

# Quantitative Profiling of Signaling Pathways Using PTMScan<sup>®</sup> Direct

### Introduction

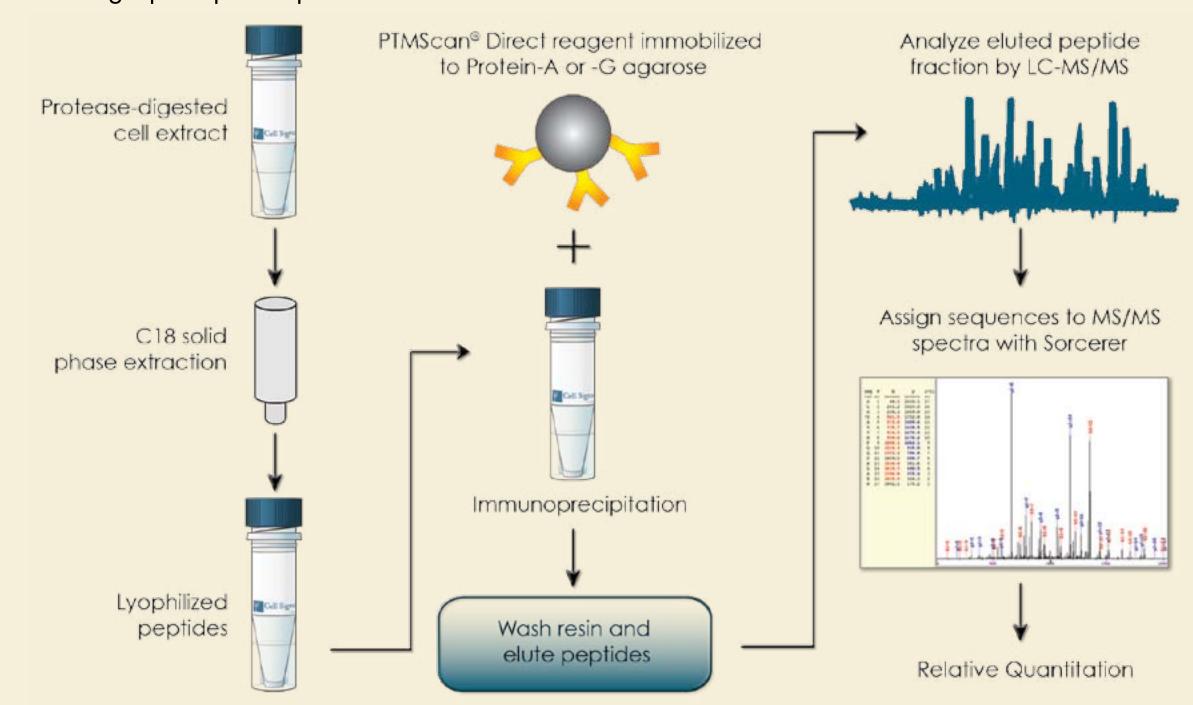
A challenge to proteomic studies of post-translational modifications (PTMs) is the ability to focus on proteins of interest rather than randomly identified proteins typified by data dependent analysis. Methods that target specific proteins and pathways known to be critical to cellular signaling are therefore desirable. Cell Signaling Technology (CST) has developed an immunoaffinity-based LC-MS/MS method, called PTMScan<sup>®</sup> Direct, for identification and quantification of hundreds of peptides from particular signaling pathways or protein types. Cell lines, tissues, or xenografts can be used as starting material. Quantification may be performed using isotopic labeling methods such as SILAC or label-free quantification of chromatographic peak areas. PTMScan Direct is a powerful new method that combines the specificity of antibody-based methods with the sensitivity and high duty cycle of LC-MS/MS analysis allowing thousands of measurements.

### PTMScan<sup>®</sup> Direct Reagents

- **1.** Multipathway Reagent
- 2. Serine/Threonine Kinases Reagent
- **3.** Tyrosine Kinases Reagent
- **4.** Akt / PI3K Pathway Reagent
- 5. DNA Damage / Cell Cycle Reagent
- 6. Apoptosis / Autophagy Reagent

### PTMScan<sup>®</sup> Direct Method

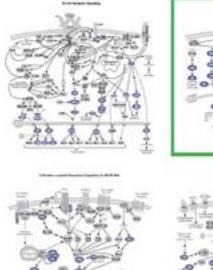
Figure 1: PTMScan Direct work flow diagram. PTMScan Direct is a published method (Stokes et al, 2012) adapted from the original PhosphoScan<sup>®</sup> method developed at Cell Signaling Technology (Rush et al, 2005, Patent # 20030044848). Cell lines, tissues, xenografts, or other biological starting materials are lysed under denaturing conditions, digested with trypsin, and desalted over C18 columns. Target peptides are immunoprecipitated using the appropriate PTMScan Direct reagent. Immunoprecipitated peptides are separated on a reversed-phase Magic C18 AQ column and data-dependent MS methods are performed with an LTQ-Orbitrap Velos or Elite mass spectrometer. MS/MS spectra are evaluated using SEQUEST 3G and the SORCERER 2 v4.0 platform from Sage-N Research (Lundgren et al., 2009). Quantification is performed using chromatographic peak apex areas or intensities.



TECHNOLOGY<sup>®</sup>

Cell Signaling

## Multipathway Reagent Selected Targets



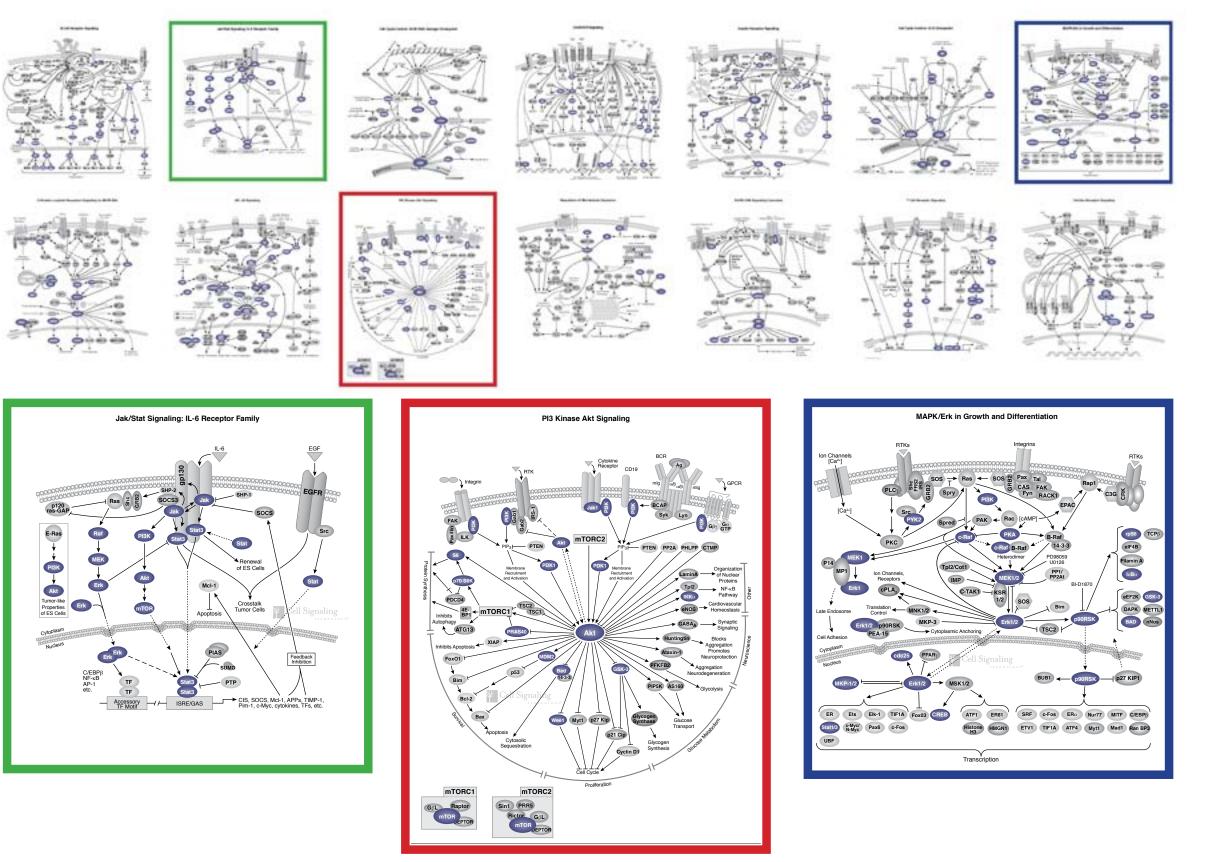
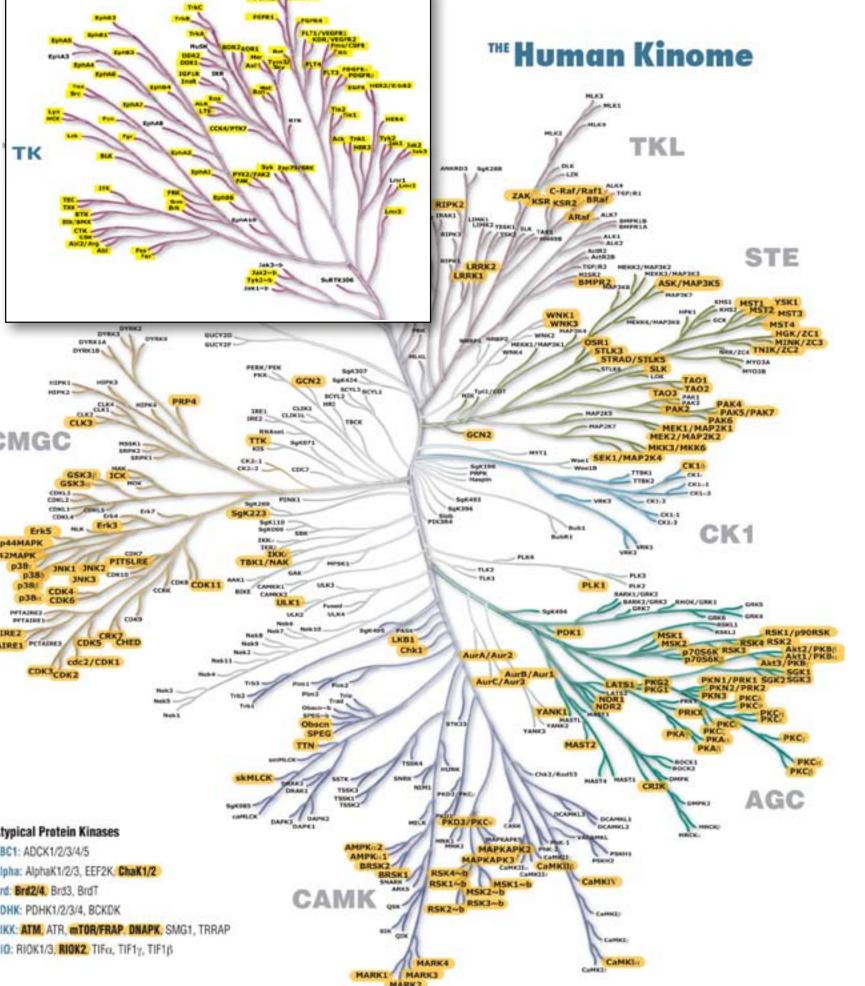
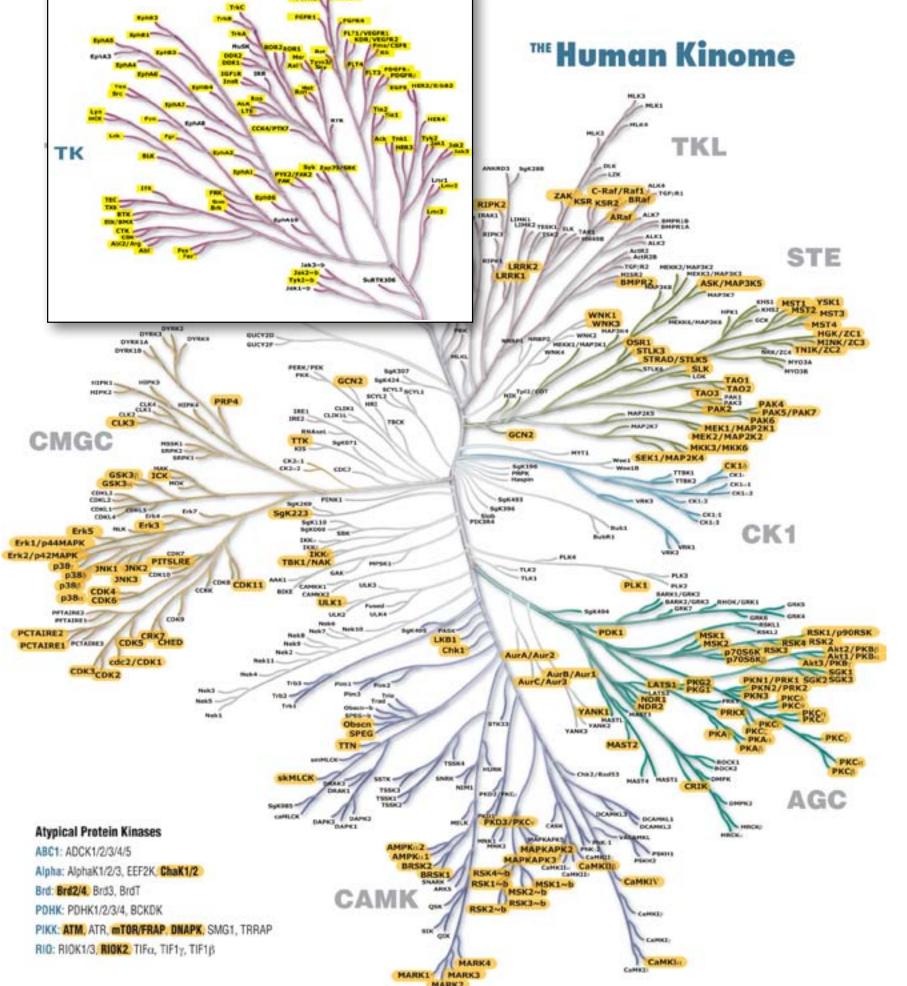


Figure 3: PTMScan Direct Ser/Thr Kinases Reagent and Tyr Kinases Reagent targets mapped onto the human kinome tree. Yellow highlighting indicates kinases identified using the reagents. Tyr Kinases Reagent coverage is shown in the inset ("TK").





### Figure 2: Coverage of selected pathways using the PTMScan Direct Multipathway

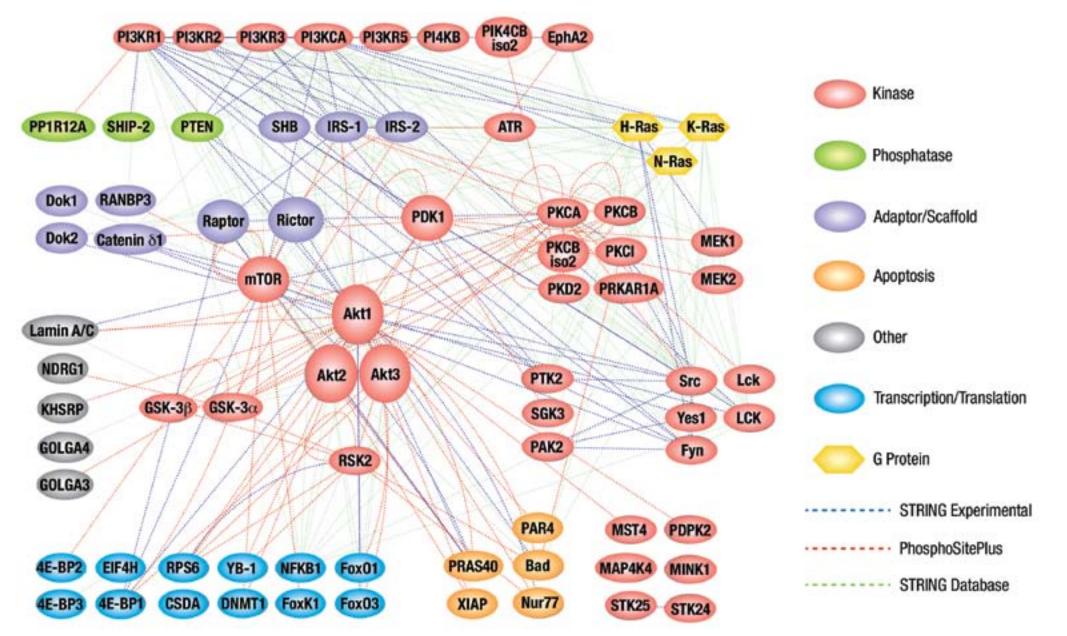
**Reagent.** Peptides from purple proteins are identified using the Multipathway Reagent.

### Ser/Thr Kinases and Tyr Kinases Reagent Targets

## Akt / PI3K Pathway Reagent

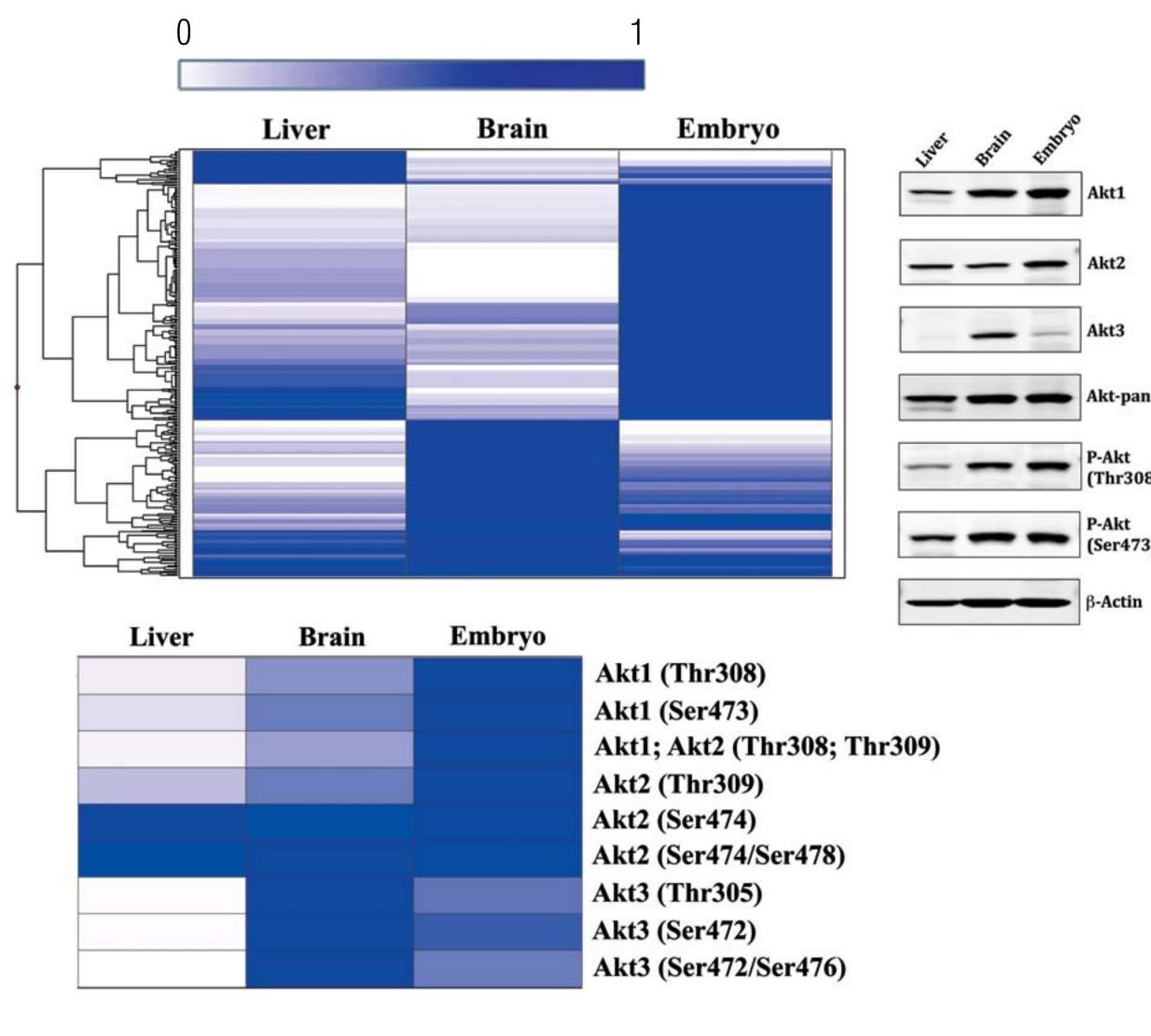
### **INTERACTION MAP**

**Figure 4: PTMScan Direct: Akt/PI3K Pathway Reagent interaction map.** Interactions were derived from STRING using high confidence scores (>0.700) from experimental, database, and textmining lines of evidence and the substrate search function in PhosphoSitePlus<sup>®</sup>. Node colors/ shapes denote different protein classes. Edge color denotes interaction type.



### **MOUSE TISSUE STUDY**

Figure 5: Akt/PI3K Pathway Reagent mouse tissue study. Heat map of relative intensities for mouse liver, brain, and embryo. Each row represents a different peptide ion and each column represents a different tissue. The max intensity across the three tissues was set to 1 and the other two intensities normalized to the max. **Blue** indicates higher intensity.



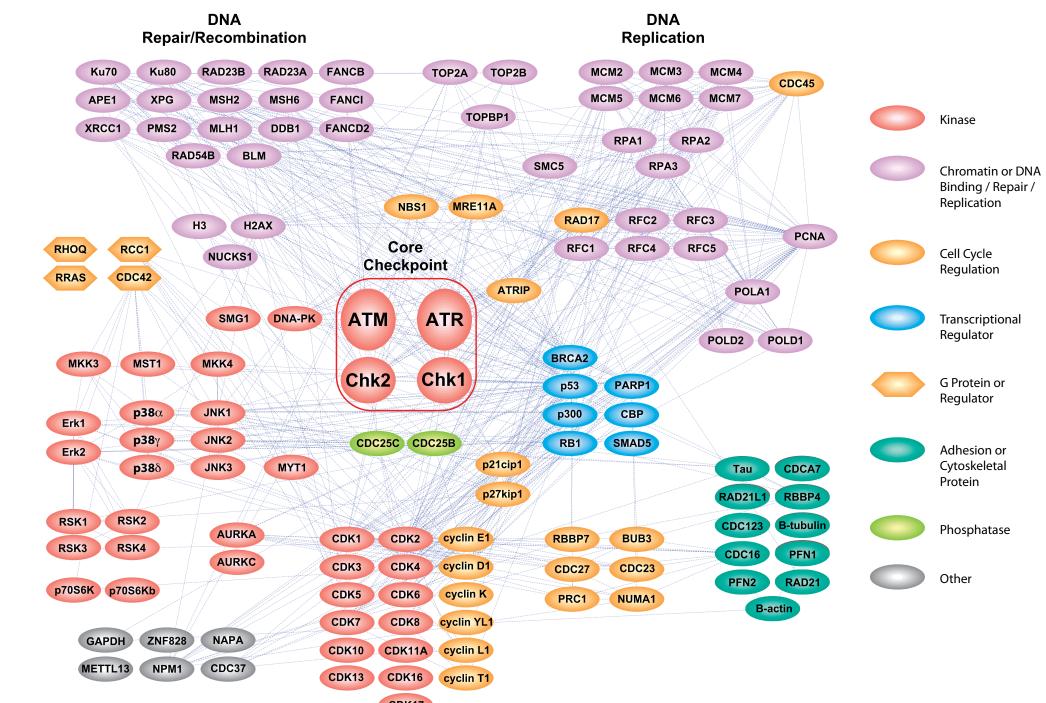
Matthew P. Stokes, Jeffrey C. Silva, Charles L. Farnsworth, Xiaoying Jia, Roberto Polakiewicz, Michael J Comb Cell Signaling Technology, Inc., Danvers, MA

## : DNA Damage/Cell Cycle Reagent :

#### **INTERACTION MAP**

Figure 6: PTMScan Direct: DNA Damage / Cell Cycle Reagent interaction map. Interactions were derived from STRING using using high confidence scores (>0.700) from

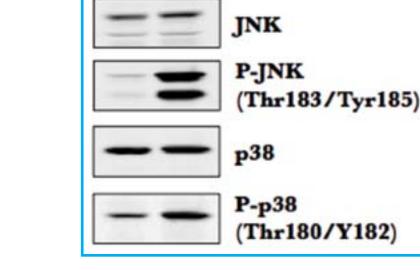
experimental, database, and textmining lines of evidence. Node colors/shapes denote different protein classes.



#### HeLa -/+ UV DAMAGE

Figure 7: DNA Damage/Cell Cycle Reagent: HeLa cell UV damage study. HeLa cells were untreated or treated with 500 mJ/cm<sup>2</sup> UV light and harvested at 2 hr post treatment. Fold changes were calculated from chromatographic peak heights/areas. Green cells indicate peptides more abundant with UV damage, **Red** cells indicate peptides less abundant. Selected DNA damage response proteins are shown in detail with accompanying western blots.

	Fold Change UV : Control	Protein Name	Site	Peptide	Max Inter
i.	1.1	ATM		QAGIIQALQNLGLCHILSVYLK	299,73
	5.5	ATM	§1981	SLAFEEGS*QSTTISSLSEK	20,664,
	-1.6	ATR		YIVLHFGR	46,40
	-1.2	ATR	§428	M#ESM#EIIEEIQCQTQQENLSSNSI	43,72
	-1.3	Chk1	-	FFHQLM#AGVVYLHGIGITHR	48,71
	3.2	Chk1	§317	YSSS*QPEPR	698,57
	4.1	Chk1	§345	LVQGISFS*QPTCPDHM#LLNSQLLG	
	-1.1	Chk2		FTTEEALRHPWLQDEDM#K	7,865,5
	9.1	Chk2	§379	ILGETS*LM#R	1,010,0
	260.2	JNK1; JNK3	§183, §185; §221, §223	TAGTSFM#MT*PY*VVTR	8,543,0
	98.0	JNK2	§183, §185	TACTNFMM#T*PY*VVTR	15,427,
	4.7	p38-alpha	§180, §182	HTDDEMT*GY*VATR	18,211,
	4.1	p38-gamma	§183, §185	QADSEM#T*GY*VVTR	1,421,7
				<ul> <li>UV</li> <li>ATM</li> <li>P-ATM (Ser198</li> <li>ATR</li> <li>P-ATR (Ser428)</li> </ul>	
		ilete il ut		– UV јмк	

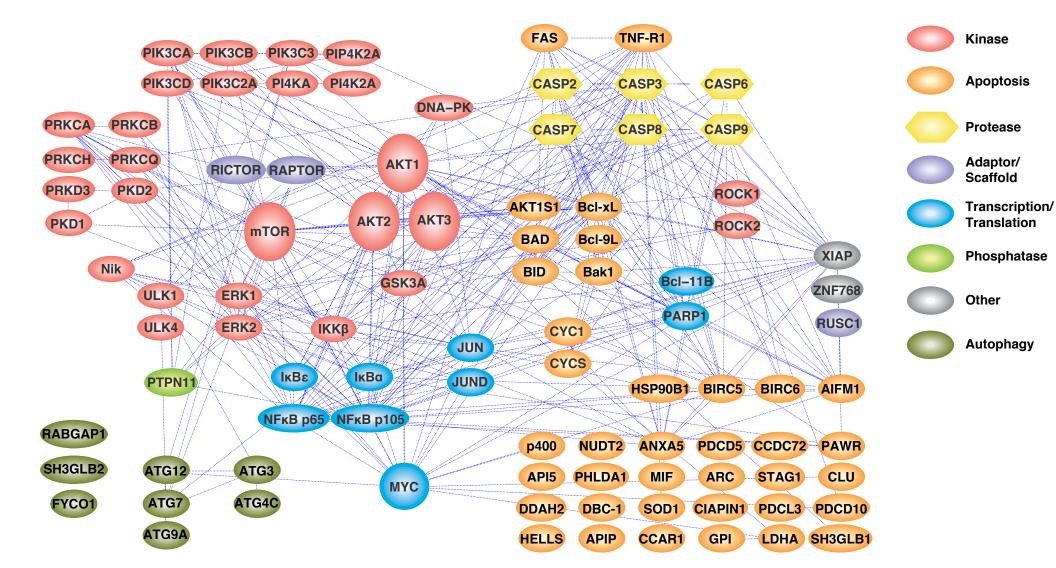


### Apoptosis / Autophagy Reagent

### INTERACTION MAP

### Figure 8: PTMScan Direct: Apoptosis / Autophagy Reagent interaction map.

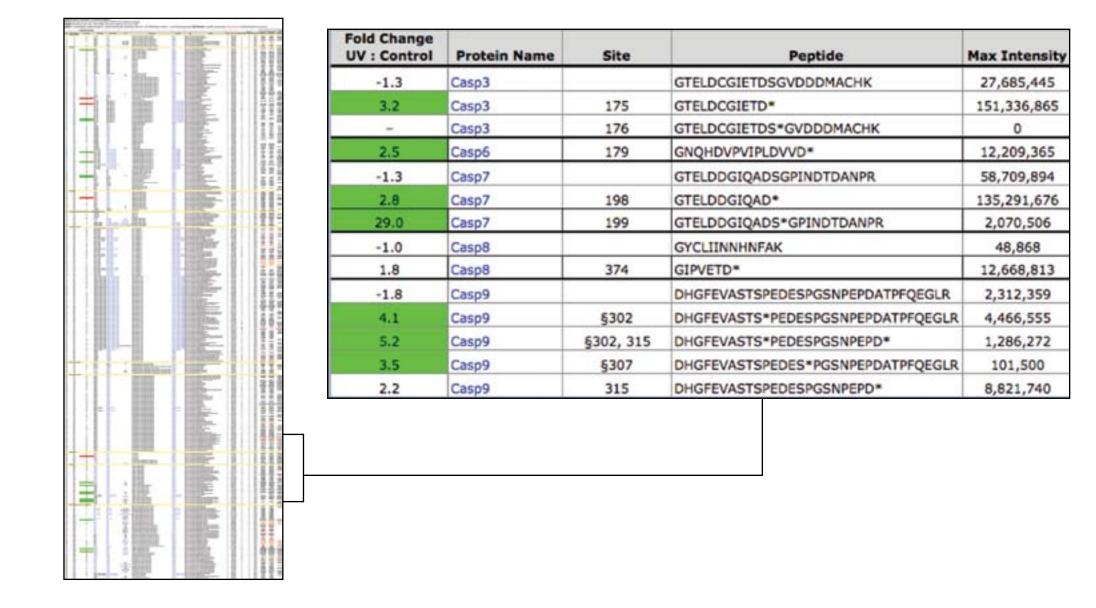
Interactions were derived from STRING using using high confidence scores (>0.700) from experimental, database, and textmining lines of evidence. Node colors/shapes denote different protein classes



#### HeLa -/+ UV DAMAGE

#### Figure 9: Apoptosis / Autophagy Reagent: HeLa cell UV damage study.

Quantitative data for caspase peptides identified with the Apoptosis/Autophagy Reagent used to probe UV damaged HeLa cells (500 mJ/cm<sup>2</sup>, harvest 2 hr post treatment). Green indicates ncreased abundance relative to control.



### Summary

PTMScan Direct is a novel method that allows identification and quantification of hundreds of peptides from selected protein types or signaling pathways. This allows focus on proteins of interest, bypassing the random sampling of peptides that occurs in traditional data-dependent proteomic analysis. PTMScan Direct is widely applicable in drug development and discovery, as well as in any application where monitoring of known signaling pathways is desired.

### References

- **1.** Rush J. et. al. (2005) *Nat. Biotechnol.* 23: 94-101
- **2.** Lundgren, et. al. (2009) *Curr. Protoc. Bioinformatics*. 13: Unit 13-3
- **3.** Stokes M.P. et. al. (2012) *Mol. Cell Proteomics.* Epub Feb 9th

## **Contact Information**

Matthew P. Stokes, Cell Signaling Technology, Inc. Email: mstokes@cellsignal.com **PTMScan Services Department** Email: ptmscan@cellsignal.com • web: www.cellsignal.com/services

