Applications of Mass Spectrometry Targeted Assays for Quantitative Analysis of Cancer Signaling Proteins

Penny Jensen¹; Bhavin Patel¹; Leigh Foster¹; Aaron Gajadhar²; Sebastien Gallien³; Jonathan R. Krieger⁴; Jiefei Tong⁵; Michael F. Moran^{4,5,6}; Rosa Viner²; Andreas Huhmer²; Kay Opperman¹; John Rogers¹ ¹Thermo Fisher Scientific, Rockford, IL; ²Thermo Fisher Scientific, San Jose, CA; Thermo Fisher Scientific, PMSC, Cambridge, MA; ⁴SPARC Biocentre, The Hospital for Sick Children, Toronto, Canada; ⁵Program in Cell Biology, The Hospital for Sick Children, Toronto, Canada; ⁶Department of Molecular Genetics, University of Toronto, Canada

ABSTRACT

Purpose: The AKT/mTOR pathway plays a central role in tumor progression and drug resistance. Quantitative measurement of alterations in the expression of pathway proteins and post-translational modifications (PTM) is necessary for understanding cancer biology. Highly accurate monitoring of these pathway proteins has not been achieved, due to poor reproducibility, unreliable quantitation, and lack of standardized methods and reagents. To overcome these challenges, the novel Thermo Scientific[™] SureQuant[™] pathway panels have been applied, which utilize an optimized multiplex immunoprecipitation to targeted mass spectrometry (mIP-tMS) workflow. SureQuant assays can quantitate multiple proteins, PTMs and interacting partners, which creates new possibilities for a broad range of applications, including cancer, drug development, and research into precision medicine.

Methods: The SureQuant total and phospho pathway panels contain two modules: 1) The IP-MS Sample Prep Module includes reagents necessary to immunoenrich AKT pathway, RAS, or TP53 proteins, and perform MS sample preparation in one day 2) The Absolute or Relative Quantitation Modules include a Thermo Scientific™ Pierce™ LC-MS/MS System Suitability Standard, AQUA Ultimate Heavy and/or AQUA Ultimate Light Peptides, and verified MS instrument and data analysis methods. Serum-starved, inhibitor-treated (LY294002/NVP-BEZ235/Rapamycin) HCT116, A549, and MCF7 cells were stimulated with hIGF-1. SureQuant AKT pathway panels (total and phospho) were used to determine the absolute concentration of target peptides using targeted MS analysis. The panels were benchmarked against Western blotting using three unstimulated, hIGF-1 stimulated or inhibited cell lysates, as well as several tissue/xenograft lysates.

Results: Previously, we verified antibodies and target peptides to AKT and RAS pathways using an optimized mIP-tMS workflow. From the standard curve, all target peptides were monitored with <20% CV, 3 orders of magnitude dynamic range, linearity (R²) >0.97, and accuracy of 80-120% in a complex matrix. Using the SureQuant pathway panels, absolute quantitation of 37 target peptides in unknown samples was achieved with <20% CV across multiple cancer cell lines. The SureQuant pathway analysis workflow allowed absolute quantitation of target peptides from positive control lysate with <15% individual operator %CV and <20% combined %CV using PRM analysis. Kit performance was evaluated through analysis of abundance levels between three different cancer cell lines, A549, HCT116, and MCF7, using the SureQuant AKT Total and Phospho assay showed preferences for certain inhibitors in specific cell lines treated with hIGF-1. The PI3K inhibitor LY294002 functioned the best in HCT116 cells whereas the dual PI3K/Rapamycin inhibitor NVP-BEZ235 worked predominantly in A549 cells. Analysis by mass spectrometry allowed for more accurate and informative data with the determination of fmol levels of protein expression and capability to discriminate between isoforms of many proteins that are unable to procure with western blot analysis. Absolute quantitation of 12 phosphorylated AKT pathway targets was obtained from five patient derived lung tumor xenograft samples. Additionally, all 12 total and 12 phospho AKT pathway targets were quantitated from three different tissue lysates, Lung, Large Intestine, and Breast tumor.

INTRODUCTION

Multiplex Immunoprecipitation to Mass spectrometry (IP-MS) kits from Thermo Fisher Scientific are developed for simultaneous enrichment and quantitation of total abundance and phosphorylation levels of multiple proteins from the AKT/mTOR Signaling Pathway. The immunoenriched, digested samples are spiked with heavy peptide internal standards, which can then be processed using discovery MS (DDA) and targeted MS (PRM) methods for analysis.

Figure 1. SureQuant Multiplex IP to Targeted MS Modules & Kits

P/MS sample prep module	Absolute quantitation module	AKT pathway: total targets	AKT pathway: phospho targets
10 reactions (samples) Biotinylated antibodies mix Positive control lysate Streptavidin magnetic beads Wash buffers (A and B) Trypsin Sample prep buffer Reduction/alkylation reagents 10% TFA Low-binding tubes	 10 reactions (samples) AQUA heavy peptide mix AQUA light peptide mix 6 protein digest matrix 7 x 5 system suitability standard Peptide Diluent Instrument method with Skyline doc Low-binding tubes Relative quantitation module	GSK3a GSK3b PRAS40 (AKT1S1) mTOR	AKT1/AKT2 pSer473/pSer474 PTEN pSer380 IRS1 pSer312 IGF1R pTyr1162/1163 GSK3a pSer21
	 10 reactions (samples) AQUA heavy peptide mix 7 x 5 system suitability standard Peptide Diluent Instrument method with Skyline doc Low-binding tubes 		GSK3b pSer9 PRAS40 pThr246 mTOR pSer2448 P70S6K pThr389 TSC2 pSer939

Figure 3. AKT Pathway

Figure 2. Experimental Workflow for Multiplex IP-MS Assays





charcoal stripped FBS for 24 hours before stimulation with 100 ng/mL of IGF for 15 minutes. Breast, Lung, and Large Intestine Tumor and Normal Adjacent Tissue were purchased from BioIVT. Cells and tissue samples were lysed with Phosphatase inhibitor cocktail (PN#78440).

Multiplex Immunoprecipitation to MS Sample Preparation and MS Quantitation: The SureQuant IP and MS Sample Preparation Modules for AKT Pathway (PN# A40081, A40086, A40091), were used to immunoenrich relevant protein targets. The SureQuant Absolute Quantitation Modules for AKT Pathway (PN# A40083, A40093) was used to generate calibration curves and determine concentrations of target peptides from unknown samples.

curves or unknown samples. IP-enriched and trypsin digested samples were then desalted on-line using the Thermo using the Thermo Scientific™ Dionex™ UltiMate™ 3000 RSLCnano System and Thermo Scientific™ Q Exactive™ HF Hybrid Quadrupole-Orbitrap[™] Mass Spectrometer. Verified instrument acquisition methods were used, as well as inclusion lists relevant to each Absolute Quantitation Module.

sequence coverage, unique peptides, areas/intensities of identified peptides, and PTMs. For targeted MS data analysis, Skyline software (University of Washington) were used to measure limit of quantitation (LOQ) from the calibration curve and target analyte concentration from unknown samples.

RESULTS

Figure 5. Pierce[™] LC-MS/MS System Suitability Standard (7 x 5 Mixture)



magnitude dynamic range and linearity $(R^2) \ge 0.9800$

Figure 6. Determination of Quantitation Limits and Linearity of AQUA Light Peptides for SureQuant Pathway Kits



All 30 AKT pathway target peptides were monitored with linear guantitation $(R^2 \ge 0.9800)$ and 2-3 orders of magnitude (LLOQ ≤ 0.5 fmol on column)



- and DMEM Media, respectively, with 10% FBS/1xPenStrep to ~70-80% confluency. Cells were serum starved with 0.1% Thermo Scientific[™] Pierce [™] IP Lysis buffer (PN#87788) supplemented with 1X Thermo Scientific[™] Halt Protease and
- Liquid Chromatography and Mass Spectrometry: Pierce LC-MS/MS System Suitability Standard (7 x 5 Mixture) (PN# A40010) was used to assess dynamic range and sensitivity (LLOQ) of the nanoLC-MS system prior to running calibration Scientific[™] Acclaim[™] PepMap[™] 100 C18 Trap Column (PN#164564) followed by seperation using a Thermo Scientific[™] EASY-Spray[™] C18 column (PN#ES800). For discovery MS and targeted PRM-MS analysis, the samples were analyzed
- MS Data Analysis: Discovery MS data were analyzed with Thermo Scientific[™] Proteome Discoverer[™] to assess percent

AKT SureQuant Pathway Mass Spec Assay kits allowed absolute quantitation of target peptides from positive control lysate with <15% individual CV and <20% combined CV using PRM analysis.

Figure 8. Quantitation of Total and Phospho AKT Pathway Proteins using Different Positive Control Lysate Amounts



Total and phospho AKT pathway PRM analysis allowed absolute quantitation of all target peptides from 50 µg to 1000 µg of positive control lysate. Overall linear correlation between lysate amount and target quantitation was observed.

Figure 9. Absolute or Relative Quantitation of AKT/mTOR Signaling Pathway Proteins Using SureQuant



AKT/mTOR pathway proteins were enriched through multiplex immunoprecipitation using the SureQuant AKT Pathway or AKT Phospho Pathway Mass Spec Assay kit. Analyses were performed on a Q Exactive HF Orbitrap mass spectrometer using directed discovery (DDA) and targeted MS (PRM) acquisition methods. DDA and PRM data were analyzed in Proteome Discoverer and Skyline software, respectively. PRM analysis using the calibration curve allowed absolute quantitation of each target peptide from positive control lysate. Western Blot data showed differential expression for phosphorylated AKT pathway proteins with hIGF-1 stimulation and inhibitor treatments



Calculated concentration (fmol) values were used to summarize data where an increase in phosphorylation was designated by at least a 40% increase compared to untreated. Inhibition was designated if below 40% hIGF-1 treated value. All pathway phosphorylated proteins showed increase in abundance with hIGF-1 treatment across all three cell lines, whereas inhibitors functioned differently among three cell lines:

- The PI3K inhibitor (LY294002) was more effective than the mTOR inhibitor (Rapamycin) with effectiveness HCT116 > MCF7 > A549
- Combined LY294002 + Rapamycin treatment was most effective in HCT116 cells with effectiveness HCT116 > A549 > MCF7 Dual inhibitor (NVP-BEZ235) worked best in A549 cells compared to LY294002 and Rapamycin together with
- effectiveness A549 > HCT116 > MCF7

Figure 11. Quantitation of Total and Phospho AKT Pathway Proteins in Lung, Breast, and Large Intestine Normal and Tumor Tissue Samples



SureQuant AKT Total and Phospho Pathway kits allowed accurate quantitation from lung, breast, and large intestine normal and tumor tissue lysates. Skyline Software was used to generate standard curves and calculate absolute concentrations of unknown samples.



Accurate quantitation was obtained from five patient derived lung tumor xenograft samples using the SureQuant AKT Phospho Pathway kit. Patient derived xenograft samples were lysed in 2% SDS before multiplex IP enrichment followed by Targeted MS (PRM) analysis using the Q Exactive HF-X and EASY-nLC 1200 system. A 50cm EASY-Spray column (ES803) in direct-injection setup was used for all LC-MS data acquisition and Skyline Software was used to generate standard curves (data not shown) and calculate absolute concentrations of unknown samples.

CONCLUSIONS

- Pierce™ LC-MS/MS System Suitability Standard (7 x 5 mixture) achieves appropriate linearity and dynamic range to assess system performance prior to acquisition of unknown samples.
- SureQuant Multiplex IP-MS and Absolute Quantitation Modules for AKT pathway proteins allowed simultaneo quantitation of multiple total and phospho AKT pathway proteins in treated cell lines and tumor samples with high accuracy and precision (CV <20%).
- Analysis of abundance levels between three different cancer cell lines using the SureQuant AKT Total and Phospho kits revealed preferences for certain inhibitors in specific cell lines, with PI3K inhibitor LY294002 demonstrating highest efficacy in HCT116 cells, whereas the dual PI3K/Rapamycin inhibitor NVP-BEZ235 was most effective in A549 cells. Orthogonal evaluation between PRM assays and Western Blot analysis of three cancer cell lines treated with hIGF-1 and various inhibitors, showed similar trends in the protein expression changes, albeit the level of precision and dynamic range achieved with the SureQuant kit is difficult or impossible to achieve with Western Blot analysis.
- SureQuant AKT pathway kits are amenable to diverse sample sources and allowed identification of target proteins from cell lysate, tissues and patient derived xenograft tissue samples.
- SureQuant AKT pathway kits are amenable to diverse sample sources and allowed identification of target proteins from cell lysate, tissues and patient derived xenograft tissue samples.

REFERENCES

- Logue JS, Morrison DK. Complexity in the signaling network: insights from the use of targeted inhibitors in cancer therapy. Genes Dev. 2012 Apr 1; 26(7):641-50.
- Carr SA, Abbatiello SE, Ackermann BL et al. Targeted Peptide Measurements in Biology and Medicine: Best Practices for Mass Spectrometry-based Assay
- Development Using a Fit-for-Purpose Approach. Mol Cell Proteomics. 2014 Mar; 13(3):907-17. Ackermann BL. Understanding the role of immunoaffinity-based mass spectrometry methods for clinical applications. Clin Chem. 2012 Dec; 58(12):1620

For Research Use Only. Not for use in diagnostic procedures.

TRADEMARKS ANDS LEGAL INFORMATION

© 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of

PO65554-EN0519S



Mendoza MC, Er EE, Blenis J. The Ras-ERK and PI3K-mTOR pathways: cross-talk and compensation. Trends Biochem Sci. 2011 Jun;36(6):320-8

