

Beyond TurboTMT: Phi-SDM Super-resolution Methods for Next-generation Highly-multiplexed Quantitative Ultrasensitive and Single-Cell Proteomics via TMTPro Complement Ion Deconvolution

Hendrik Wesseling¹; Konstantin Aizikov²; Dmitry Grinfeld²; Arne Kreutzmann²; Daniel Marc Mourad²; Oliver Lange²; Alexander A. Makarov²; Erik Hett¹; David H. Perlman¹
¹**MERCK** Cambridge Exploratory Science Center, Cambridge, MA
²**ThermoFisher** Scientific (Bremen) GmbH, Bremen, Germany

Introduction

Single-Cell Proteomics has emerged as an extremely promising new field capable of unlocking the mechanisms behind the vast cellular heterogeneity that defines biology, disease, and therapeutic response. However, enormous challenges exist, including how to maximize both quantitative signal and experimental throughput. We utilize the multiplexed method (e.g., SCoPE-MS¹, SCoPE2^{2,3}) in which individual cells' proteomes are labeled, pooled, and combined with abundant carrier material using tandem mass tag (TMT) reagents, improving both peptide identification and cell throughput. Still, entire proteomes must be gleaned through short single-shot nanoLC-MS runs that are plagued by limited ion current. Under these circumstances, relatively long trap fill-times are necessary to obtain sufficient peptide signal, and, optimally, both identification and accurate quantification need to be achieved from single MS2 fragmentation spectra for sufficient throughput. Unfortunately, accurate quantitation of TMT reporter ions in MS² remains challenging, due to the co-isolation artifacts they suffer.

TMT complement (TMTc) ions, consisting of the peptide plus TMT balancer moieties, have the potential to provide quantitative read-outs in the MS² spectra that are essentially free of co-isolation artifacts, so long as the ions can be sufficiently resolved. Specifically, in order to deconvolute the nearly isobaric ¹⁵N and ¹³C TMTc species of the TMTProTM products so as to exploit these reagents' full multiplexing power, ultra-high spectral resolution is required.

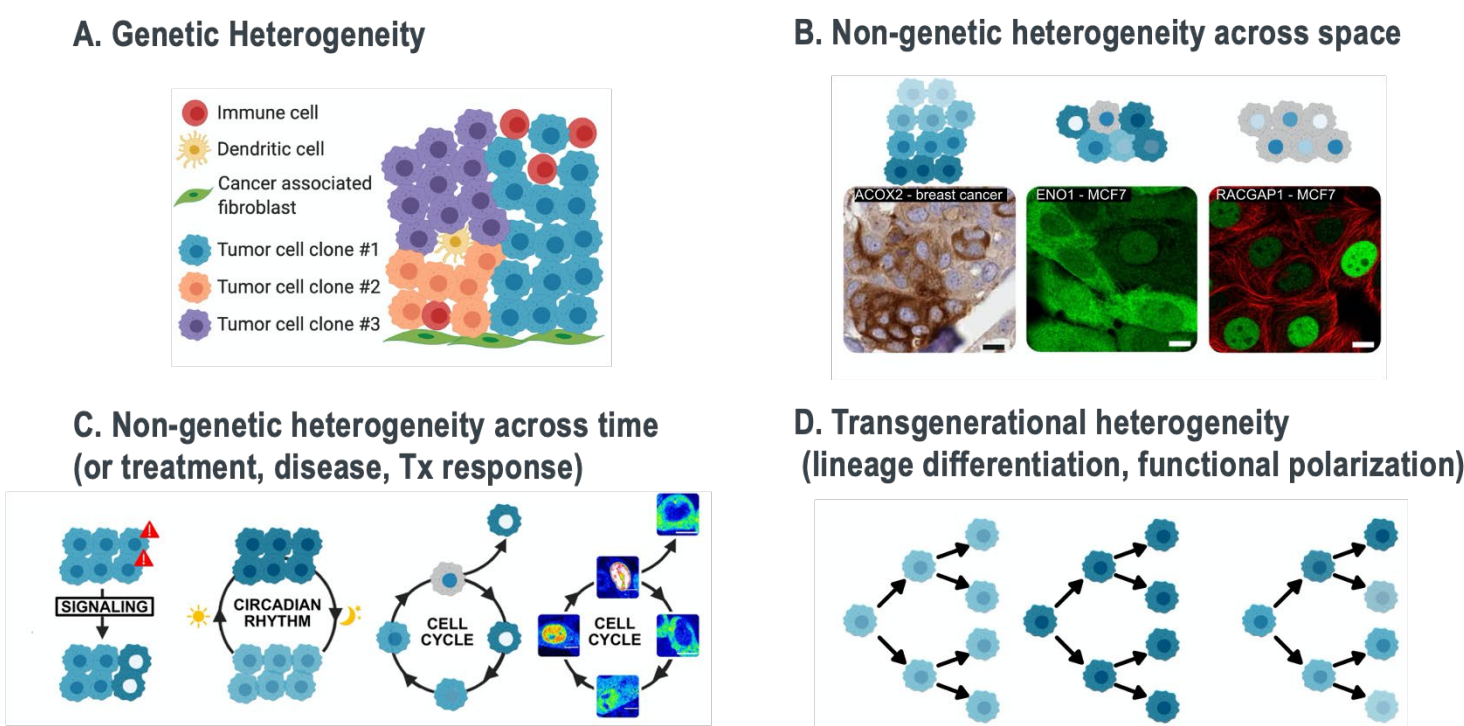
Recently, Phi-SDM signal processing, which is capable of resolving TMT reporter ions from transients 2-3-times shorter than required by conventional eFT processing^{1,5}, was introduced and commercialized as the TurboTMTTM feature in the Orbitrap EclipseTM and ExplorisTM MS instrument lines. Phi-SDM is similarly capable of rendering 2-3-times higher resolution spectra than eFT for any given transient length. Here, we have explored whether Phi-SDM directed not at the TMT reporter ions, as in TurboTMT, but rather at the high-mass TMTc ions might provide sufficiently high-resolution measurement to enable their full deconvolution.

¹Budnik, B., Levy, E., Harmange, G. et al. Mass spectrometry of single mammalian cells quantifies proteome heterogeneity during cell differentiation. *Genome Biol* 19, 161 (2018)
²Specht H, Emmott E, Petelski A, Huffman RG, Perlman DH, Serra M, Kharchenko P, Koller A, and Slavov N. Single-cell proteomic and transcriptomic analysis of macrophage heterogeneity using SCoPE2. *Genome Biol* volume 22, 50 (2021)
³Petelski A, Emmott E, Leduc A, Huffman RG, Specht H, Perlman DH. Multiplexed single-cell proteomics using SCoPE2. *Nature Protocols* (in press) (2021)
⁴Grinfeld D, Aizikov K, Kreutzmann A, Damoc E, and Makarov A. Phase-Constrained Spectrum Deconvolution for Fourier Transform Mass Spectrometry. *Anal. Chem.* 89, 2, 1202–1211. (2017)
⁵Kelstrup CD, Aizikov K, Bath TS, Kreutzman A, Grinfeld D, Lange Q, Mourad D, Makarov A, Olsen JV. Limits for Resolving Isobaric Tandem Mass Tag Reporter Ions Using Phase-Constrained Spectrum Deconvolution. *J Proteome Res* 17 (11), 4008-4016. (2018)

Methods

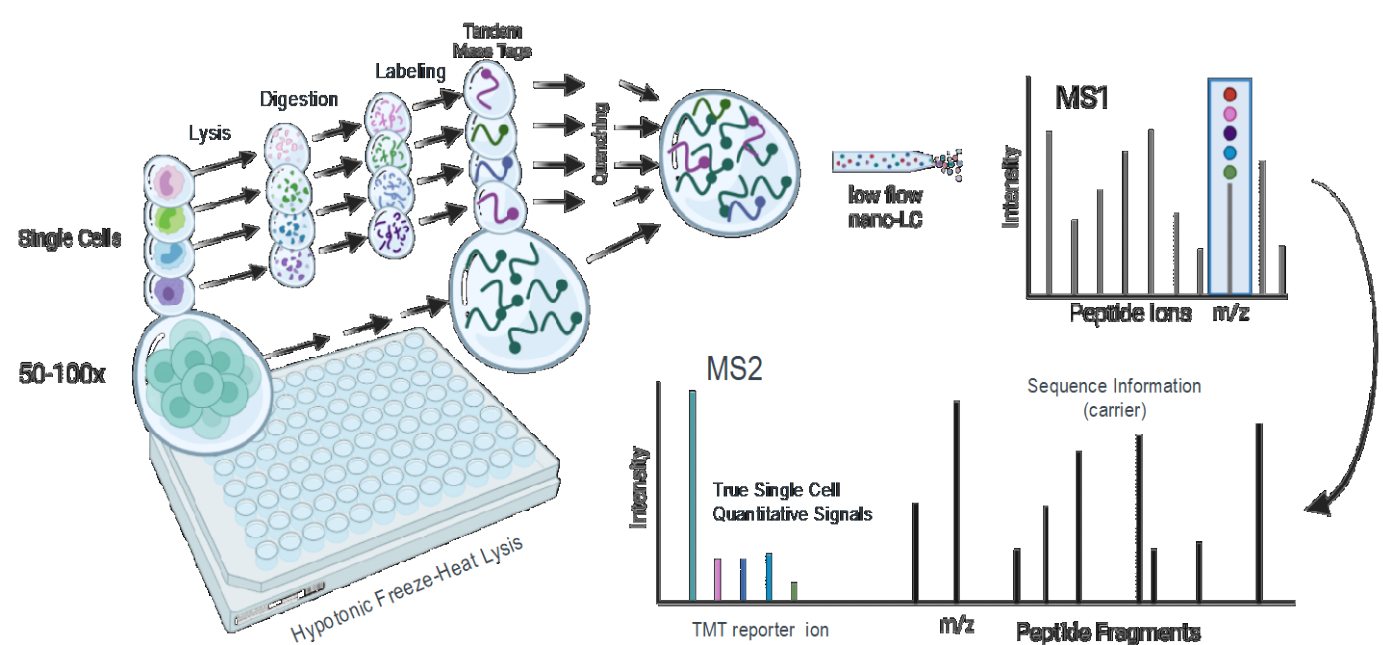
Single-Cell Proteomics dilution standards were made using commercial purified peptide standards (MassPREPTM Peptide Mixture, Waters; PierceTM HeLa Digest/PRTC Standard, Thermo ScientificTM; Mass Spec-Compatible Human Protein Extract, Promega) labeled with TMTProTM reagents (Thermo Scientific), according to the SCoPE-MS paradigm. Samples were injected on the single-cell level and analyzed on a Q ExactiveTM HF-X mass spectrometer (Thermo Scientific) outfitted with an EvosepTM Plus nano-flow LC device (Evosep), or an Orbitrap ExplorisTM 480 mass spectrometer outfitted with a VanquishTM Neo UHPLC system (Thermo Scientific). Research-grade versions of Tune instrument control software were employed that enabled high-resolution phase-constrained spectral deconvolution (Phi-SDM) across specified ranges, including TMTc ion regions. Data was acquired at various instrument resolution settings (as indicated), using either eFT or Phi-SDM processing, with ion fill-times adjusted to the maximum afforded for complete parallelization with transient acquisition. Spectra were visualized using XcaliburTM software (Thermo Scientific).

Cellular Heterogeneity is Fundamental to Human Biology



Mammalian cells are defined by their phenotypic and proteomic heterogeneity. This heterogeneity can be classified into at least 4 separate categories: (A) Genetic heterogeneity, (B) Non-genetic heterogeneity across space, (C) Non-genetic heterogeneity across time, and (D) Transgenerational heterogeneity.

The SCoPE-MS Method For Multiplexed Single-Cell Proteomics

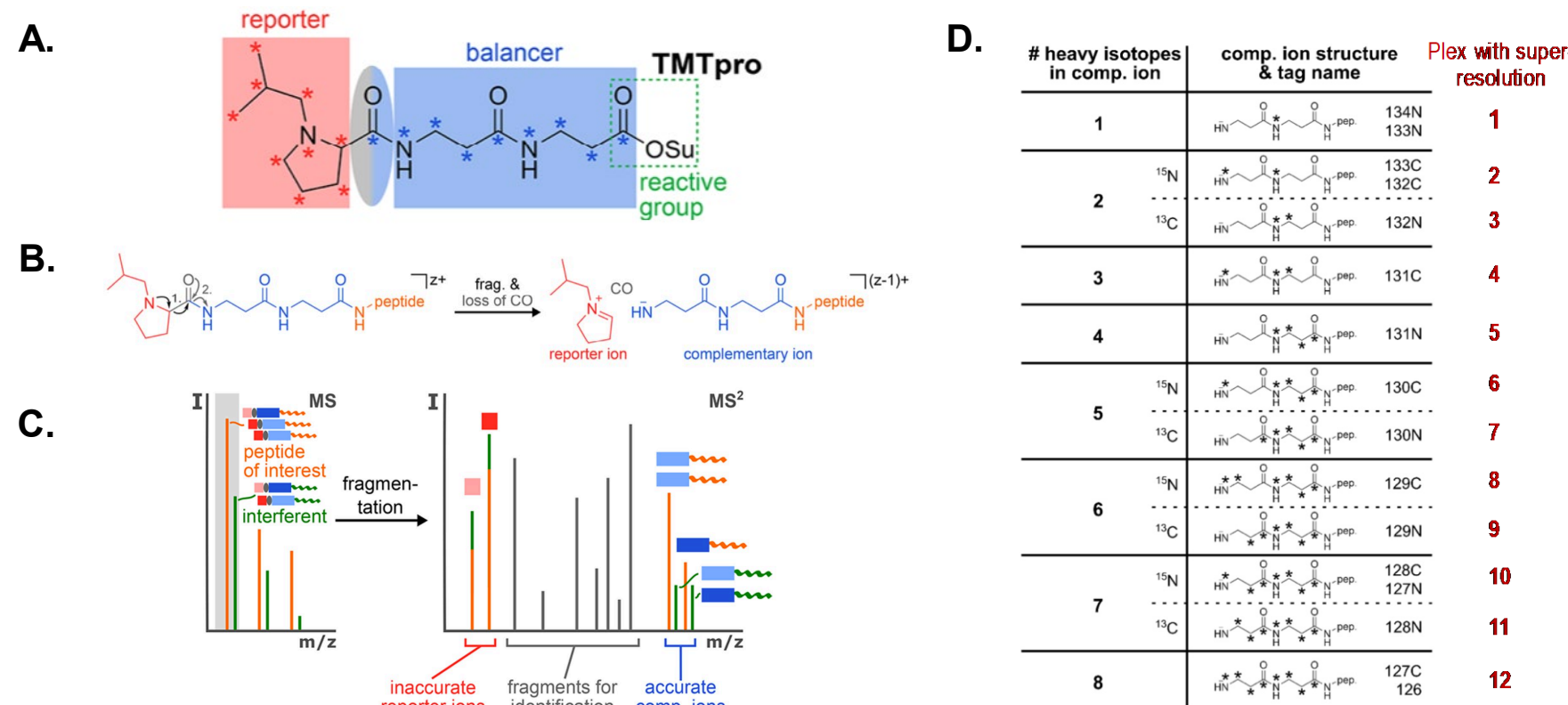


In the SCoPE-MS method for Single-Cell Proteomics 1-3, single mammalian cells are subject to a mass-spec compatible lysis free of harsh detergents or chaotropes, followed by digestion, TMT labeling, and pooling with an abundant (e.g., 20-200 cell-equivalent) TMT-labeled carrier material. Upon LC-MS/MS of the pooled sample, fragmentation spectra contain abundant peptide backbone ion signals lent to the sample from the carrier, which facilitate identification, and also minute, authentic single-cell TMT reporter ion signals that read out relative peptide abundance in each single cell.

Challenging Features of Single-Cell Proteomic LC-MS

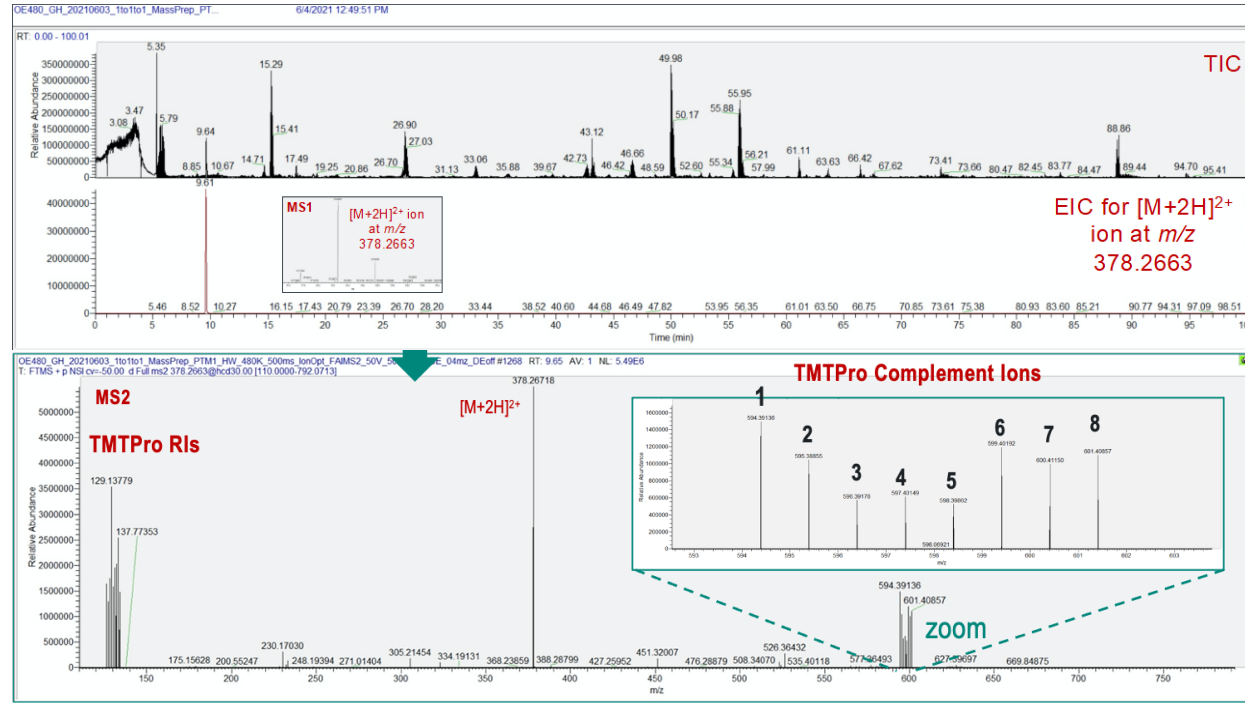
- Single shot LC-MS
 - "All your eggs in one basket": demands highest chromatographic resolution and extreme frugality to achieve maximal analytical depth.
- Signal starved
 - Majority of ions require long fill-times (e.g., 256ms, 512ms, etc.) for each MS2 scan.
 - Coordinates well with longer transient lengths (e.g., 120K res, 240K res settings) for high resolution Orbitrap measurement.
- Duty-cycle constraints
 - Due to time pressure and signal-intensity constraints, identification and quantification must be accomplished in MS². MS³-based methods provide diminishing returns.
- Need to maximize cell throughput
 - Necessitates maximal multiplexing, short LC runs, etc.
- TMT reporter ion species suffer MS² co-isolation artifacts
 - Caused by the presence of isobaric peptides, TMT-quench products (1+ and multiply charged), and TMT-labeled protein material (semi-intact trypsin, etc.)
 - Long fill times increase opportunity for co-isolation

TMT Complement Ions for Accurate Multiplexed Quantitation



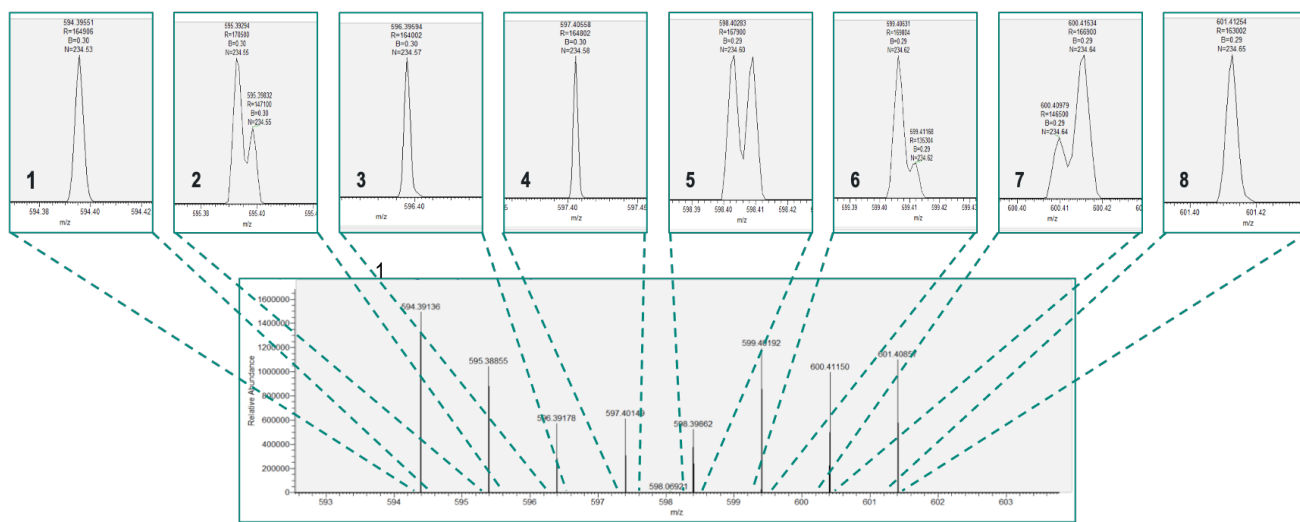
Quantitation via TMTc. (A) The structure of the TMTPro family of labels consisting of the reporter moieties joined to the balancer regions. (B) During gas-phase fragmentation and TMT-RI formation, the TMTc ion-species is simultaneously formed from the TMTPro balancer moiety fused to the parent peptide, thus it has a peptide-specific, plus TMT-tag specific mass. (C) When a TMT-labeled peptide of interest (red, MS) is co-isolated along with an interferent (green), it produces a fragmentation spectrum (MS²) in which the TMT-RLs from both the peptide and interferent are coincident, while their TMTc ion species retain distinct masses and are distinguishable. (D) The TMTPro 16-plex family forms 8 nominal mass TMTProC species, which can be further separated into 12 distinct species with ultra high-resolution methods.

Detection of TMTPro Complement Ions in labeled Peptide STD



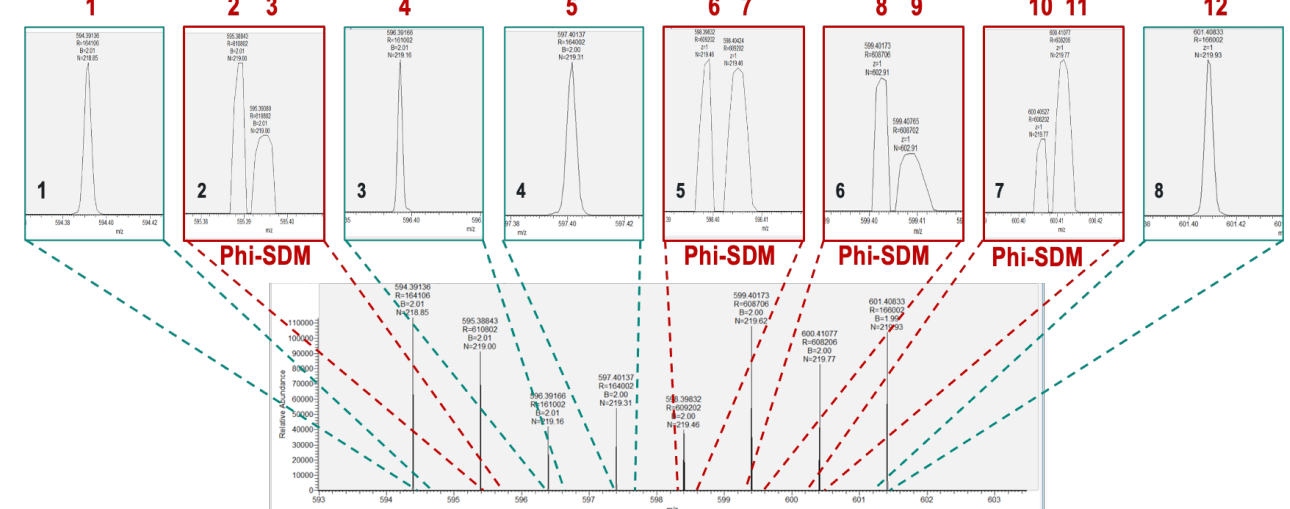
LC-MS/MS of TMTPro labeled peptide standards yields MS² spectra displaying both low-mass TMTPro-RI species, and also prominent TMTProC ion clusters in the higher mass region. Zooming in on these TMTProC peaks show that they consist of the predicted 8 nominal mass species.

Insufficient Resolution of TMTProC Ions at 240k eFT Setting



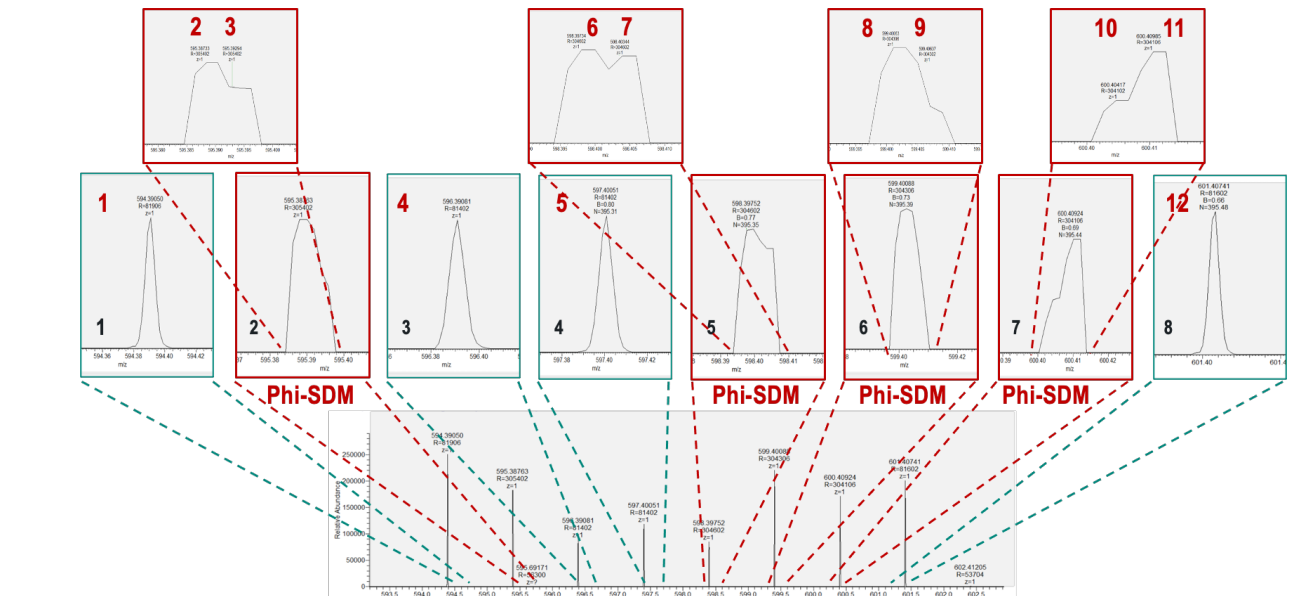
At the maximum resolution setting of the commercial Q Exactive HF-X instrument (240K using eFT) we are unable to completely resolve the 12 TMTProC species from the 8 nominal mass peaks. Zoomed in regions around each ion are shown above the 8-peak TMTProC cluster spectrum.

PhiSDM TMTProC Ion Deconvolution from a 512ms Transient



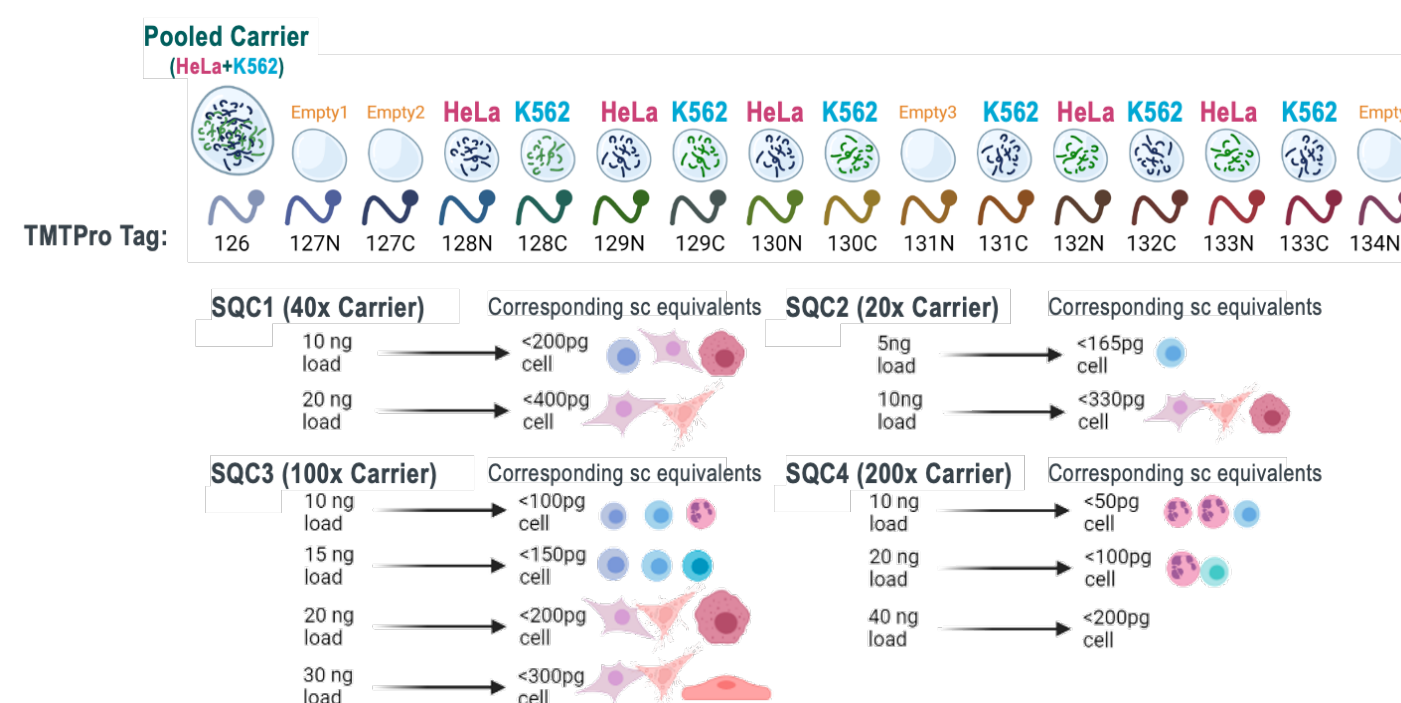
Enabling phi-SDM processing over the TMTProC region on the same Q Exactive HF-X instrument using the same transient acquisition settings as above (240K), allows us to completely resolve all 12 TMTProC species. Estimates of the empirical resolution at the peaks subjected to phi-SDM are >600K, while the resolution at the neighboring peaks subjected to eFT is ~160K

Sufficient Peak Calling with PhiSDM of a 256ms Transient



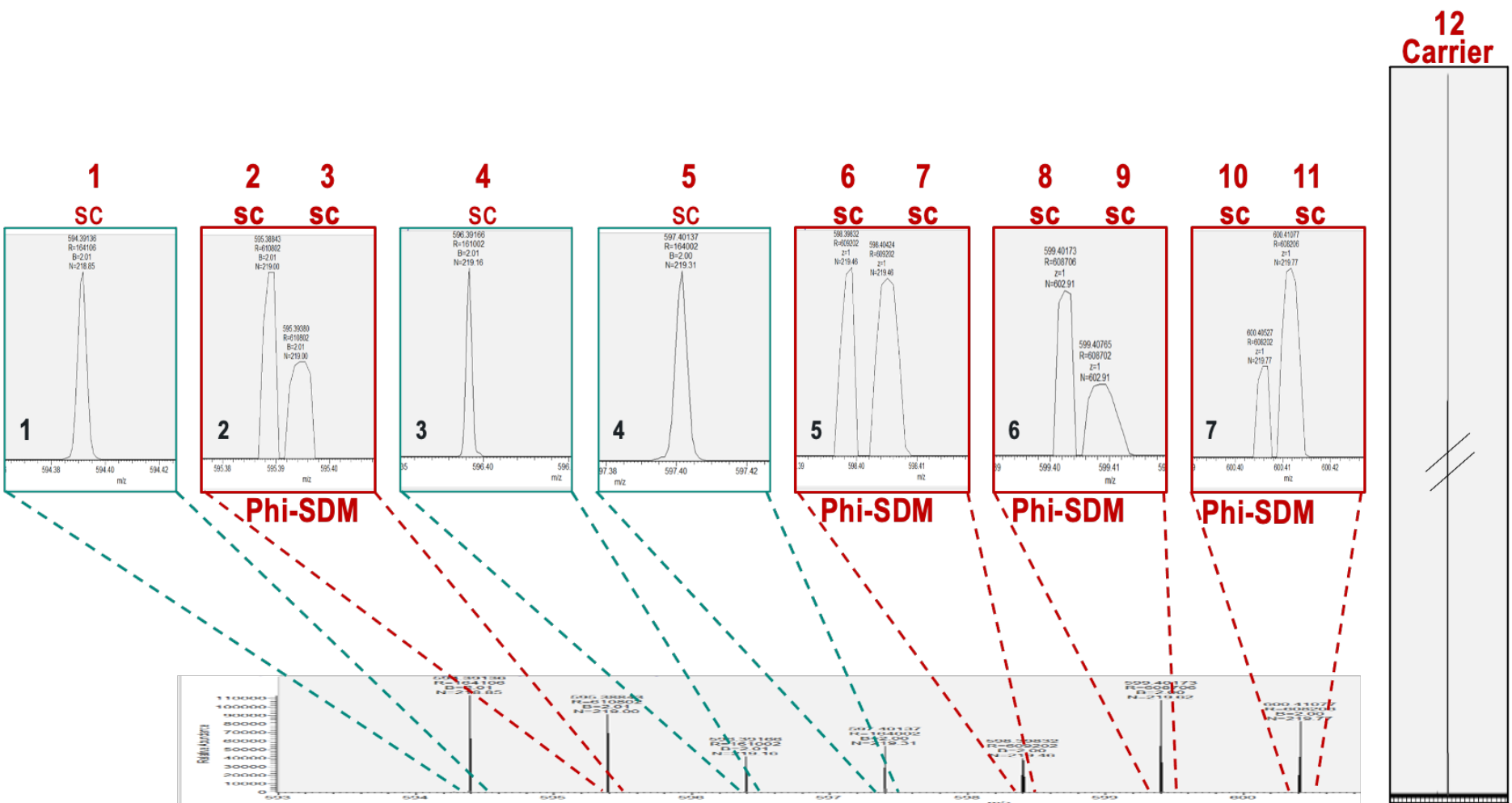
Reducing the transient length by half (nominally a 120K setting) and performing phi-SDM over the TMTProC regions as above, yields incompletely resolved peaks with partial splitting, at least as rendered by Xcalibur. However, zooming into these peaks shows that the phi-SDM algorithm was indeed able to identify and call out both of the peaks in each case.

Design of SCoPE-style Single Cell Standards



A set of SCoPE-style Single-Cell Proteomics QC standards (SQC) was designed to contain alternating peptides from 2 different cell types (HeLa and K562) in the single-cell TMT channels, along with a pooled carrier from the 2 cell types, and empty channels (to assess carryover). Different variants of the standards (SQC1-SQC4) were formulated to contain different ratios of carrier peptides to single cell channel peptides, from 20x to 200x, such that with a 10ng SQC sample load we could model the proteomic yield of different sizes of cells in the single cell channels.

Resolution and Quantification of Single Cell STD TMTProC Ions



We are able to resolve and quantify each of the 12 TMTProC ions from our Single-Cell Proteomic SQC samples using phi-SDM processing across the TMTProC region on an Exploris 480 instrument. Notably the 126-labeled carrier (Peak 12) is well separated from the lower-mass single-cell channel signals and does not interfere with them. Therefore, with this design, 10 single cells plus 1 empty control well could be incorporated into each multiplexed Single-Cell Proteomics set.

Summary

Here we demonstrate that Phi-SDM applied across mass ranges spanning TMTc ions present in the tandem mass spectra of TMTPro-labeled peptides and single-cell standards deliver effective resolution values across these regions that surpass any achievable on current commercially available Orbitrap platforms. Moreover, we demonstrate baseline resolution of the nearly isobaric ¹⁵N and ¹³C TMTc species (delta mass ca. 6.3 mDa) at transient lengths of ca. 512ms and 256ms, corresponding to the nominal 240K and 120K eFT resolution settings, which synchronize well with maximum ion-fill times commonly used during single-cell analyses. This novel application of Phi-SDM beyond TurboTMT is a demonstration of the potential of this super-resolution method and brings us closer to an operational workflow for TMTPro complement ion-based quantification in highly-multiplexed ultrasensitive and single-cell proteomics.