# Mapping deregulated pathways downstream of RTKs in cell line models

# Abstract

In virtually all epithelial tumors, growth factor receptor activity is dereg- method of TMT labeling coupled with serial peptide immunoprecipitaulated by activated mutations, genomic amplification, and autocrine tion to profile drug-treated and untreated RTK pathway-addicted loops. The dependence of tumor cell survival upon the driving oncogene cancer cell lines. We also examined the ratio of protein levels in treated has been called "oncogene addiction" and demonstrates the acute cell lines. Quantitative analysis was performed to look at the effects of sensitivity of cancer cells to inhibition of the pathways driving their drugs on RTK-dependent cell lines: MKN45 (Met) with proliferation, growth and survival. Major questions regarding the sensi- Crizotinib, H1703 (PDGFRlpha) with Gleevec, and H3255 (EGFR) with tivity of solid tumors to targeted kinase inhibitors are why some tumors Iressa. Phosphotyrosine, acetyl-lysine, methyl arginine, ATM/ATR respond while others do not and why tumors which respond still are substrate (s/tQ), and AGC/CAMK/STE family kinase motif were all able to develop resistance through the activation of other pathways. selected for use in peptide immunoprecipitation which allowed us to Robust, but not durable, response to receptor tyrosine kinase inhibitors characterize and quantify a broad PTM space. In this study we were able (RTKIs) highlights the need to inhibit pathway activity at multiple levels. to identify and quantify hundreds of modification sites for phosphorylation,

In this study, we identify extensive signaling networks downstream sensitive, including all PTM types, were used for pathway analysis of of receptor tyroine kinases (RTKs) across multiple spaces, including signaling downstream of each RTK disease driver. phosphorylation, acetylation, and methylation. We used the established

arginine methylation, and lysine acetylation. Those sites found to be drug-

# Materials and Methods

Cells were cultured in DMEM medium or RPMI 1640 medium supple mented with 10% fetal bovine serum and processed in the appropriate amount of a 8 M urea lysis buffer. Cells were starved overnight and then treated with the following drugs for 1, 3, 6, and 24 hrs: MKN45 (Met) with Crizotinib, H1703 (PDGFR $\alpha$ ) with Gleevec<sup>®</sup>, and H3255 (EGFR) with Iressa<sup>®</sup>. Samples were then reduced and alkylated before being digested with trypsin. After digestion, samples were purified over Sep-Pak<sup>®</sup> C18 columns and eluted peptides were lyophilized. After peptide quantitation 2 mg of each sample was labeled using TMT reagents with 100 µg used for total proteome analysis and the rest used to profile PTM spaces. Six plex were organized the following way:

Labeled samples were then combined and subjected to sequential immunoprecipitations using motif antibodies for PTMs and in parallel basic reverse phase (bRP) chromatography was performed to profile the full proteome. Samples were purified over Sep-Pak<sup>®</sup> C18 columns before being processed by LC-MS/MS/MS analysis. Data generated from the MS analysis were processed through several modules in CORE, ending with quantitated data.

Channel

Sample	TMT Channel	Sample	TMT Channel	Sample	T
H3255 Control	126	H1703 Control	126	MKN 45 Control	12
H3255 Control	127	H1703 1 hr Gleevec	127	MKN 45 Control	12
H3255 1 hr Iressa	128	H1703 3 hr Gleevec	128	MKN 45 1 hr Criz.	12
H3255 3 hr Iressa	129	H1703 6 hr Gleevec	129	MKN 45 3 hr Criz.	12
H3255 6 hr Iressa	130	H1703 24 hr Gleevec	130	MKN 45 6 hr Criz.	13
H3255 24 hr Iressa	131	H1703 24 hr Gleevec	131	MKN 45 24 hr Criz.	13

Ab Target	Clone #	Motif
pY-1000	D1G10/D2D1	Phospho Tyr
AGC/PSD	D3E5/D8D9/D4E2/D8B11	AGC/CAMK/STE kinase acitivation loop
Akt substrate	100B7	RXX(s/t)
AMPK	D72H3/D78G9	LXRXXT(S)
ATM/ATR substrate	D23/D69	(s/t)Q
ATM/ATR substrate	D14/D86	(s/t)Q
Ac-K	D10G3/11D/16E	Acetyl lysine
K-Me	D4P3J, D3Z9J, D8R1C	K-Me
R-Me	Me-R4-100	R-Me

Figure 1. Motif antibodies chosen for PhosphoScan<sup>®</sup> profile (MS/MS).





Figure 2. Full proteome. Samples are harvested in 8M urea, homogenized if necessary, reduced, and alkylated before being digested. Samples were then fractionated over an ACCELA® HPLC using a 4.6X150 mm ZORBAX® 300 Extend-C18 (5 µm particle size with 300 A-pore). 96 fractions were collected over the 60-minute gradient. The fractions were then combined to give a total of 12 and these fractions were again purified over a C18 micro column. Each purified fraction was then subjected to MS/MS/MS analysis. Data processing was performed using CORE. (Protein Sieve, Protein Assembler, Protein Quant.)



Figure 3. Experimental flowchart. Cell lysates were prepared from RTK inhibitor treated and untreated cell lines, proteins were digested with trypsin, peptides were TMT-labeled, and samples were mixed for 6-plex analysis. The peptide mixtures were serially fractionated through immunoprecipitation with antibodies specific for phosphotyrosine, ATM/ATR substrate motif, AGC/CAMK/STE family kinase motif, acetyl-lysine, and methyl-arginine, though any motif-recognizing antibody could be used with this protocol. The enriched peptides from each immunoprecipitation were analyzed by LC-MS/MS using an LTQ<sup>®</sup> Orbitrap<sup>®</sup> Elite, with quantification of the TMT labels enabled by HCD fragmentation.





Figure 5. Number of sites affected by RTK inhibitors across spaces

#### Klarisa Rikova, Benjamin Hall, Anthony Possemato, Keri Mroszczyk, Kimberly A. Lee, Jian Min Ren, Ailan Guo, Daniel Mulhern, Yi Wang, Sean A. Beausoleil, Michael J. Comb

Figure 6. Pathway diagrams for each cell line were generated using Cytoscape®. These diagrams include all proteins with modification sites increasing or decreasing at least 2x upon relevant drug treatment. Proteins with PTM site peptides decreasing upon treatment (dark green protein names) are distinguished from those with sites increasing upon drug treatment (red protein names). Proteins identified in different PTM classes (phosphotyrosine, ATM/ATR substrate phosphorylation, AGC/CAMK/STE kinase family motif phosphorylation, lysine acetylation, and arginine methylation) are distinguished by different colored protein symbols. Shapes of protein symbols signify protein functionality (cytoskeletal, adaptor, kinase, adhesion protein, etc.). First order interacting partners with the driving RTK are presented in bright yellow.

#### Summary

- Intelligent fractionation through sequential immunoprecipitation can be used to probe numerous PTM spaces, allowing in-depth pathway mapping of valuable samples
- Deep data sets generated through enrichment of different types of modified peptides allow elucidation of novel biological relationships

### Contact Information

Klarisa Rikova Cell Signaling Technology, Inc. Tel: (978) 867-2330 Email: krikova@cellsignal.com

#### www.cellsignal.com