

Efficient Enrichment of Endogenous SUMOylated Peptides from Alpha-lytic Protease Digest Using K-ε-GG Remnant Immuno-affinity Purification

Hongbo Gu¹, Xiaoying Jia¹, Jianmin Ren¹, Elizabeth Komives², Matthew Stokes¹
Cell Signaling Technology, Inc., Danvers MA USA 01923
Department of Chemistry & Biochemistry, UCSD, La Jolla, CA USA 92033

INTRODUCTION

The family of Small Ubiquitin-like Modifier (SUMO) proteins in humans includes four distinct genes: SUMO1, SUMO2/3 and SUMO4. SUMOylation regulates the function of various proteins by reversible, covalent, isopeptide-bond attachment between the C-terminus of SUMO and a free ε-amine group of the target protein's lysine. The function of SUMO conjugation is less well-understood, but SUMOylated proteins are involved in gene expression, DNA repair, nuclear import, heat shock, cell motility, and lipid metabolism. SUMO targets are generally low-abundance proteins, and the amount of the modification at steady-state is also low. In recent years several groups have developed methods for analysis of SUMOylated proteins. However, many of these involve overexpression of mutant SUMO sequences. Therefore, methods that allow proteome-level identification of endogenous SUMOylation sites are needed.

We recently described the novel application of wild-type α-lytic protease (WaLP) to proteome digestion for shotgun proteomics. Although relatively relaxed specificity was observed, WaLP prefers to cleave after threonine residues and never cleaves after arginine (Fig. 1b). We show here that WaLP cleaves at the C-terminal TGG sequence (in SUMO 2/3 and 4) leaving a SUMO-remnant KGG at the position of SUMO attachment. The resulting KGG-containing peptides can then be identified using methods already developed for Ub-profiling (Fig 1d). The method allows identification of SUMO attachment sites under completely native conditions using the Ub-profiling workflow by simply substituting WaLP for trypsin. The same sample can be subjected to analysis of both Ub-attachment and SUMO-attachment by digesting the sample with either trypsin or WaLP respectively.

METHODS

Hela cells were heat shocked and lysed in urea buffer. Protein digestion was performed in 2M urea by alpha-lytic protease (WaLP) that preferentially cut at the C-terminus of T, S, A and V resulting a GlyGly remnant left on the lysine residue previously carrying SUMO-1 or SUMO-2/3. Digested peptides were subject to immunoaffinity purification (IAP) using the K-ε-GG remnant antibody and LC-MS/MS analyzed for identification and quantification of SUMOylation sites. To validate the KGG sites identified by the method, specific sumo-proteases were used to remove endogenous SUMO tails from modified proteins, which were then digested by WaLP and trypsin, respectively. Label-free quantification was applied to determine the quantitative changes of identified KGG peptides from WaLP and trypsin digest in parallel.

CONCLUSIONS

There were about 1,100 non-redundant KGG peptides identified from 5mg Hela lysate digested with WaLP and enriched with KGG antibody IAP including known SUMOylated proteins including TRIM28, TRIM33, SUMO1 and SUMO-2/3. By using specific SUMO-proteases (SENP 1 and 2), both SUMO-1 and SUMO-2/3 were efficiently removed *in-vitro*, as confirmed by western blots; while the total ubiquitination level was not affected by SUMO-proteases. We compared the quantitative changes of identified KGG peptides from WaLP and trypsin digest with and without SUMO-protease treatment. 612 out of 724 WaLP-digested KGG peptides (SUMOylated sites) decreased in abundance, while only 16 out of 9,031 trypsin digested peptides (Ubiquitinated sites) decreased with SUMO protease treatment. The results validated that the KGG sites identified by our method truly come from SUMOylation instead of ubiquitination.

Protein function classification placed proteins involved in 1) transcriptional regulation; 2) translation, 3) RNA processing, and 4) DNA binding and repair as top protein groups modified by SUMOylation that indicated a potential correlation between SUMOylation and epigenetics. More importantly, more than 2/3 of the KGG sites identified by our method are novel SUMOylation sites, providing a rich resource for the field of SUMOylation.

REFERENCES

Meyer, J. G. et al. Expanding proteome coverage with orthogonal-specificity α-lytic proteases. Mol Cell Proteomics 13, 823-835, doi:10.1074/mcp.M113.034710 (2014).

Udeshi, N. D., Mertins, P., Svinkina, T. & Carr, S. A. Large-scale identification of ubiquitination sites by mass spectrometry. Nat Protoc 8, 1950-1960 (2013).

Becker, J. et al. Detecting endogenous SUMO targets in mammalian cells and tissues. Nat Struct Mol Biol, 525-531, doi:10.1038/nsmb.2526 (2013).

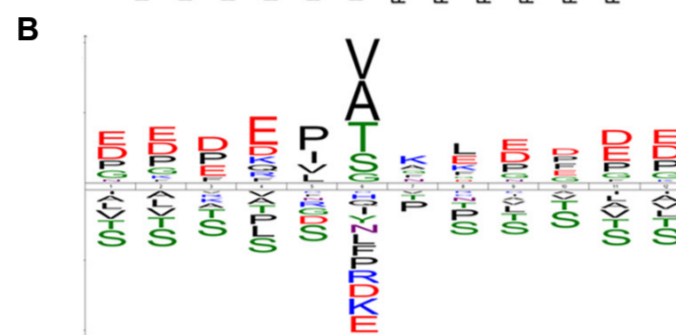
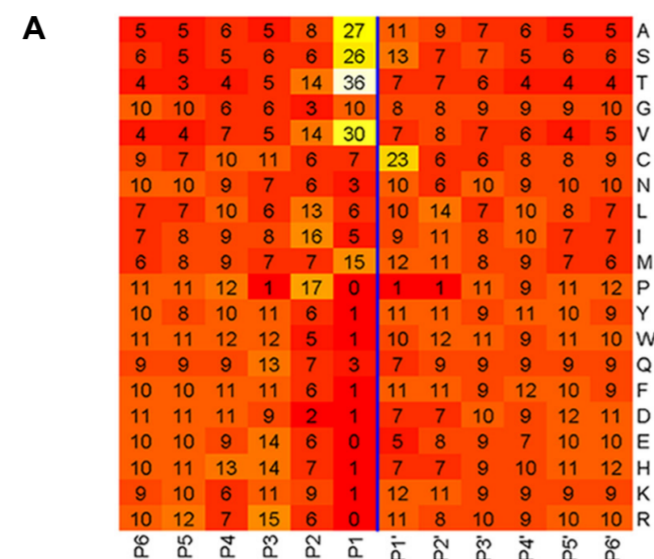


Figure 1: Heat map summarizing the analysis of the occurrence of each amino acid at each position in the identified peptide from the WaLP digest in reference 1. (A), WaLP peptides, counts normalized for the occurrence of each amino acid at each position. (B), iceLogo depicting the enrichment and depletion of specific amino acids at each position in the WaLP peptides, with residues colored according to property (acidic, red; basic, blue; hydrophobic, black; small/neutral, green). WaLP yields peptides with the following P1 (C-terminal) residues: A, 20%; V, 20%; S, 16%; T, 16%; G, 8%; L, 6%. The cleavage site is marked by a vertical line in the heat map

A

sp P0CG48 UBC_HUMAN	QKESTLHLVLRRLRGG
sp P63165 SUMO1_HUMAN	EEEDVIEVYQEQTGG
sp P61956 SUMO2_HUMAN	EEDTIDVFQQQTGG
sp P55854 SUMO3_HUMAN	EEDTIDVFQQQTGG
sp Q6EEV6 SUMO4_HUMAN	EEDTIDVFQQPTGG

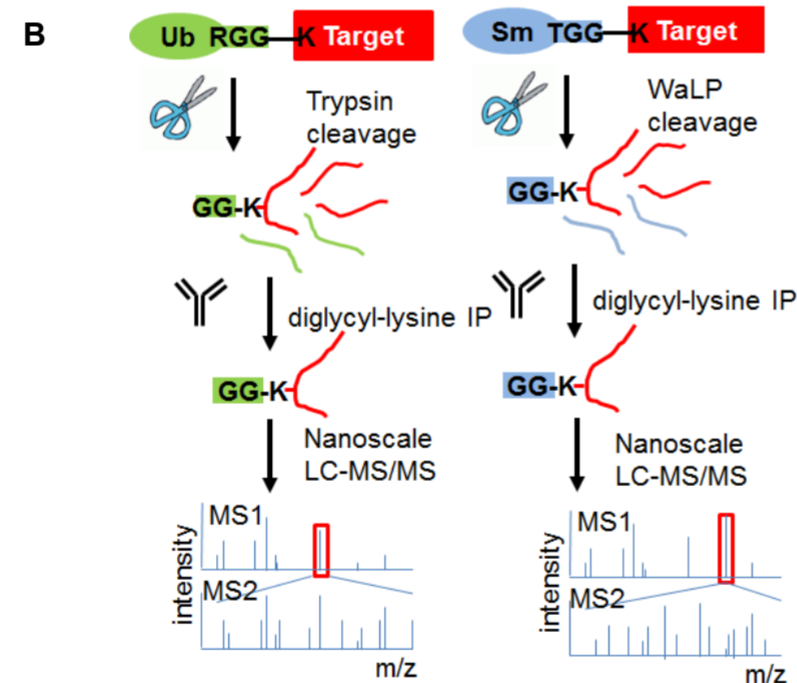


Figure 2: A Strategy for mapping endogenous SUMO1/2/3 attachment sites. (A), C-terminal sequence alignment of mature human SUMO 1-4 and Ubiquitin. (B), Cartoon showing ubiquitin-remnant and SUMO-remnant mapping strategy. Proteins modified by ubiquitin digested with trypsin and modified by SUMO digested with WaLP leave a diglycyl attached to the ε-amine of the lysine where ubiquitin or SUMO was attached. An antibody specific for the K-GG remnant is used to enrich peptides from ubiquitinated sites and SUMOylated sites that are then identified by nano LC-MS/MS.

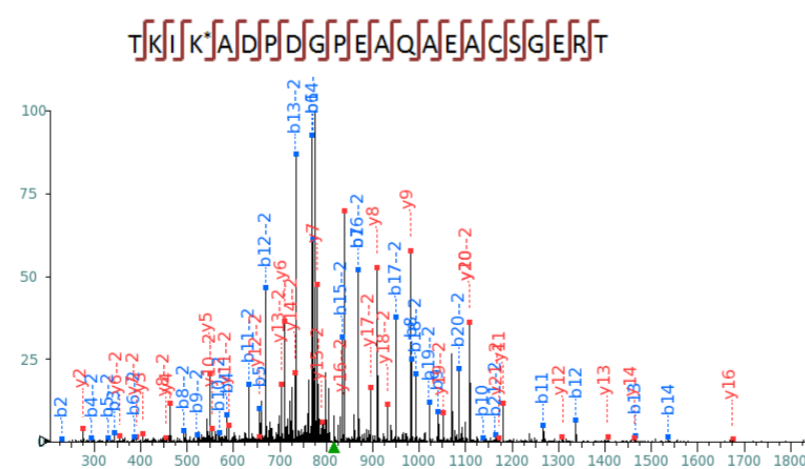
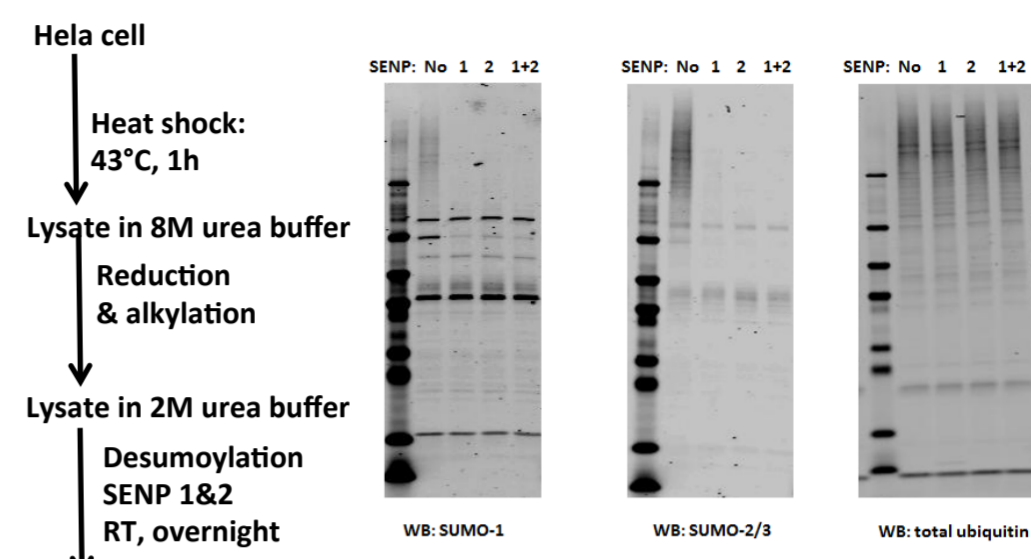


Figure 3: A known SUMOylation site (K5 of Human NHP2, Q9NX24) was confidently identified from WaLP digest of Hela cells by K-GG antibody enrichment and LC-MS/MS. Green triangle indicates the m/z of the precursor peptide.



Western blot for SUMO1, SUMO 2/3 and Ubiquitin

Figure 4: A strategy to validate identified SUMOylation sites. Endogenous SUMO attachment was removed by SUMO proteases SENP 1 and SENP 2. Endogenous ubiquitin levels were not affected by SENP 1 and SENP 2 treatment.

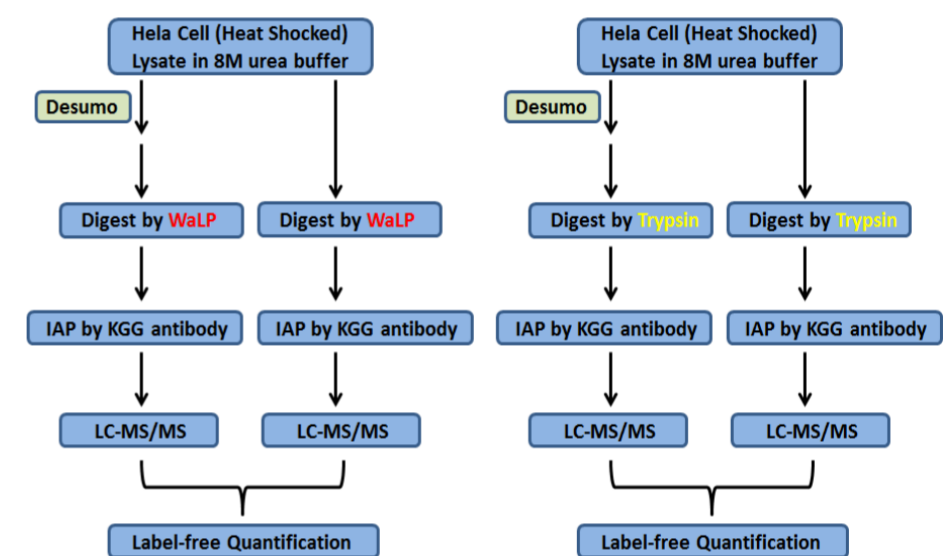


Figure 5: A parallel quantitative analysis of K-GG peptides enriched from WaLP and Trypsin digest of Hela cells. Label-free quantification was applied to quantify peptides from cells treated with SENP 1+2 vs. untreated.

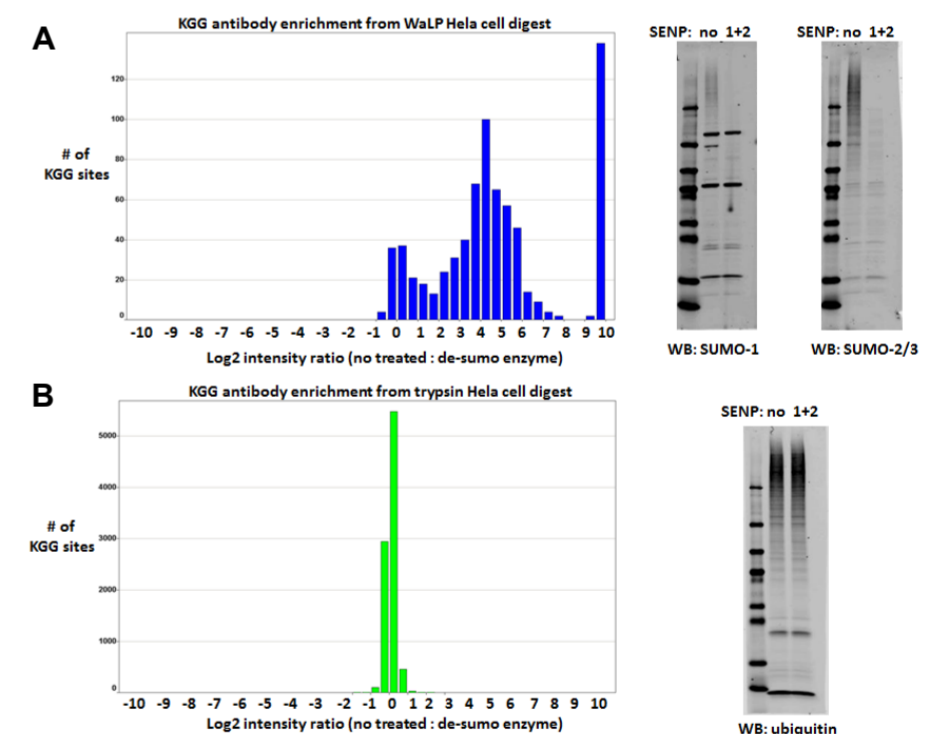


Figure 6: (A), Bar graphs of results from quantification of K-GG peptides after treatment with SENP 1+2 showing the dramatic reduction of K-GG-sites observed after WaLP digestion. Importantly, for over 400 sites, SENP 1+2 treatment resulted in no detectable K-GG modification upon WaLP digestion. These sites were assigned as a log₂ value of 10. (B), No significant reduction in intensity of K-GG sites was observed after tryptic digestion. Both quantitative results are consistent with western blots.

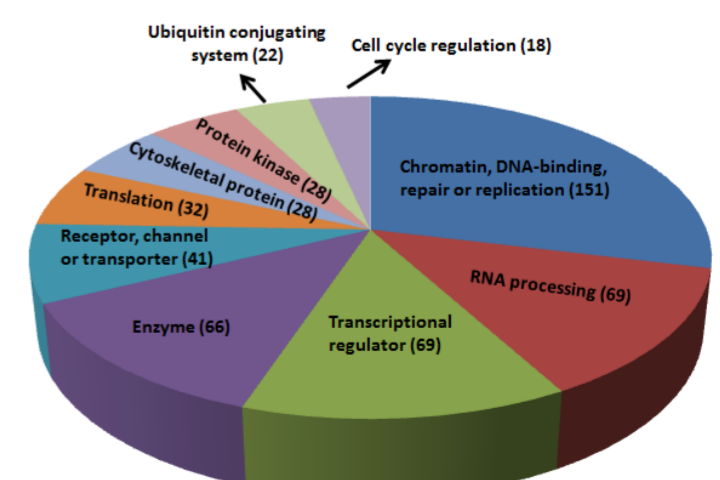


Figure 7: Classification of identified SUMOylated proteins.