# **Complex Peptide Mixture Fractionation via Parallel Isoelectric Focusing on a digital ProteomeChip for Direct LC-MS/MS Analysis**

### Overview

ores

**Purpose:** To develop an LC-MS/MS compatible workflow for peptide analysis

### Introduction

Protein Forest Inc. has developed the digital ProteomeChip (dPC) that traps peptides in acrylamide gel features based on charge, with a pH resolution of less than 0.1 units and run times less than one hour. The dPC is composed of 41 gel plugs each at a specific pH, creating a discrete 2.0 pH gradient across the dPC with a resolution of 0.05 pH units. This workflow employs buffers and conditions that have been designed to be MS compatible, eliminating the need for post-separation sample clean-up and dramatically reducing artifacts introduced by the separation process.

### Methods

### **Sample preparation for dPC fractionation**

Frozen Human A431 cell lysates were reduced with tributylphosphine and alkylated with iodoacetamide. Proteins were precipitated in methanol/chloroform and dissolved in 50 mM ammonium bicarbonate. Trypsin was added and the sample was allowed to digest overnight at 37°C. After proteolysis, peptides were reconstituted with dPC running buffer. Chips were run for a maximum of 45 min, briefly rinsed with water and gel plugs ejected into microcentrifuge tubes.

#### **Sample preparation for LC-MS/MS analysis**

Gel plugs were pooled in groups of 5, corresponding to 0.25 pH units, and extracted with 200 µL of 0.2% formic acid in 50% acetonitrile at 37°C for 1 hour. Extracts were collected in fresh microcentrifuge tubes and concentrated to a volume of 50  $\mu$ L. Samples were then loaded onto a Thermo Scientific Micro AS, equipped with a 25 µL syringe, a 2.4  $\mu$ L sample needle, a 200  $\mu$ L buffer tubing and a 10  $\mu$ L sample loop. The wash solvent used was 0.1 % formic acid in 30% acetonitrile. The pump was a Surveyor MS pump plus equipped with a static split at a ratio of 800:1 at 250 µL/min at 95% A (0.1 % formic acid in water), 5% B (0.1% formic acid in acetonitrile) with a 75  $\mu m \times 15$  cm column packed with Michrom Bioresources Magic 5  $\mu m C_{18}$  media. The Thermo LTQ<sup>™</sup> linear ion trap mass spectrometer was used in a "Top 8" configuration.

#### **Data Analysis**

Raw files from the LTQ mass spectrometer were searched using SEQUEST® through Thermo Scientific Bioworks 3.3.1 SR1 software package against the Refseq Human database, with static carbamidomethyl modified cysteines and differentially modified oxidized methionines. SEQUEST data was exported and analysed further by the Protein Forest Mass Spectrometry Research Analysis Tool (MSRAT<sup>®</sup>)

Dayana Argoti<sup>1</sup>, William Kuhlman<sup>1</sup>, James R. Dasch<sup>1</sup>, Stephen Haralampu<sup>1</sup> <sup>1</sup>Protein Forest, Inc., Lexington, MA 02421;



Figure 1. Top. Fractionation by parallel isoelectric focusing on the dPC. Bottom. Peptides and proteins are charged when in an environment that is either above or below their isoelectric point (pI) allowing them to migrate in an electric field until they encounter a gel that is at or near their pI and lose both charge and mobility.

## Workflow

Complex dPC Peptide Peptide ID Peptides Protein Separation LCMS & concentrate Sample ~ 45 min Search ~90 min dPC Fractionato Data 1) Harvest plug into Eppendorf Analysis tube or 96 well plate: **2**) Add 100 µl 50% ACN, 0.2% formic acid, incubate 1 hr @ b) Transfer fluid to fresh tube, SpeedVac to  $\sim 50 \ \mu$ l 4) Sample is ready for LC-MS

**Figure 2.** Overview of integrated workflow for maximum peptide recovery

5 10	15 20	25	30	35
Red paperson	Yellow peptide			
and and a second se				
5 10	15 20	25	30	35
Rednermon	Vellow peptide			

**Figure 3.** Comparison of two separate dPC runs on the 4.2-6.2 chip, each loaded with 50  $\mu$ g aliquot of an A431 cell lysate, 25  $\mu$ g of red and 10  $\mu$ g of yellow flourescent peptides.



Figure 4. Focusing reproducibility between two separate dPC runs across the current working range. Each dPC was loaded with a 50 µg aliquot of an A431 tryptic digest. Grey to black boxes represent pools of plugs in which the same peptides were found in each of the two chips. The number of peptides observed is represented by the degree of intensity towards black.



**Figure 5.** Number of proteins identified in each pH range and reproducibility observed between two technical replicates at each pH range.

### Advantages

•Reproducible peptide focusing with overall RSD <20%

# Conclusions

3.5-8.0

### References

Immobilized pH Gradient Isoelectric Focusing as a First-Dimension (1) Separation in Shotgun Proteomics; Journal of Biomolecular Techniques 2005,16, 181–189. G. Zilberstein et al, Parallel isoelectric focusing chip; Proteomics 2004, 4, (2) 2533–2540.

Acknowledgements Protein Forest would like to extend our thanks to David Sarracino at Thermo Fisher/ Scientific BRIMS center for his input and suggestions.



Figure 6. LC-MS/MS base peak chromatograms for 5-plug pH fraction 5.95-6.20 from two 4.20-6.20 dPC technical replicates each loaded with a 50 µg aliquot of an A431 tryptic digest. Illustrating the typical run to run degree of reproducibility between both dPC fractionation and LC-MS/MS analysis.

Minimized sample handling prior to LC-MS/MS analysis

•Fractionation to MS ready samples in 3 hours

•Only three simple steps to LC-MS/MS analysis

•Harvest plug  $\rightarrow$  Elute peptides  $\rightarrow$  Concentrate

•More than 3,500 peptides were found using three dPCs with a combined pH range of

•Approximately 1,600 unique proteins were reproducibly identified across the entire working pH range

•Future work will further optimize the workflow for targeted quantitaive assays and extend pH range to pH = 9.0