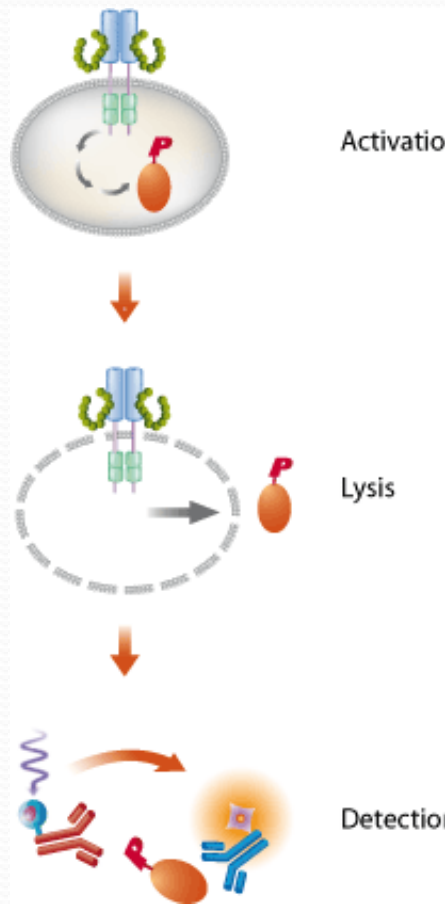
1. Seed cells in a 96 well plate. Stimulate cells with ligands. Fix, permeabilize, and block cells.

2. Add primary antibodies (rabbit anti-phospho-Akt (S473) and mouse anti-total Akt).

3. Add secondary antibodies (HRP-conjugated anti-rabbit IgG and AP-conjugated anti-mouse IgG).

4. Add fluorogenic substrates F1 and F2 and measure fluorescence.

~\$4.70/well



Activation 5-30 min

30 min

Freeze Lysates

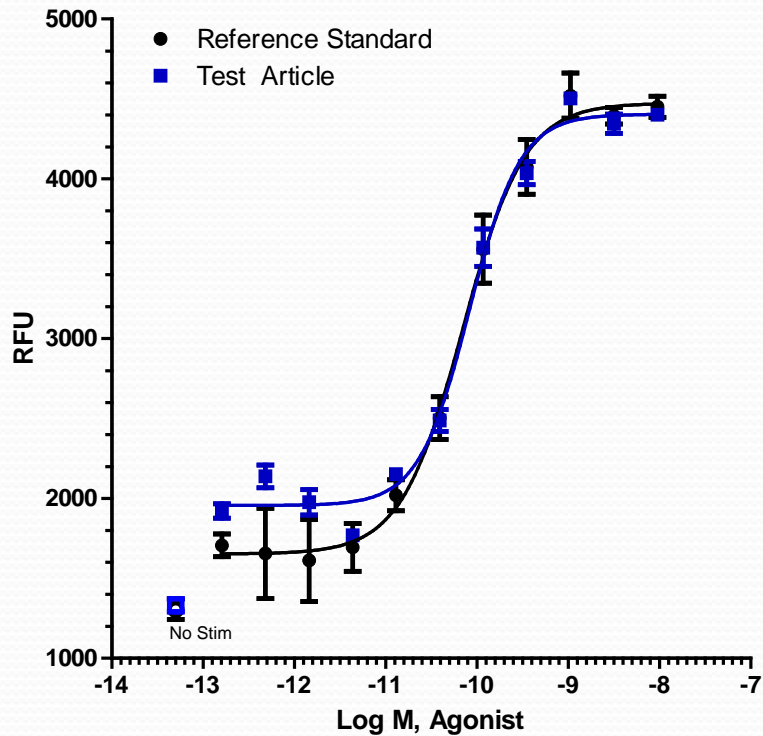
~3h

Detection

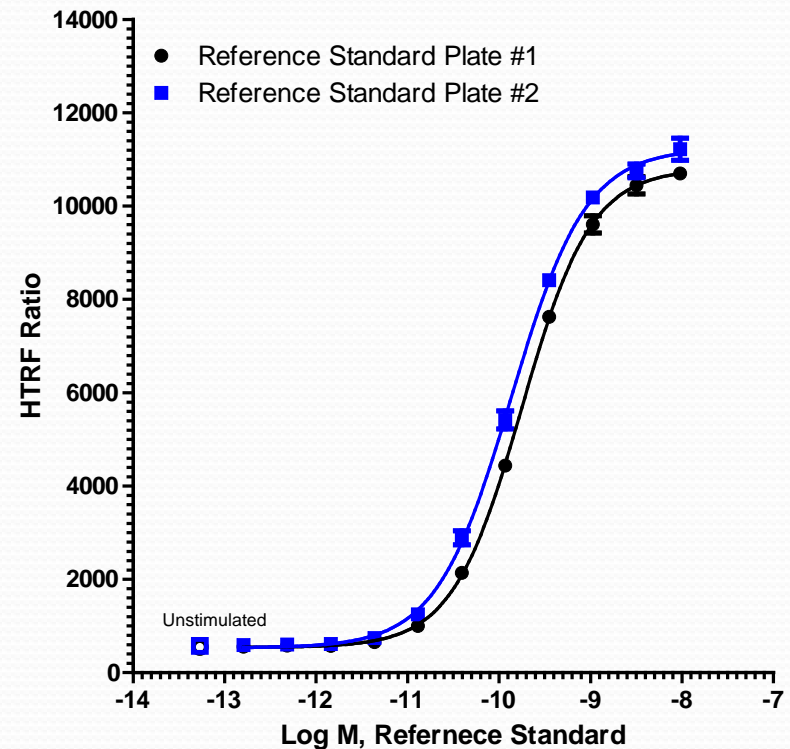
~\$0.70/well

# Comparison of Assay Performance:

## pAkt-Ser<sup>473</sup> Cell Based

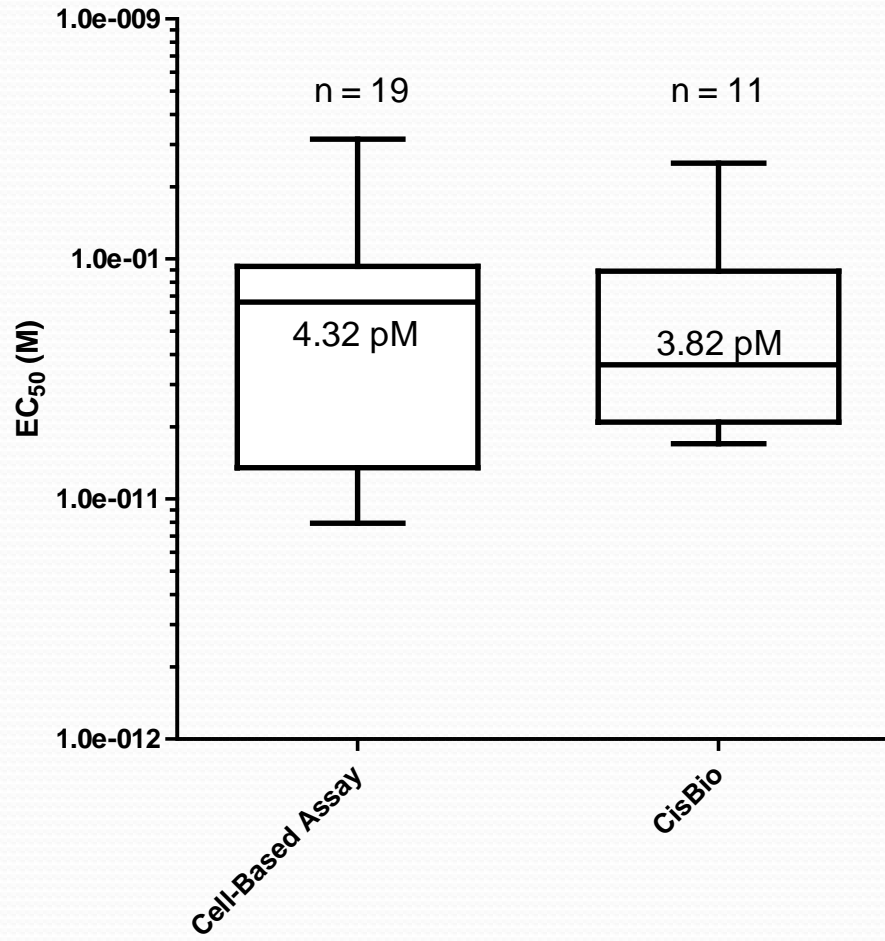


## pAkt-Ser<sup>473</sup> Cisbio



- Both Assays went through standard optimization parameters
  - Cell number
  - Stimulation time
  - Lysis volume (Cisbio)

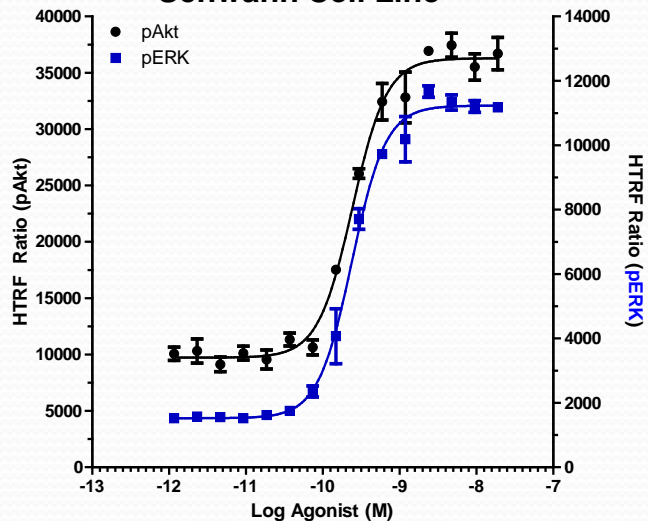
# Comparison of Assay Performance: pAkt-Ser<sup>473</sup>



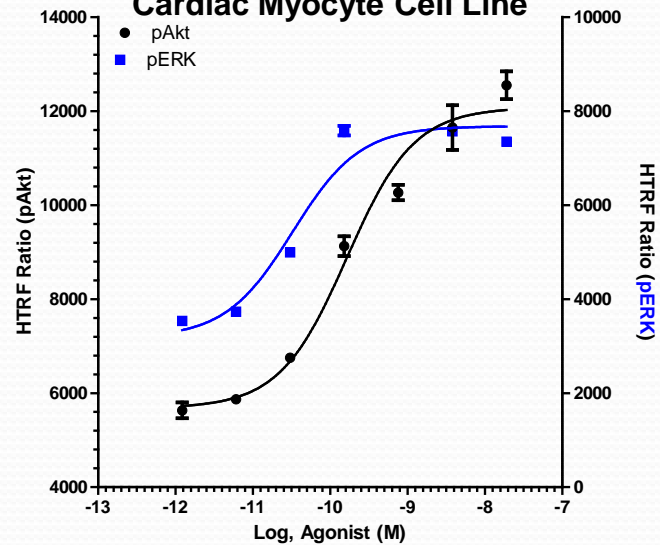
# Activity Assays:

## Comparison of pAkt-Ser<sup>473</sup> and Cellul'ERK in Multiple Cell Lines

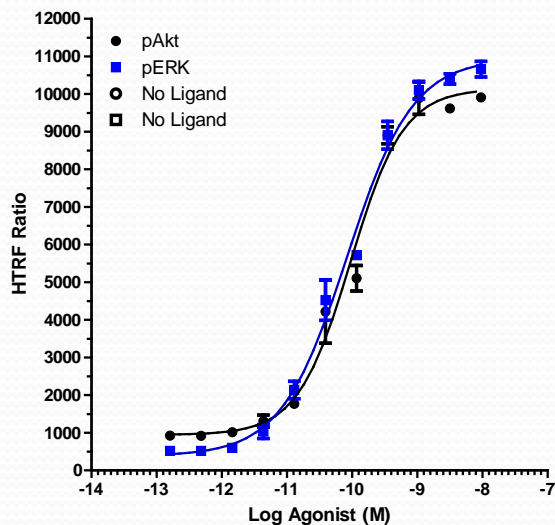
### Schwann Cell Line



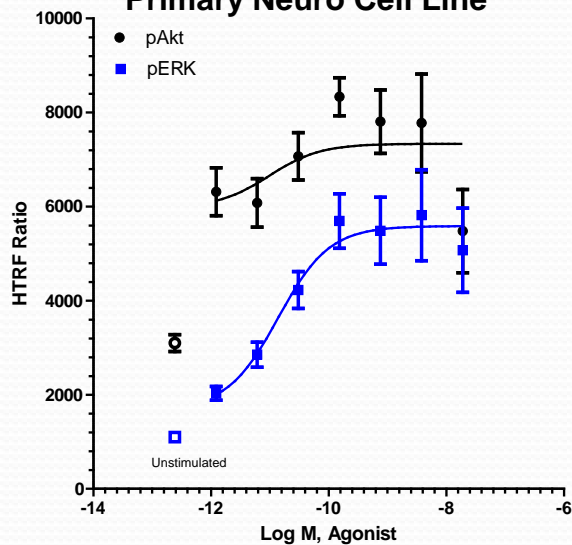
### Cardiac Myocyte Cell Line



### Breast Cancer Cell Line



### Primary Neuro Cell Line



# Summary

- HTRF is a robust methodology with several key benefits:
  - Time savings
    - Day-to-day
    - Assay development
  - Cost savings
    - Reagent
    - Personnel time
  - Equal if not better assay performance than standard assay formats
  - Flexible conditions
- Other uses:
  - Screening for novel molecules
  - Multiplexed assay
    - Using two acceptor assay systems
  - Moving assays into qualification/GXP studies
  - ATA Assay???