

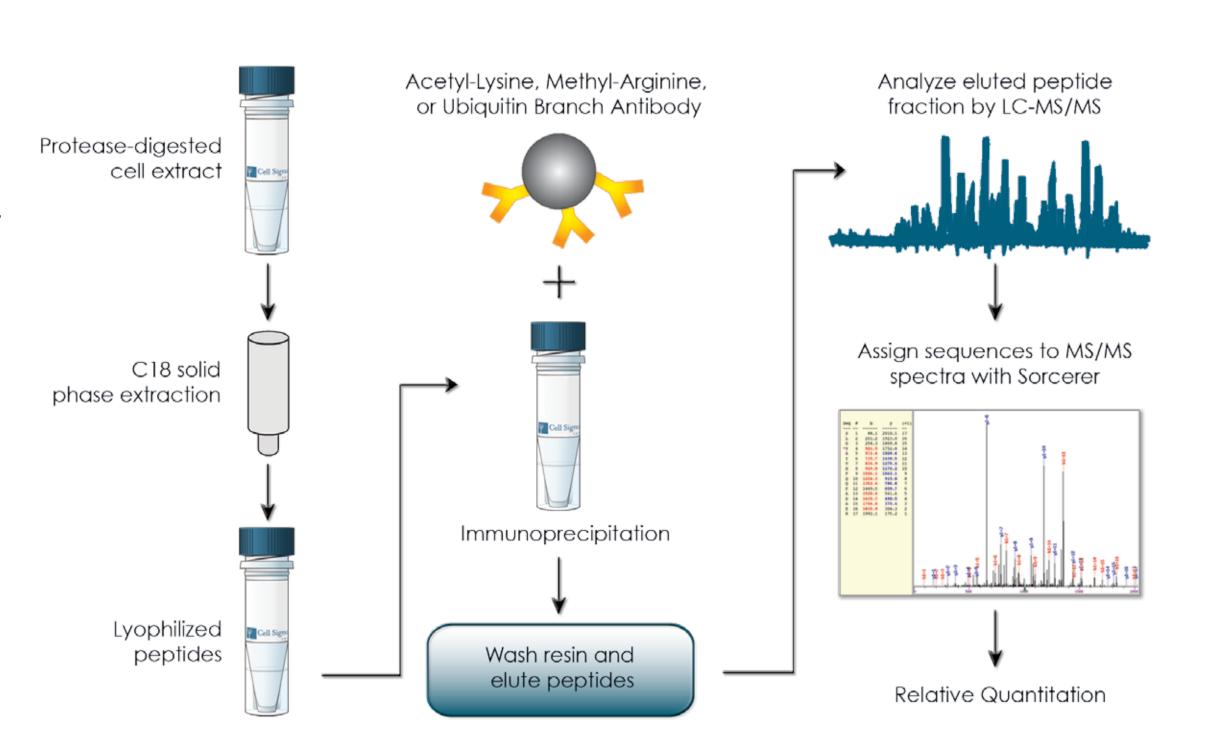
# Network and Pathway Analysis of Post-translationally Modified Peptides Identified by Immunoaffinity-based Proteomics

Jeffrey C. Silva<sup>1</sup>; Stuart Tugendreich<sup>2</sup>; Hongbo Gu<sup>1</sup>; Charles L. Farnsworth<sup>1</sup>; Kimberly A. Lee<sup>1</sup>; Xiaoying Jia<sup>1</sup>; Jian Min Ren<sup>1</sup>; Matthew P. Stokes<sup>1</sup> <sup>1</sup>Cell Signaling Technology, Danvers, MA USA 01923; <sup>2</sup>Ingenuity Systems, Redwood City, CA 94063

#### Introduction

Post-translational modification of proteins provides critical regulation of protein activity, stability, and localization in nearly all aspects of cellular signaling. We have developed antibodies that recognize specific post-translational modifications (phosphorylation, ubiquitination, acetylation, methylation), consensus kinase substrate motifs, or peptides from proteins that reside in the same signaling pathway or pathways. The antibodies are used to immunoaffinity purify post-translationally modified peptides from cell lines, tissues, xenografts, or other biological materials which are then identified using LC-MS/MS, enabling identification of thousands of post-translationally modified peptides in a single LC-MS/MS run. Ingenuity® Systems pathway analysis software IPA® has been applied to these large datasets in order to facilitate interpretation of the LC-MS/MS data and provide a biological context to the results.

PTMScan® analysis (Rush et al. 2005, Lee et al. 2011, Kim et al. 2011) was performed on HCT-116 human colon carcinoma and mouse liver tissue peptides with the Acetyl-Lysine, Mono-Methyl-Arginine, or Ubiquitin Branch Antibodies from Cell Signaling Technology. Samples were run on LTQ-Orbitrap VELOS<sup>™</sup> or Elite LC-MS/MS instruments and searches were performed using Sorcerer<sup>™</sup> software (Lundgren et al. 2009). Results from each study were combined and imported into the Ingenuity® Systems IPA® software package. Core analyses were confidence predicted relationships.



motif antibodies for enrichment of acetylated (AcetylScan®), methylated (MethylScan<sup>™</sup>), and ubiquitinated (UbiScan<sup>®</sup>) peptides.

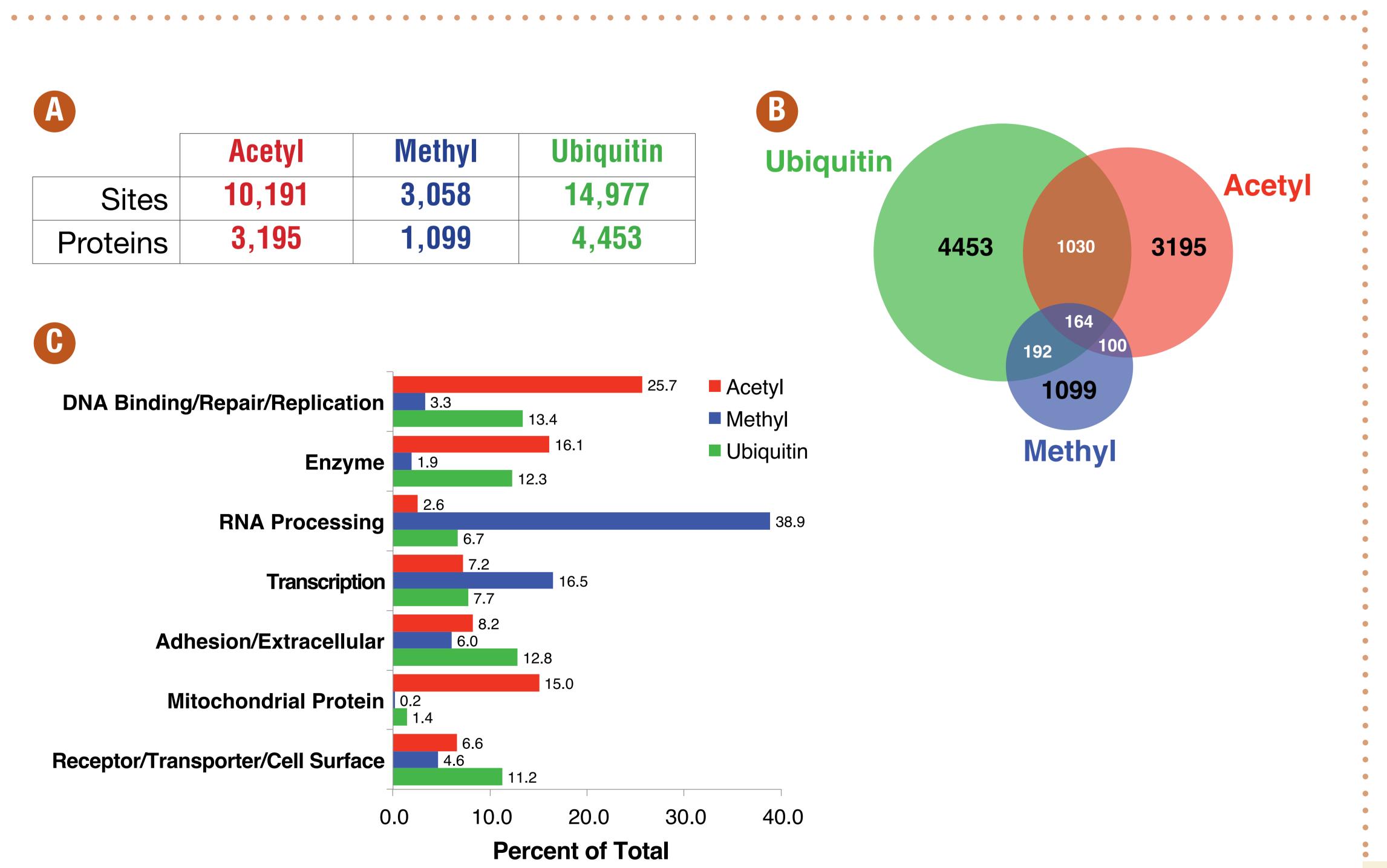
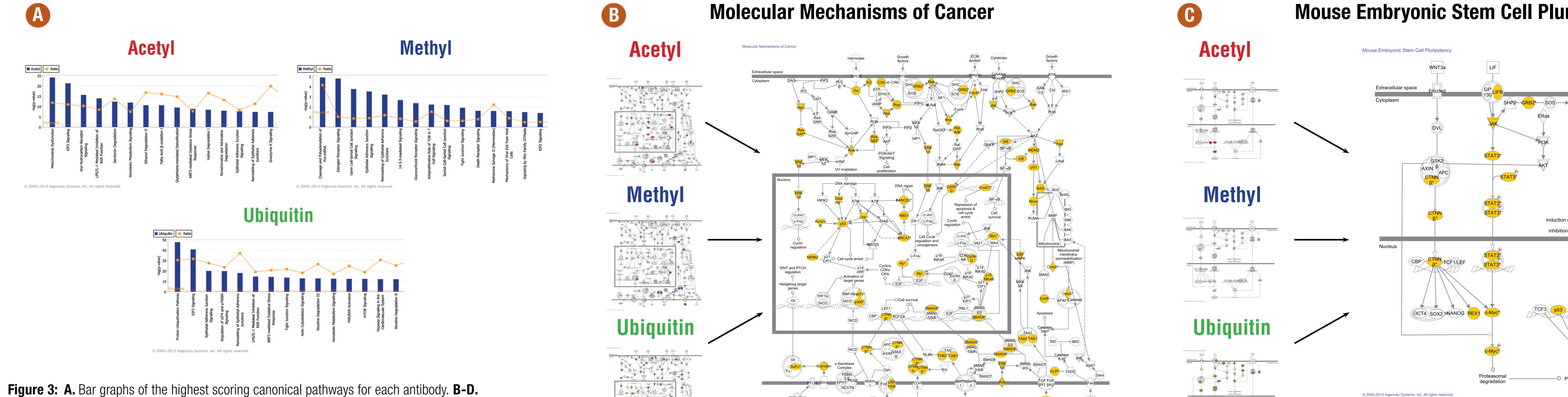
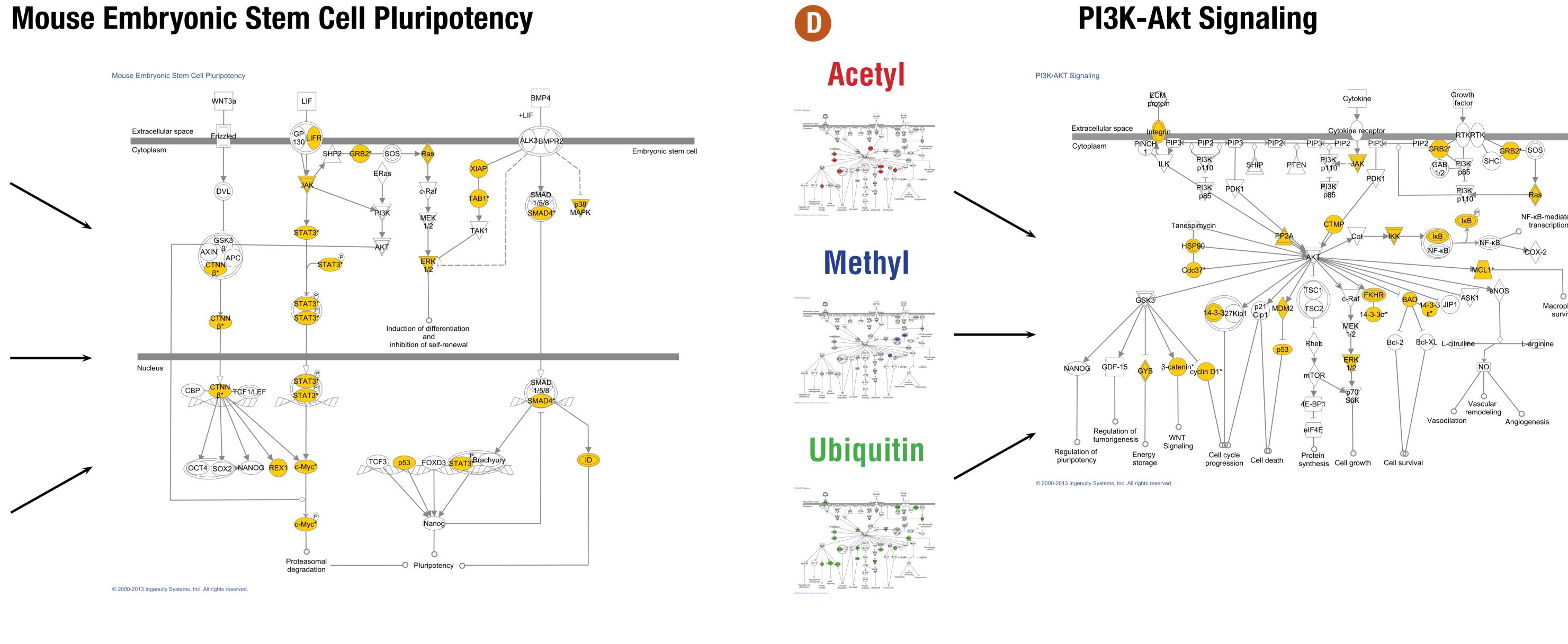


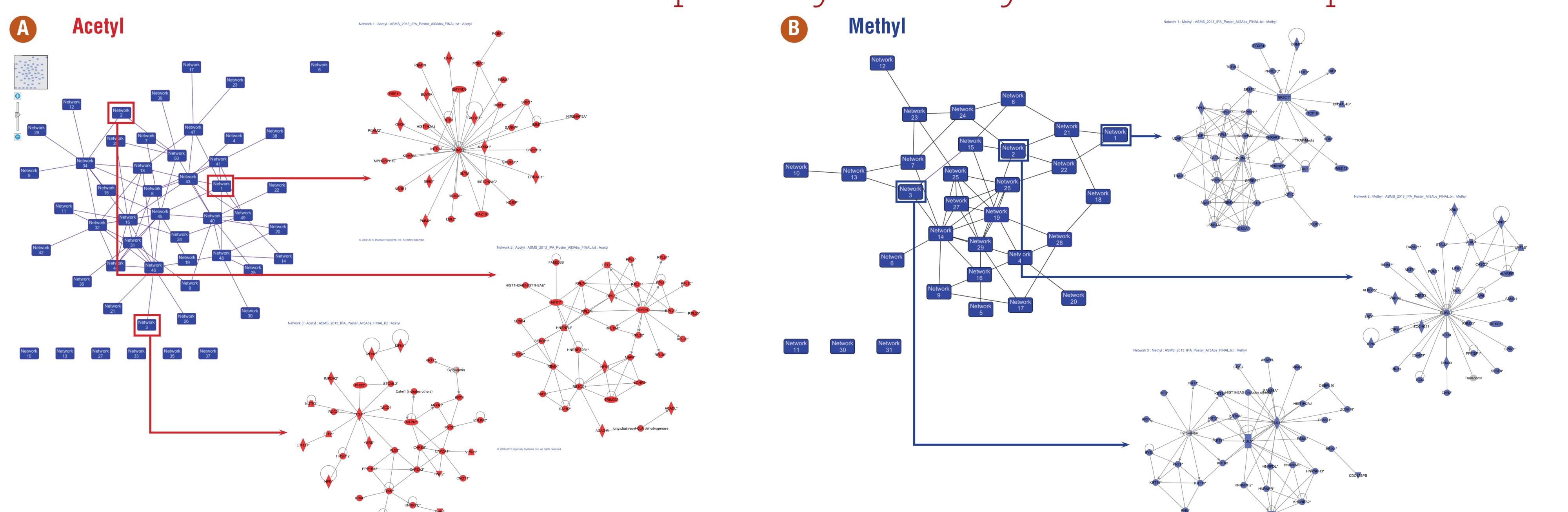
Figure 2: PTMScan® results using Acetyl-lysine, Mono-methyl-Arginine, and Ubiquitin Branch motif antibodies. A. Number of nonredundant sites/proteins identified for each antibody. B. Area proportional Venn diagram showing the number of proteins identified using each antibody (black numbers) and overlap between antibodies (white numbers). C. Top seven most highly represented protein classes identified with each antibody. Protein type information from the PhosphoSitePlus® database.

# IPA® Canonical Pathways: Mapping identified proteins onto pre-defined signaling networks





## IPA®-Generated Networks: De novo pathway assembly from identified proteins



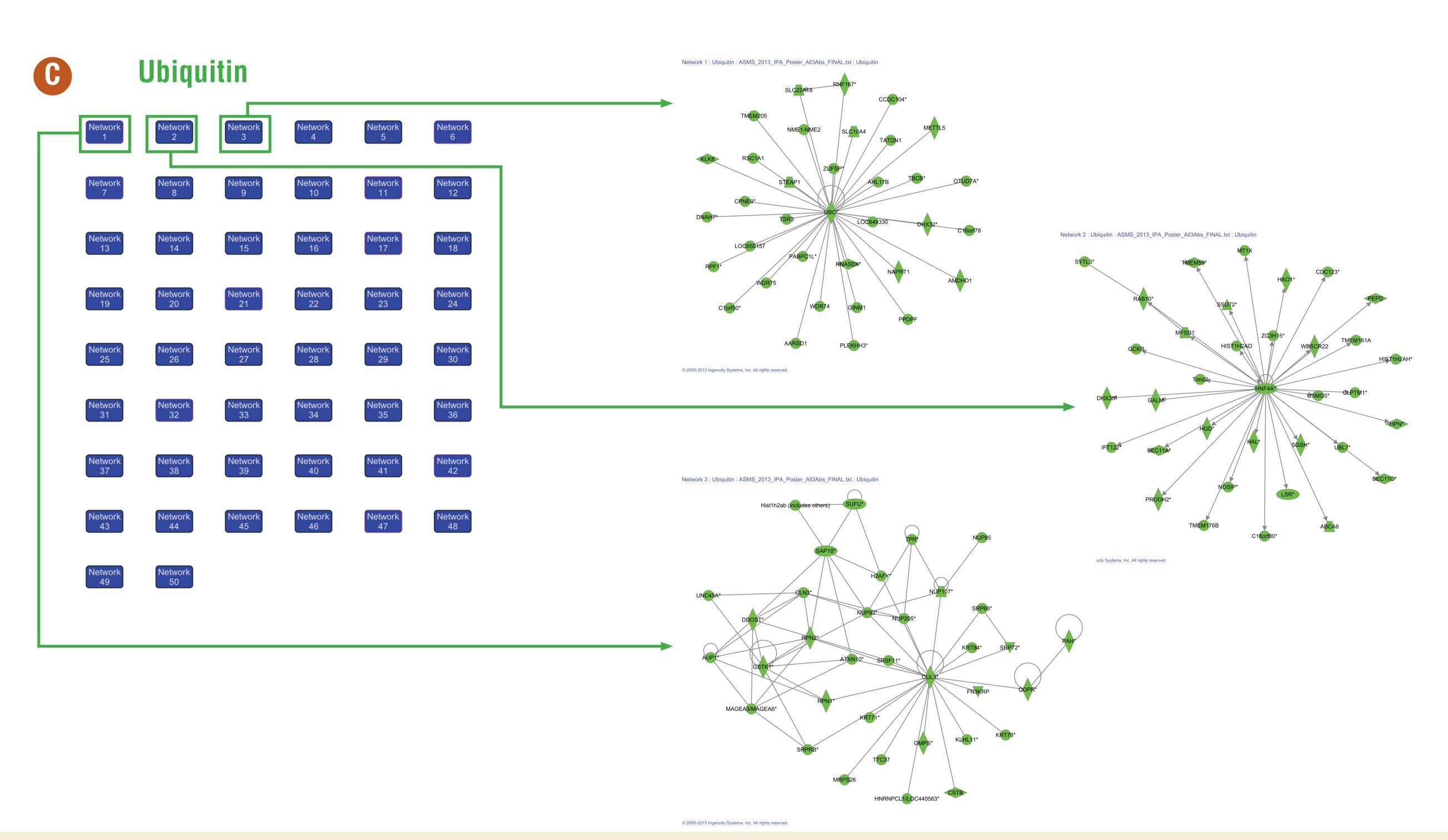


Figure 4: The top three IPA®-generated networks are shown for each antibody (A = Acetyl, B = Methyl, C = Ubiquitin). The overlap of the networks created for each antibody are shown with the selected networks displayed in detail.

#### Conclusion

Selected canonical IPA<sup>®</sup> Signaling Pathways are color-coded for each antibody (Acetyl = Red,

Methyl = **Blue**, Ubiquitin = **Green**). A composite pathway is shown for each (Yellow highlights).

Ingenuity® Systems IPA® software provides in-depth analysis of the complimentary biological networks assembled using the Acetyl-Lysine, Mono-Methyl-Arginine, and Ubiquitin-Branch motif antibodies in PTMScan®.

### References

**1.** Rush, J. et. al. (2005) *Nat. Biotechnol.* 23, 94–101. 2. Lee, K. A. et. al. (2011) *J. Biol. Chem.* 286(48):41530–41538.

**3.** Kim, W. et. al. (2011) *Mol. Cell.* 2011 Oct 21;44(2):325–340. 4. Lundgren, D. H. et. al. (2009) Curr. Protoc. Bioinformatics Chapter 13, Unit 13 13.