

Enhanced Detection of Hemoglobin Variants in Clinical Research by DBS and High-Resolution Accurate Mass (HRAM) MS

Yvonne E. Song, Kerry Hassell, Ed Goucher, Jingshu Guo, and Tanis Correa
Thermo Fisher Scientific, 355 River Oaks Parkway, San Jose, CA, 95134

Abstract

Purpose: to develop the top-down approach for hemoglobin variant detection in clinical research using dried blood spots and Orbitrap Exploris 240 mass spectrometer.

Methods: A 3.2 mm disc was punched and put into a 96-well plate for direct protein extraction and precipitation. LC-MS analysis was performed by Thermo Scientific™ Vanquish™ Flex UHPLC system with Thermo Scientific™ MAbPac™ RP column (1 x 100 mm, 4 μm, Part No. 303183) coupled to Thermo Scientific™ Orbitrap Exploris™ 240 mass spectrometer with BioPharma Option installed. Sequence search was performed using ProSighPD™ with a modified database and quantitation was performed using Thermo Scientific™ TraceFinder™ software 5.1.

Results: Identification of variants applied top-down analysis to characterize the structural mutation while quantitation was performed by extracting isotopic m/z of precursors to quantify ratios between target chains or variants. Minimal sample preparation (< 1 hr) in 96 well plate format, easily transferrable to automation using a robotic system. Confident data processing and review was successfully developed using automated software tools, ProSightPD™ and TraceFinder™.

Introduction

Hemoglobinopathies and thalassemia are the most common genetically determined disorders. They are caused by pathogenic variants in genes that control the production of hemoglobin subunits. Common techniques used in Hb analysis are electrophoretic and chromatographic assays. However, there are challenges to differentiate isomers (i.e., between Hb C and Hb E), very similar masses (i.e., < 1 Da difference between normal beta and Hb C, Hb E, and Hb D-Punjab), or new variants due to their lack of resolution. Recently, top-down approach has emerged to target intact protein biomarkers owing to high-resolution accurate-mass (HRAM) mass spectrometry such as Orbitrap. Here, we present the top-down analysis for enhanced detection of various hemoglobin variants using Orbitrap Exploris mass spectrometry for clinical research.

Materials and methods

Sample Preparation

Normal hemoglobin (Sigma-Aldrich, P/N H7379) and hemoglobin S (HbS, Sigma-Aldrich, P/N H0392) were used. Different concentrations of HbS calibrator stock were spiked into 2-EDTA stabilized normal human whole blood (BioIVT, Westbury, NY), which generated final concentrations ranging from 0, 0.50, 1.0, 2.5, 5.0, 10, 20, 50, and 100 mg/mL. A 3.2 mm disc was punched and put into a 96-well plate for direct protein extraction and precipitation as shown in Figure 1.

Test Method(s)

LC-MS analysis was performed by Thermo Scientific™ Vanquish™ Flex UHPLC system with Thermo Scientific™ MAbPac™ RP column (1 x 100 mm, 4 μm, Part No. 303183) coupled to Thermo Scientific™ Orbitrap Exploris™ 240 mass spectrometer with BioPharma Option installed. Conditions are described in Table 1.

Data Analysis

Sequence search was performed using ProSighPD™ with a modified database. The modified database manager and workflow are shown in Figure 2. Protein Quantitation was performed using TraceFinder™ software 5.1.

Figure 1. Experimental workflow

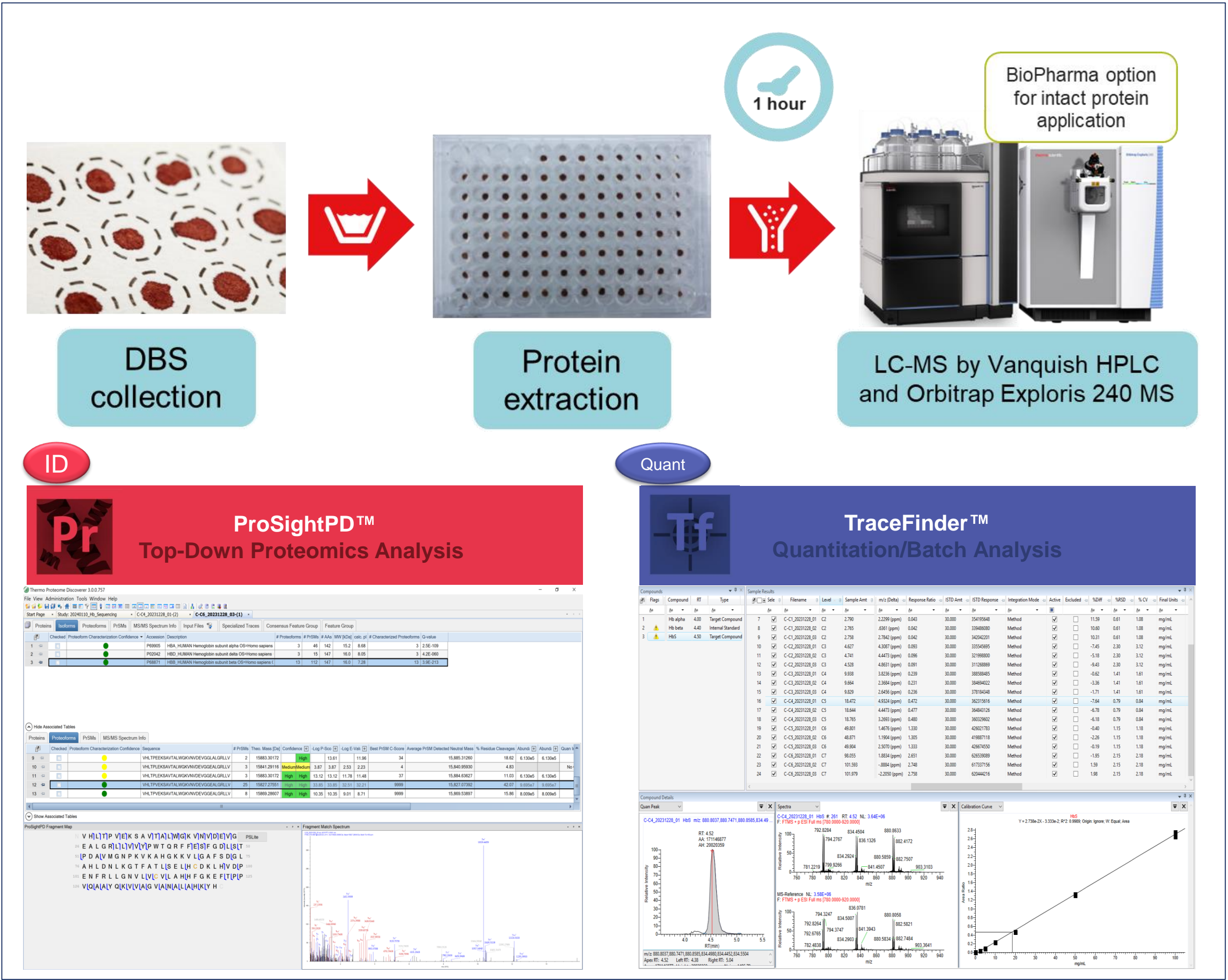


Table 1A. LC conditions

LC gradient			
Time (min)	% A	% B	Curve
0.0	75	25	5
0.5	75	25	5
0.7	68	32	5
8.5	57	43	5
9.2	10	90	5
9.7	10	90	5
9.8	75	25	5
12	75	25	5
Separation conditions			
Mobile phase A	0.1 % formic acid+ 0.02 % trifluoroacetic acid in water		
Mobile phase B	0.1 % formic acid + 0.02 % trifluoroacetic acid in 10: 20: 70 water: isopropanol: acetonitrile (v/v/v)		
Flow rate	0.12 mL/min		
Column temperature	80 °C		
Injection volume	2 μL		

Table 1B. MS conditions

MS global parameters	
Application mode	Intact protein
Pressure mode	Low pressure
Source type	Heated electrospray ionization (H-ESI)
Polarity	Positive
Spray voltage (V)	3800
Sheath gas (Arb)	35
Aux gas (Arb)	7
Sweep gas (Arb)	1
Ion Transfer tube temp (°C)	325
Vaporizer temp (°C)	150
Divert Valve A	0.0 min: position 1-6 (waste) 0.5 min: position 1-2 (MS) 9.5 min: position 1-6 (waste)
MS scan parameters	
Orbitrap resolution	120,000
Scan range (m/z)	500-3000 for intact mass confirmation 780-920 for targeted identification and quantitation
RF lens (%)	60
AGC target	Standard
Maximum injection time mode	Custom
Maximum injection time (ms)	250
Microscans	3
Intensity	1.00E+04
Charge state properties	18
ddMS2 settings	
Isolation window (m/z)	2.5
HCD collision energies (V)	20, 22, 26, 28, 30
Orbitrap resolution	90000
Scan range (m/z)	200-2000
AGC target	Custom
Normalized AGC target(%)	100
Maximum injection time (ms)	250

Results

Figure 2. ProSightPD workflow and database manager

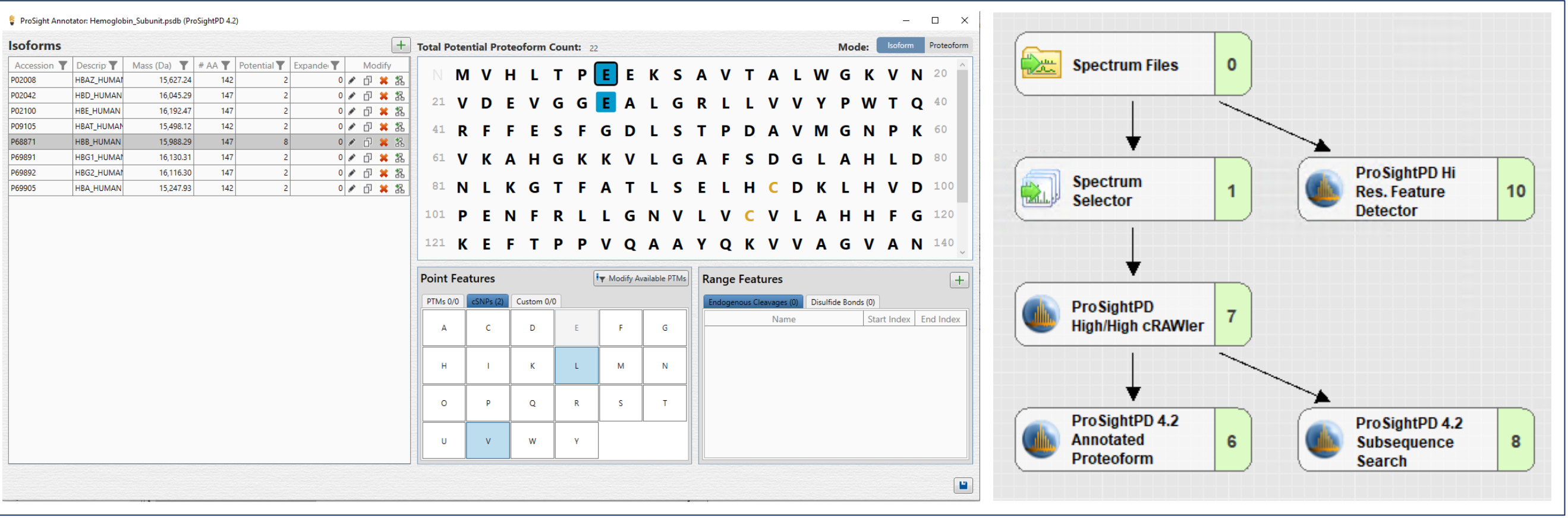


Figure 3A. Chain/variant identification with ProSightPD

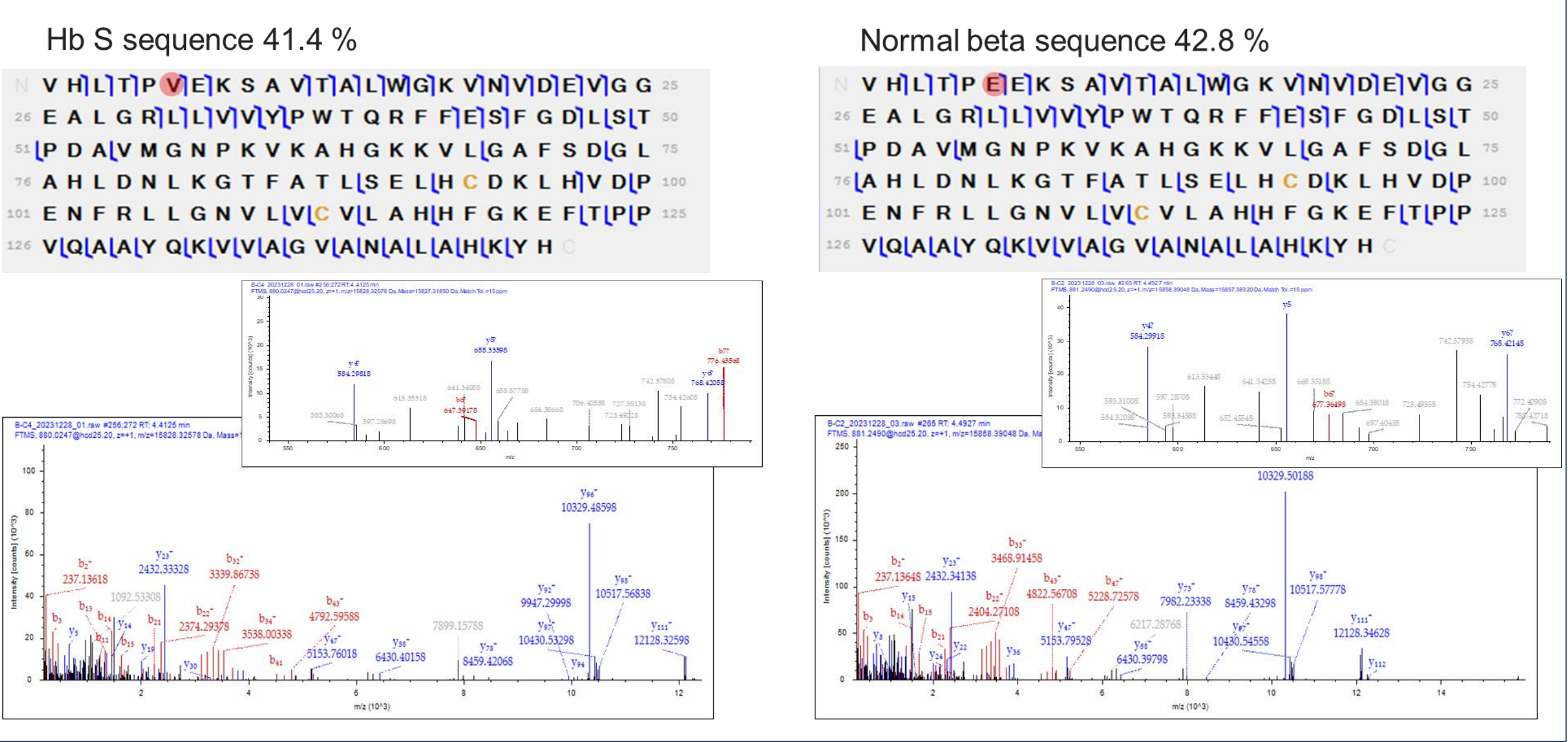


Figure 3B. Chain/variant identification with ProSightPD

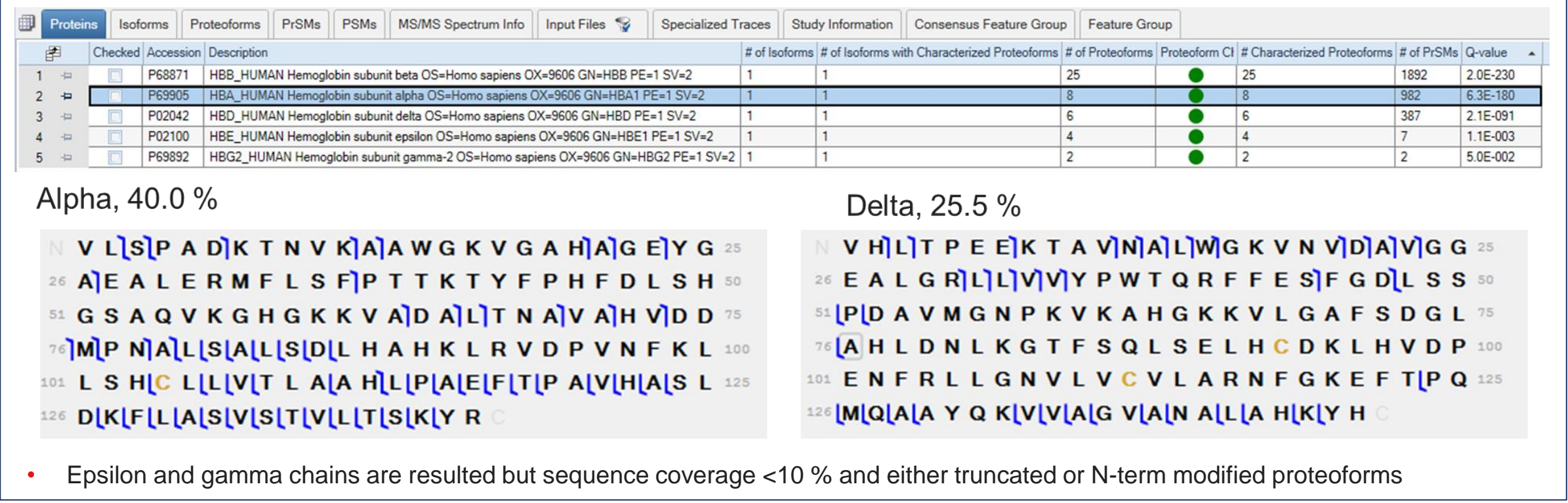


Figure 4. MS spectrum and chromatograms of 4 chains/variant

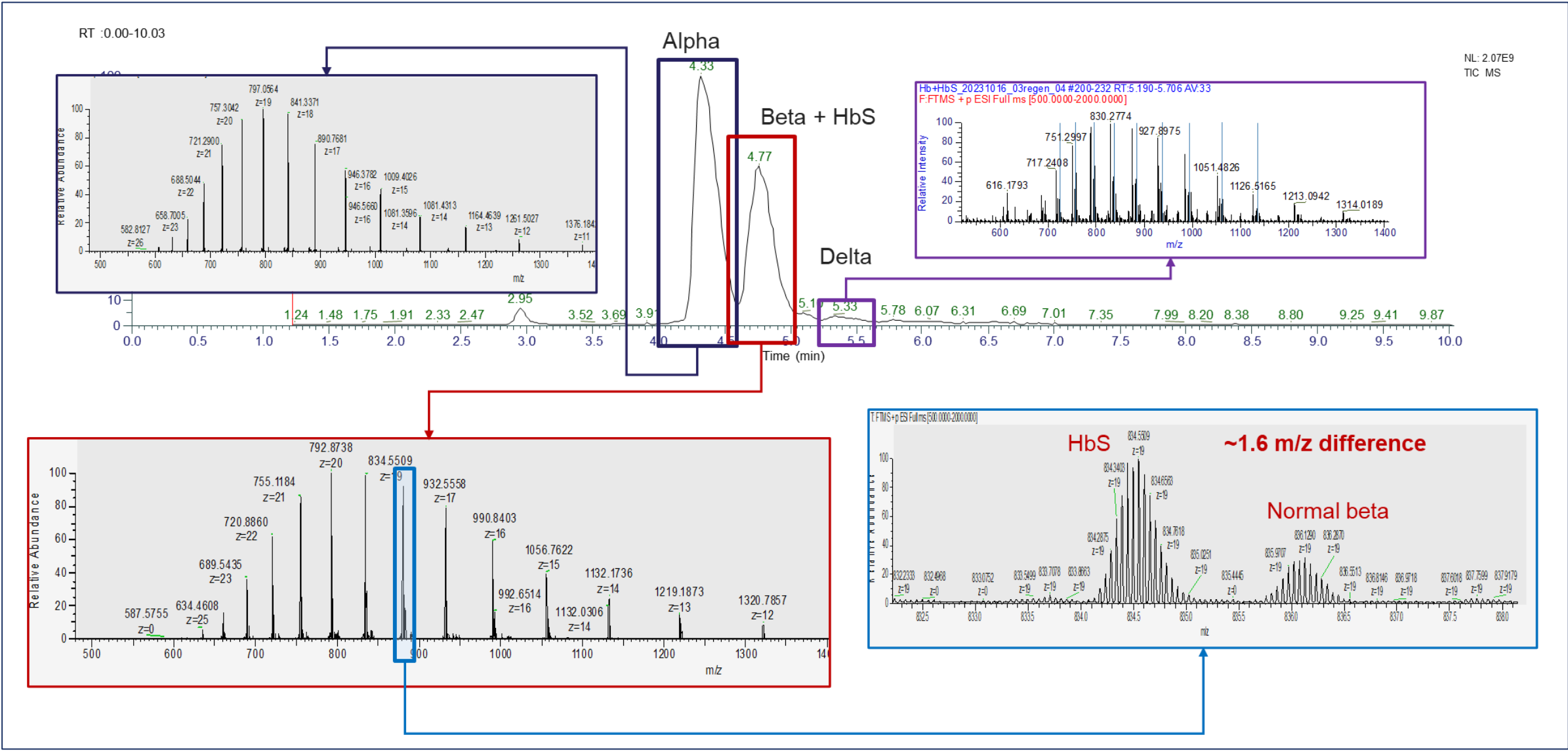
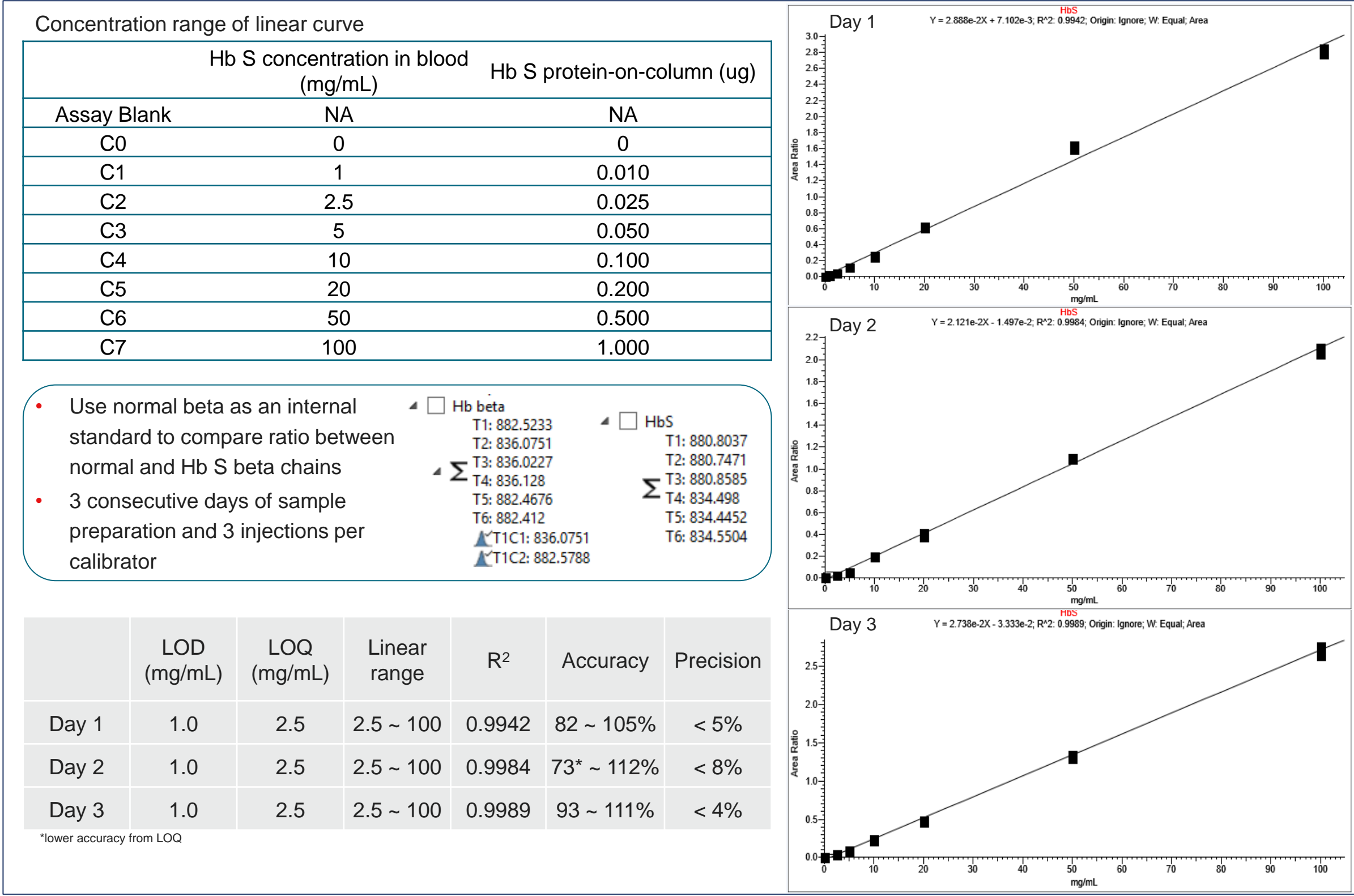


Figure 5. Quantitation strategy and summary



Conclusions

- Direct protein extraction from dried blood spots for protein identification and quantitation
- Identification: using top-down analysis for variant sequencing to characterize the structural mutation
- Quantitation: using intact protein quantitation by extracting isotopic m/z of precursors to quantify ratios between target chains or variants
- Minimal sample preparation (< 1 hr) in 96 well plate format, easily transferrable to automation using a robotic system
- Confident data processing and review using automated software tools, ProSightPD and TraceFinder

Trademarks/licensing

© 2024 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 others.