

# Development of a Multiplex Screening Panel for Akt Signaling Pathway Biomarkers in Cell and Tissue Lysate Models

Thomas Miller, Karen Tressler, Jill Dunty, Paula Eason, Leonid Dzantiev, Sripriya Ranganathan, Laura Schaefer, David Stewart, Pankaj Oberoi, and Jacob N. Wohlstatter

Meso Scale Discovery® , Rockville, Maryland, USA

## 1 Abstract

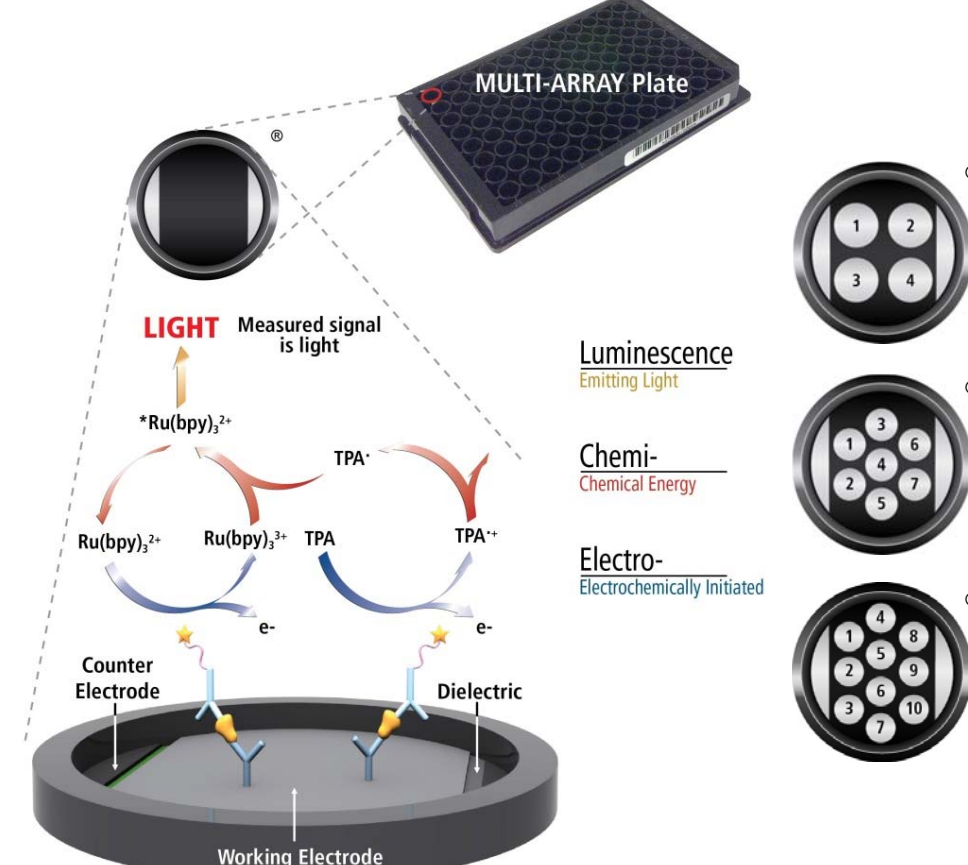
Major challenges face the large-scale implementation of intracellular tumor biomarkers for clinical diagnosis and therapeutic development. These challenges include non-quantitative results, insufficient assay sensitivity, and lack of multiplexing capability. For many studies, sample volumes are limited (e.g., tumor lysates); however, it is essential to extract as much biomarker information as possible from a given sample while maintaining the quality and consistency necessary to sustain ongoing studies. To address these challenges, MSD® has developed multiplex panels for assaying cell signaling biomarkers using fit-for-purpose methods that emphasize optimal multiplex combinations and rigorous development of critical reagents. Here we report the development and verification of a multiplex screening panel of immunoassays for simultaneous measurement of total and phosphorylated analytes of the Akt signaling pathway using MSD technology. Markers in the panel include phosphorylated and total forms of GSK3B, p70S6K, FOXO3a, PTEN, Akt, S6RP, PRAS40, and ERK1.

We demonstrate the ability of these assays to measure analyte levels in multiple tumor-derived cell lines and human tissue samples (normal and tumor) with excellent sensitivity and performance. Most analytes can be quantified using no more than 10 μg of sample. By using recombinant proteins to calibrate some of the assays, we were able to quantify analytes in cultured cell lysates, including MCF-7 and Jurkat T-cells. Lack of analyte specificity is a well-known issue when multiplexing intracellular signaling analytes, especially those in a common signaling pathway such as the Akt pathway. Through rigorous optimization, we achieved non-specific binding less than or equal to 2% for all analytes in this panel. Spike recovery and dilution linearity were matrix tolerant in tumor cell line and tissue lysates. In addition, the assays exhibited a dynamic range between 3 and 4 logs. This characterization allowed for quantification of analytes at different abundance levels without using multiple dilutions.

In conclusion, MSD MULTI-SPOT® Akt signaling assay panels have been developed and verified for measurement of intracellular tumor biomarkers of relevance to clinical diagnosis, therapeutic development, and treatment of various cancers. The accuracy, reliability, ease of use, and high-throughput features of these multiplex assays make them ideally suited for use in large-scale clinical studies.

## 2 Methods

MSD's electrochemiluminescence detection technology uses SULFO-TAG™ labels that emit light upon electrochemical stimulation initiated at the electrode surfaces of MULTI-ARRAY® and MULTI-SPOT® microplates. The multiplex panels were developed as two 8-plex panels designed for optimal performance. The assays use a common Total-analyte specific capture antibody and an assay specific total or phospho-specific detection antibody. Assay controls were prepared by diluting blended Jurkat and MCF7 cell lysates to 400 μg/mL, 100 μg/mL, and 25 μg/mL in lysis buffer supplemented with protease and phosphatase inhibitors. We performed a complete analytical validation of the kits.



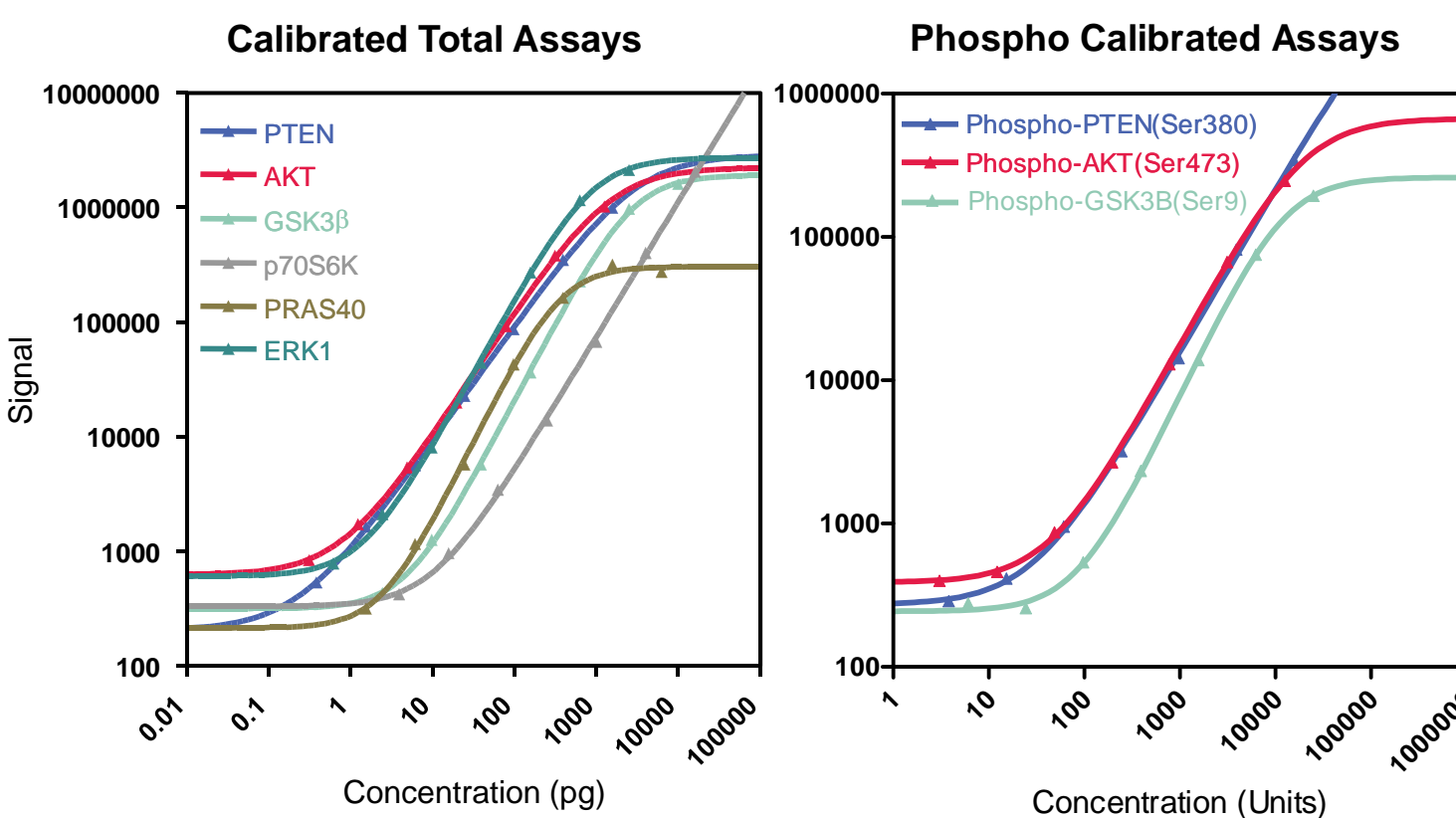
## Electrochemiluminescence Technology

- Minimal non-specific background and strong responses to analyte yield high signal-to-background ratios.
- The stimulation mechanism (electricity) is decoupled from the response (light signal), minimizing matrix interference.
- Only labels bound near the electrode surface are excited, enabling non-washed assays.
- Labels are stable, non-radioactive, and directly conjugated to biological molecules.
- Emission at ~620 nm eliminates problems with color quenching.
- Multiple rounds of label excitation and emission enhance light levels and improve sensitivity.
- Carbon electrode surface has 10X greater binding capacity than polystyrene wells.
- Surface coatings can be customized.

## 3 Protocol and Calibration Curves

The panels use a simple 4-step protocol that can be completed in 3-3.5 hours. Representative calibration curves are shown below for each of the 9 Calibrated assays. While only calibrators for the 9 specified assays were available at the time this study was completed, MSD is continuing to develop calibrators for the remaining assays.

- Add 150 μL Blocker A. Incubate for 1 hour at room temperature (RT).
- Wash with TBS-T. Add 25 μL of calibrator, control, or diluted sample. Incubate for 1 hour at RT.
- Wash with TBS-T. Add 25 μL of detection antibody. Incubate for 1 hour at RT.
- Wash with TBS-T. Add 150 μL of 1X Read Buffer T. Read on MSD imager.



## 4 Sensitivity

The lower limit of detection (LLOD) is a calculated concentration corresponding to the signal 2.5 standard deviations above the background (zero calibrator).

GSK3B			p70S6K			PTEN			
Conc. (pg)	Average Signal	%CV	Conc. (pg)	Average Signal	%CV	Conc. (pg)	Average Signal	%CV	
Cat-1	10000.00	1596397	9.6	16000.00	2059368	0.2	1562.50	988802	8.3
Cat-2	2500.00	971158	14.5	4000.00	401055	5.0	390.63	348848	6.8
Cat-3	625.00	225626	1.6	1000.00	67571	6.9	97.68	86694	0.7
Cat-4	156.25	36455	8.4	250.00	13887	3.6	24.41	22713	0.8
Cat-5	39.06	5728	8.5	62.50	3445	5.5	6.10	5632	9.0
Cat-6	9.77	1265	11.9	15.63	966	4.1	1.53	1641	1.2
Cat-7	2.44	432	12.3	3.91	422	8.9	0.38	536	10.4
Cat-8	0	144	6.9	0	234	10.0	0	198	0.0
	<b>1.25</b>	<b>LLOD</b>		<b>2.50</b>	<b>LLOD</b>		<b>0.09</b>	<b>LLOD</b>	

AKT			PRAS40			ERK1			
Conc. (pg)	Average Signal	%CV	Conc. (pg)	Average Signal	%CV	Conc. (pg)	Average Signal	%CV	
Cat-1	1250.00	1024694	3.9	6250.00	274312	26.8	2500.00	2116114	0.7
Cat-2	312.50	380428	4.6	1562.50	319782	6.4	625.00	1155337	4.1
Cat-3	78.13	92211	0.1	390.63	163216	2.3	156.25	270121	0.7
Cat-4	19.53	19866	3.1	97.66	42839	2.1	39.06	44137	4.6
Cat-5	4.88	5411	2.8	24.41	5702	16.3	9.77	8059	0.9
Cat-6	1.22	1727	5.4	6.10	1164	1.2	2.44	2101	2.8
Cat-7	0.31	838	2.6	1.53	317	8.5	0.61	785	4.1
Cat-8	0	531	0.5	0	154	12.9	0	426	8.5
	<b>0.10</b>	<b>LLOD</b>		<b>1.02</b>	<b>LLOD</b>		<b>0.27</b>	<b>LLOD</b>	

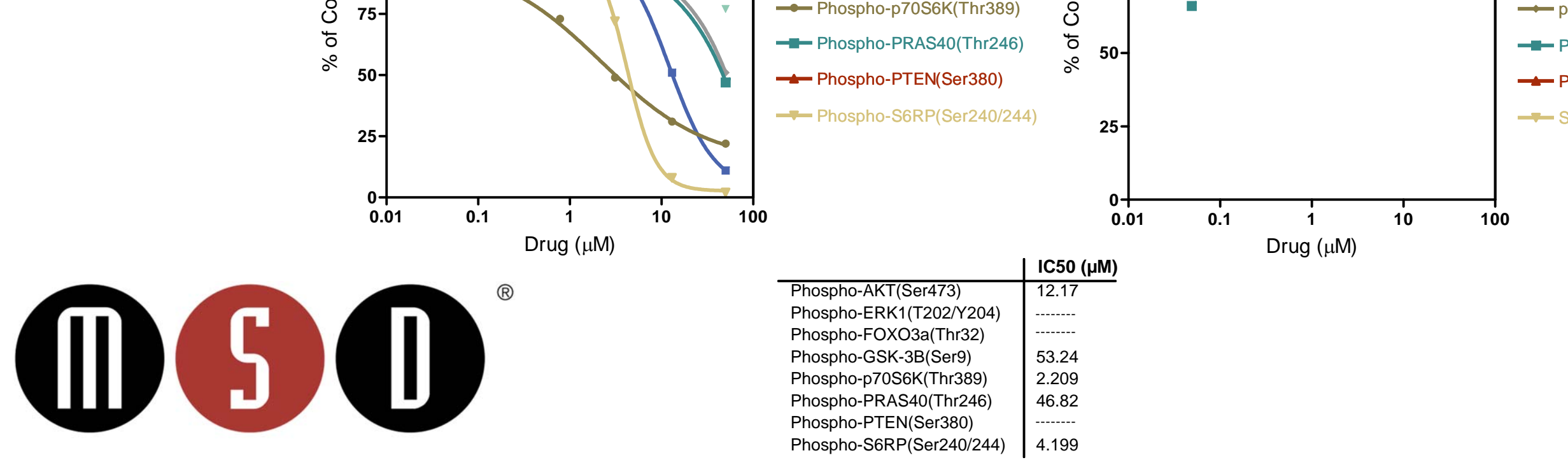
Phospho-GSK3B(Ser9)			Phospho-PTEN(Ser380)			Phospho-AKT(Ser473)			
Conc. (Units)	Average Signal	%CV	Conc. (Units)	Average Signal	%CV	Conc. (Units)	Average Signal	%CV	
Cat-1	100000	208121	10.0	15625	314200	4.9	12500	150284	7.1
Cat-2	25000	61834	6.1	3906	67487	4.7	3125	35749	5.6
Cat-3	6250	8955	3.3	977	11238	3.4	781	5835	2.9
Cat-4	1563	1319	5.2	244	2958	3.0	195	1453	4.1
Cat-5	391	348	8.4	61	806	6.2	49	754	2.7
Cat-6	98	199	7.0	15	405	4.9	12	537	6.0
Cat-7	24	162	28.6	4	329	4.5	3	508	7.5
Cat-8	6	152	21.4	1.0	377	17.1	0.8	473	7.0
	<b>277</b>	<b>LLOD</b>		<b>16.1</b>	<b>LLOD</b>		<b>29</b>	<b>LLOD</b>	

## 5 Specificity and Reproducibility

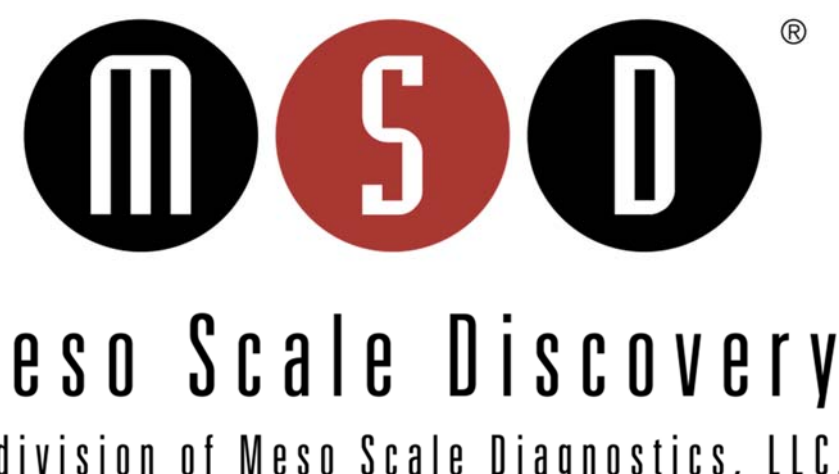
Each assay in the 16-plex panel was evaluated for cross-reactivity with other analytes on the panel as well as related analytes that are not on the panel. The cross-reactivity for all assays was <2%. To evaluate reproducibility, three levels of controls spanning the quantifiable range of the panels were prepared and measured over multiple runs. The %CV for each assay of the 16-plex assays is shown below. The controls were prepared for single-use, and 7 runs were conducted.

CTL-1	Non-Specific Interaction	Total Detect								
		PTEN	Akt	GSK3b	FOXO3a	PRAS40	ERK1	p70S6K	S6RP	
Capture	PTEN	100.00%	0.00%	0.02%	0.01%	0.05%	0.02%	0.01%	<0.1%	
	Akt	0.06%	100.00%	0.15%	0.01%	0.05%	0.07%	0.02%	<0.1%	
	GSK3b	0.65%	<0.1%	100.00%	0.02%	0.00%	<0.1%	0.14%	0.73%	
	FOXO3a	0.08%	0.02%	0.01%	100.00%	0.04%	0.01%	0.49%	0.31%	
	PRAS40	0.09%	<0.1%	0.00%	0.42%	100.00%	0.00%	0.00%	0.02%	
	ERK1	0.03%	0.04%	0.01%	0.01%	0.04%	100.00%	0.01%	0.08%	
	p70S6K	<0.1%	<0.1%	0.00%	0.00%	<0.1%	0.01%	100.00%	<0.1%	
	S6RP	0.08%	0.04%	0.01%	0.05%	0.00%	0.01%	0.32%	100.00%	
CTL-1	Non-Specific Interaction	Phospho Detect								
		PTEN	100.00%	0.00%	0.02%	0.36%	0.03%	0.14%	0.00%	0.04%
		Akt	0.05%	100.00%	0.15%	0.05%	0.00%	0.27%	0.00%	0.03%
		GSK3b	0.65%	<0.1%	100.00%	0.02%	0.00%	<0.1%	0.17%	0.00%
		FOXO3a	1.02%	0.02%	0.00%	100.00%	0.00%	0.16%	0.00%	1.21%
		PRAS40	1.34%	<0.1%	<0.1%	0.05%	100.00%	0.01%	0.00%	0.01%
		ERK1	0.14%	1.16%	0.00%	0.09%	0.02%	100.00%	0.00%	0.09%
		p70S6K	0.02%	0.02%	0.01%	0.01%	0.00%	<0.1%	100.00%	0.01%
S6RP	1.28%	0.04%	0.04%	1.91%	0.01%	<0.1%	0.00%	100.00%		

Assay	Sample ID	AVG Intra-run Signal	CV	AVG Intra-run Conc CV
GSK3B	Chi-1	6.6%	5.7%	23.0%
	Chi-2	5.5%	4.4%	6.2%
	Chi-3	5.1%	4.2%	4.1%
p70S6K	Chi-1	2.6%	2.4%	11.7%
	Chi-2	1.8%	1.5%	6.0%
	Chi-3	3.7%	1.5%	4.4%
FOXO3a	Chi-1	2.8%	N/A	N/A
	Chi-2	4.1%	N/A	N/A
	Chi-3	2.0%	N/A	N/A
PTEN	Chi-1	0.8%	0.8%	5.2%
	Chi-2	4.6%	4.6%	3.5%
	Chi-3	4.4%	4.6%	3.4%
AKT	Chi-1	4.4%	3.9%	2.6%
	Chi-2	7.2%	6.3%	3.6%
	Chi-3	4.6%	4.3%	3.3%
S6RP	Chi-1	1.3%	N/A	N/A
	Chi-2	1.7%	N/A	N/A
	Chi-3	4.4%	N/A	N/A
PRAS40	Chi-1	3.9%	3.5%	N/A
	Chi-2	6.0%	4.9%	N/A
	Chi-3	2.2%	1.8%	N/A
ERK1	Chi-1	4.8%	4.5%	N/A
	Chi-2	3.8%	3.1%	N/A
	Chi-3	2.2%	1.8%	N/A



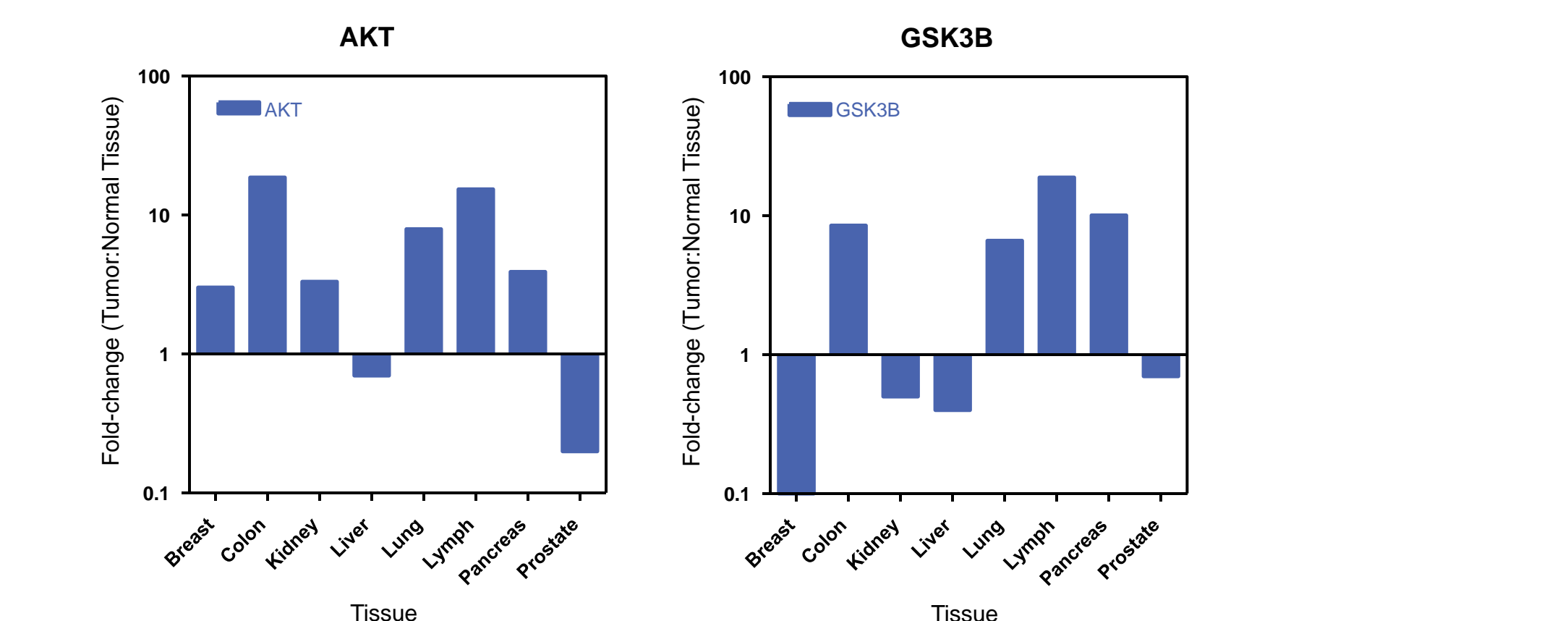
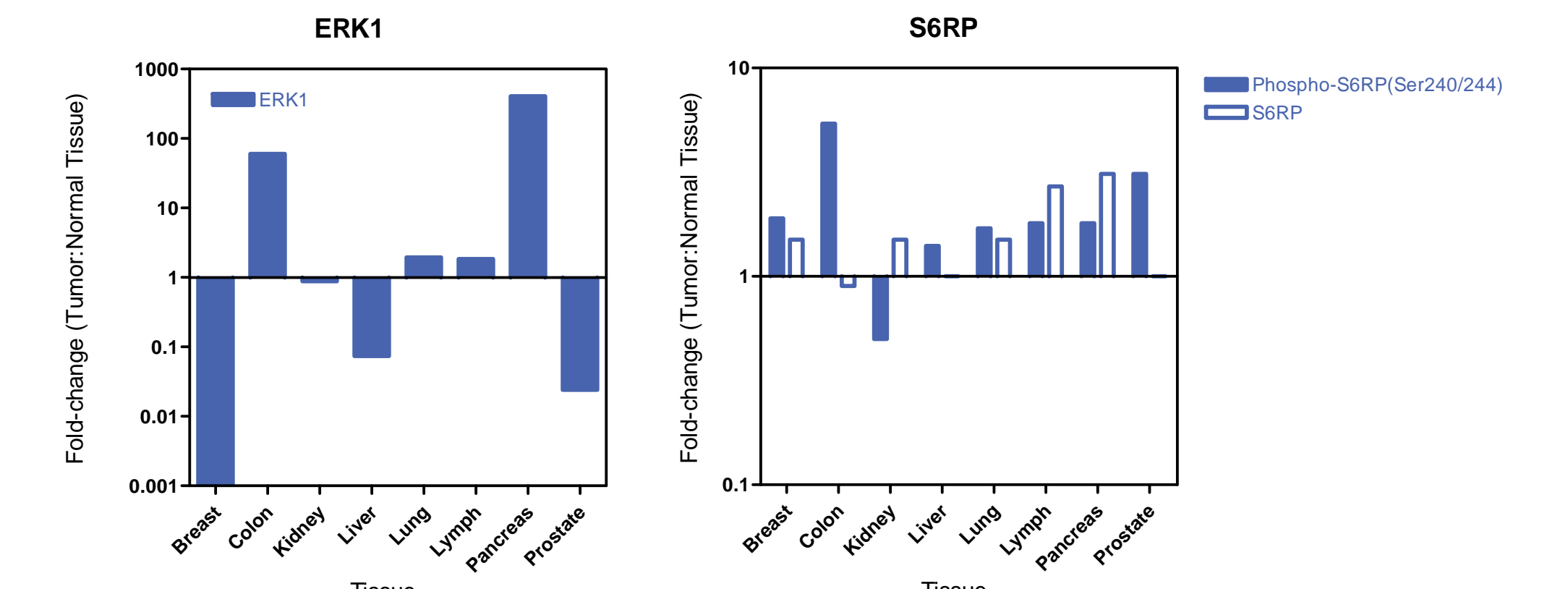
IC50 (μM)	
Phospho-AKT(Ser473)	12.17
Phospho-ERK1(T202/Y204)	-----
Phospho-FOXO3a(Thr32)	-----
Phospho-GSK-3B(Ser9)	53.24
Phospho-p70S6K(Thr389)	2.209
Phospho-PRAS40(Thr246)	46.82
Phospho-PTEN(Ser380)	-----
Phospho-S6RP(Ser240/244)	4.199



## 8 Quantitation in Human Tissue Samples

Total and phospho-analytes were measured in normal and tumor human tissue lysates (obtained from Novus Biologicals) from 5 μg of sample. Analyte levels varied across tissue types and normal/tumor status indicated by calculated concentration for calibrated assays and mean signal for non-calibrated assays in the table. ND indicates not detectable. Optimization of lysis conditions may improve analyte detection. Samples were not collected with optimal concentrations of phosphatase inhibitors. Graphs below are representative of analyte changes across tissues expressed as fold change in tumor to normal.

		Phospho-AKT (Ser473)	Phospho-ERK1 (T202/Y204)	Phospho-GSK3B (Ser9)	p70S6K	Phospho-p70S6K (Thr389)	Phospho-PTEN (Ser380)	FOXO3a	Phospho-FOXO3a (Thr32)	Phospho-S6RP (Ser240/244)					
											AKT (pg)	ERK1 (pg)	GSK3B (pg)	(Units)	(pg)
Normal	Breast	1.05	ND	587.12	ND	13.00	ND	3.07	ND	ND	ND	4731	1103		
	Colon	ND	ND	5.10	ND	3.04	ND	2.06	ND	ND	ND	4740	1893		
	Kidney	0.04	ND	98.48	ND	2.79	ND	1.60	ND	ND	ND	4819	1758		
	Liver	0.28	4.46	103.31	ND	1.83	47.73	1.43	ND	0.06	ND	4938	1874		
	Lung	0.08	ND	82.09	ND	0.74	ND	1.60	ND	0.07	ND	2632	1233		
	Lymph	0.21	ND	173.26	ND	0.46	ND	1.45	ND	0.08	ND	2861	1072		
	Pancreas	0.01	1.22	0.08	ND	0.34	ND	1.05	ND	0.05	ND	3105	1178		
	Prostate	2.29	3.56	335.18	ND	9.67	17.45	1286	0.36	ND	4856	ND	7702	973	
Tumor	Breast	4.46	3.74	0.29	ND	1.66	ND	4.58	ND	0.43	ND	9235.00	ND	7072	2130
	Colon	10.25	18.27	308.82	5476.00	21.80	257.33	8.98	764.00	2.75	ND	5357.00	ND	4074	10283
	Kidney	1.41	ND	86.87	ND	1.18	ND	2.01	ND	0.11	27.98	ND	ND	7153	909
	Liver	0.06	ND	7.37	ND	0.48	ND	1.73	ND	0.07	ND	ND	ND	4708	2687
	Lung	4.63	ND	157.14	3569.00	6.64	ND	2.29	ND	0.17	ND	3852	2050		
	Lymph	11.98	21.31	317.60	ND	13.95	66.71	21.38	1495.00	0.80	14.18	7801.00	1333.00	7658	1930
	Pancreas	1.68	5.53	255.20	ND	7.19	161.03	2.66	ND	0.19	ND	ND	ND	9594	2142
	Prostate	0.05	ND	8.37	ND	6.68	ND	2.12	ND	0.10	ND	4007	1868.00	7550	2975



## 9 Conclusion

A multiplex screening panel has been developed and validated for use in quantitating Akt signaling analytes in a variety of human samples. The assays in the panel have sufficient sensitivity, reproducibility, accuracy, and sample performance to allow testing of multiple sample types in cancer biology applications. Consistent results were demonstrated for calibration curves as well as for control samples over multiple runs on different days by different operators. All of the assays in the panel can be run using two fit-for-purpose-designed 8-plex panels that use a simple three-step protocol. Biomarker screening can be performed with as little as 5 μg of sample. Evaluation of various disease samples using this panel demonstrates that it can be an effective tool for cancer biomarker and drug discovery.

