



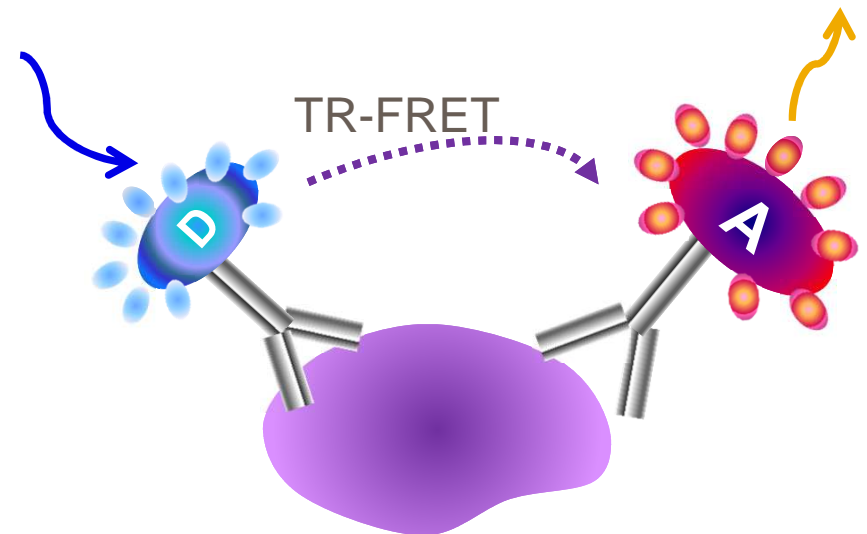
## Cellular High Throughput Screening assays using HTRF detection of cytokines

HTRF in Drug Discovery Symposium

April 2013

Lorena Kallal

GlaxoSmithKline, Collegeville, PA USA

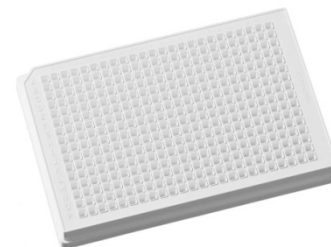


# GlaxoSmithKline, Upper Providence (Collegeville), Pennsylvania, USA

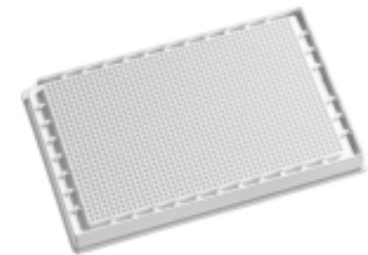


- Assay development and screening environment
- 384 and 1536 well cell based assays

**384 well**



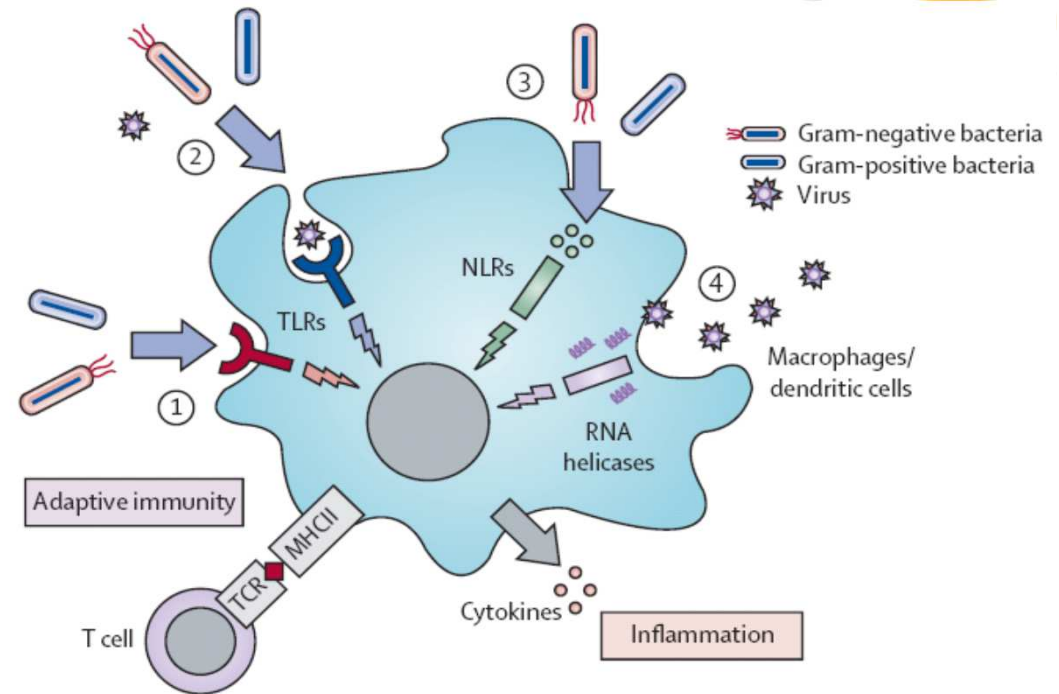
**1536 well**



# Pattern Recognition Receptors: immunity

NOD, Toll, TLR, NLRPs, etc.

Activated by pathogenic & endogenous danger signals



Stimulate:

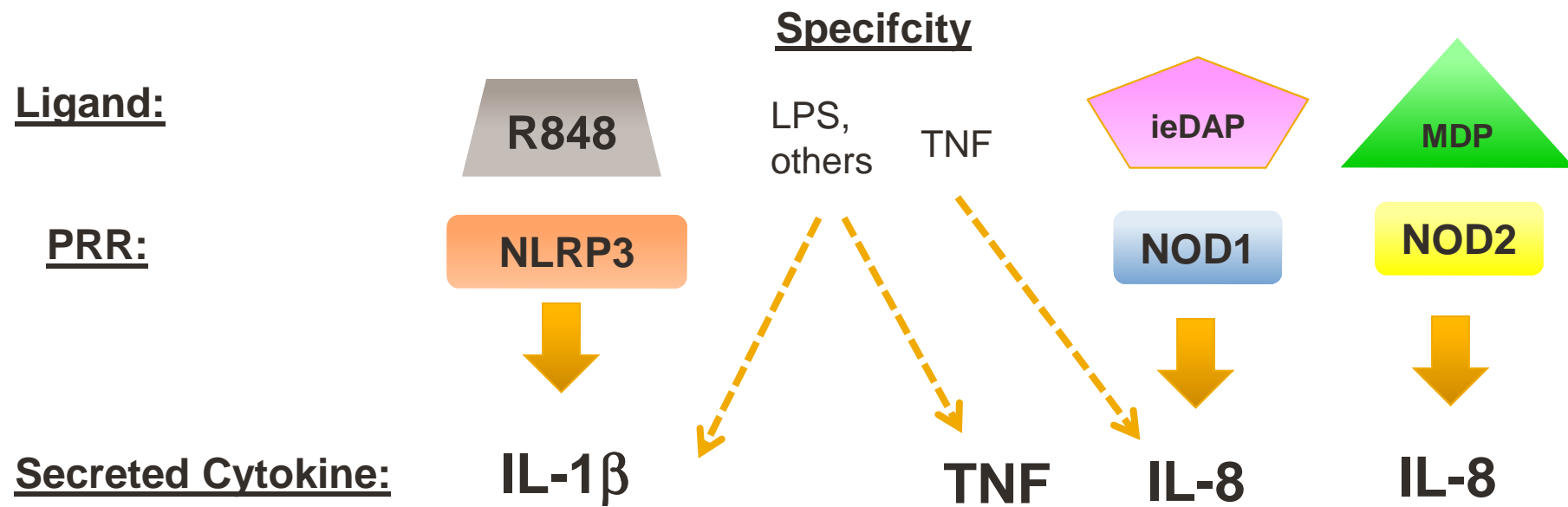
–**Cytokine production**

–inflammatory signal transduction cascades

Therapeutics:  
inflammatory diseases

# HTRF assays for the detection of cytokines

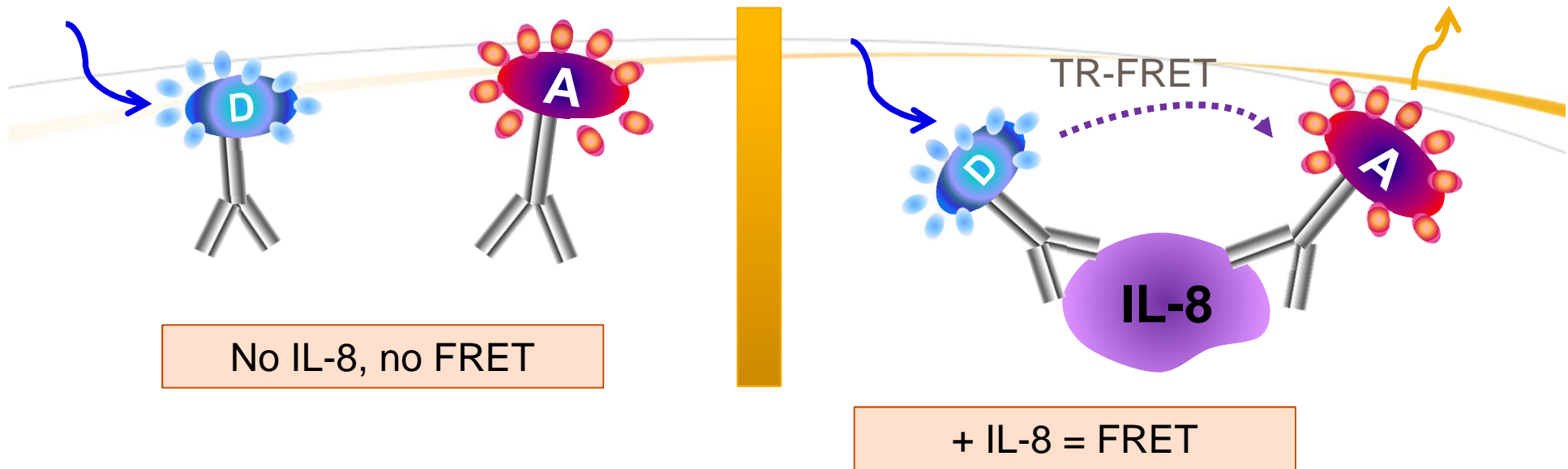
- Interested in NOD1, NOD2, and NLRP3 signaling
- Pathway specificity: different activators, different cytokines



**Cell system:** PMA differentiated THP1 cells:

Engineered HEK cells, stably expressing NOD1 or NOD2:

## Assay development strategy: NOD1 and NOD2 HTSs:



- Goal: Establish 1536-well homogeneous cell based assays
- Also: Will assay-ready frozen cells work for the assay?

## Assay development for high throughput screen (HTS):

HTS: ~2 M compounds; @10  $\mu$ M

1536 well plates

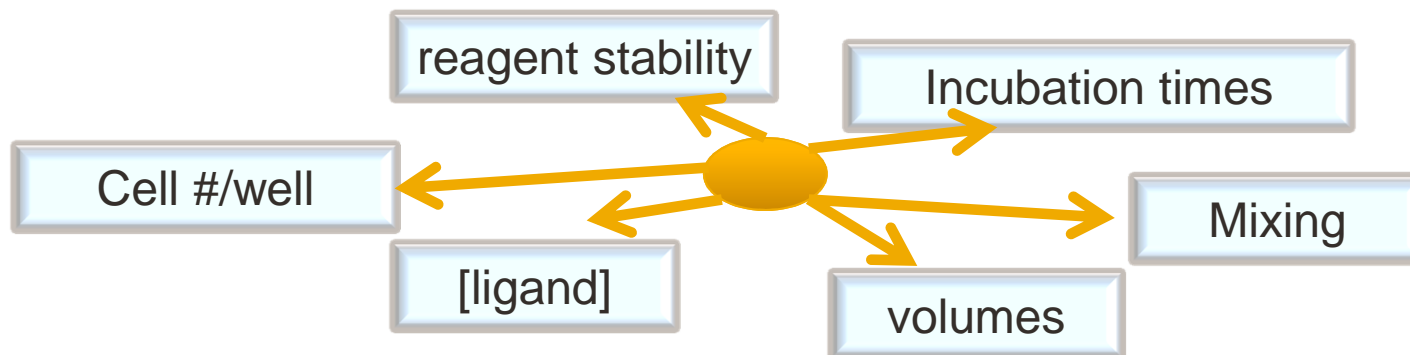
XC50 curves; primary assay + specificity assay

384 well plates

Other specificity assays and downstream assays

→ tox, off target...

- Test many conditions during assay development: will share only examples

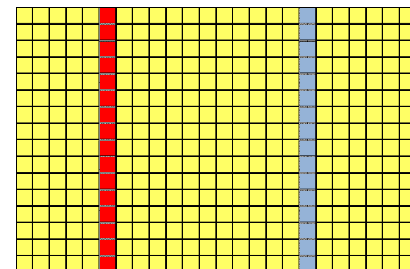


### ■ Statistics

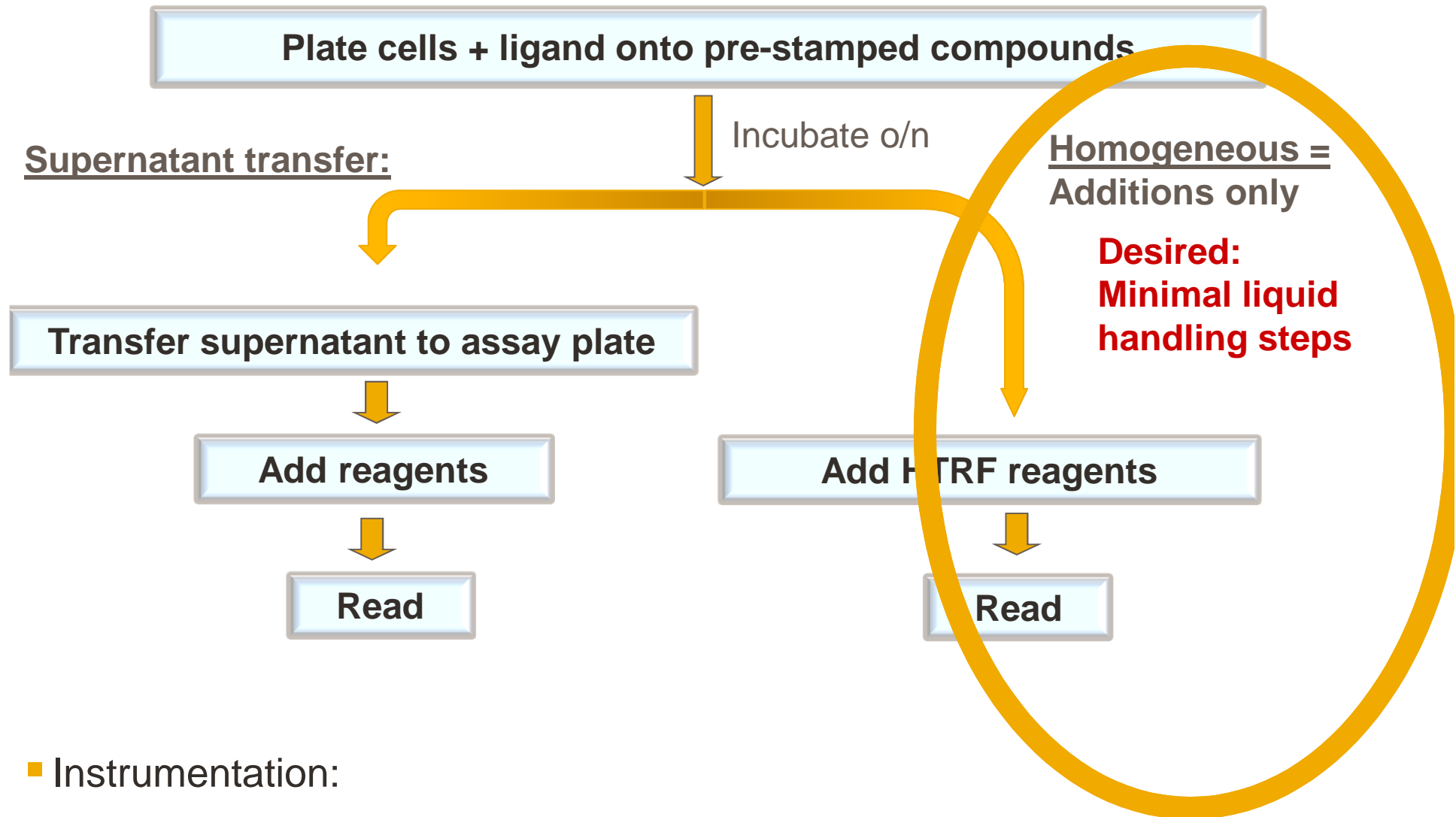
—Signal / background (**S/B**)

—**Z'** gives “variability”: want  $\geq 0.4$ ; 1 = perfect

Control columns



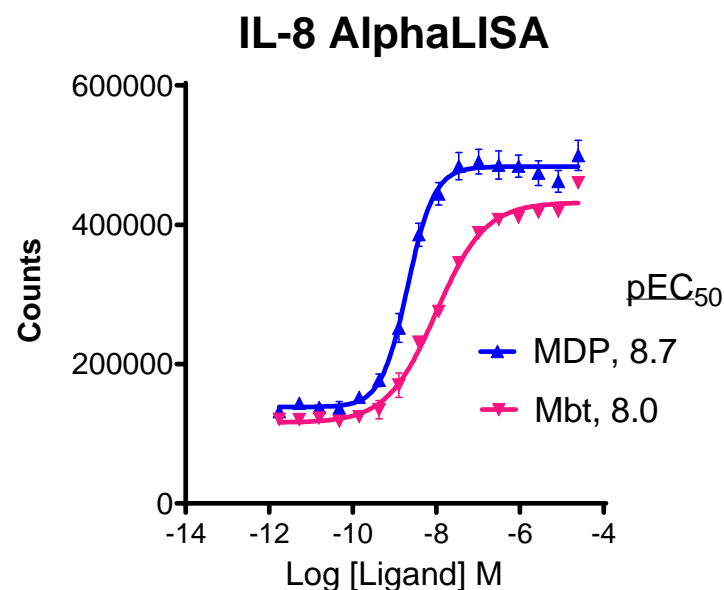
Flow chart: supernatant transfer vs homogeneous assays:



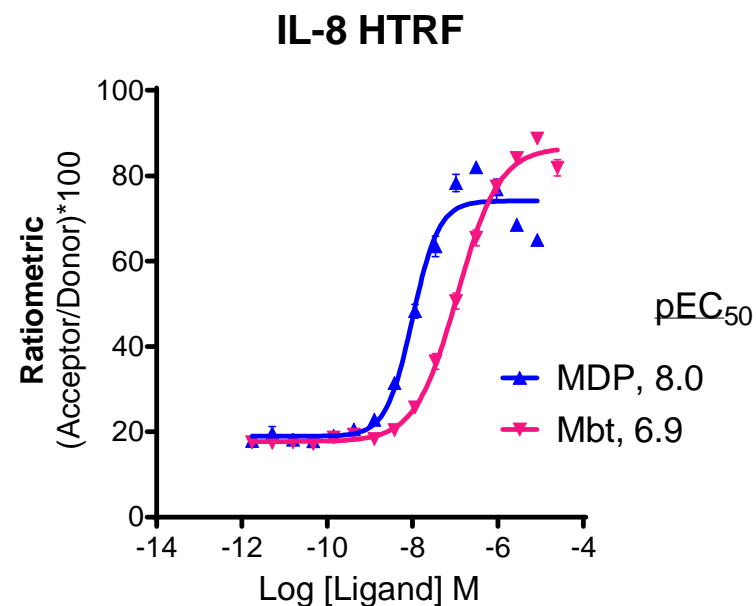
■ Instrumentation:

- Envision: PMT-based, sensitive, ~ 7 – 35 min per plate
- Viewlux: CCD camera based, less sensitive, ~ 2 - 4 min per plate); ☺

## NOD2 IL-8: AlphaLISA (Perkin Elmer) vs HTRF comparison



Ligand	S/B	Z'
MDP	2.3	0.45
MBt	2.2	0.33

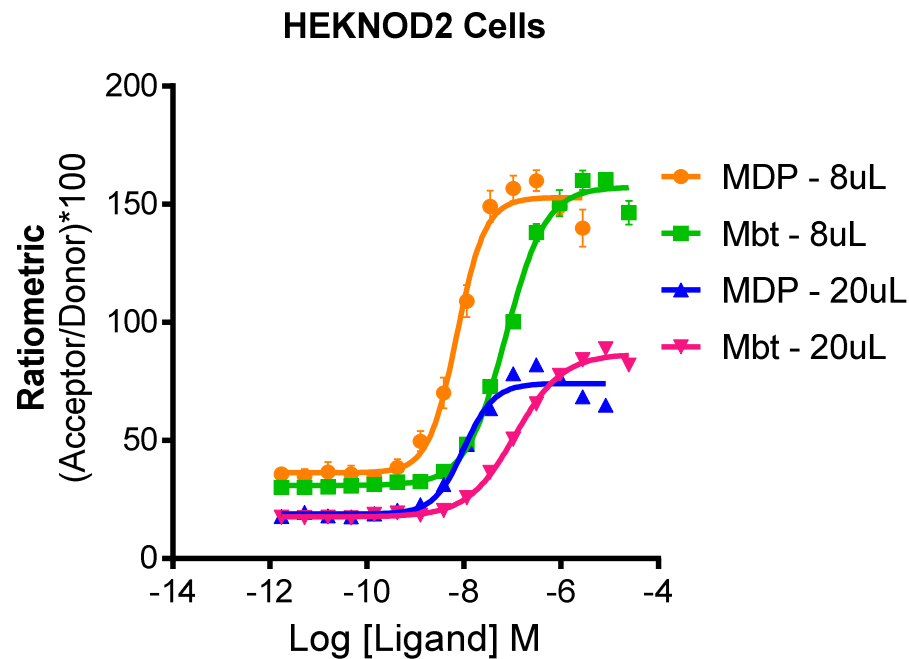


Ligand	S/B	Z'
MDP	2.9	0.56
MBt	2.9	0.57

- AlphaLISA :higher pEC<sub>50</sub>s, lower Z'
  - more steps, reagents less stable
  - Envision: longer read time (7 vs. 2 min on a Viewlux)
- HTRF assay was chosen for screening**

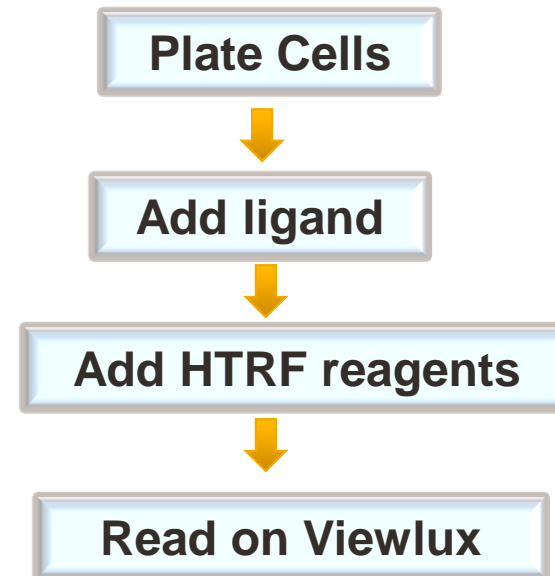


## NOD2 HTRF Assay: miniaturization: 8 $\mu$ L vs. 20 $\mu$ L (low vol 384):



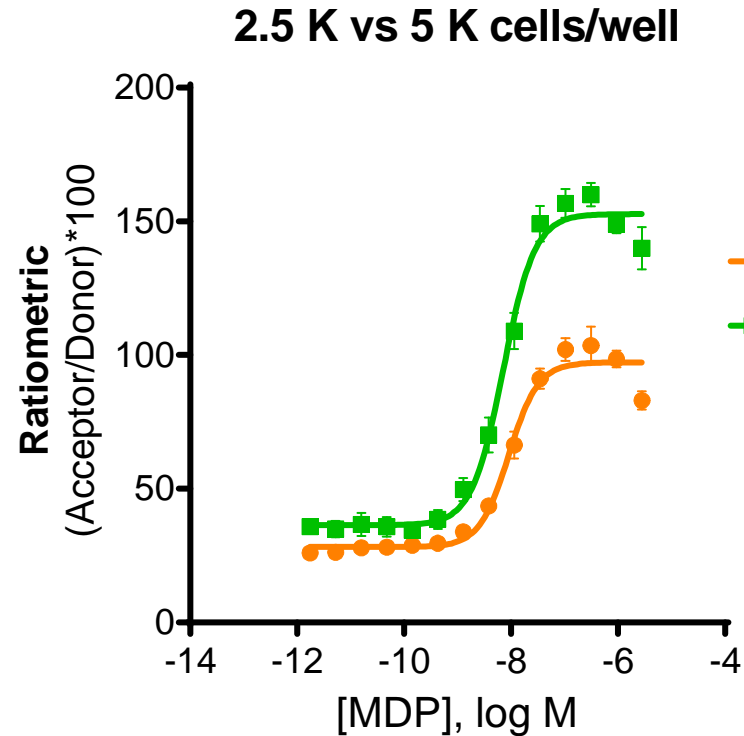
ligand	volume	S/B	Z' (16/16)
MDP	8	3.3	0.58
MDP	8	3.4	0.55
Mbt	20	2.9	0.56
Mbt	20	2.9	0.57

### Assay flow chart: "homogeneous"



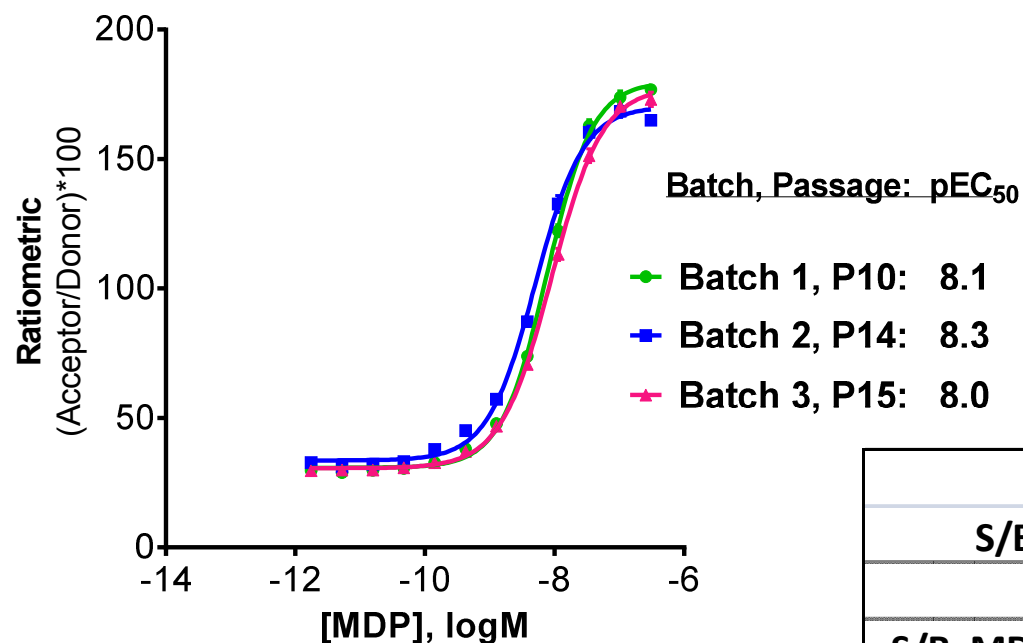
- S/B higher in 8  $\mu$ L than 20  $\mu$ L assay
- EC<sub>50</sub>s and Z' comparable → low vol assay OK → good for 1536

## NOD2 Cell Density (#cells/well) optimization :



- 5 K per well gave slightly higher S/B and Z' values than 2.5 K

## NOD2: Assay-Ready Frozen cells were acceptable:

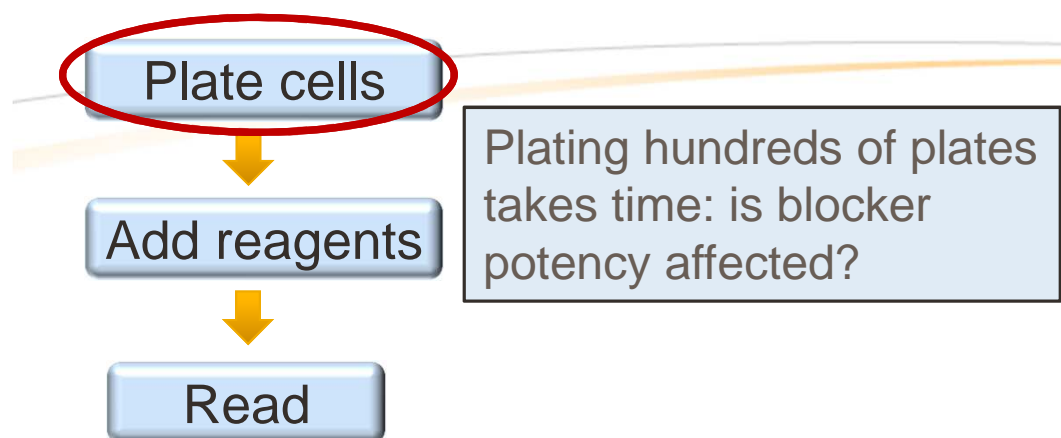


### S/B and Z' statistics:

	Batch 1	Batch 2	Batch 3
S/B: MDP / DMSO	4.4	4.3	4.2
Z'	0.60	0.75	0.59
S/B: MDP / MDP + blocker 1	4.6	4.3	4.5
Z'	0.64	0.74	0.62
S/B: MDP / MDP + blocker 2	6.3	6	5.6
Z'	0.66	0.78	0.64

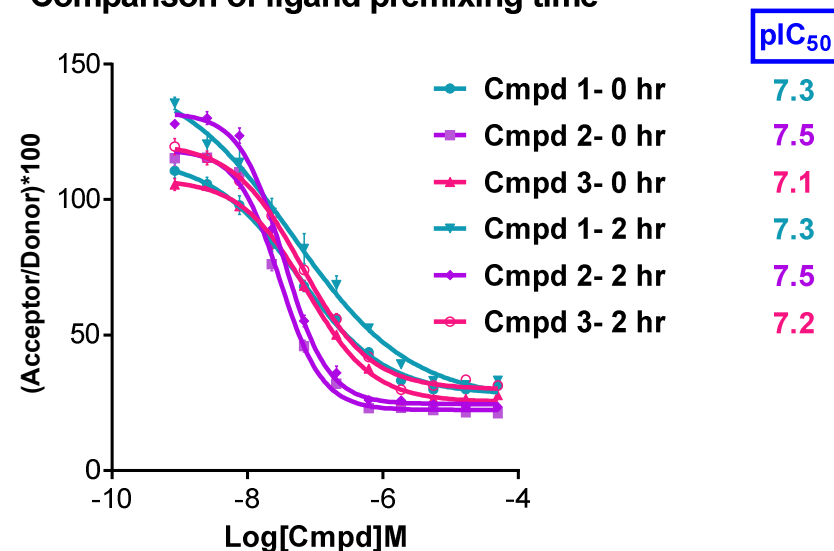
- Similar and consistent responses were obtained with assay-ready frozen cells

## Addressing cell plating lag time:



30nM EC80	Plating time			
	0 hr		2hr	
	S/B	Z'	S/B	Z'
DMSO	4.0	0.72	4.5	0.73
Blocker 1	3.9	0.71	4.1	0.74
Blocker 2	4.8	0.74	4.8	0.76
Blocker 3	3.8	0.70	3.4	0.70

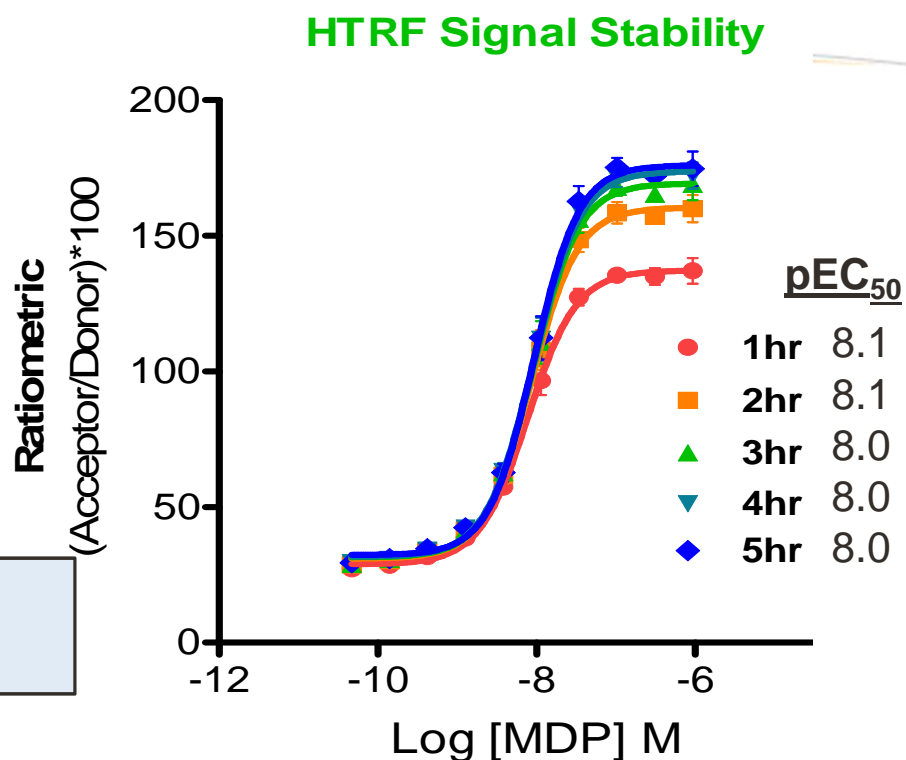
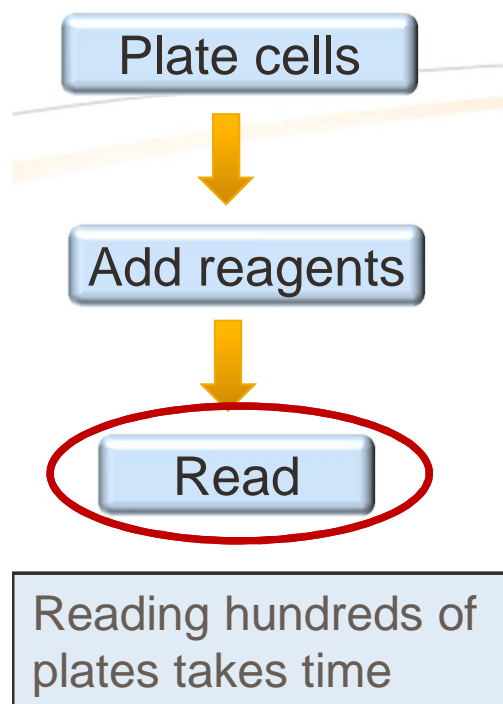
## Comparison of ligand premixing time



Low volume 384

- Blocker potencies did not change with cell pre-exposure to ligand for 2 hr

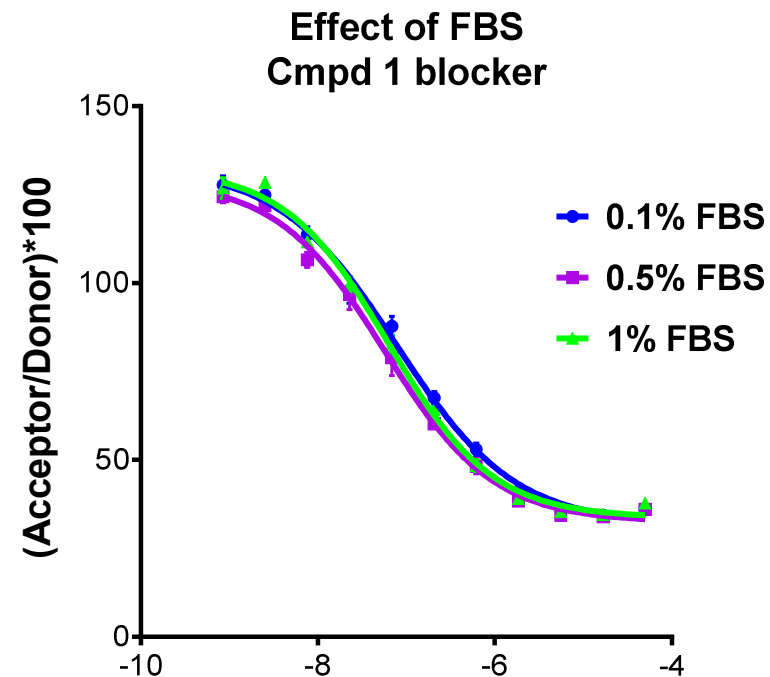
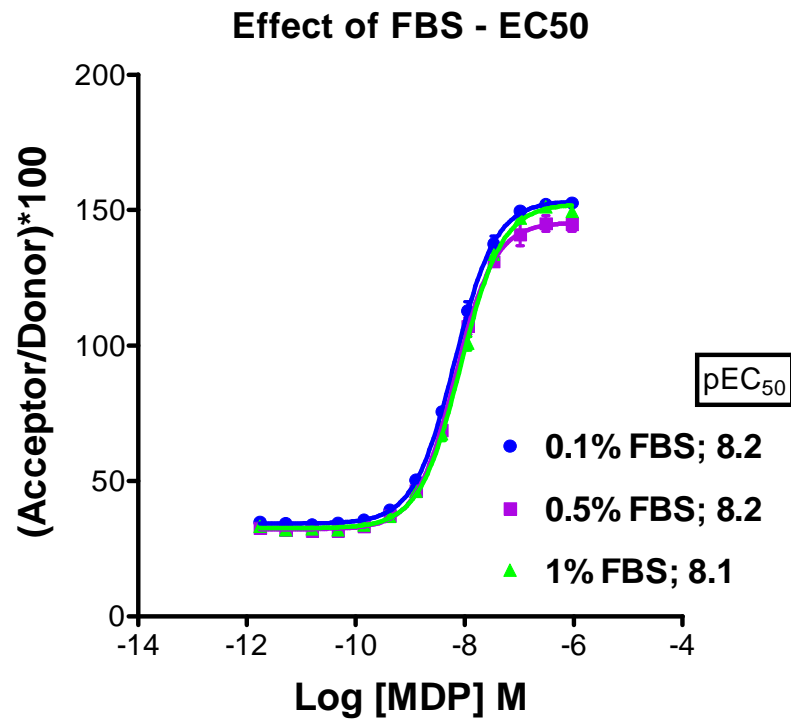
## Addressing plate reading lag time:



TIME (hr)	Z'
1	0.66
1	0.55
2	0.66
2	0.56
3	0.64
3	0.55
4	0.63
4	0.55
5	0.64
5	0.55

- The HTRF signal was stable 1-5 hrs after reagent addition

## Low serum had no effect on the NOD2 IL-8 assay:

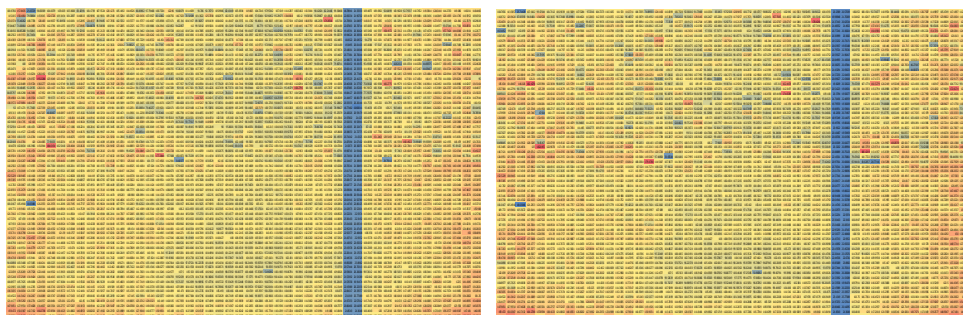


FBS:	0.1%	0.5%	1%
S/B	3.5	3.4	3.5
Z'	0.63	0.62	0.67

- Lowering the serum: no effect on MDP potency or compound blocker pIC<sub>50</sub>

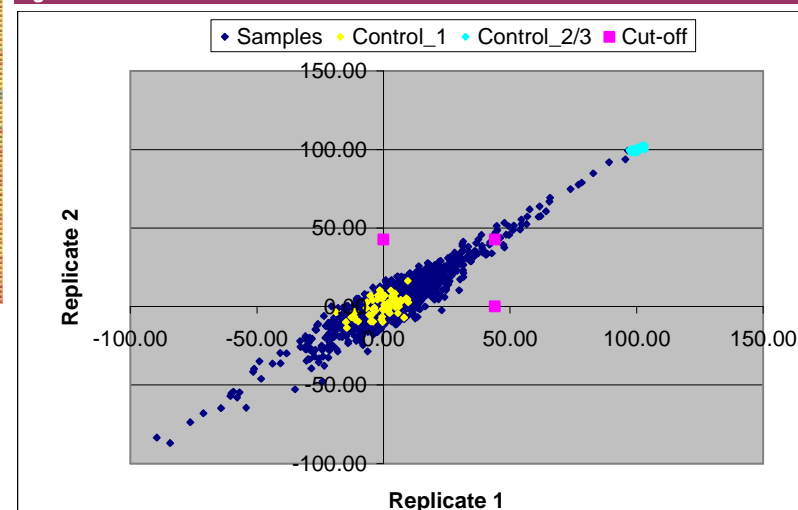
## NOD2 assay statistics: small compound set testing: “robustness”:

2 replicates



	rep 1	DMSO 1	rep 2	DMSO 2
<b>S/B</b>	4.4	4.2	4.5	4.2
<b>Z'</b>	0.77	0.72	0.79	0.72

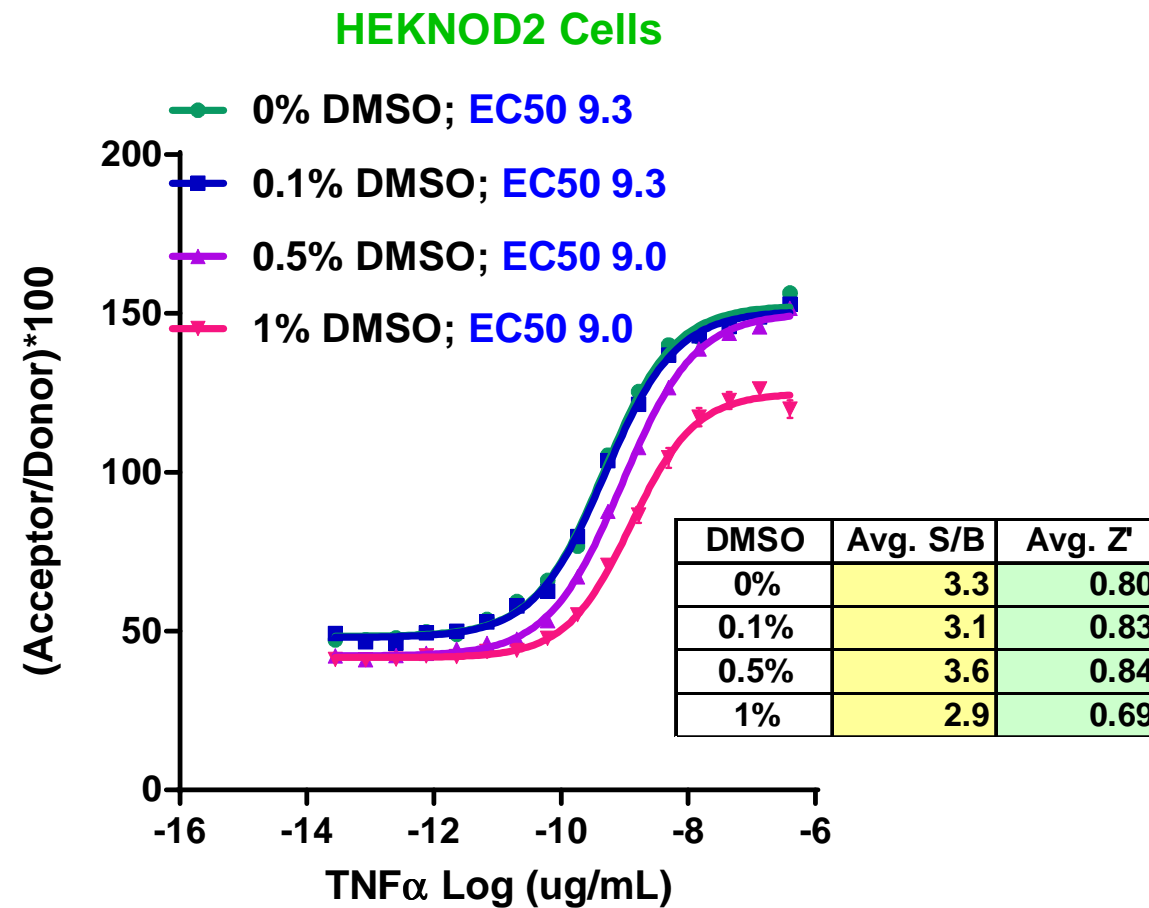
Figure 4. Correlation Plot



Replicate	Number of samples	% Inh Cut-off	N hits	HR(%)	Correlation Coeff.
<b>1</b>	1408	44	34	2.4	0.93
<b>2</b>		43	34	2.4	

- Random compound set of 1408 compounds used to assess statistics
  - Excellent for a cell based assay (Z', hit rate, correlation)

## NOD2 Specificity Assay; TNF $\alpha$ Stimulation of IL-8:



- TNF $\alpha$ -driven assay = similar responses/stats as the MDP NOD ligand



## NOD1,2 IL-8 HTRF Homogenous Protocol (low vol. 384 or 1536):

Plate 5  $\mu$ L NOD1 or NOD2 cells + ligand onto pre-dispensed compounds



Incubate O/N @ 37°C

Add 3  $\mu$ L HTRF reagents



Incubate @ RT for 2 hours

Read on Viewlux

- Simple protocol, amenable for HTS, screening 2M compounds

## Summary of NOD1, NOD2 work:

- HTRF assays for IL-8 performed better than the AlphaLISA assay (Z')
- HTRF assays performed well in 1536 well format: excellent cellular assay Z'
- Assay-ready frozen cells for HTSs = no need for continuous culture

**HTS 1.9 M**



**4000 XC<sub>50</sub>s**



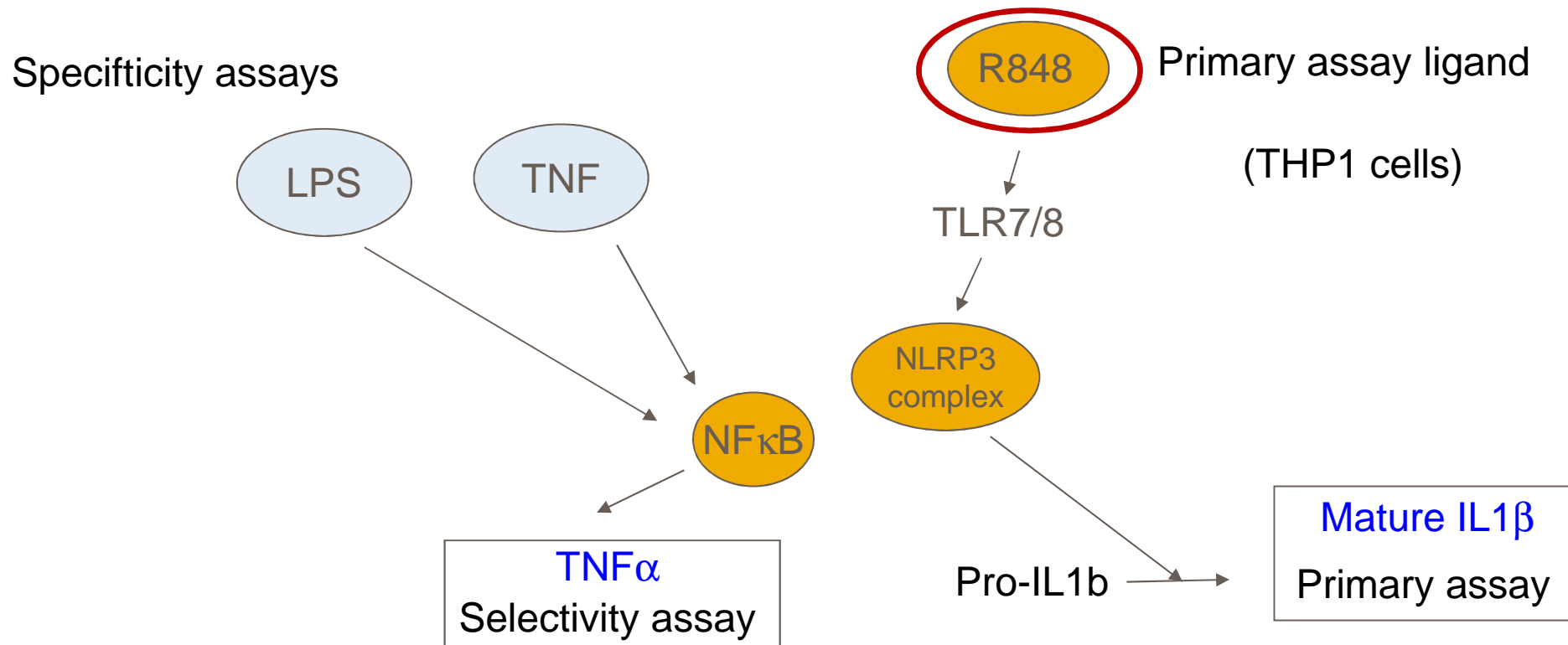
**Other specificity assays**



**= 1 NOD2 specific cmpd**

GSK Paper submitted

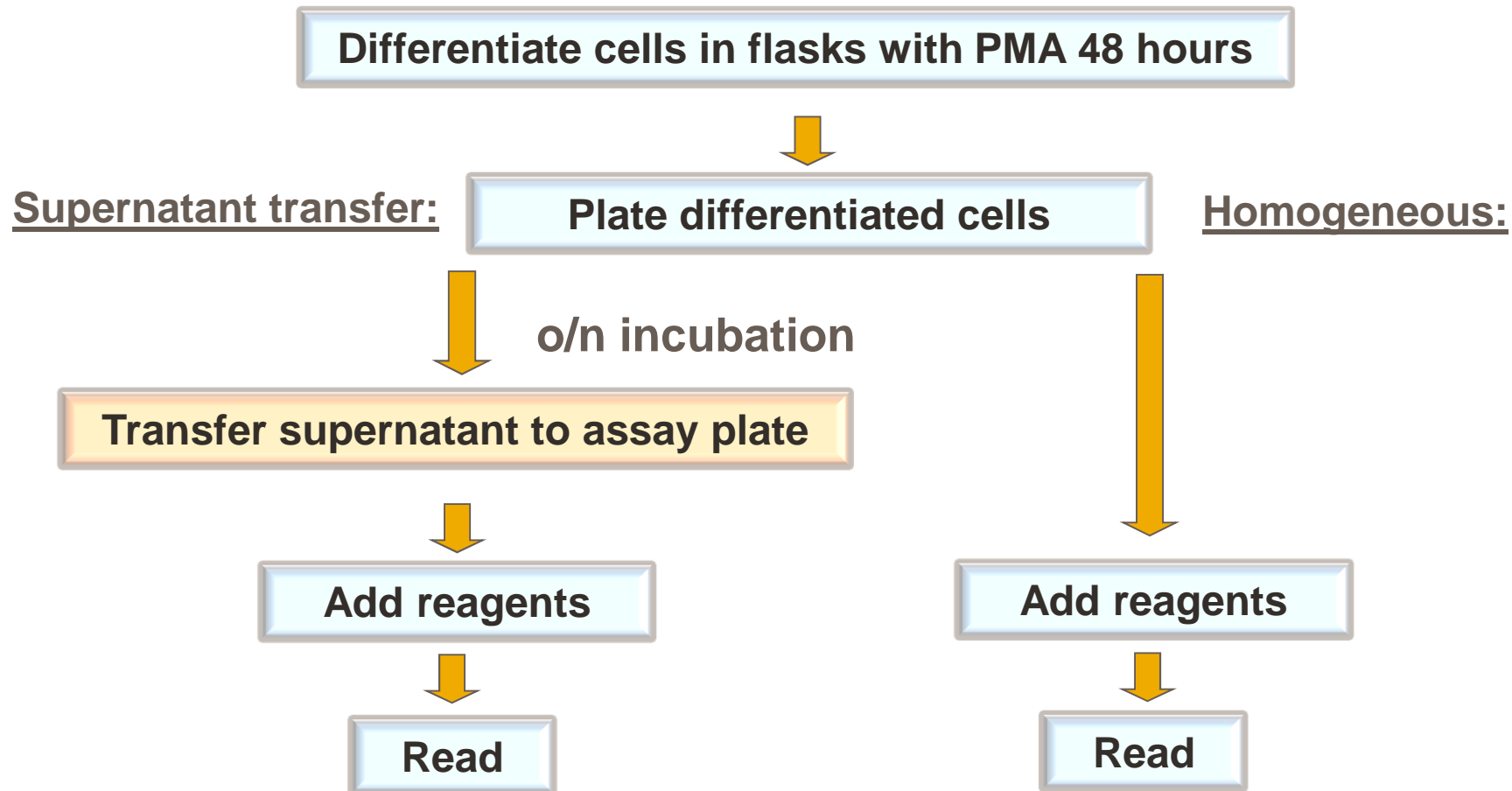
## NLRP3 primary and specificity cellular assays:



## IL-1 $\beta$ and TNF $\alpha$ assays:

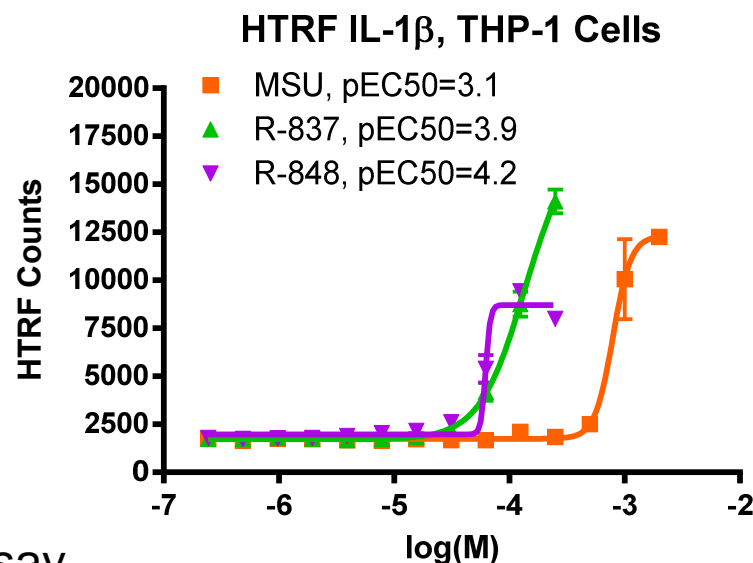
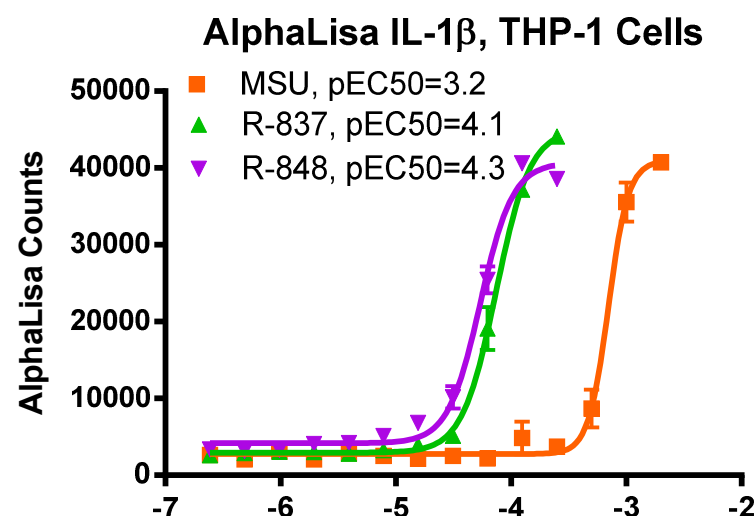
- HTRF (Cisbio) and AlphaLISA (Perkin Elmer) kits were available for TNF $\alpha$
- HTRF for IL-1 $\beta$ , yes; **AlphaLISA: no. Only pre-market test reagents (PE)**

## Flow chart for NLRP3 assays: supernatant vs homogeneous assays:



- Time on Envision instrument to read a plate: ~7 – 35 min
- Time on Viewlux instrument to read a plate: ~2 – 4 min

# IL-1 $\beta$ : supernatant assay: both AlphaLISA and HTRF OK:



- assay in the supernatant transfer assay

## Single concentration stats:

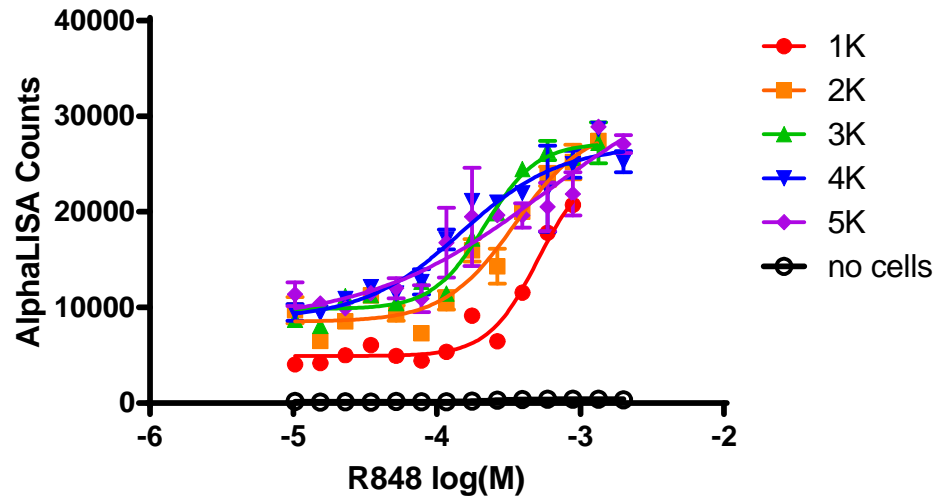
AlphaLISA (n=2)				
Cells	Ligand	pEC <sub>50</sub>	Z'	S/B
THP-1	MSU	3.4	0.7	14
	R837	3.7	0.6	16
	R848	4.2	0.7	14

HTRF (n=2)				
Cells	Ligand	pEC <sub>50</sub>	Z'	S/B
THP-1	MSU	3.2	0.8	10
	R837	3.5	0.7	10
	R848	4.0	0.7	7

- Both kits OK for IL-1 $\beta$  in a supernatant transfer assay

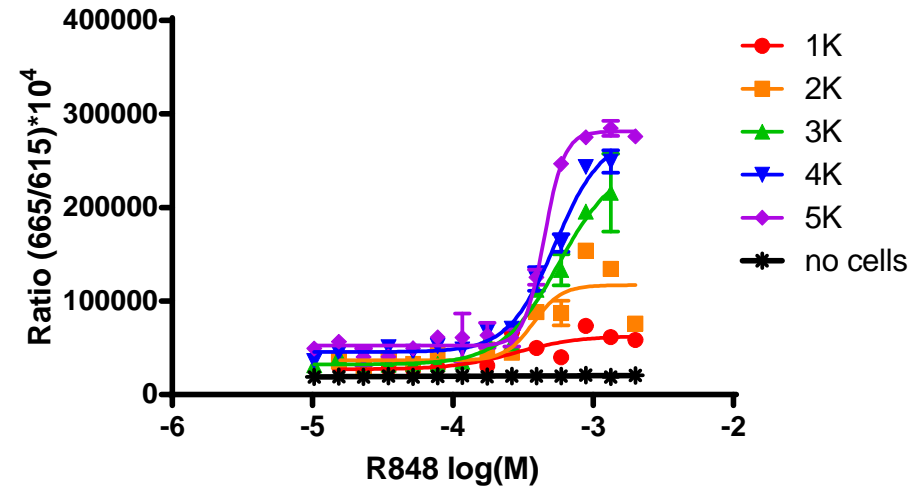
# IL-1 $\beta$ homogeneous: HTRF performed better:

IL-1 $\beta$  AlphaLISA



A-LISA	
Cells/well	S/B
1K	5
2K	5
3K	4
4K	3
5K	3

IL-1 $\beta$  HTRF

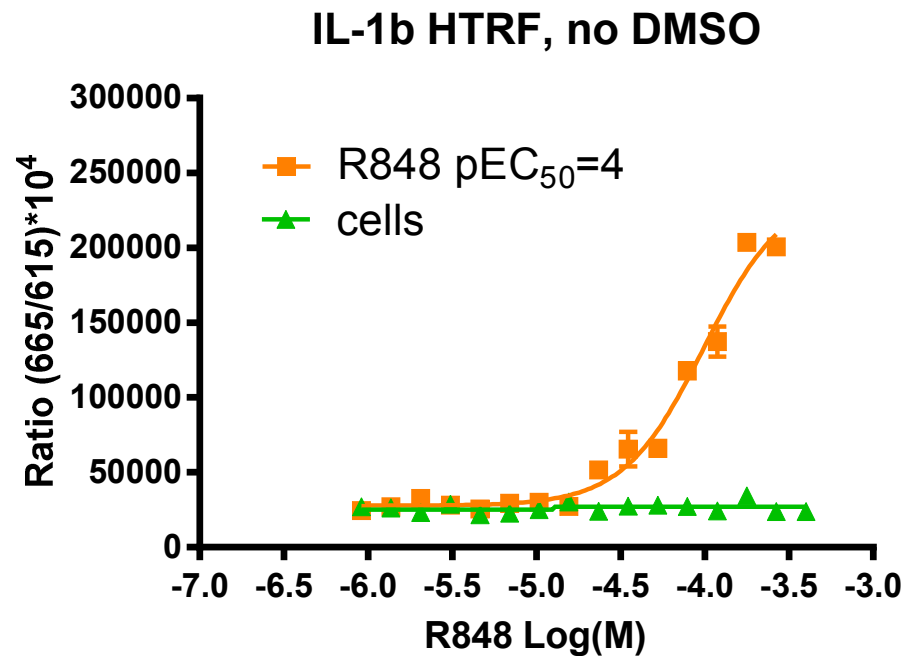


HTRF	
Cells/well	S/B
1K	2
2K	3
3K	8
4K	6
5K	5

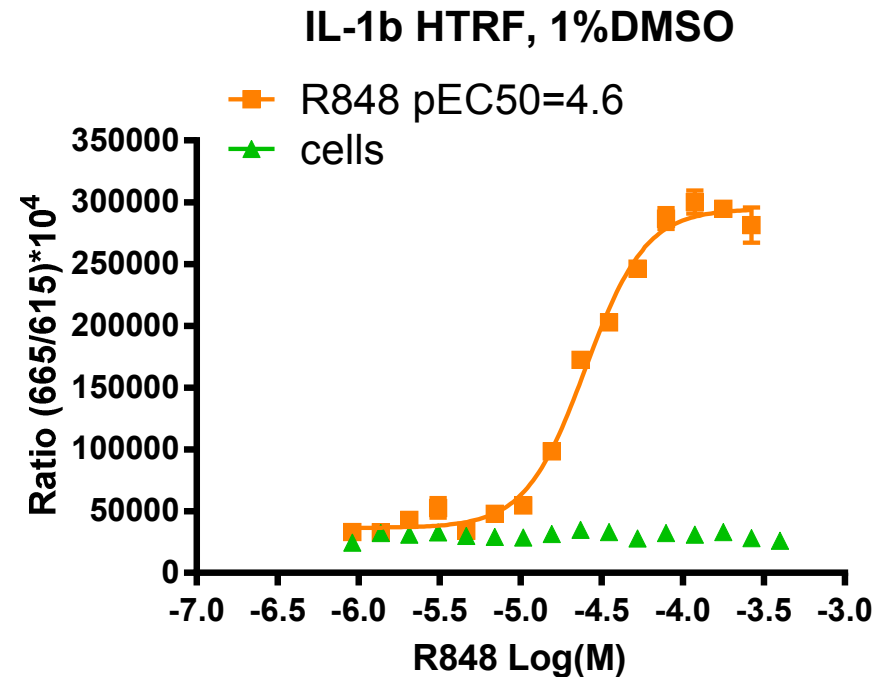
- IL-1 $\beta$  AlphaLISA did not perform as well as HTRF in homogeneous format

# IL-1 $\beta$ : DMSO helped the R848 ligand:

No additional DMSO



1% DMSO



- DMSO increased the R848 potency by approximately 0.5 log

## Homogeneous HTRF assay protocol for Profiling and Screening:

Differentiate cells in cell stacks in PMA, 48 hr



Plate cells + ligand (5 uL) onto 50 nL cmpds



16-18hr @ 37°C

Add 3 uL mixed HTRF reagents



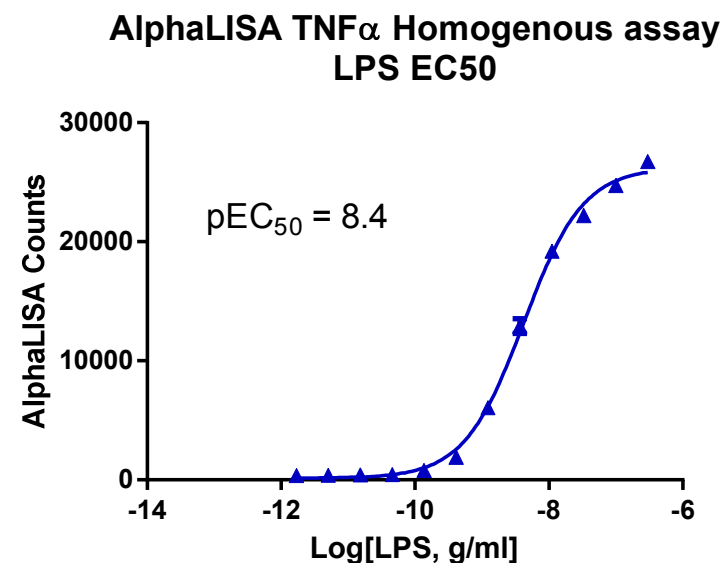
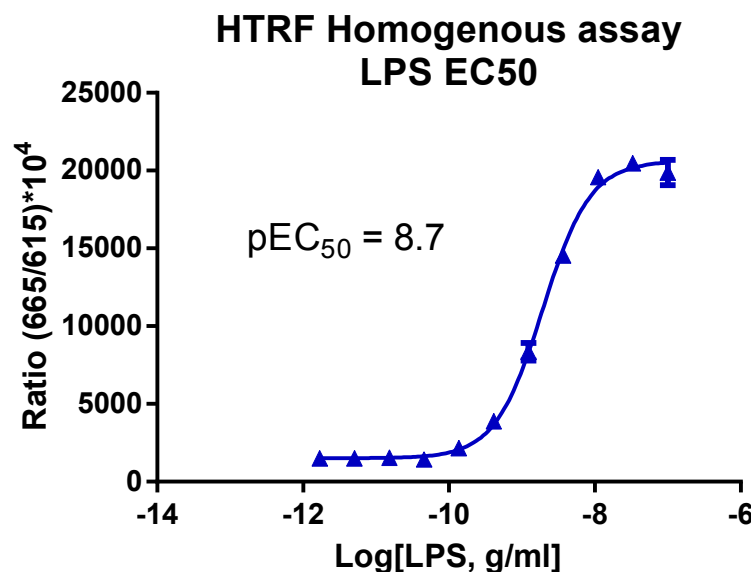
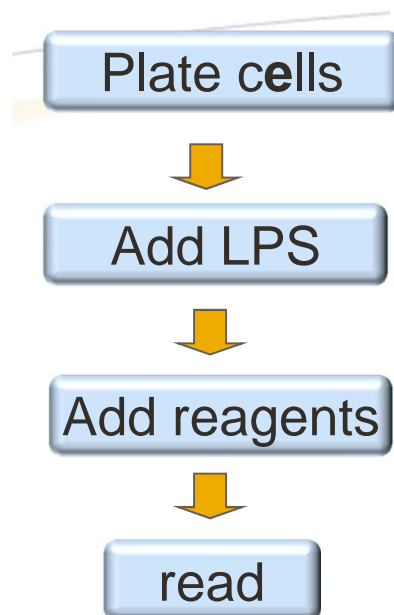
2hr @ RT

Read on Envision or Viewlux

- Cellular model: PMA differentiated THP1 cells
- With much effort, frozen cells did not respond well: fresh cells for HTS



## NLRP3 Specificity assay; TNF $\alpha$ AlphaLISA vs. HTRF



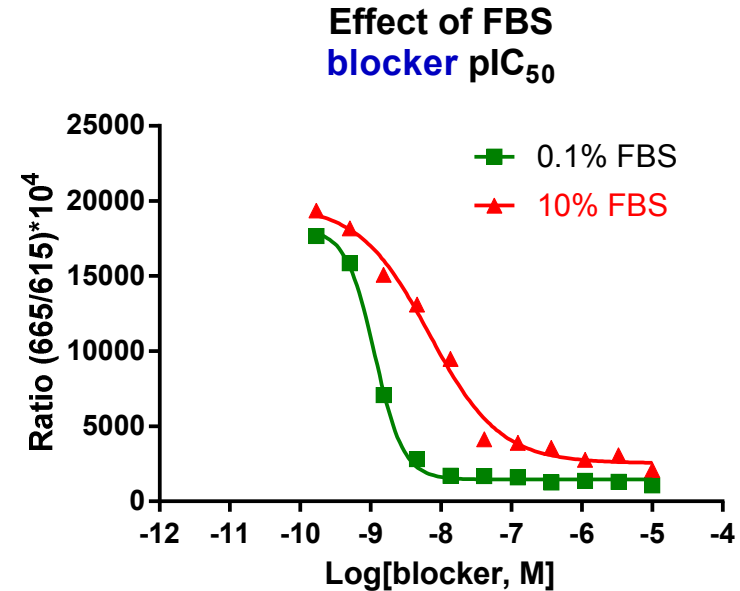
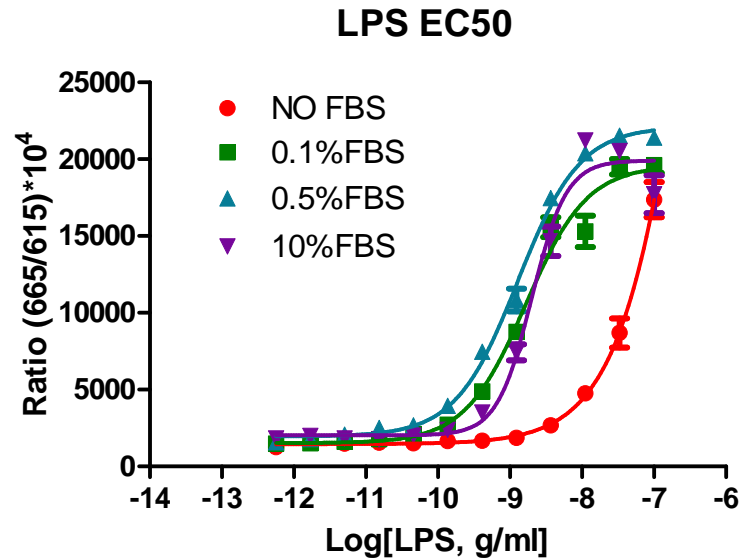
	Z'	S/B
AlphaLISA	0.88	56
HTRF	0.69	9

- For TNF $\alpha$ , the AlphaLISA performed better than HTRF
  - Higher Z', S/B
- HTRF statistics were acceptable
  - Used HTRF to match the primary screening assay

# FBS effect on blocker potency:TNF example:

Ligand:

Inhibitor:



- LPS ligand response: required at least 0.1% FBS
- Compound blocker lost potency in 10% FBS
  - common observation: reason to use reduced FBS in cellular assays

## Summary:

- HTRF assays for cytokines performed well in homogeneous format
  - 384 or 1536 well formats, excellent statistics
  - Have run numerous full HTS campaigns using HTRF for cellular assays
- Comparisons to AlphaLISA; in some cases performed better
  - In the assays shown, HTRF was better than AlphaLISA for IL-1 $\beta$
  - AlphaLISA better than HTRF for TNF $\alpha$
- Ease of use was better for HTRF than for AlphaLISA in our hands
  - HTRF reagents less light sensitive than AlphaLISA
  - Envision read times were longer than Viewlux
    - AlphaLISA cannot be read on a Viewlux
- Not all cells could be adapted to an “assay-ready” frozen cell format

## Currently:

- Using HTRF for the detection of an intracellular nuclear protein
- Using dual acceptor approach (one donor fluor, 2 acceptor fluors)

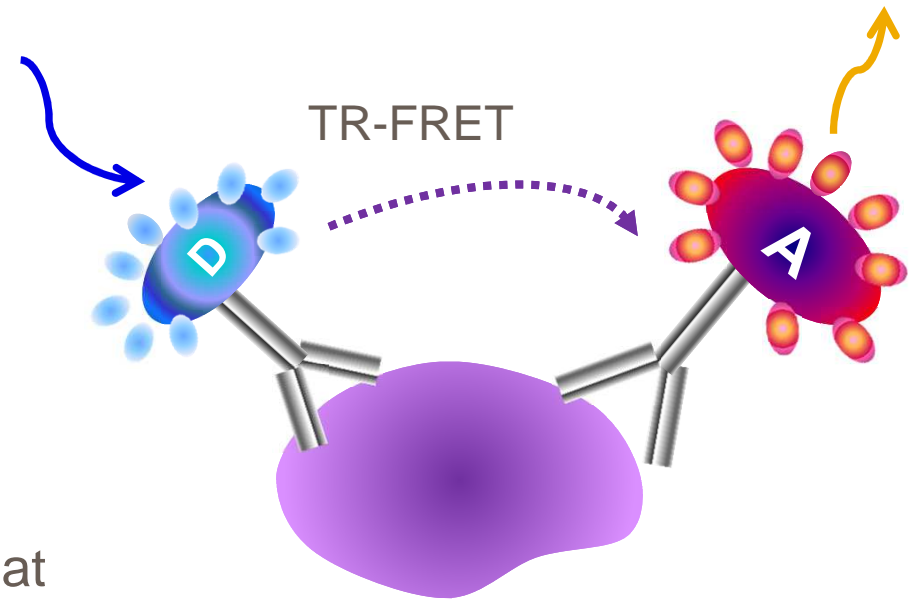
About **HTRF** .....

**H**as **T**he **R**ight homogeneous **F**ormat

**H**elps you **T**o **R**un assays **F**ast

Cisbio **H**elps **T**est new **R**eagents to **F**ind new assays with custom labeling

Scientists are **H**appy **T**o **R**un these **F**RET assays..... (OK maybe not....)



# Acknowledgements

## Assay dev and screening:

- Tushar Chordia
- Anna Waszkiewicz
- Maggie Truong
- Jonathan Basilla
- Xin Zeng
- Jim Fornwald
- Monica Montoute
- Zining Wu
- Derek Poore
- Hu Li
- Guofeng Zhang
- Lisa Payne
- Tom Sweitzer
- Subhas Chakravorty
- Elizabeth Davenport
- Bob Ames

## Biology and Chemistry

- David Rickard
- John Emery
- Viera Kasparcova
- Chuck Hanning
- Pamela Haile
- Jong Yu
- Mike Cook
- Clark Sehon
- Patrick Eidam
- Robert Marquis
- Joshi Ramanjulu
- Carol Anne Ogden
- Peter Gough
- John Bertin