

Quantitative profiling of signaling pathways using immunoaffinity purification and LC-MS/MS

Matthew P. Stokes, Jian Min Ren; Kimberly A. Lee; Xiaoying Jia; Jeffrey C. Silva

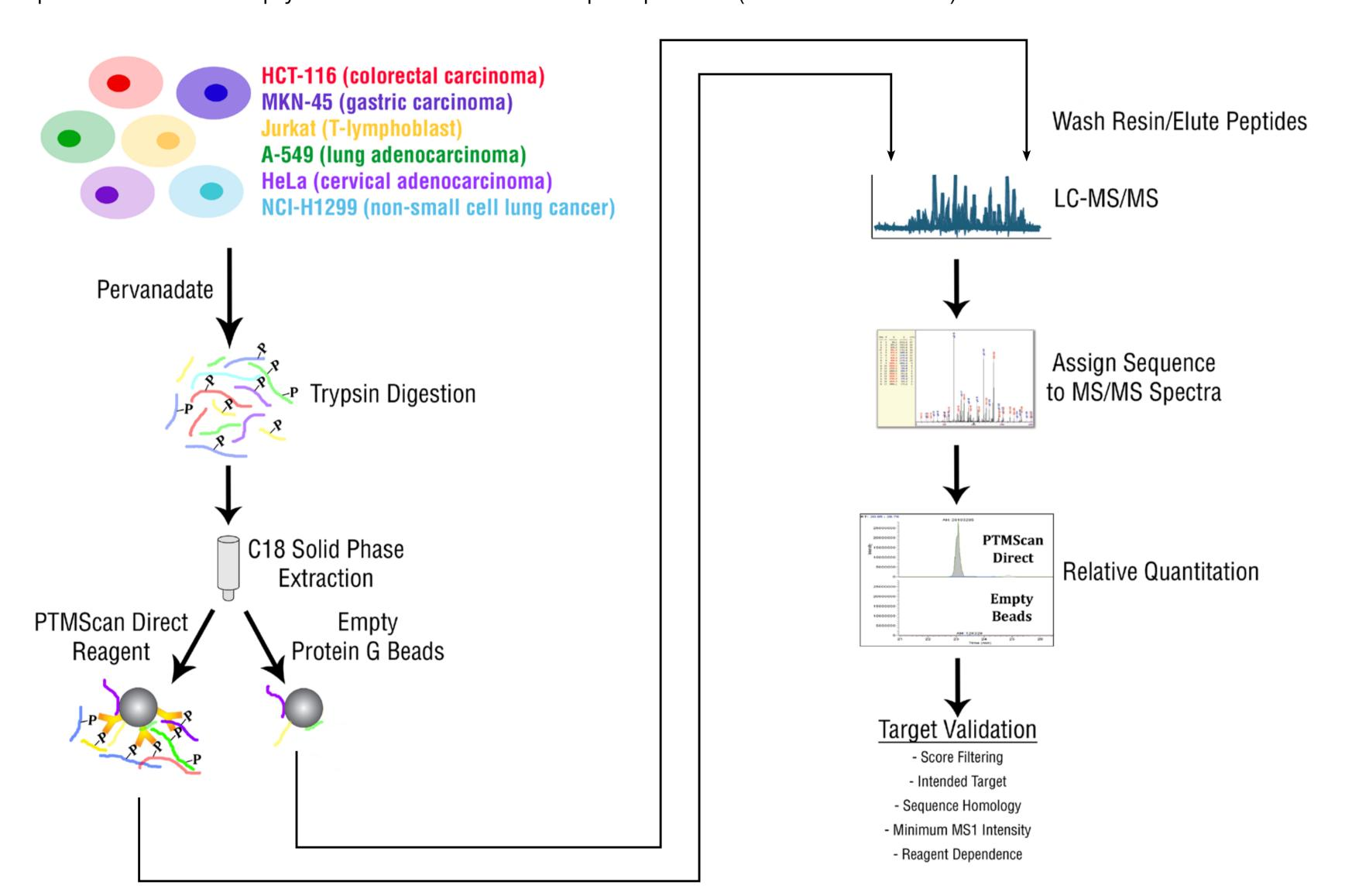
Cell Signaling Technology, Danvers MA USA 01923

Introduction

Proteomic analysis of post-translationally modified peptides has traditionally employed methods that broadly sample the proteome but are unfocused with respect to the sites identified. We have developed a novel immunoaffinity method for identification and quantitation of post-translationally modified peptides from proteins that reside in the same signaling pathway or pathways, allowing a global view of pathway activation from a single LC-MS/MS run. Six different antibody reagents have been prepared that focus on diverse signaling areas: Ser/Thr Kinase activity, Tyr Kinase activity, PI3K/Akt signaling, Cell Cycle/DNA Damage signaling, Apoptotic/Autophagolytic pathways, and a Multipathway reagent for detection of critical signaling proteins across many different pathways. Reagents were validated using human and mouse samples with a variety of treatments (inhibitors, growth factors, etc.). This technology is broadly applicable to any experimental system in which quantitative profiling of specific critical signaling molecules is desirable.

PTMScan® Direct Method

Figure 1: PTMScan® Direct Reagent validation strategy. PTMScan® Direct is a published method (Stokes et al. 2012) adapted from the original PhosphoScan® method (Rush et al. 2005) developed at Cell Signaling Technology. PTMScan Direct Reagents are validated using mixtures of pervanadate-treated human cancer cell lines digested with trypsin. Peptides are desalted over C18 columns and immunoprecipitated with either the PTMScan Direct reagent or empty Protein G beads. Immunoprecipitated peptide mixtures are analyzed by LC-MS/MS and relative quantitation is performed. Validated peptides meet several strict criteria: they are targeted by the reagent or are homologous to a target, they must pass score filtering (Lundgren et al. 2009) and signal intensity thresholds, and they must be present in higher abundance in the PTMScan Direct reagent immunoprecipitation than the empty Protein G beads immunoprecipitation (Stokes et al. 2012).



PTMScan® Direct

- 1. Multipathway (updated)
- 2. Ser/Thr Kinases
- 3. Tyr Kinases
- 4. Akt/PI3K
- 5. DNA Damage/Cell Cycle
- 6. Apoptosis/Autophagy

Interaction maps were generated from IPA® pathway analysis or the STRING database (string.embl.de) using high confidence scores (>0.700) from experimental, database, and text mining lines of evidence. Interactions were also defined from the substrate search page of PhosphoSitePlus® (www.phosphosite.org). Node colors and shapes denote different protein classes. Edge color denotes interaction type.

Multipathway

Monitoring more than 1,000 phosphorylation sites on over 400 critical signaling proteins.

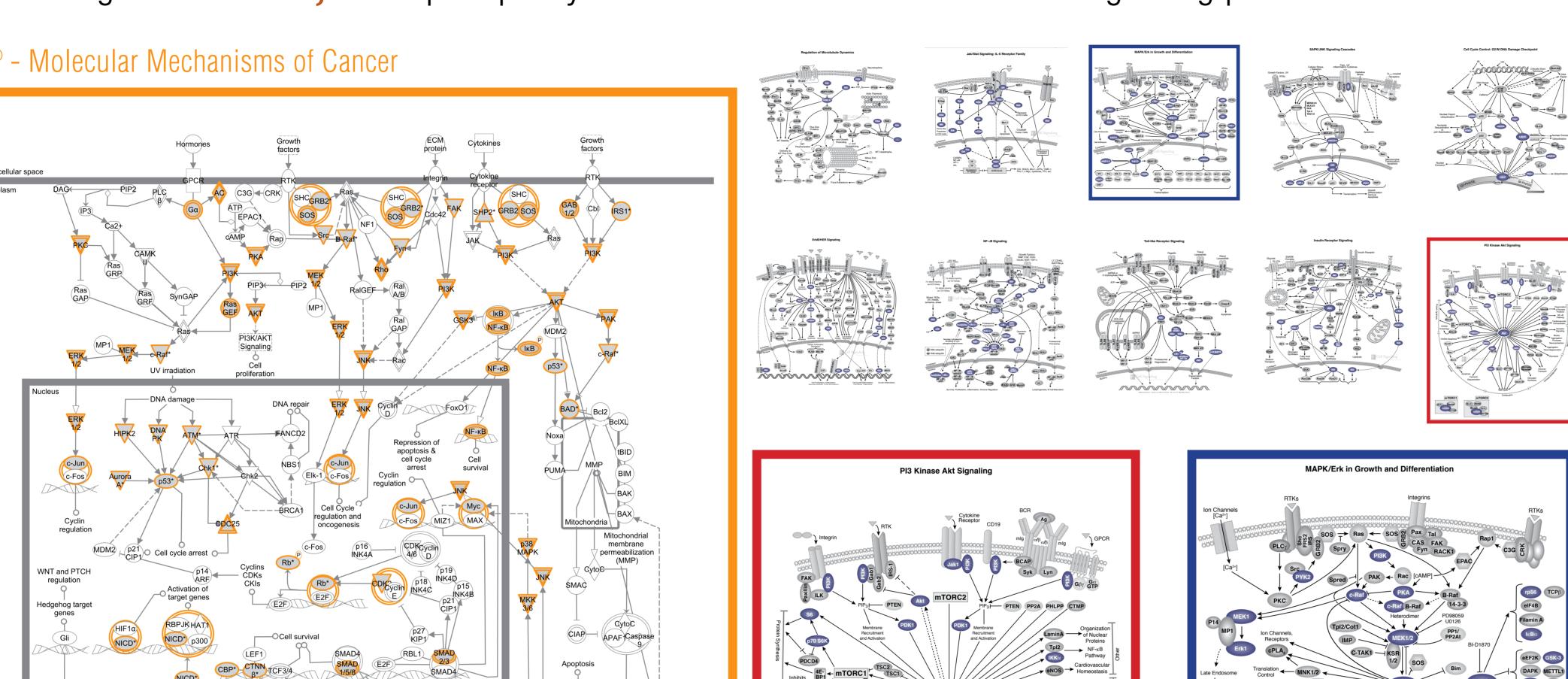


Figure 2: Multipathway coverage of selected signaling pathways.

Ser/Thr Kinases and Tyr Kinases

Profiling sites of activation and inhibition on protein kinases.

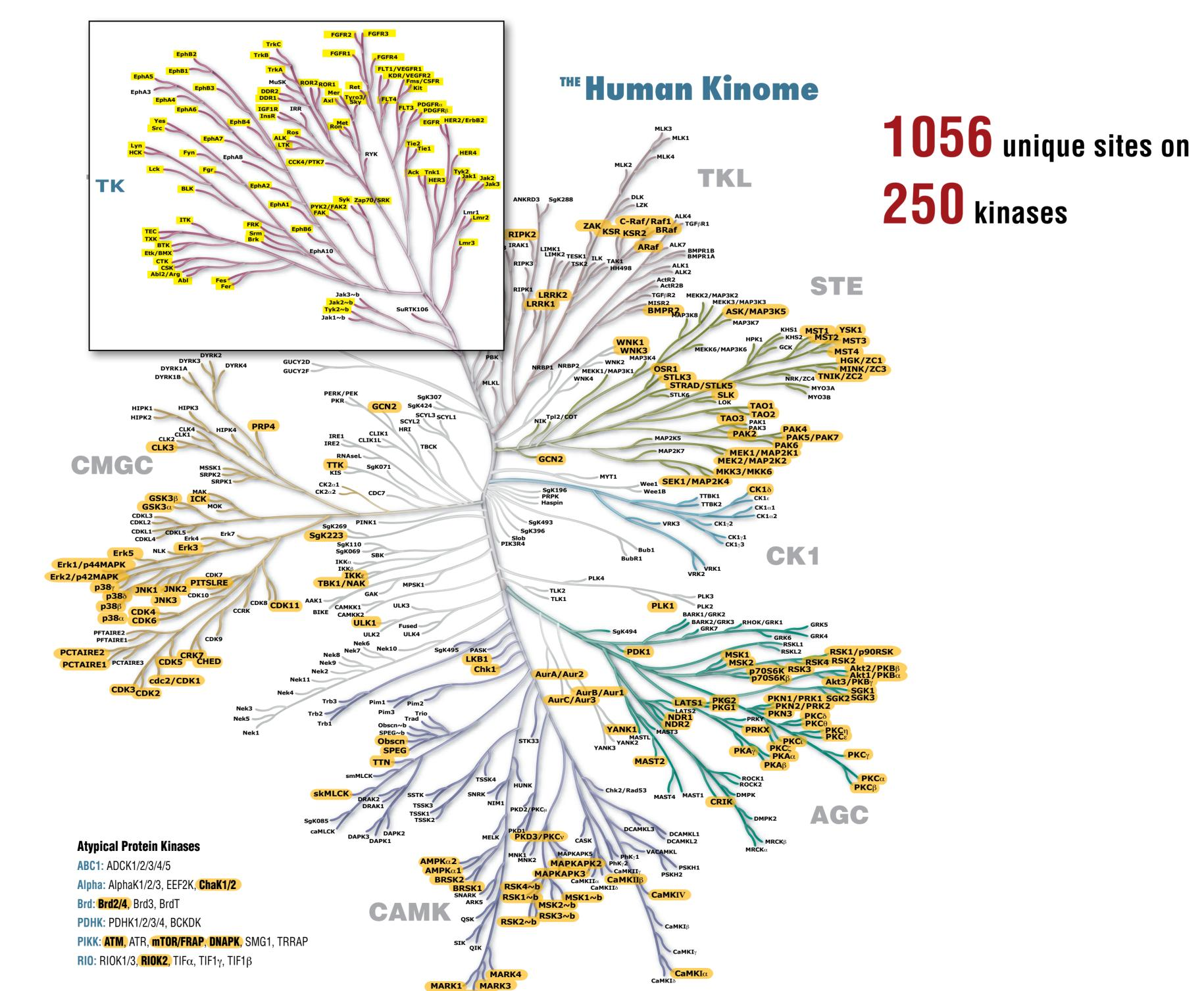


Figure 3: Ser/Thr Kinase and Tyr Kinase targets mapped onto the human kinome tree. Yellow highlighting indicates kinases for which peptides are identified using the reagents. Tyr Kinase coverage is shown in the inset ("TK").

: Akt/PI3K Pathway

In-depth analysis of PI3K/Akt and associated signaling pathways.

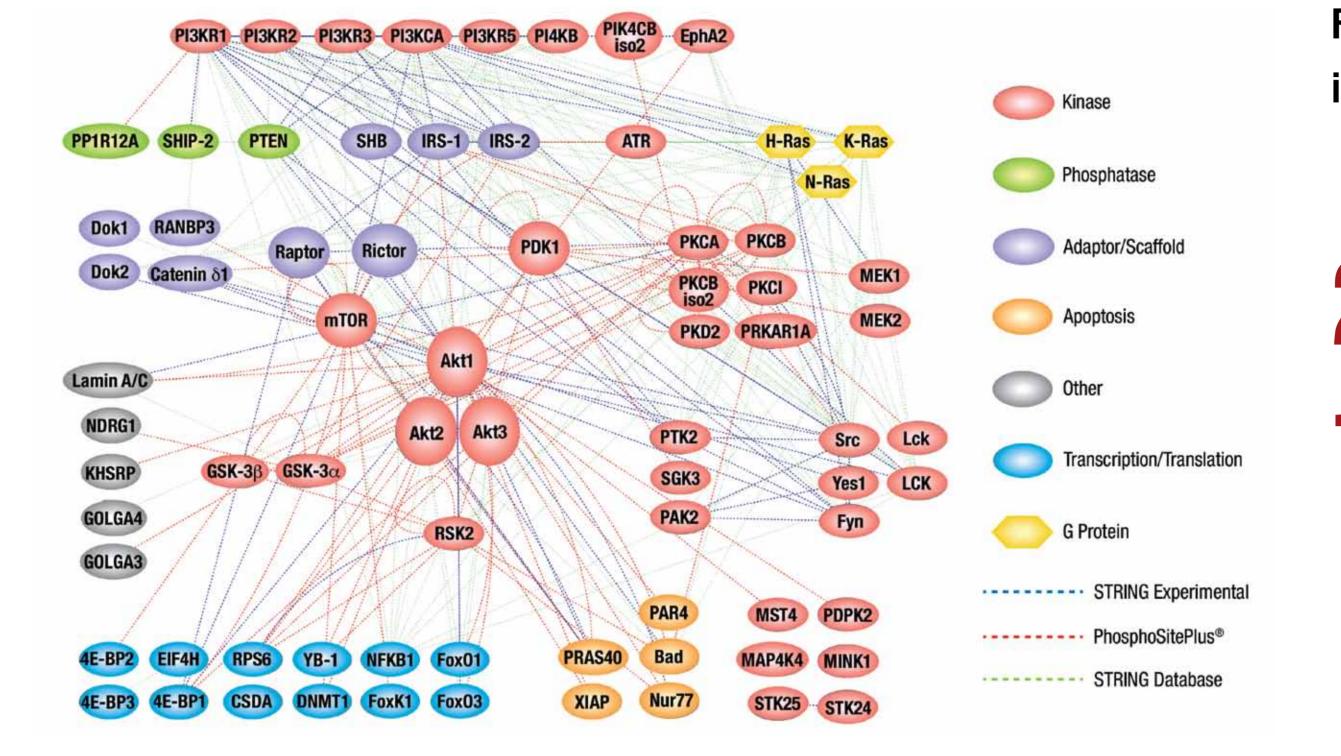


Figure 4: Akt/PI3K Pathway - interaction map of validated targets.

296 unique sites on105 proteins

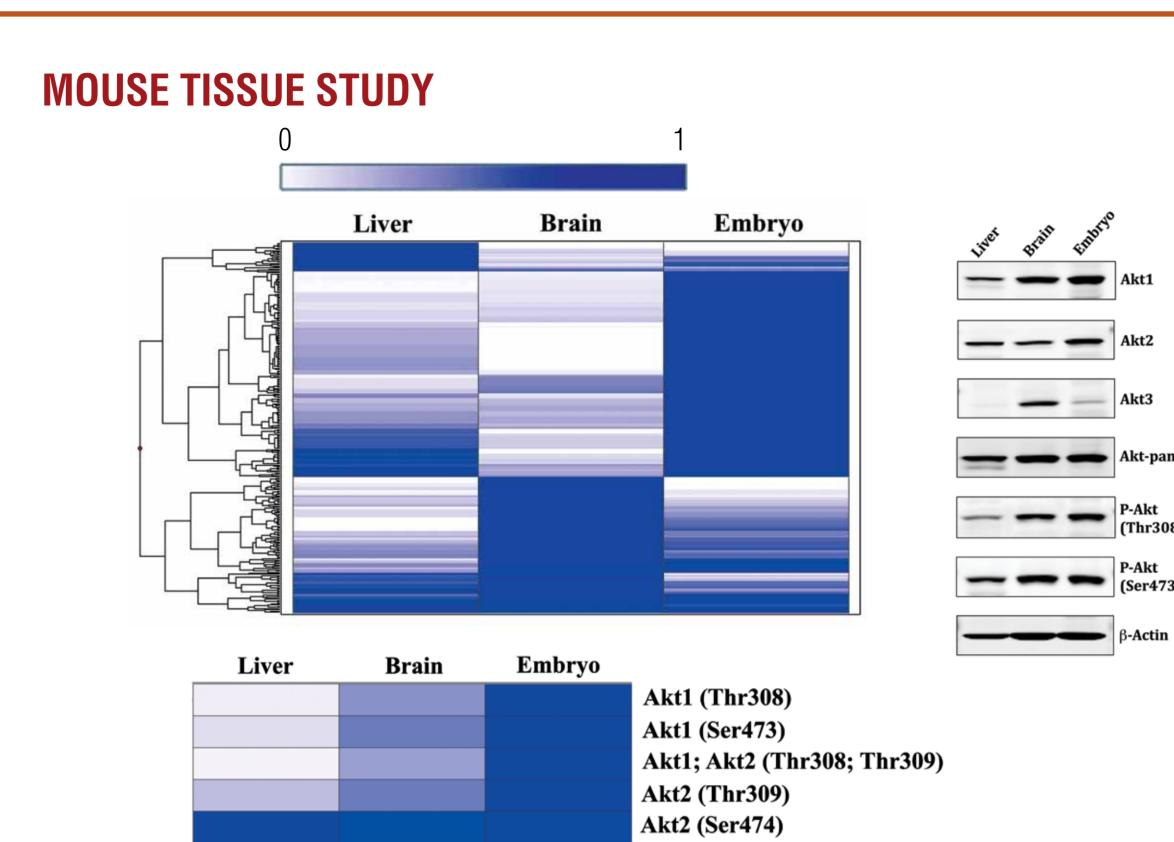


Figure 5: Akt/PI3K Pathway mouse tissue profiling. Hierarchical clustering of relative intensities for mouse liver, brain, and embryo. Each row represents a different validated Akt/PI3K pathway peptide. The maximum intensity across the three tissues was set to 1 and the other two intensities normalized to the max. Blue indicates higher intensity. Selected kinase peptides are shown in detail with accompanying western blots.

DNA Damage/Cell Cycle

A comprehensive view of the DNA damage response.

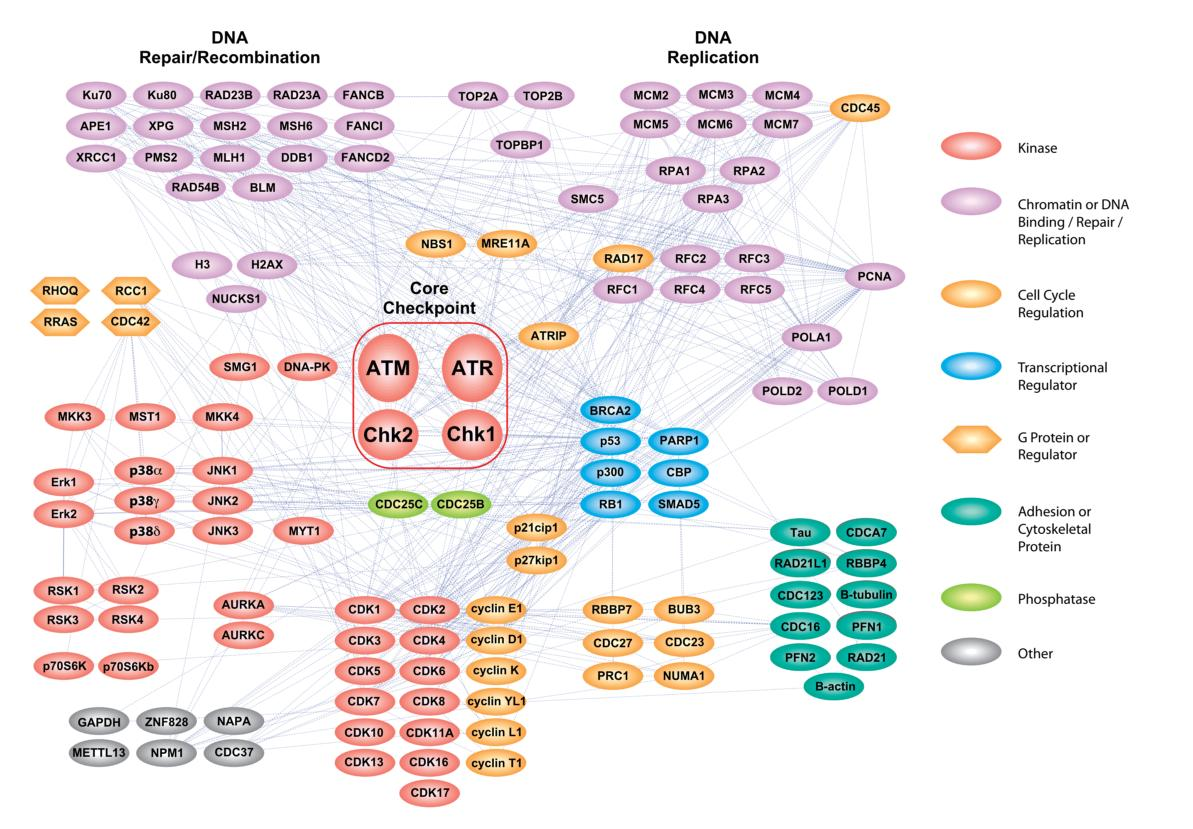


Figure 6: DNA Damage/Cell Cycle - interaction map of validated targets.

263 unique sites on137 proteins

-1.2 ATR §428 M#ESM#EIIEEIQCQTQQENLSSNSI 43,724

-1.3 Chk1 - FFHQLM#AGVVYLHGIGITHR 48,715

3.2 Chk1 §317 YSSS*QPEPR 698,572

4.1 Chk1 §345 LVQGISFS*QPTCPDHM#LLNSQLLG 1,840,956

-1.1 Chk2 - FTTEEALRHPWLQDEDM#K 7,865,595

9.1 Chk2 §379 ILGETS*LM#R 1,010,005

260.2 JNK1; JNK3 §183, §185; §221, §223 TAGTSFM#MT*PY*VVTR 8,543,090

98.0 JNK2 §183, §185 TACTNFMM#T*PY*VVTR 15,427,858

4.7 p38-alpha §180, §182 HTDDEMT*GY*VATR 18,211,944

4.1 p38-gamma §183, §185 QADSEM#T*GY*VVTR 1,421,705

Figure 7: DNA Damage/Cell Cycle profiling of response to

UV damage. Hela cells were untreated or treated with 500 mJ/cm²

UV light and harvested at 2 hr post treatment. Fold changes were calculated from chromatographic peak heights/areas. Green cells

ndicate peptides more abundant with UV damage; Red cells indicate

P-ATR (Ser428)

P-ATR (Ser428)

peptides less abundant. Selected DNA damage response proteins are shown in detail with accompanying western blots.

hoptosis/Autophagy

onitoring activity of apoptotic proteins and autophagolytic pathway

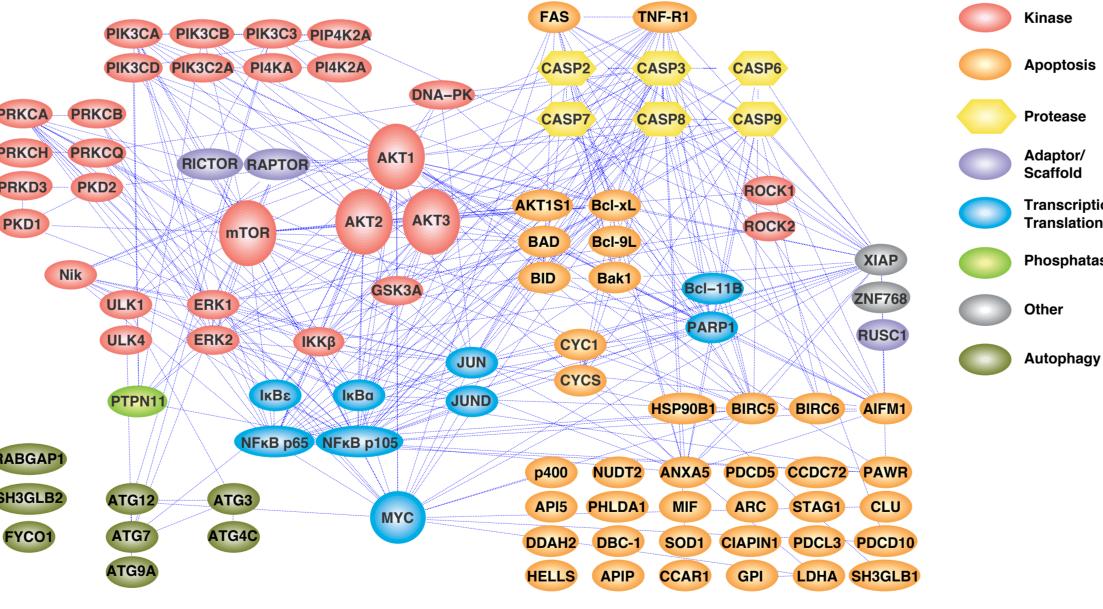


Figure 8: Apoptosis/Autophagy - interaction map of validated targets.

175 unique sites on100 proteins

HeLa -/+ UV DAMAGE

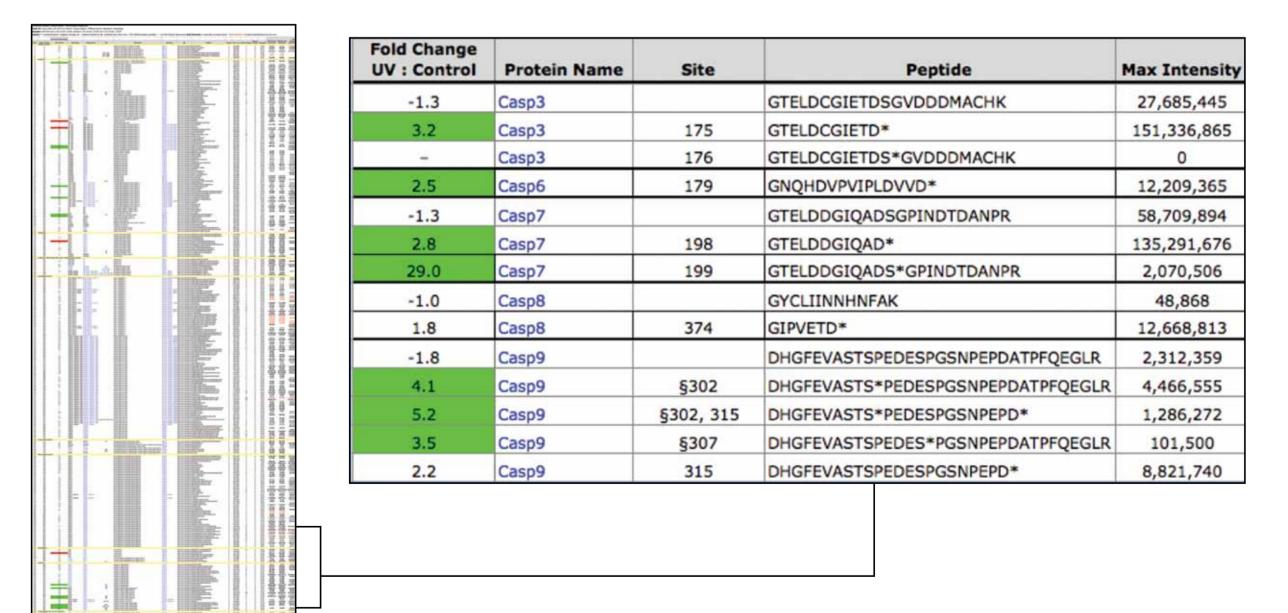


Figure 9: Apoptosis/Autophagy profiling of response to UV damage. Hela cells were untreated or treated with 500 mJ/cm² UV light a harvested at 2 hr post treatment. Fold changes were calculated from chromatographic peak heights/areas. Green cells indicate peptides morabundant with UV damage; Red cells indicate peptides less abundant.

Conclusion

PTMScan® Direct is a recently published method that allows identification and quantification of hundreds of peptides from selected protein types or signaling pathways. This approach allows focus on proteins of interest instead of 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

Stokes, M.P. et al. (2012) *Mol. Cell Proteomics* 11, 187–201.
 Rush, J. et al. (2005) *Nat. Biotechnol.* 23, 94–101.
 Lundgren, D.H. et al. (2009) *Curr. Protoc. Bioinformatics* 13, 13.3.1–13.3.21.
 Stokes, M.P. et al. (2012) *Int. J. Mol. Sci.* 14, 286–307.

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