

Mass Spectrometry Immunoassay Coupled with Peptide Enrichment to Detect Thyroglobulin by Capillary Flow LC/MS/MS in Clinical Research

Kerry M. Hassell ¹, Adrian Sanchez-Woehler ². ¹Thermo Fisher Scientific, Somerset, NJ ²Thermo Fisher Scientific, Bannockburn, IL

ABSTRACT

Purpose: To develop an assay to capture Thyroglobulin in biological matrix and quantitate.

Methods: Using immuno-capture sample preparation and capillary chromatography detect Thyroglobulin with a highly sensitive mass spectrometer.

Results: Thyroglobulin was quantitated down to 0.1 ng/mL in plasma with analytical reproducibility, accuracy and precision.

INTRODUCTION

As one of the most common endocrine cancers, thyroid cancers the measurement of serum thyroglobulin (Tg) is important for diagnostics and follow-up treatment. Immunoassays have been the standard technique for quantitation, however a high amount of false negative results occur. Anti-Tg auto antibodies found endogenously block the binding epitope leading to these results. A user friendly, automated technique with low dead volume and improved performance for washes over beads was utilized to determine if this method would have utility for clinical research. Sample preparation techniques of using mass spectrometric immunoassay (MSIA) with stable isotope standards and capture by anti-peptide antibodies (SISCAPA) in conjunction with capillary flow LC/MS on a triple quadrupole provides high analytical selectivity and specificity at low levels of Tg found in serum.

MATERIALS AND METHODS

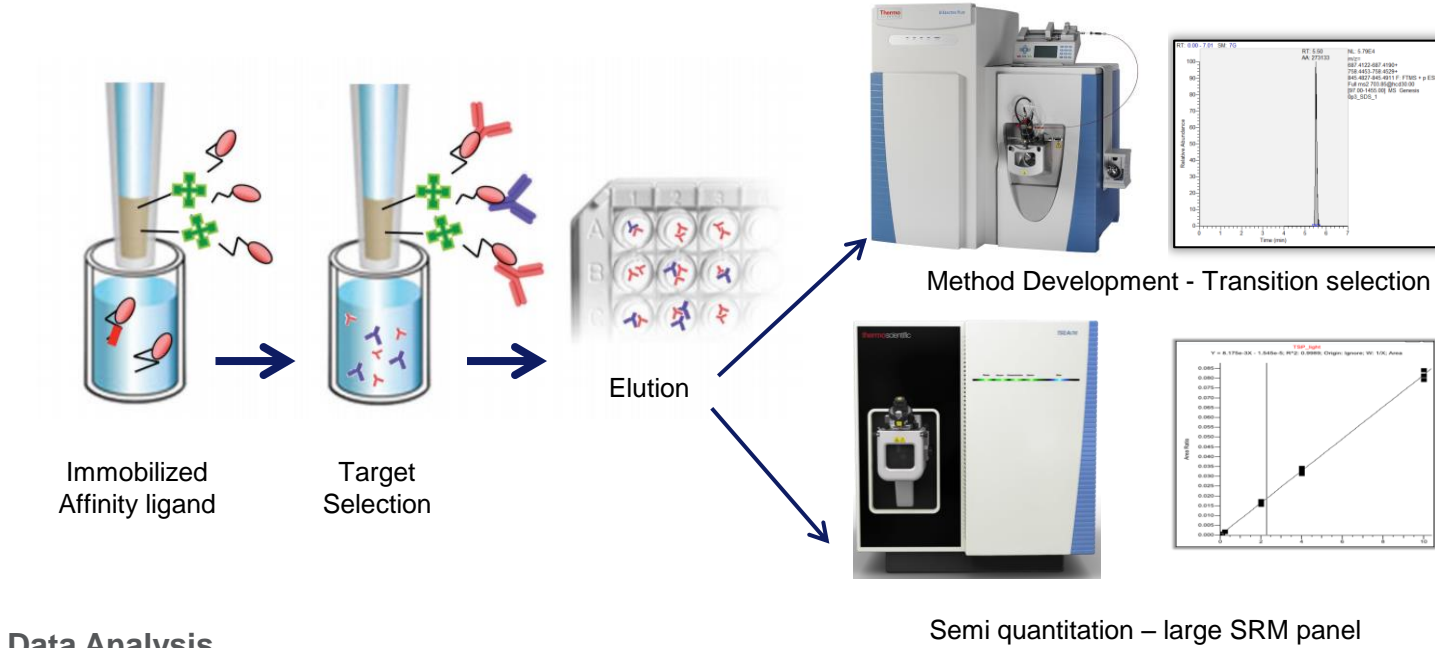
Sample Preparation

Human thyroglobulin was spiked into human heparin plasma and reduced and alkylated, followed by digestion with trypsin. Immuno-capture was then executed with a Thermo Scientific™ Finnpipe™ Novus i pipeteter with Thermo Scientific™ MSIA™ D.A.R.T.™S™ derivitized with anti-TG FSP antibody (SISCAPA). Elution of the peptide of interest was placed into a 96 well plate and injected into the LC/MS. Workflow of this process is showed in Figure 1.

LC/MS Parameters

Method development was performed on a Thermo Scientific™ Dionex™ UltiMate™ 3000 LC and Thermo Scientific™ Q Exactive™ mass spectrometer. The chromatography was performed with a Thermo Scientific™ ProSwift™ RP4H, 500um x 10cm column and run with a 20 minute long method. The quantitation experiments will be verified with a Thermo Scientific™ UltiMate™ 3000 RSLCnano™ LC system equipped with a capillary flow selector. Chromatographic separation was performed using a 150 um x 15 cm Thermo Scientific™ Acclaim™ PepMap™ column packed with 2.2 um C18 using a 30 minute gradient. LC/MS parameters are found in Figure 2.

Figure 1. Workflow of MSIA immuno-capture of Tg peptide and detection through LC/MS.



Data Analysis

Quantitation data was processed in Thermo Scientific™ TraceFinder™ software and retention times were scheduled using Skyline software.

Figure 2. LC/MS parameters for quantitation on the Thermo Scientific™ TSQ Altis™ triple quadrupole MS, a) gradient from liquid chromatography and b) SRM parameters for FSP peptide.

| a) | No. | Time (min) | Flow Rate (µL/min) | Temp (°C) | Column |
|----|-----|------------|--------------------|-----------|--------|
| | | | | | |
| A | 1 | 0.00 | 0.00 | 10 | 1 |
| | 2 | 0.00 | 0.00 | 10 | 1 |
| | 3 | 0.00 | 0.00 | 10 | 1 |
| | 4 | 0.00 | 0.00 | 10 | 1 |
| | 5 | 0.00 | 0.00 | 10 | 1 |
| B | 1 | 0.00 | 0.00 | 10 | 1 |
| | 2 | 0.00 | 0.00 | 10 | 1 |
| | 3 | 0.00 | 0.00 | 10 | 1 |
| | 4 | 0.00 | 0.00 | 10 | 1 |
| | 5 | 0.00 | 0.00 | 10 | 1 |

RESULTS

Method Development

Sample conditions and liquid chromatography parameters were developed on the Q Exactive MS platform and UltiMate 3000 RSLC nano system. Transitions of FSP peptide are shown in the high resolution data in Figure 3. In order to get the peptide to stick to the column, the gradient had to begin at 5% mobile phase B. Chromatography of the peptide of interest can be seen comparing to retention time standard PRTC in Figure 4. In addition, in order for the peptide to elute off the MSIA tips higher percent organic was needed, but good peak shape and retention was achieved with 0.4% TFA and 20% ACN as a reconstitution solvent. In this solution conditions a higher injection can be used without fronting of the chromatography, demonstration of this is shown in Figure 5.

Figure 3. XIC of FSP peptide on the Q Exactive MS and the bottom panel is the MS/MS spectrum.

