

Reference Yeast and Human Whole Cell Protein Extracts for Mass Spectrometry Instrument Performance Monitoring and Method Development



Sergei Saveliev, Ethan Strauss, Mike Rosenblatt and Marjeta Urh

Promega Corporation, Madison, WI sergei.saveliev@promega.com

1. Introduction

To monitor mass spec instrument performance for proteomics applications and optimize protein mass spec sample preparation, a reference protein material of high complexity is required. Whole cell protein extracts provide this desired sample complexity. However, to be compatible with mass spec applications, such extracts should meet a number of design requirements:

- ✓ compatibility with LC/MS (free of detergents, etc.)
- ✓ high protein integrity (minimal level of protein degradation and non-biological PTMs)
- ✓ compatibility with common sample preparation methods such as proteolysis, PTM enrichment and mass-tag labeling
- ✓ Lot-to-lot reproducibility

Here we describe whole cell protein extracts from yeast and human cells that meet the above criteria. Two extract formats have been developed:

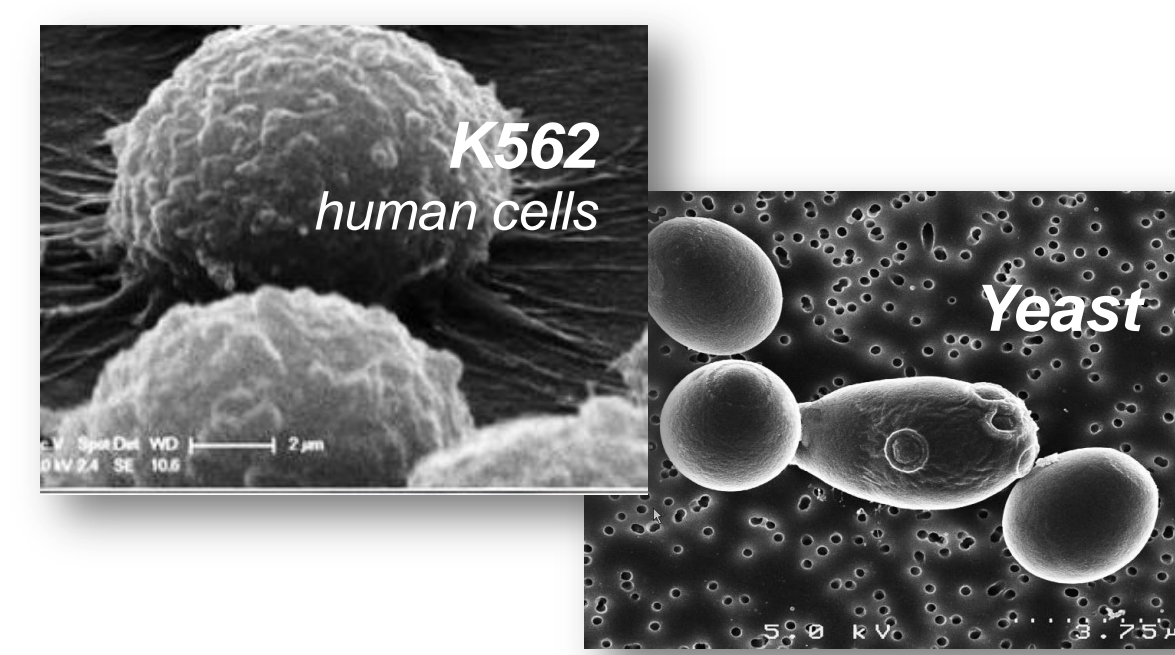
- Pre-digested extracts (peptides); primary use - instrument performance monitoring
- Intact protein extracts; primary use – optimization of sample preparation

2. The extract features

Source

Human extract - from K562 cells
Complex human proteome with a large dynamic range.

Yeast extract - from *Saccharomyces cerevisiae*
~6,600 ORFs; all proteins quantified previously.



Extract preparation

Cell lysis

- The process is optimized to allow for minimal protein fragmentation.
- Instant inactivation of endogenous proteases to assure virtual lack of nonspecific cleavages.
- Only the reagents compatible with LC and MS are used.

Sample preparation and digestion

- Strict control over the procedural steps assures that non-biological PTMs are maintained at lowest level.
- Proteins are exhaustively digested (less than 8% of missed tryptic cleavages).

Peptide clean-up and lyophilization

- Non-peptide material is removed with solid phase extraction (C₁₈).
- Peptides are dispensed in 100 µg aliquots with high accuracy and lyophilized for long term stability.

Mass spec analysis report for the extracts

Extract	Human	Yeast
Identified proteins	2120	1637
Total spectra	26440	21264
Spectrum match	65%	52%
Unique peptides	16375	12079
Deamidation spectra	4.9%	6.5%
Oxidation spectra	0.6%	1.3%
Carbamylated spectra	0.2%	0.6%
Nonspecific cleavages	1%	2%
Missed tryptic cleavages	4.9%	4.8%

Q Exactive (Thermo), 2h gradient
1 µg extract injection

Strict control over all the steps of preparation protocol assures that the extracts meet the critical requirements of reference materials for mass spectrometry.

3. Pre-digested extracts

Reference material for instrument performance monitoring

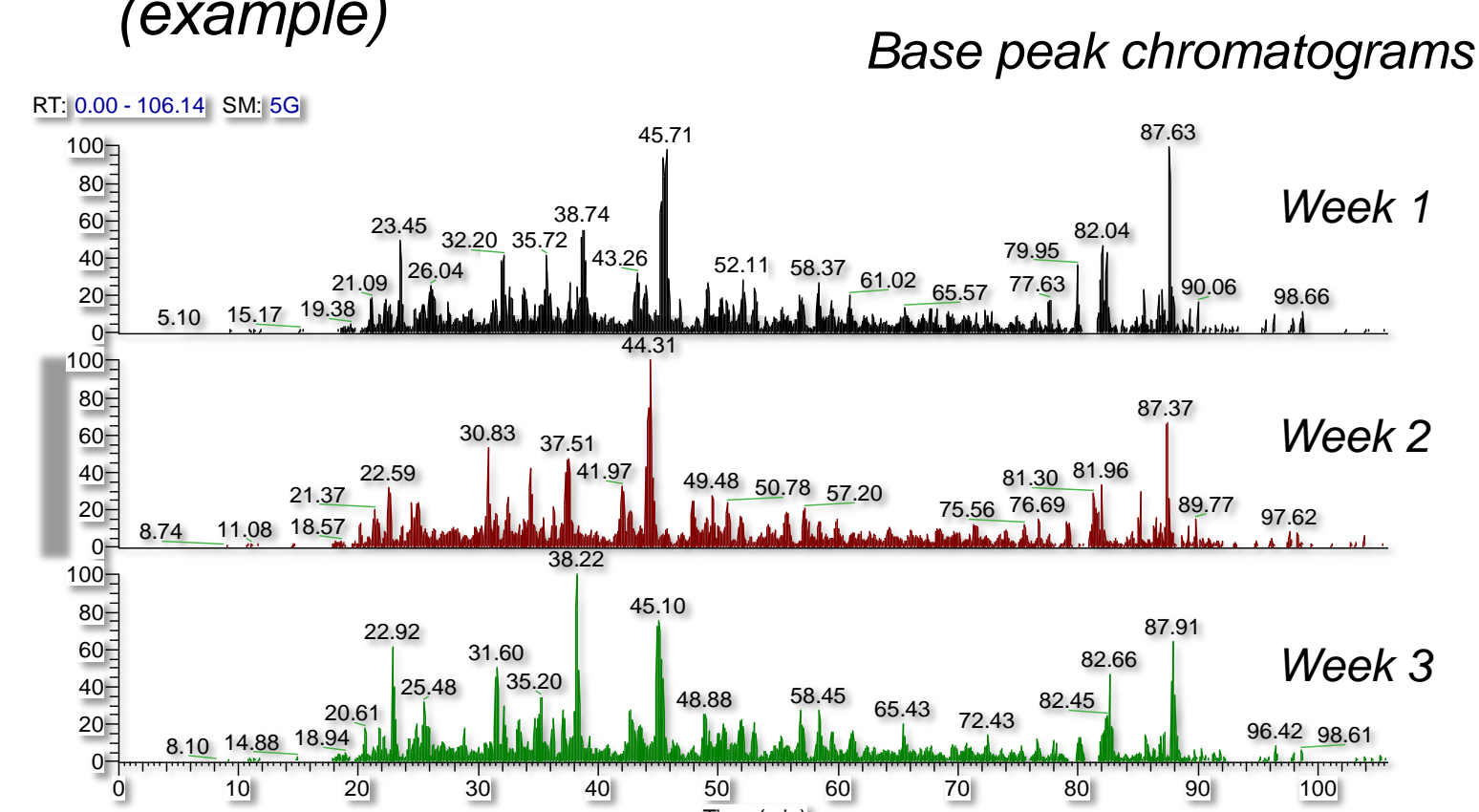
Features

- ✓ Full compatibility with LC and MS
- ✓ Ready for analysis - no need to reduce, alkylate, digest, or clean-up the sample
- ✓ Efficiently digested (less than 8% missed cleavages)
- ✓ Highly pure (SPE C₁₈ clean-up)
- ✓ Select material complexity you need – relatively small yeast proteome or complex human proteome with a large dynamic range

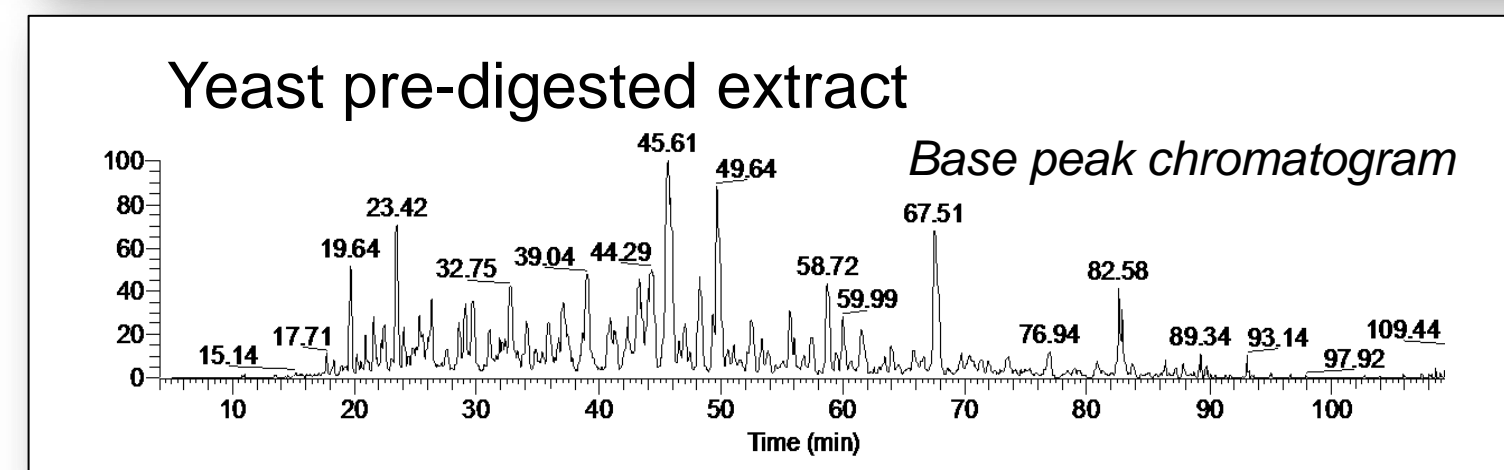
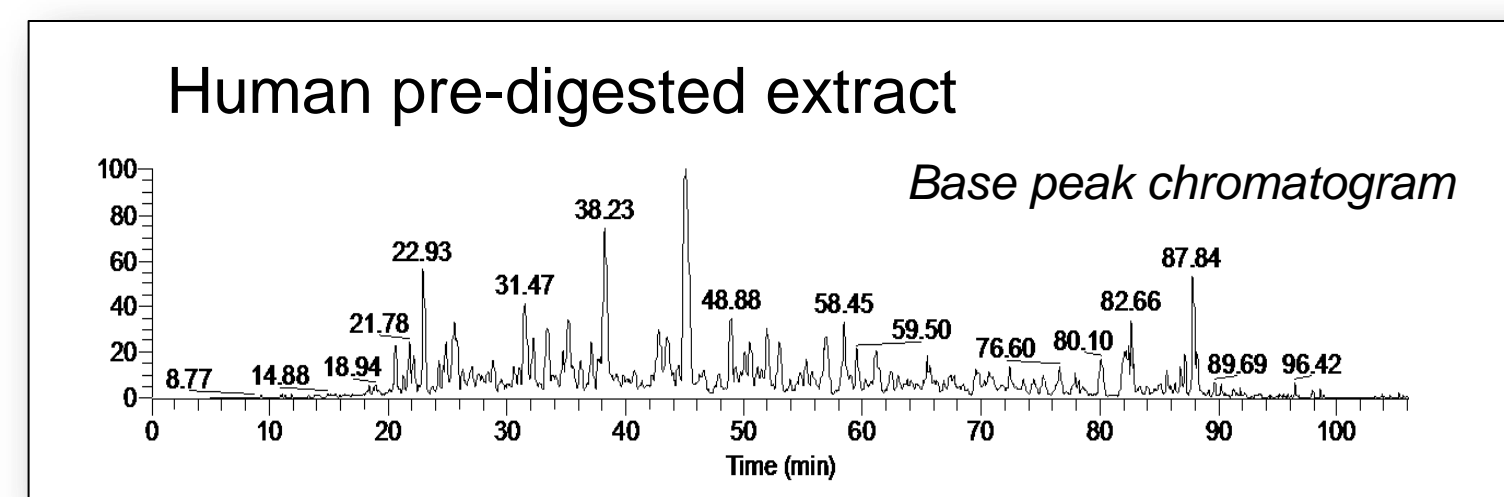
Applications

- ✓ Instrument performance monitoring
- ✓ Mass spec method optimization
- ✓ Interlaboratory instrument benchmarking

Monitoring LC/MS instrument performance (example)



Instrument performance (Waters NanoAquity HPLC system interfaced to a ThermoFisher Q Exactive) was monitored with human pre-digested extract (1 µg injection, 2h gradient).



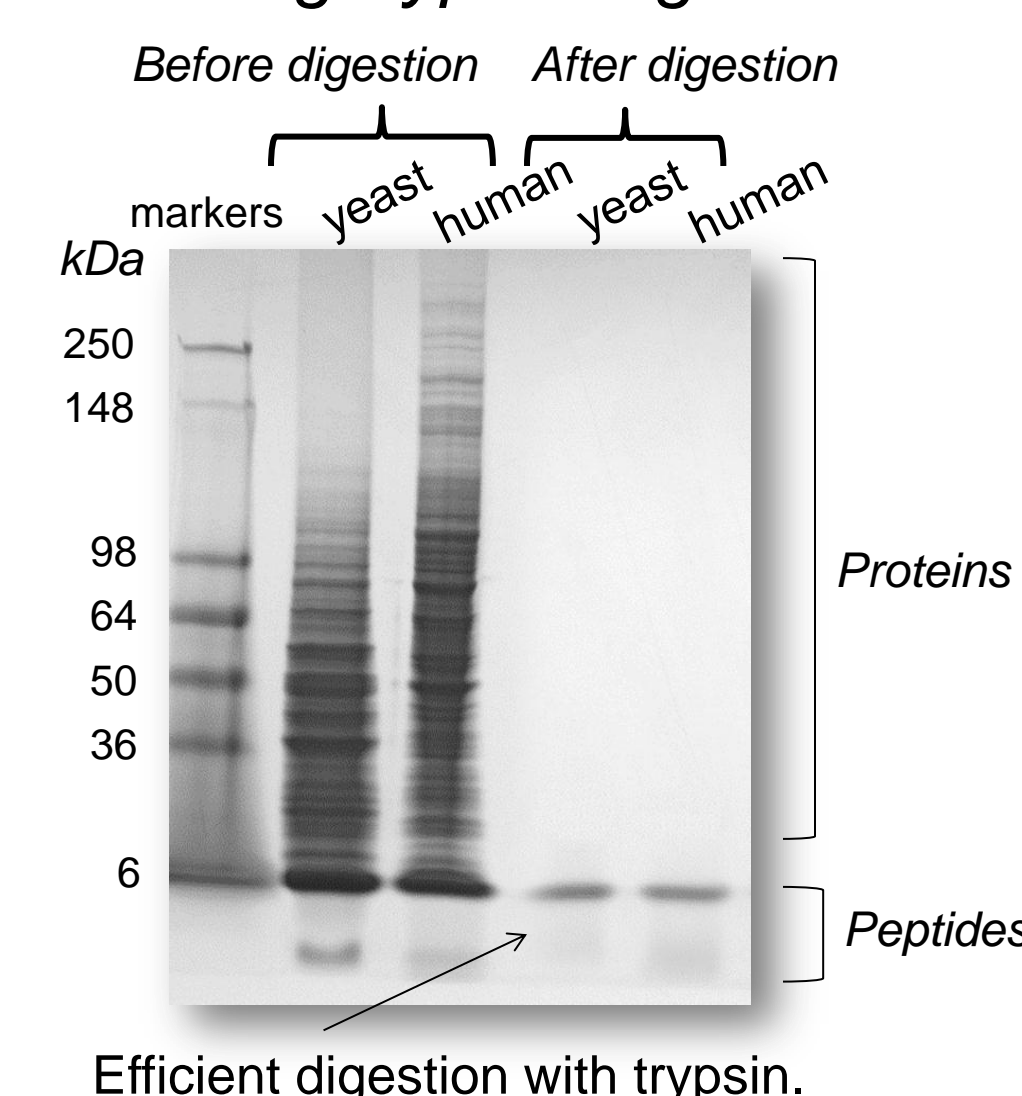
Test results by different users

Instrument, laboratory and method	Pre-digested extract	Identified unique peptides	Identified proteins
Instrument #1, Lab #1, method #1	yeast	12427	1567
	human	18040	2370
Instrument #1, Lab #2, method #2	yeast	27491	3096
	human	46018	4664
Instrument #2, Lab #3, method #3	yeast	11233	1220
	human	10344	1588
Instrument #3, Lab #4, method #4	yeast	14064	2703
	human	19531	3885

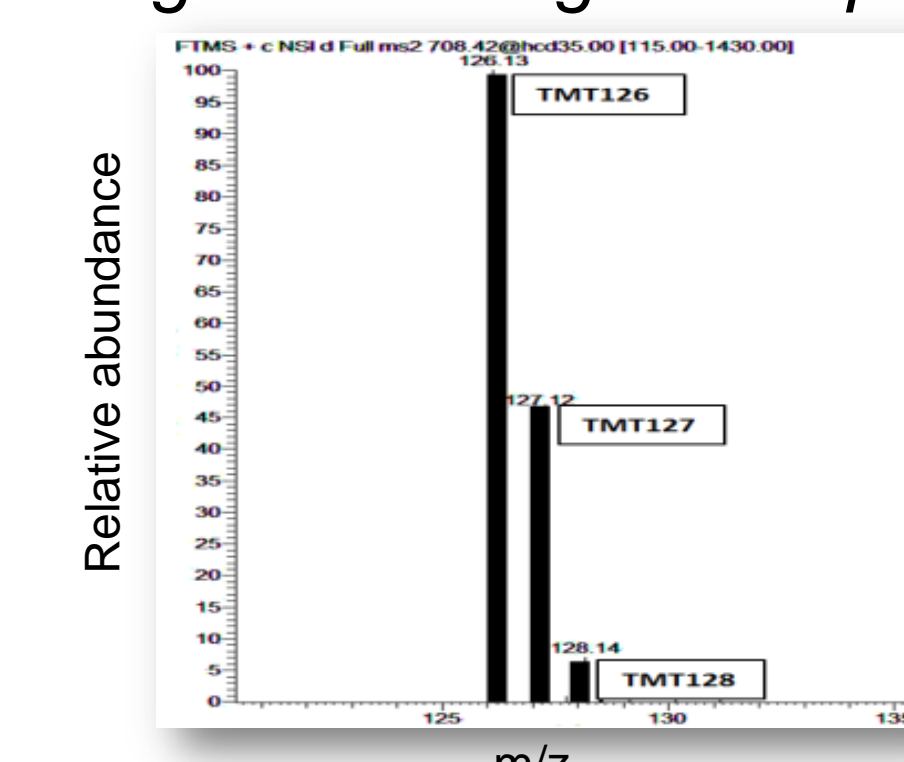
4. Intact extracts

Test material for optimization of sample preparation

Testing trypsin digestion



Optimizing isobaric tag-based protein quantitation



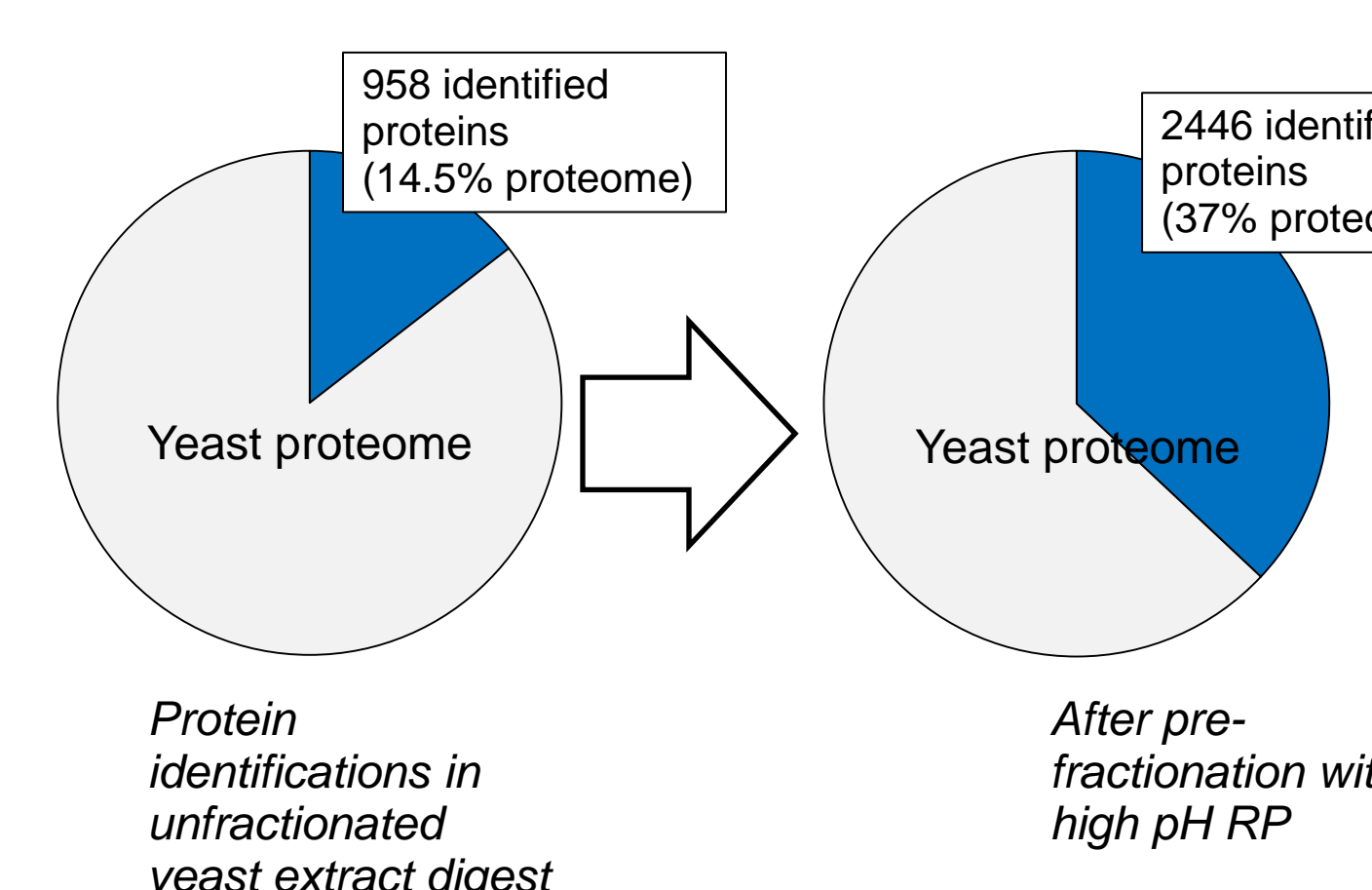
Three aliquots of the digested human extract were individually labeled with a specific TMT tag and mixed at 10:5:1 ratio. Good correlation between observed and expected protein quantitation levels was observed.

Validating phosphopeptide enrichment method

Extract	Total peptides	Phosphopeptides	Phosphopeptide enrichment
Yeast	2502	1715	68.5%
Human	1768	1537	87%

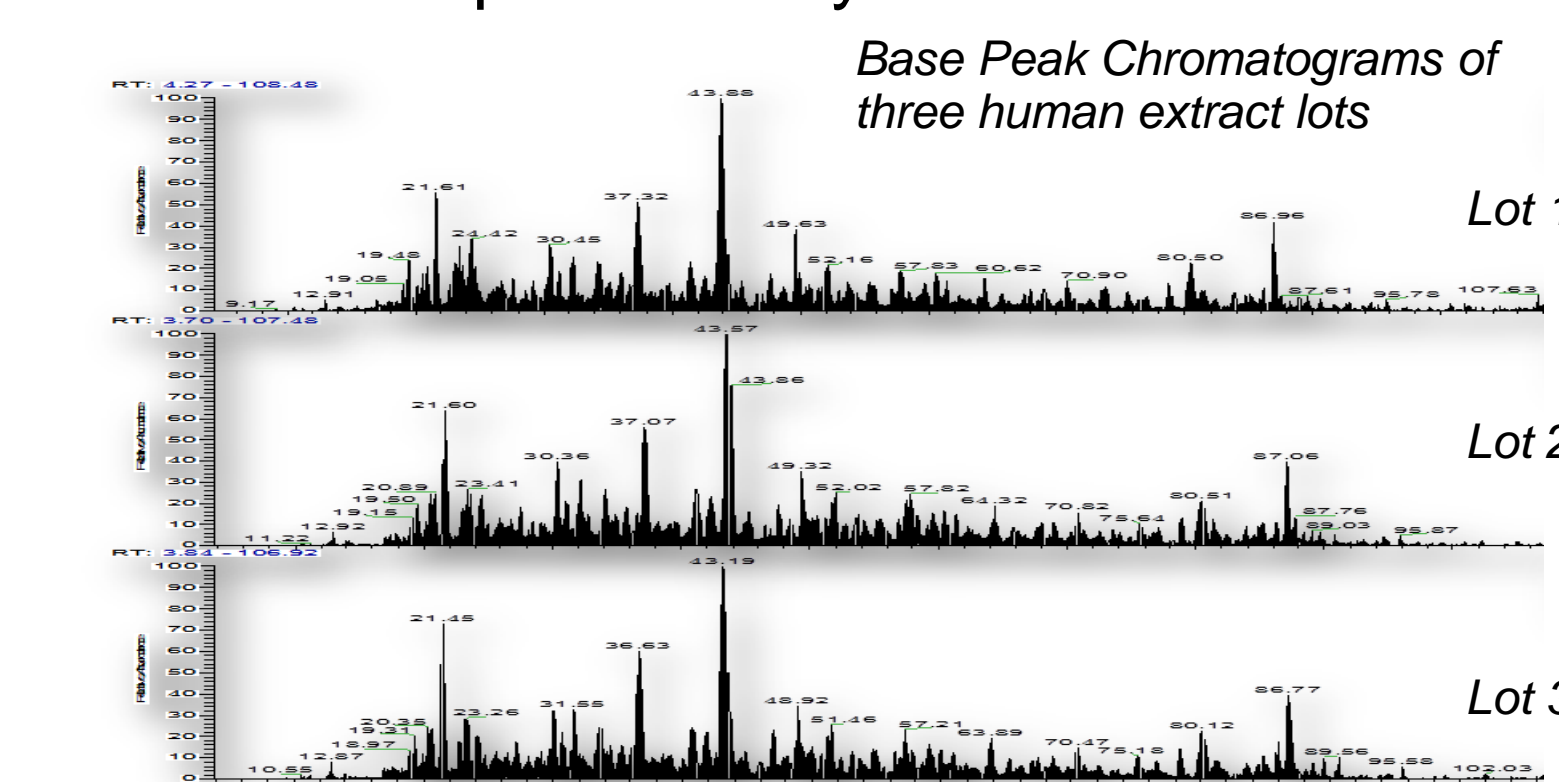
The extracts provided a good model system for establishing efficient phosphopeptide enrichment with TiO₂ resin.

Establishing a method for fractionation of a complex peptide mix



4. Lot-to-lot reproducibility

Lot-to-lot reproducibility



Waters NanoAquity HPLC system interfaced to a Thermo Fisher Q Exactive, 2h gradient, 1 µg extract injection.

Lot	#1	#2	#3
Identified proteins	2101	2076	2088
Unique peptides	25116	25821	25883
Deamidation spectra	4.9%	5.1%	5.0%
Oxidation spectra	0.6%	0.5%	0.6%
Nonspecific cleavages	1%	1%	1%
% missed tryptic cleavages	4.8%	4.8%	4.6%

Quantitation via spectral counting

Human extract Spectral counting				Yeast extract Spectral counting			
protein	lots			protein	lots		
	#1	#2	#3	Protein	#1	#2	#3
HS90B	155	157	153	YBR118W	358	375	368
ACTG	116	113	113	YGR192C	356	344	335
HS90A	134	130	129	YCR012W	369	336	304
PRKDC	126	129	116	YBR196C	153	153	146
TBB5	96	96	99	YLR249W	130	129	128
IF2P	18	17	17	YDL145C	23	22	25
ANXA2	21	23	23	YDR261C-D	31	30	28
DLDH	19	16	18	YKL035W	31	26	32
PDIA4	21	27	22	YOR375C	28	32	28

Quantitation via SRM assay
Courtesy by Dr. Koomen, Moffitt Cancer Center

Protein	Peptide	Lot	Normalized abundance*
STAT3_HUMAN	TLTDEELADWK	A	1.10
		B	0.83
		C	1.00
STAT5A_HUMAN	LSPAPGLFTSAR	A	0.92
		B	1.09
		C	1.00
	YYTPVLAK	A	1.00
		B	1.09
		C	0.93

*GAPDH normalized data

Lot-to-lot extract reproducibility with respect to protein composition and abundance

5. Conclusions

- We developed MS ready whole cell yeast and human protein extracts
- Our method allows for protein recovery with high reproducibility and minimal level of protein fragmentation or non-biological PTMs
- The extracts are provided in two ready-to-use alternative formats
 - Pre-digested extracts (tryptic peptides) serve the need of instrument performance monitoring
 - Intact protein extracts are designed for sample preparation method development and optimization
- The extracts provide reference material for comprehensive LC/MS instrument validation, performance monitoring and method development. They can also be used as a model test material for optimizing protein mass spec sample preparation.