

Efficiency assay of detergent removal columns on protein and peptide samples for mass spectrometric analysis



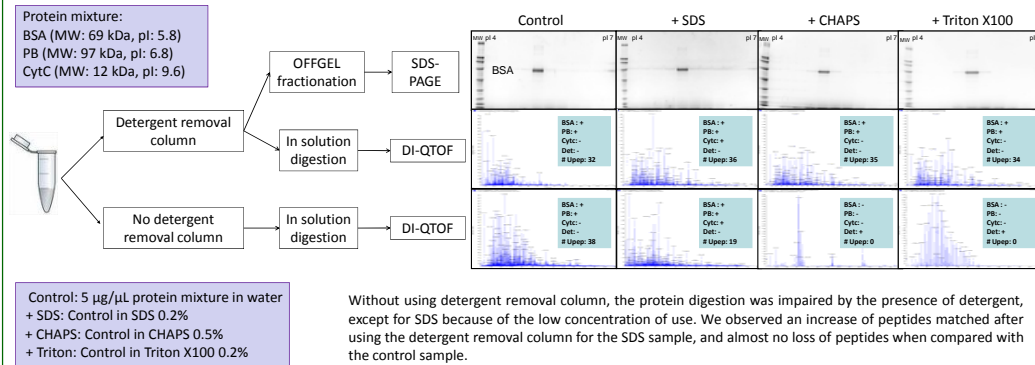
Sophie Alvarez¹, Aftab Alam², Leslie M. Hicks¹

¹Donald Danforth Plant Science Center, St Louis MO 63132, USA

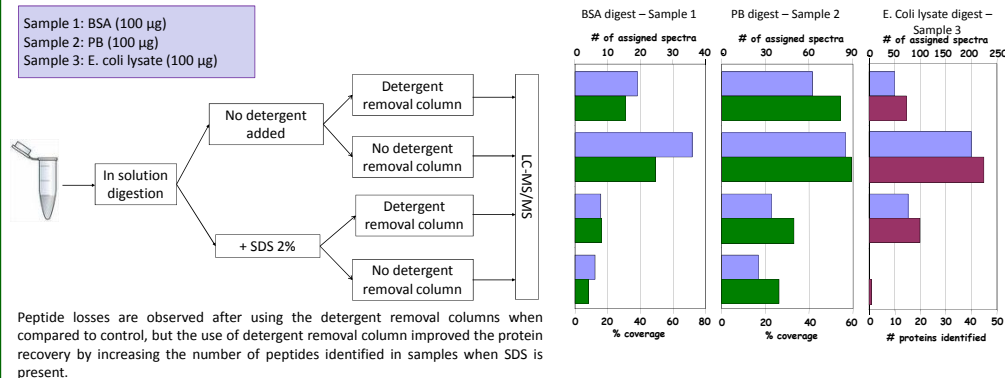
²G-Biosciences, Maryland Heights, MO 63043, USA

Introduction: Detergents are essential for protein solubility during protein extraction and sample preparation, especially when working with hydrophobic proteins (e.g. proteins tightly bound to cell membranes). However, the presence of high concentrations of detergents in protein samples can impair protease digestion of proteins and suppress peptide ionization when analyzed by mass spectrometry. New (GB-S10) detergent removal columns from G-Biosciences were tested for their efficiency to remove anionic, nonionic or zwitterionic detergents (e.g. SDS, TritonX100 or CHAPS) from protein and tryptic peptide samples with minimal sample loss for downstream analysis by mass spectrometry and other techniques.

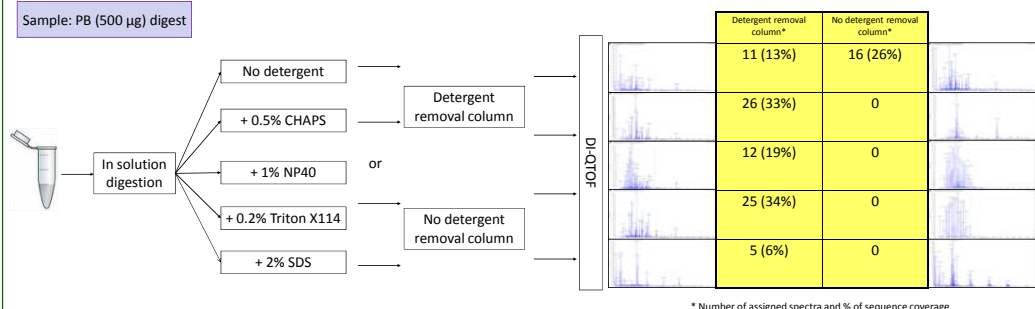
Detergent Removal from Protein Samples



Detergent Removal of High Concentrations of SDS from Simple and Complex Tryptic Peptide Samples



Detergent Removal of Various Detergents from Tryptic Peptide Samples



Summary:

Questions	0.2% SDS (anionic)	0.5% CHAPS (zwitterionic)	0.2% Triton X100 (nonionic)
Protein recovery	✓	✓	✓
Detergent removal from protein sample	✓	✓	✓
Protein electrophoresis	✓	✓	✓
Protein digestion	✓	✓	✓

Questions	2% SDS (anionic)	0.5% CHAPS (zwitterionic)	1% NP40 (nonionic)	0.2% Triton X114 (nonionic)
Peptide recovery	✓	✓	✓	✓
Detergent removal from peptide sample	✓	✓	✓	✓

Methods

OFFGEL electrophoresis and SDS-PAGE: 100 µg samples were separated on 4-7 pH strip, 12 fractions using the 3100 OFFGEL fractionator (Agilent). After applying 20 kV, fractions were collected and 1/3 was dried down and run via SDS-PAGE. The gel was stained with Sypro.

In solution digestion: Samples were reduced and alkylated before tryptic digestion.

DI-QTOF: Samples were resuspended in 5% ACN/ 0.1% FA, zipipped using C18, and infused using nanospray tips into an ABI QSTAR XL (Applied Biosystems/MDS Sciex) hybrid QTOF MS/MS mass spectrometer. TOF mass and product ion spectra were acquired using information dependent data acquisition (IDA) in Analyst QS v1.1 with the following parameters: mass ranges for TOF MS and MS/MS were m/z 300-2000 and 70-2000, respectively. Every second, a TOF MS precursor ion spectrum was accumulated, followed by three product ion spectra, each for 3 s.

LC-MS/MS: Nano-LC was performed with a nano 2D LC (Eksigent) equipped with a Dionex C18 PepMap100 column (75 µm i.d.) flowing at 200 nL/min. Peptides (5 µL injections) were resolved on a gradient from 90.5% solvent A (0.1% FA in MilliQ water) and 9.5% solvent B (0.1% FA in ACN) to 25% B in 4 minutes, then increasing to 40% B over 75 minutes, and from 40-90.5% B over the final 5 minutes. Mass spectrometer parameters are identical to those described above.

Database searching: The peptide tandem mass spectra were searched against NCBItrn using an in-house version of MASCOT v2.2 (Matrix Science Inc). The following parameters were selected: tryptic peptides with s1 missed cleavage site; precursor and MS/MS fragment ion mass tolerance of 0.8 and 0.8 Da, respectively; fixed carbamidomethylation of cysteine; and variable oxidation of methionine. Data was then compiled in Scaffold (Proteome Software). Positive identification was determined based on the following criteria: ≥ 2 peptide sequences, minimum peptide probability of 50% and minimum protein probability of 99%.