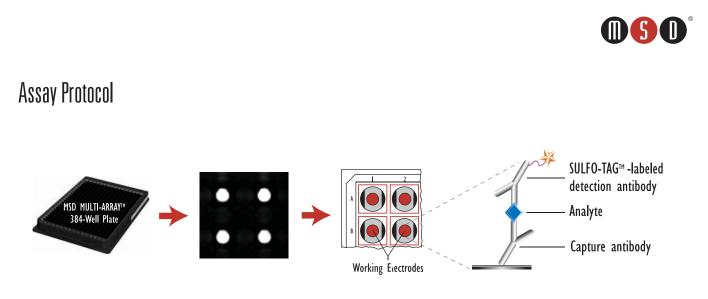


# High Throughput Assays for Biomarkers in 384-well Format

Robert M. Umek, Paula D. Eason, Sharon H. Tynan, Alan Kishbaugh, Pankaj Oberoi, and Jacob N. Wohlstadter

Monitoring biomarkers is common throughout the drug discovery process and increasingly popular in the earliest phases. Biomarker assays can be powerful readouts for cell based assays in support of focused library screening and the pursuit of structure-activity relationships. In order to maximally support these early drug discovery efforts, biomarker assays should be compatible with existing screening practices, ideally formatted in 384-well plates. We describe here a number of immunoassays for biomarkers formatted in 384-well plates. The assays are inclusive of a wide range of classes of biomarkers including intracellular phosphoproteins, the amyloid peptides of Alzheimer's Disease, and serum biomarkers. The assays are quantitative, sensitive, and retain the performance characteristics of the same assays in 96-well format. The assays are typically performed with 10 µL of sample making them well-suited to 384-well cell culture applications or the characterization of limiting amounts of fluids including cell supernatants, serum, plasma, and CSF.

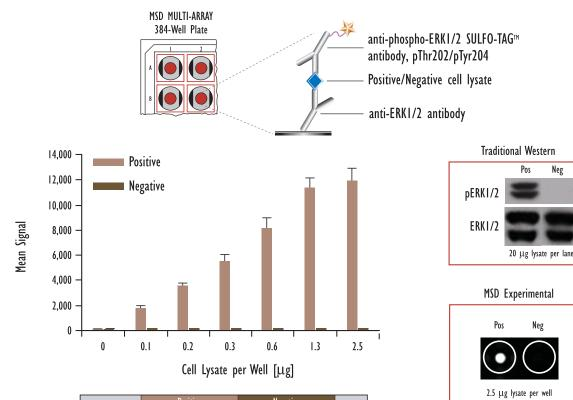


# General Protocol

- I. MSD MULTI-ARRAY 384-Well plates precoated with capture antibodies, plates are blocked and washed.
- 2. Samples are incubated in the assay plate with shaking, 10  $\mu$ L per well, and then washed.
- 3. Antibodies labeled with MSD SULFO-TAG reagent are incubated in the assay plate with shaking, 10  $\mu$ L per well, and then washed.
- 4. MSD Read Buffer T , 40  $\mu L$  per well, followed by plate analysis on an MSD SECTOR Imager instrument.



## Detection of Phosphorylated ERK1/2 (pThr202/pTyr204)



Lysates (µg)	Positive				P/N		
	Average	StdDev	%CV	Average	StdDev	%CV	171
0	116	13	- 11	108	13	12	
0.1	1,777	269	15	131	10	8	13.6
0.2	3,526	277	8	126	19	15	28.0
0.3	5,438	567	10	122		9	44.8
0.6	8,165	859	11	131	13	10	62.5
1.3	11,334	775	7	140	21	15	81.2
2.5	11,873	1,028	9	147	27	18	80.7

Jurkat cells were treated with PMA (200 nM; 15 minutes) (positive) or LY294002 (50  $\mu$ M; 2.25 hr) (negative). Whole cell lysates were added to MSD MULTI-ARRAY 384-well plates coated with anti-ERK1/2 antibody. Phosphorylated ERK1/2 was detected with anti-phospho-ERK1/2 antibody labeled with MSD SULFO-TAG reagent.



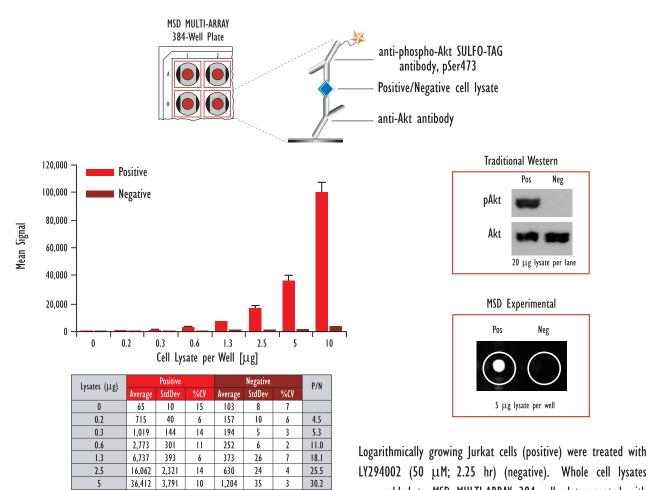
### Detection of Phosphorylated Akt (pSer473)

10

100,129 6,939

7

3,347



170

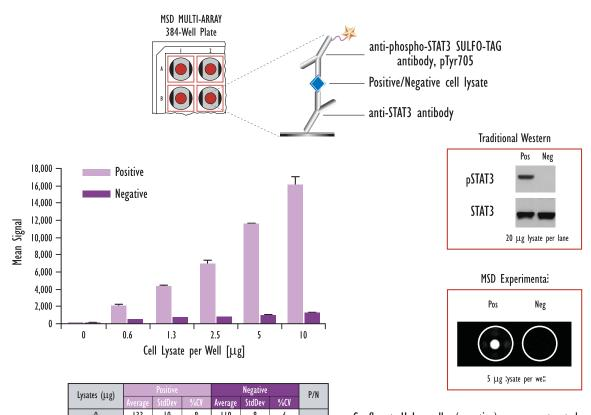
5

29.9

LY294002 (50  $\mu$ LM; 2.25 hr) (negative). Whole cell lysates were added to MSD MULTI-ARRAY 384-well plates coated with an anti-total-Akt antibody. Phosphorylated Akt was detected with anti-phospho-Akt antibody labeled with MSD SULFO-TAG reagent.



### Detection of Phosphorylated STAT3 (pTyr705)



Confluent HeLa cells (negative) were pretreated with sodium vanadate (1mM, 4h) and stimulated with Oncostatin M (40ng/mL, 5min)(positive). Whole cell lysates were added to MSD MULTI-ARRAY 384-well plates coated with an anti-total-STAT3 antibody. Phosphorylated STAT3 was detected with antiphospho-STAT3 antibody labeled with MSD SULFO-TAG reagent.

Lysates (µg)		Positive		Negative			P/N
	Average	StdDev	%CV	Average	StdDev	%СV	171
0	133	10	8	119	8	6	
0.6	2,133	164	8	491	40	8	4.3
1.3	4,395	45	1	739	63	9	5.9
2.5	6,920	460	7	834	44	5	8.3
5	11,580	124		1,052	16	2	11.0
10	16,103	982	6	1,271	74	6	12.7

# 

#### Detection of Phosphorylated c-Met (pTyr1349)

1.3

2.5

5

10

13,704

26,006

45,257

76,342

2,148

1,737

1,891

5,632

16

7

4

7

692

1,034

1,499

2,098

31

43

19

15

5

4

I

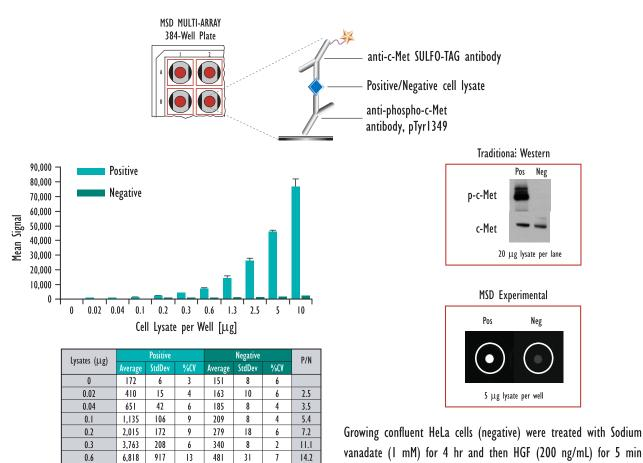
Ι

19.8

25.2

30.2

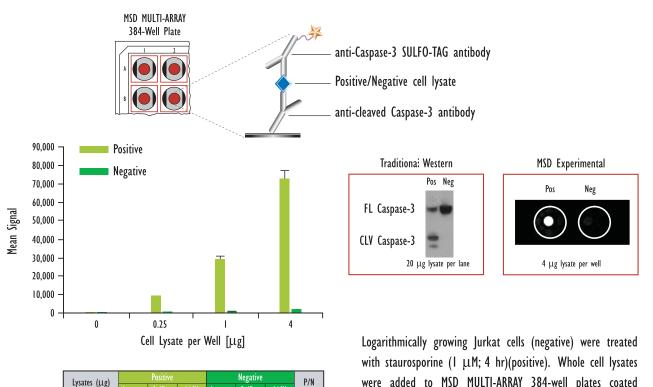
36.4



vanadate (1 mM) for 4 hr and then HGF (200 ng/mL) for 5 min (positive). Whole cell lysates were added to MSD MULTI-ARRAY 384-well plates coated with anti-phospho-c-Met antibody. Phosphorylated c-Met was detected with anti-c-Met antibody labeled with MSD SULFO-TAG reagent.

# 

#### Detection of Cleaved, Active Caspase-3

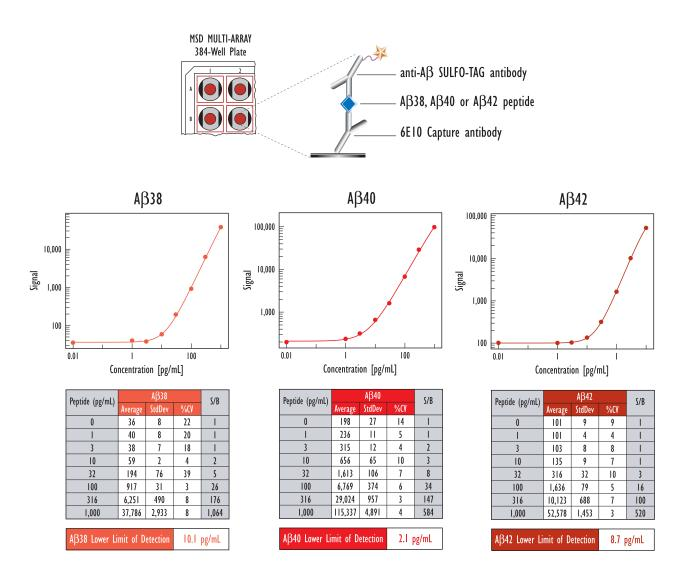


Lysates (µg)	T USILITE			negative			P/N	
Lysaics (prg)	Average	StdDev	%CV	Average	StdDev	%CV	170	
0	54	8	14.9	40	5	11.6		
0.25	8,989	417	4.6	264	42	15.8	34.1	
I	28,930	1,674	5.8	653	37	5.7	44.3	
4	72,361	4,950	6.8	1,625	81	5.0	44.5	

Logarithmically growing Jurkat cells (negative) were treated with staurosporine (1  $\mu$ M; 4 hr)(positive). Whole cell lysates were added to MSD MULTI-ARRAY 384-well plates coated with anti-cleaved Caspase-3 antibody. Cleaved, active Caspase-3 was detected with an anti-Caspase-3 antibody labeled with MSD SULFO-TAG reagent.



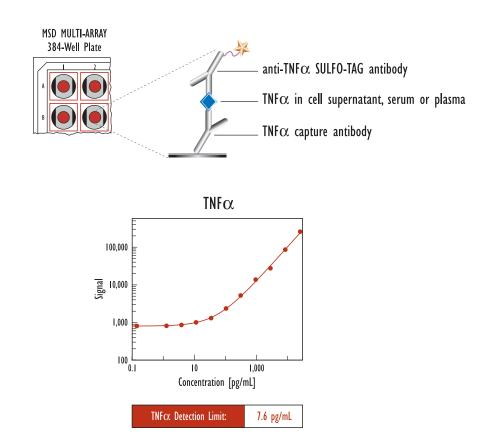
#### Ultra-Sensitive Assays for Amyloid Peptides A $\beta$ 42, A $\beta$ 40, and A $\beta$ 38



The A $\beta$  peptides are fragments of the amy:oid precursor protein (APP) formed by sequentia: c:eavage of APP by  $\beta$ -secretase and  $\gamma$ -secretase. One of the A $\beta$  peptides, A $\beta$ 42, is the major component of amy:oid p:aques, the extra-ce::u:ar protein deposits characteristica::y seen in the brains of patients with A:zheimer's Disease (AD). MSD has developed nove: antibodies to A $\beta$ 38, A $\beta$ 40, and A $\beta$ 42. U:tra-sensitive immuno-assays were developed for the amy:oid peptides in combination with the 6E:0 antibody that binds amino acids 3-8 near the N-terminus of the peptides. Here, synthetic peptides were diluted into a BSA solution and titrated in MSD MULT:-ARRAY 384-we:: p:ates coated with 6E:0 antibody. The A $\beta$  peptides were individua::y detected with MSD's peptide-specific antibodies :abe:ed with SULFO-TAG.



#### Detection of Human Cytokine TNFlpha



TNF $\alpha$  regulates a wide range of cellular and systemic effects. It is one of the most commonly studied biomarkers in drug discovery. Quantitative measurements of TNF $\alpha$  are required for cell supernatants, serum and plasma. Here, recombinant, human TNF $\alpha$  had been diluted into tissue culture medium and titrated into MSD 384-well MULTI-ARRAY plates coated with an anti-TNF $\alpha$  capture antibody. In this homogenous assay, 10  $\mu$ L of sample and 20  $\mu$ L of anti-TNF $\alpha$  detection antibody in MSD Read Buffer are added sequentially to the well. The plate is incubated 4 hours and then analyzed in an MSD Sector Imager instrument.



#### Conclusions

- We have developed numerous immuno-assays in 384-well format for a wide range of biomarkers.
- The assays retain the performance of 96-well assays but require as little as 10 microliters of sample.
- The performance of the 384-well assays for intracellular phosphoproteins makes them compatible with screening applications using 384-well tissue cultures.
- Secreted proteins can be quantified from cell culture supernatants and animal fluids including serum, plasma, CSF and others.