

Meso Scale Discovery:

**The New Kid on the Block for
HTS Protein Kinase Assays**

A Beta-Site Evaluation

David Powers
Research Scientist
Amgen Inc. Thousand Oaks, CA

Introduction

MSD Contacted Amgen Sept '02

Signed Early Access Program Feb '02

Objectives

- Does it work?

Technology, Hardware, Biology

- Will it have utility at Amgen?

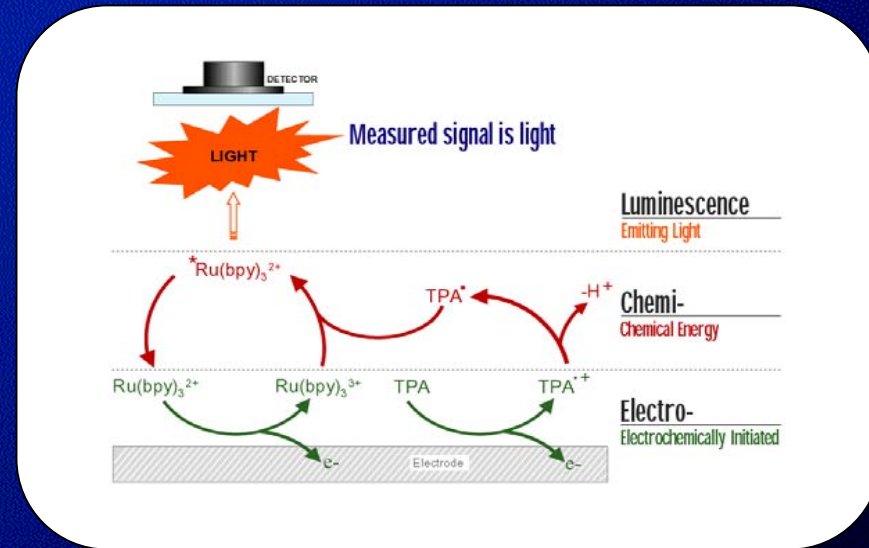
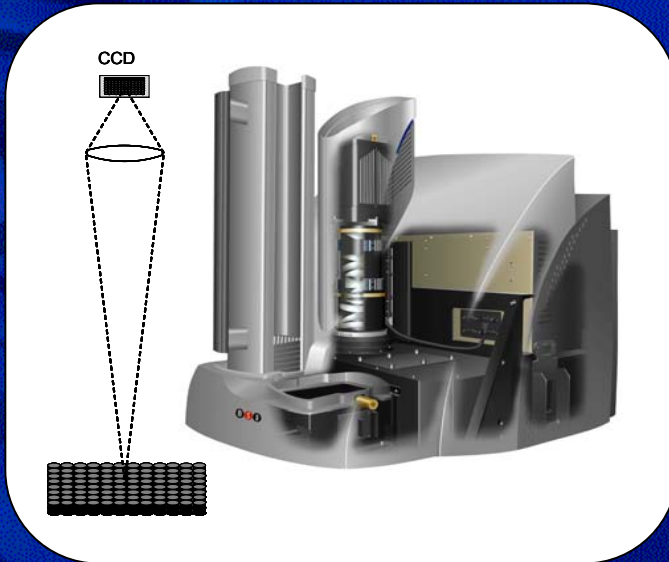
Integration it into our Existing Process?

Future?

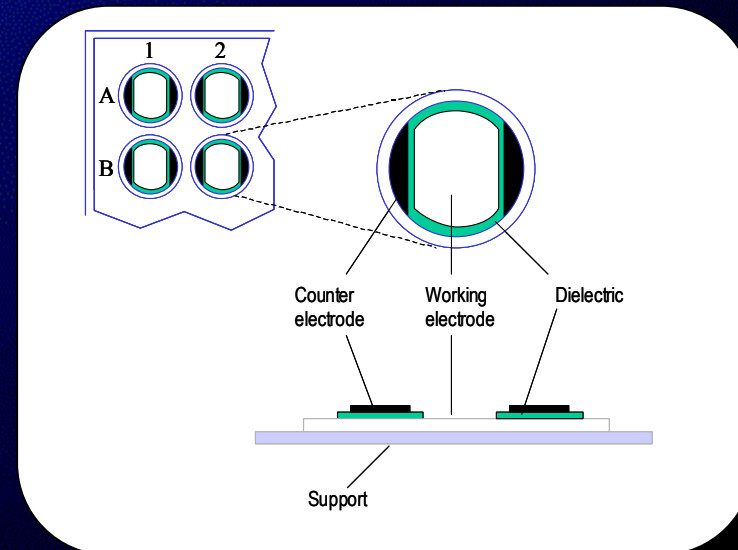
Structure

- Installation and Training
- Conduct 75,000 Multi-Array determination (Ser/Thr Kinase Screens)
- Conduct 35,000 Multi-Spot determinations (Cytokine Secretagogue)
- Comparison of Multi-Array Technology platform against TR-FRET
- Provide Feedback

Meso Scale Sector HTS



- *Plate Based*
- *Electrochemiluminescent*
- *Homogeneous*
- *Imaging Supports High Density Plates*



Hardware

Features

- CCD Imaging (1 min/plate)
- Small Footprint
- 75 plate capacity Stacker
- Stylish Design

- More than 3500 plates cycled
- 1500 plates cycled in three days
1 instrument failure (loose wire)
- No Stacker Failures
- No Communication Failures

Wish List

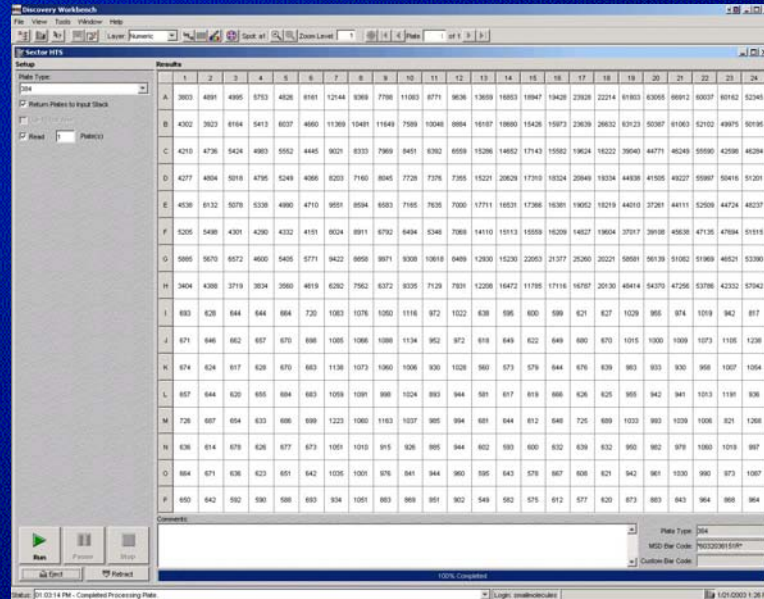
- Barcode Position Options



Software

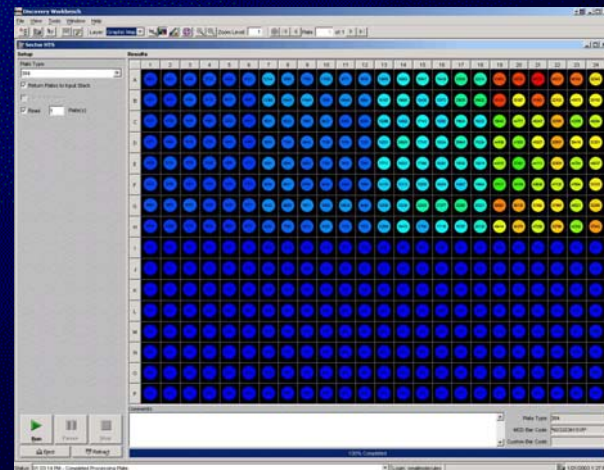
Features

- Stable Software
- Simple/Intuitive GUI
- No Frills
- Plate In/ Data Out

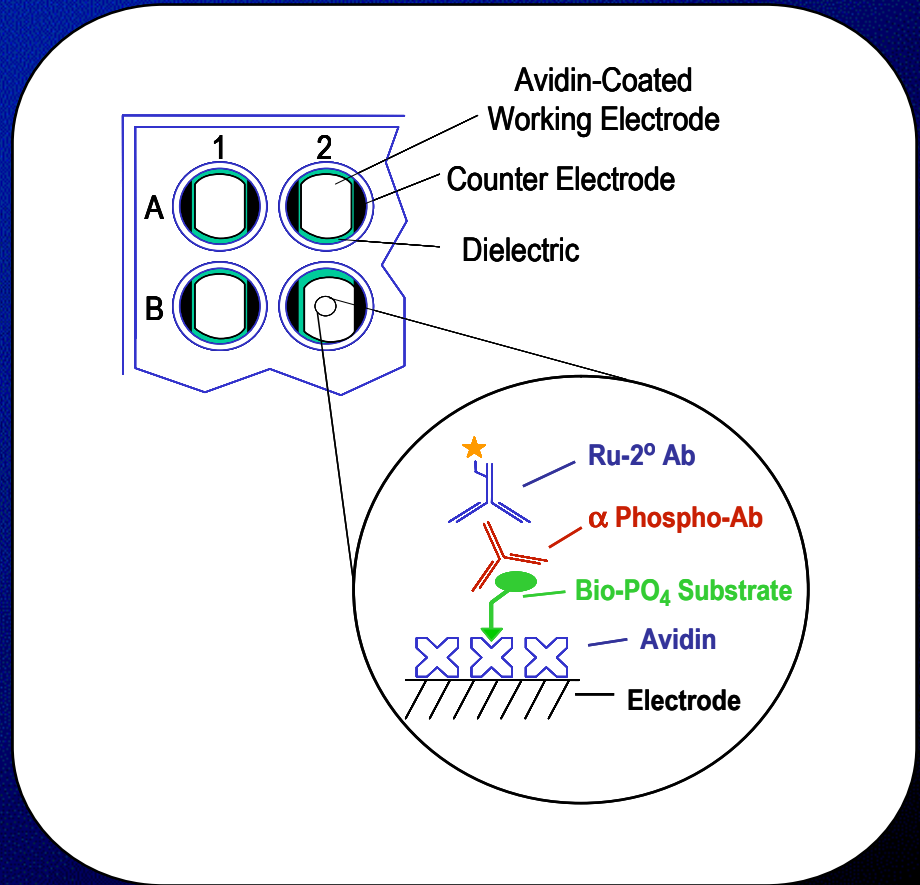
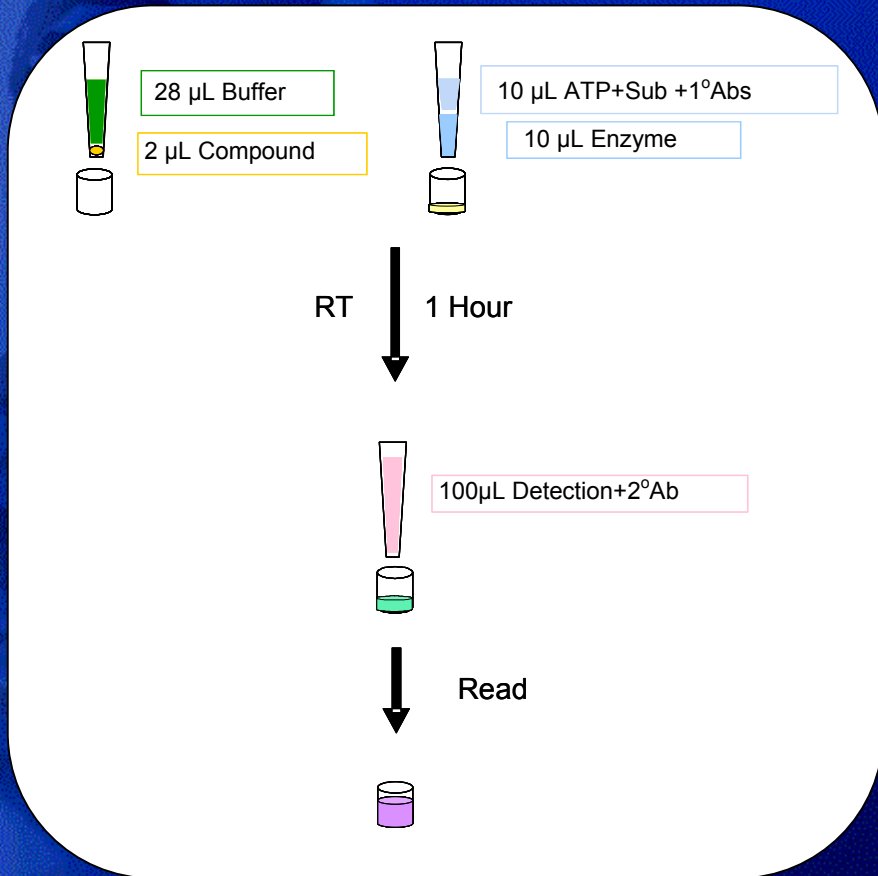


Wish List

- Assay Based Protocol
- Improved File Management



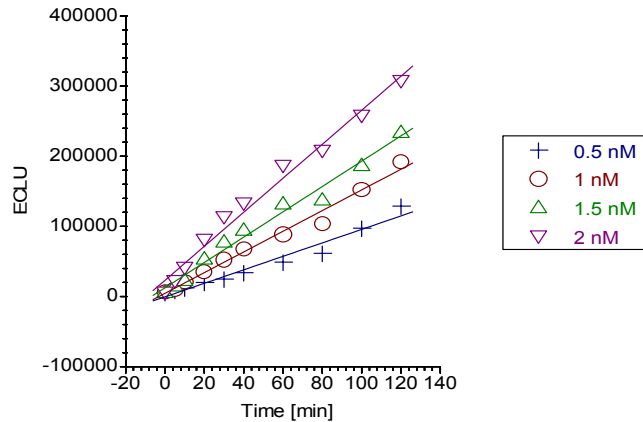
Screening Protocol



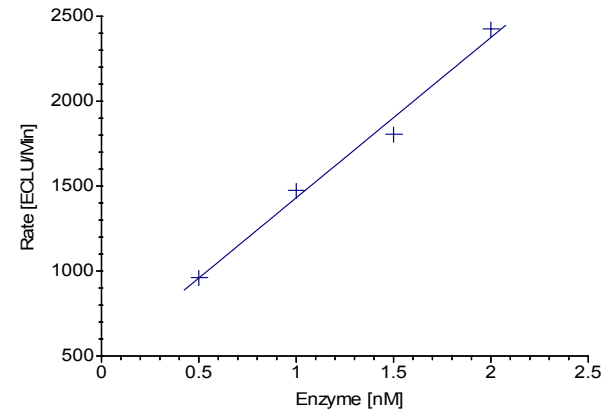
*Ser/Thr Kinase
Biotinylated Substrate
Labeled Generic 2 $^\circ$ Antibody*

Kinase Assay Development

MSD Kinase Titration/Time Course



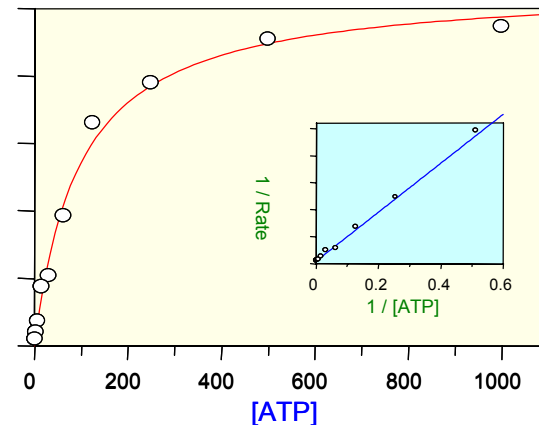
MSD Kinase Dilution



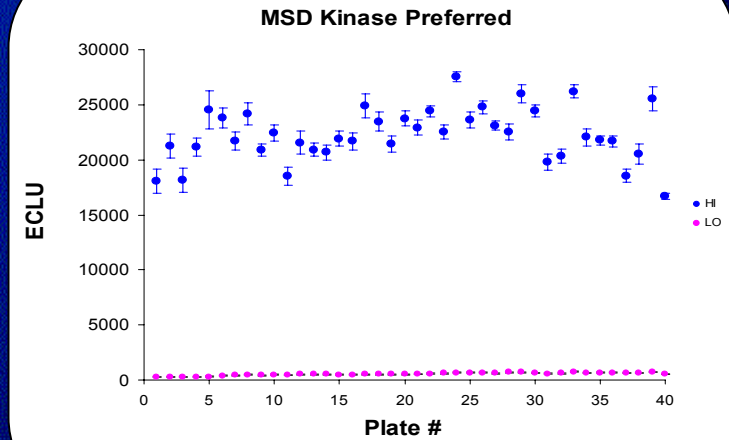
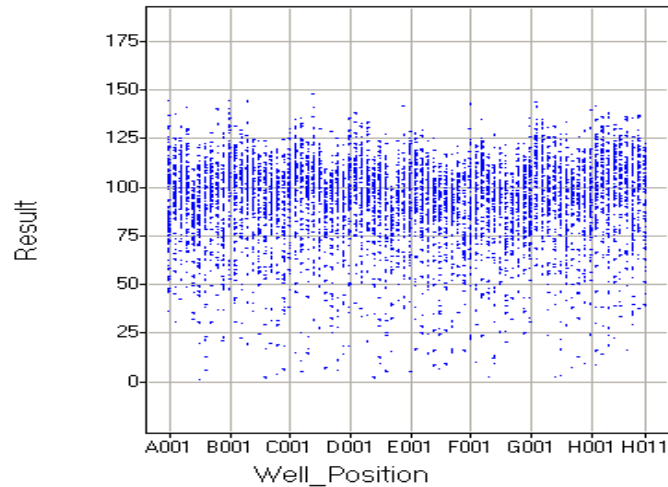
ATP $K_{m\ app}$ [μ M]		
Meso Scale	Lance	FBA
96.56	99.71	100.15

- Ser/Thr Kinase
- Transferable Assays
- Comparable To Other Platforms
- Less Enzyme (5-10 fold)

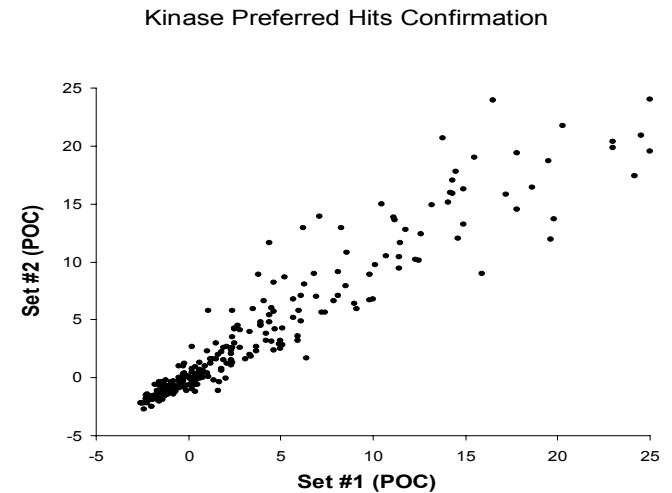
Kinase 1 MSD - K_m for ATP



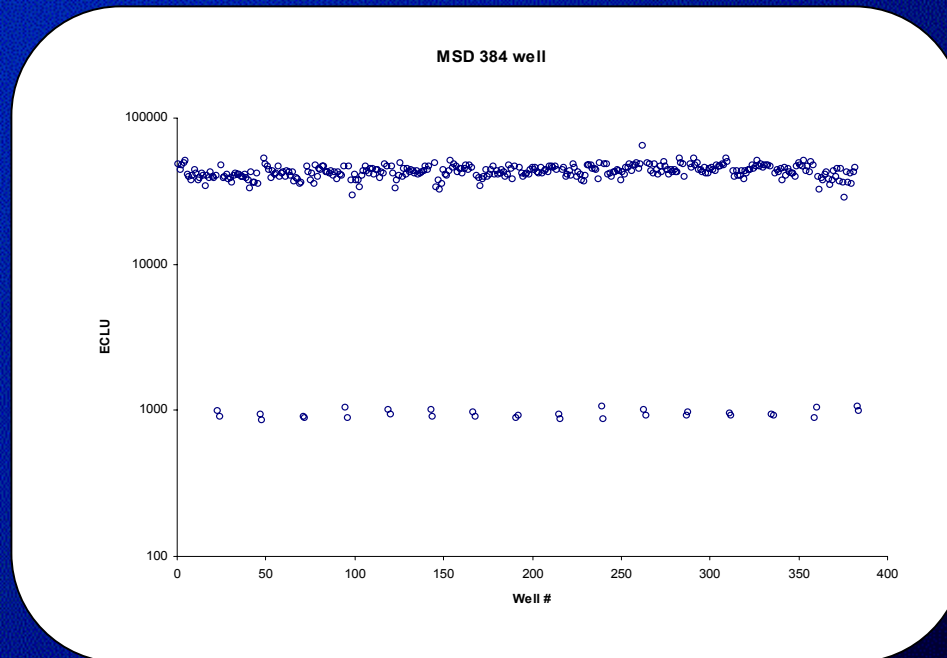
Kinase Preferred



- Preferred Collection 40 μ M
- Kinase = 2 nM
- S/B = 40
- Z' = 0.68



384 Well



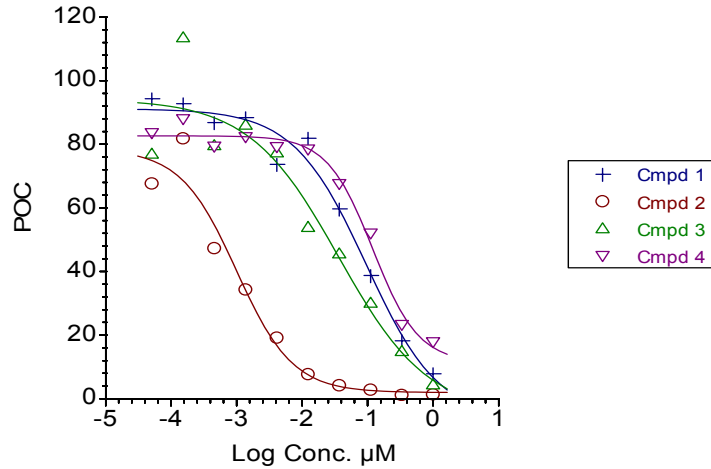
- Kinase = 5 nM
- 50 μ L Vol.
- S/B = 45
- Z' = 0.67

Read Time

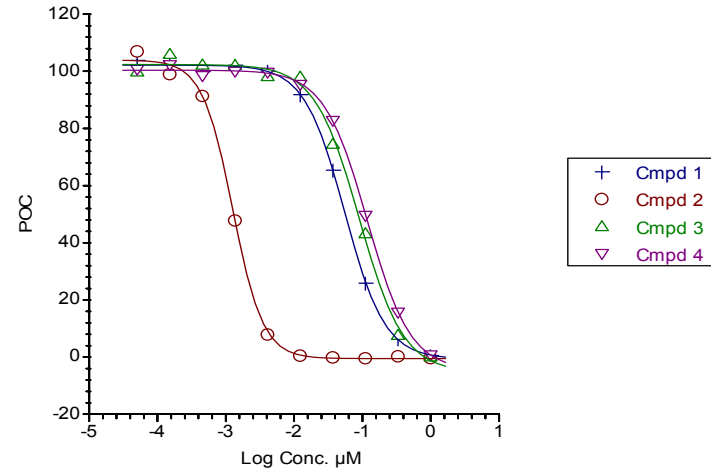
Sector HTS	Discovery
~ 1 min	>8 min

LANCE vs. MSD

MSD Dose Response



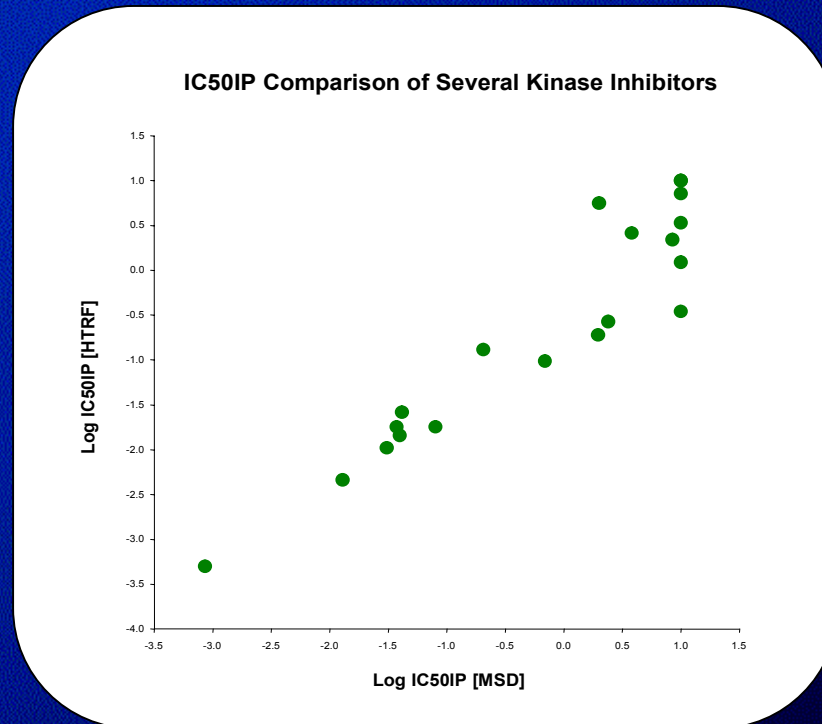
LANCE Dose Response



- Same as TR-FRET
- Intrinsic Potency
- Selectivity
- 5x less protein
- Lower 'bottom end'

	MSD [IC50 μM]	LANCE [IC50 μM]
Cmpd 1	0.0954	0.0548
Cmpd 2	0.001	0.0013
Cmpd 3	0.0362	0.0892
Cmpd 4	0.12	0.1102
S/B	37	7

LANCE vs. MSD

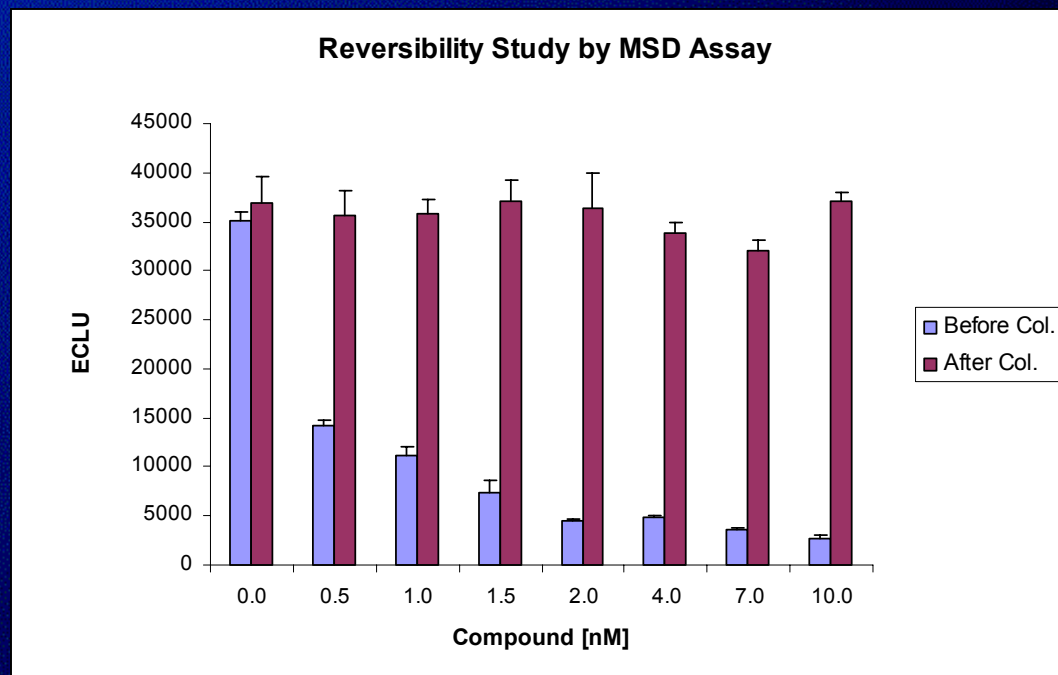


- *20 Inhibitors in side-by-side comparison*
- *IC50 similar across both platforms*
- *Some compounds appear less potent in MSD*

Mechanistic Studies

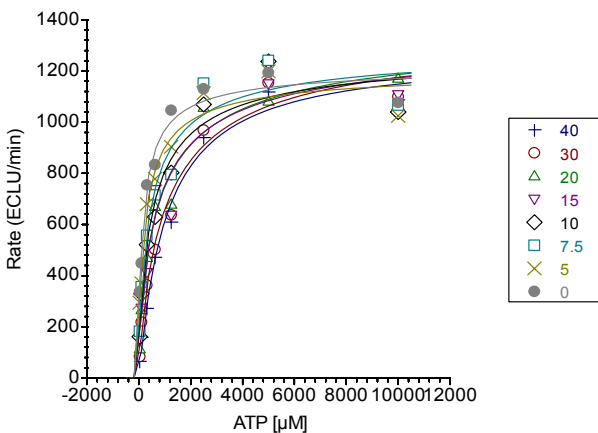


Are Inhibitors Reversible?

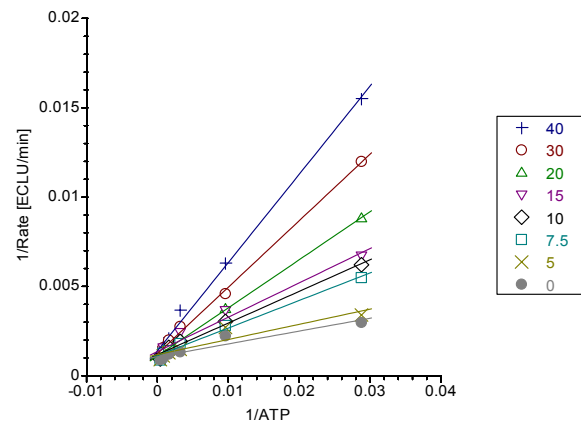


Mechanistic Studies

MSD Kinase Assay

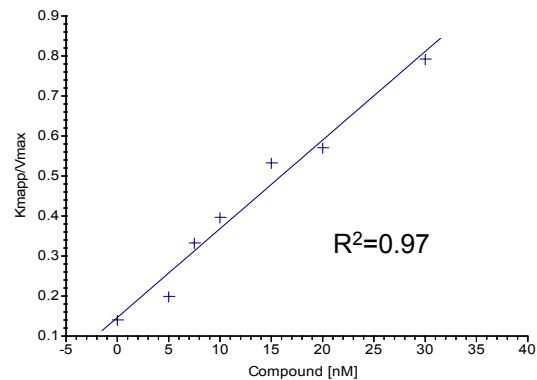


MSD Kinase



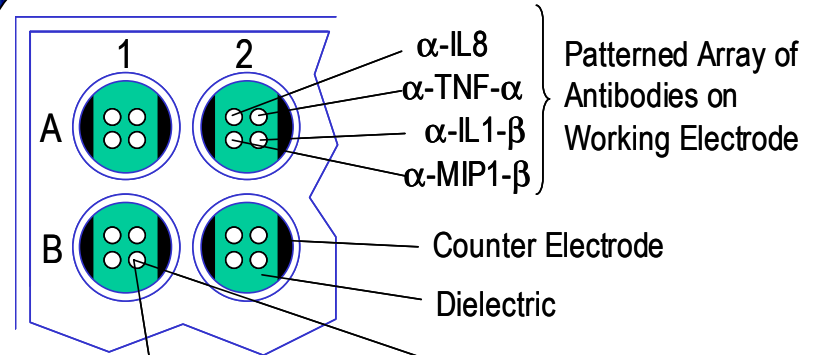
Are inhibitors ATP Competitive?

K_mapp/V_{max} vs. Compound [nM]



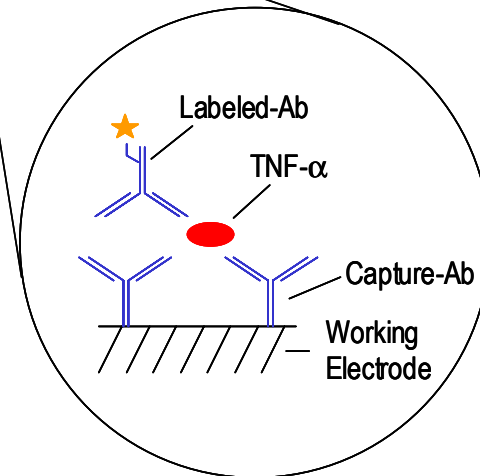
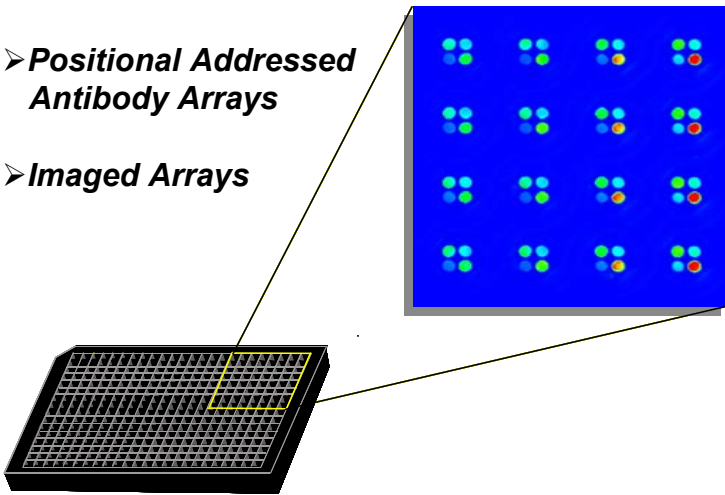
Multiplex Secretagogue Assay

- *LPS Stimulated THP-1 cells*
- *Dispense 50 μ L media*
- *Add 100 μ L Abs in MSD Buffer*
- *Incubate 2 hours*
- *Read*

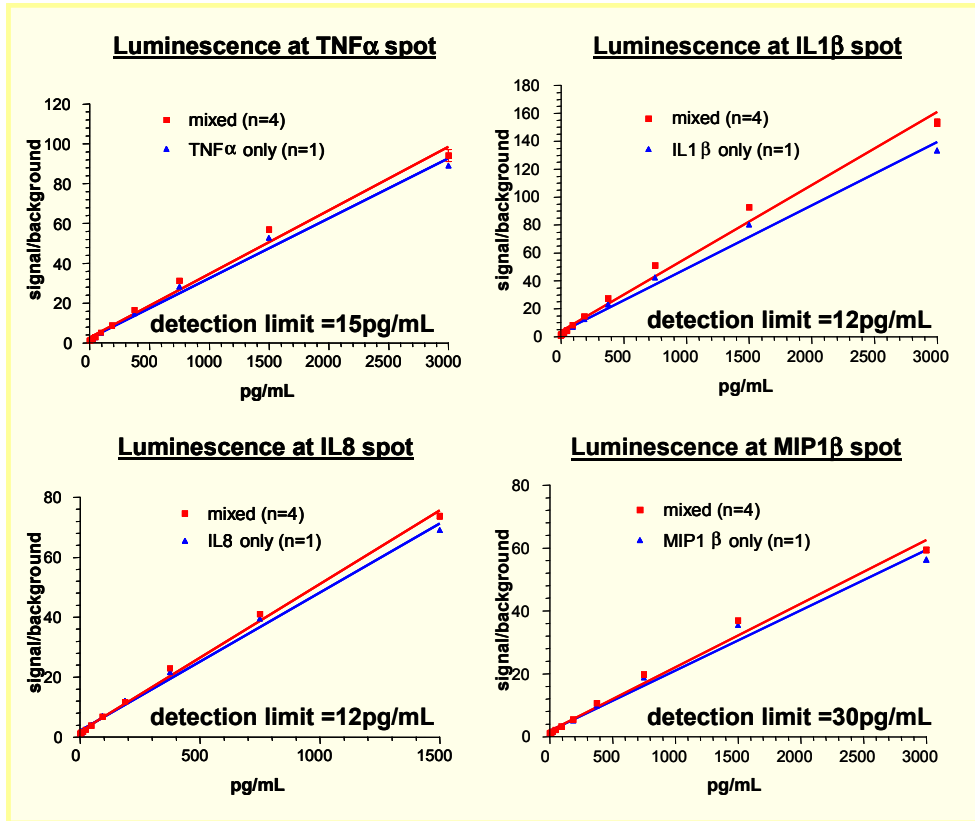


➤ **Positional Addressed Antibody Arrays**

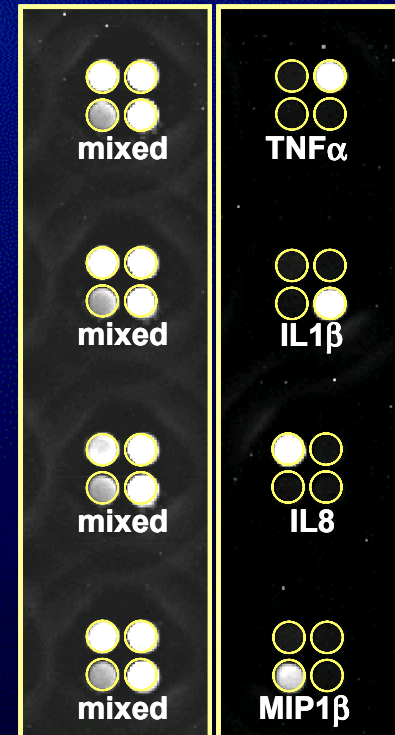
➤ **Imaged Arrays**



Standard Curves and Detection Limits



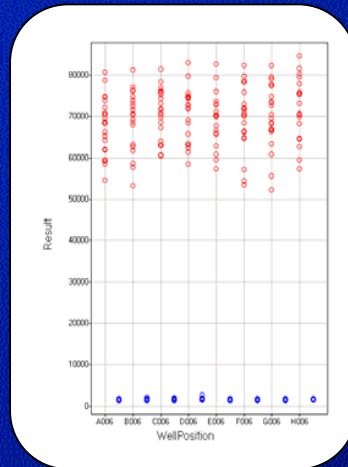
750 pg/mL cytokine:



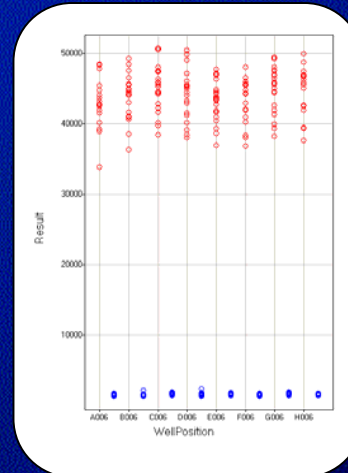
No Ab cross-reactivity or signal cross-talk

HTS Mutiplex

IL8
 $Z'=0.9$ (intra)
 $Z'=0.8$ (inter)

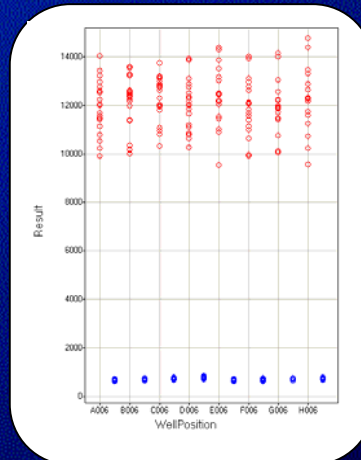


TNF α
 $Z'=0.8$ (intra)
 $Z'=0.7$ (inter)

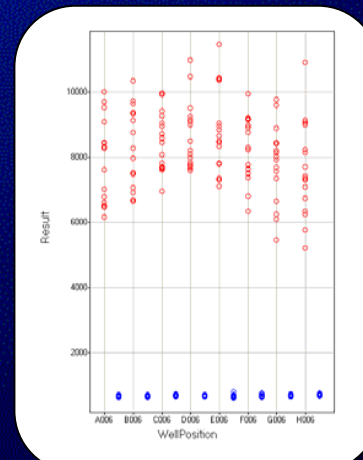


HI/LO Scatter plots
20 plates

MIP1 β
 $Z'=0.8$ (intra)
 $Z'=0.8$ (inter)



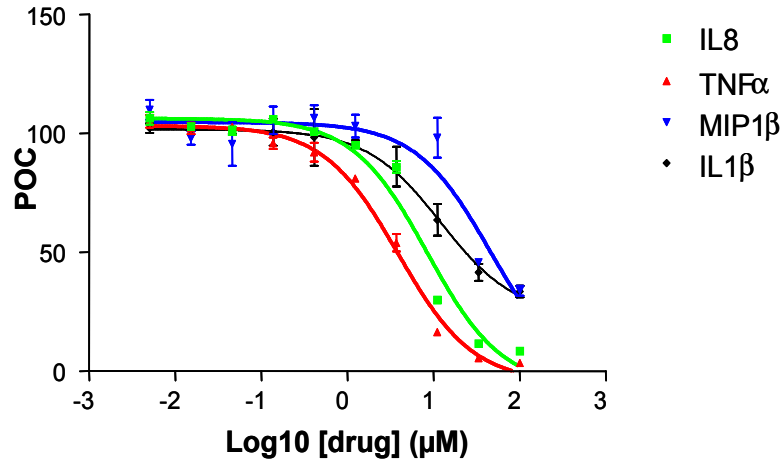
IL1 β
 $Z'=0.8$ (intra)
 $Z'=0.6$ (inter)



Total 4320 Compounds Screened @ 10 μ M

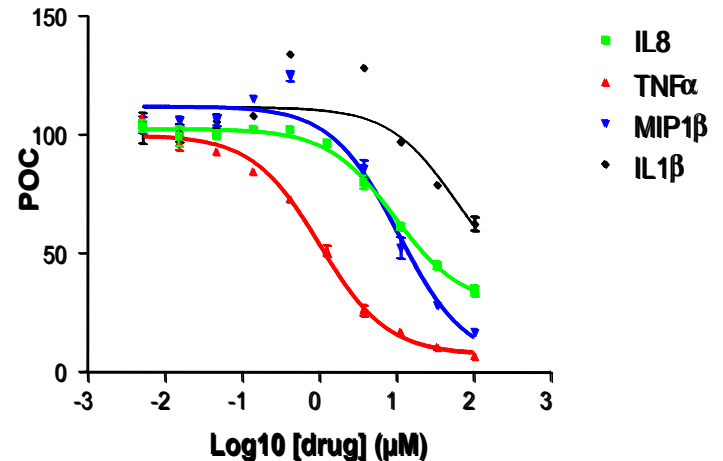
Multiplexed Dose Response

Dose-Response

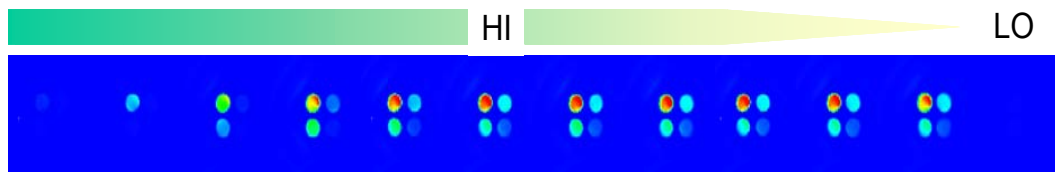


	IC50 IP	STDEV
IL8	6.76	0.30
TNFα	3.70	0.24
MIP1β	23.15	6.07
IL1β	10.60	1.36

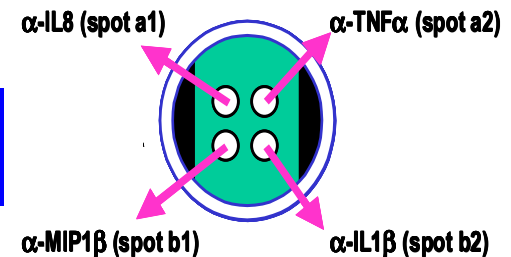
Dose-Response



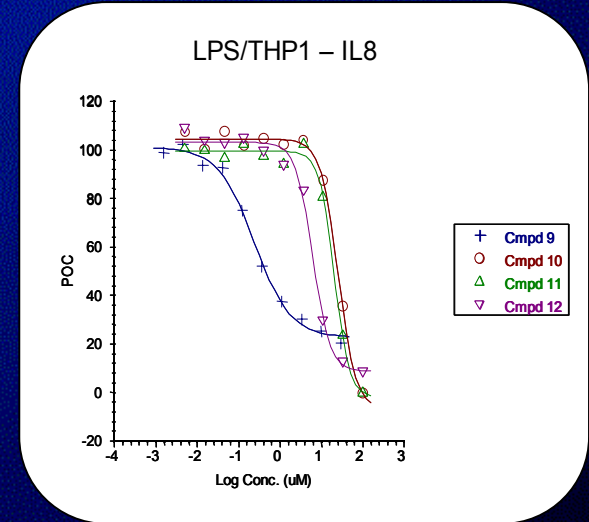
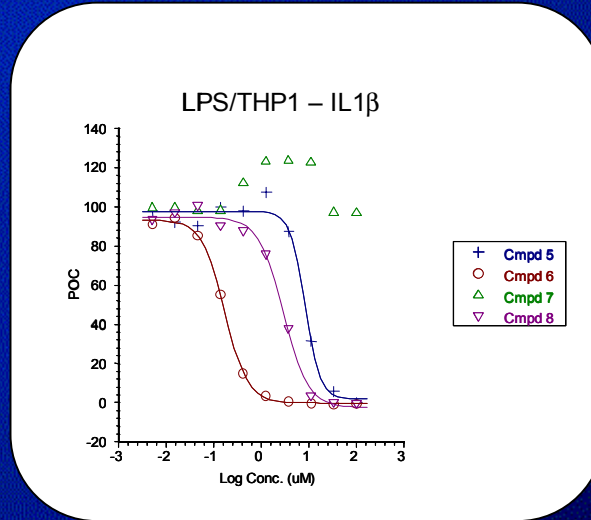
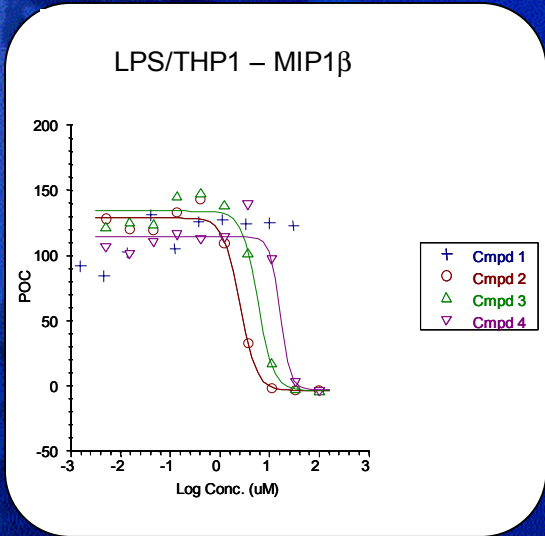
	IC50 IP	STDEV
IL8	8.32	1.64
TNFα	1.00	0.17
MIP1β	7.42	1.15
IL1β	>100	N/A



Drug concentration

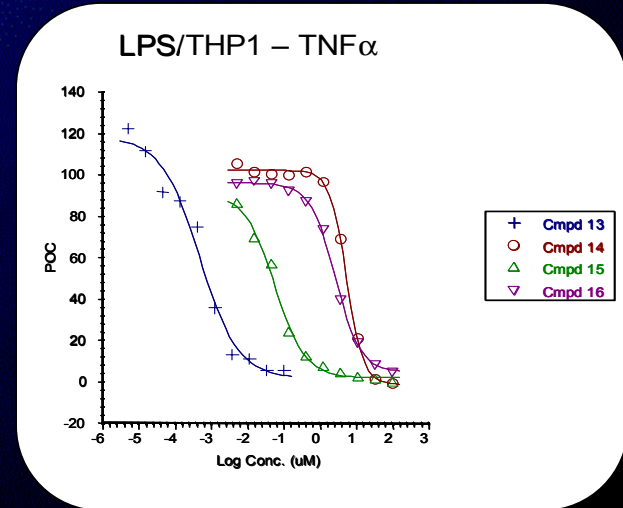
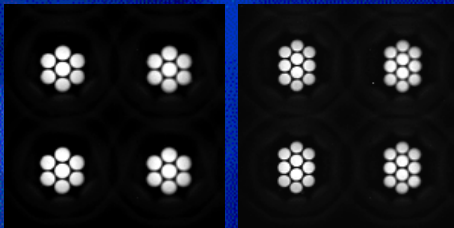


Quantitative Arrays



Enabling Capacity

Annotation of Libraries
Multiple Stimulators
Multiple Cell Lines
Multiple Species



.....towards fingerprinting

Lance vs. MSD

	Compound 1		Compound 2		Compound 3	
LANCE	TNF α	3.4 μ M	TNF α	0.0007 μ M	TNF α	0.433 μ M
MSD	TNF α	5.49 μ M	TNF α	0.0005 μ M	TNF α	0.84 μ M
	IL1 β	3.5 μ M	IL1 β	0.008 μ M	IL1 β	5.3 μ M
	IL8	6.8 μ M	IL8	0.007 μ M	IL8	3.9 μ M
	MIP1 β	>30 μ M	MIP1 β	>30 μ M	MIP1 β	>30 μ M

Comparable across platforms

Conjugation of Ab pairs limits Lance™ Application

MSD Technology Platform

Pros

- Homogeneous
- Robust Stable Signal
- Short read times 1 min/plate (high density)
- Multiplex-High Content
- 1° Antibody labeling not required

Cons

- Single Read
- Plate Based (logistics, storage)

Future Directions

- Kinetics- Solution vs. Solid Phase
- Carbon Surface Immobilization
- What biology won't work (Binning Targets)

Conclusion

MSD Technology Platform Fits into our Process

- Build assays quickly
- Identify 'hits'
- Assign Potency/Selectivity
- Determine Cellular Efficacy
- Establish *in-vivo* predictors

- Supports Higher Density
- Quantitative HTS Arrays

Identification, Stratification and Advancement of Compounds

Acknowledgments

Amgen

Yanyan Tudor

Violeta Yu

Meso Scale Discovery

Jim Wilbur

Kent Johnson

Charles Clinton

George Sigal