

# Profiling and Characterization of Polyphenol Polymers from Cinnamon Using an Ion Trap Mass Spectrometer

Min He<sup>1</sup>, Peter Wang<sup>2</sup>, Ying Xiang<sup>3</sup>, Ying Qi<sup>3</sup>, Howard Sun<sup>3</sup>, and Julian Phillips<sup>1</sup>

<sup>1</sup>Thermo Fisher Scientific, San Jose, CA, USA; <sup>2</sup>Thermo Fisher Scientific, Shanghai, China;

<sup>3</sup>Shanghai R&D, Nu Sin Enterprises, Shanghai, China

**Goal:** To develop a rapid and rugged LC/MS<sup>n</sup> method for high throughput characterization of the active components in cinnamon extract.

## Introduction

Type 2 diabetes is the most common metabolic disease worldwide. Although the prevention and control of it remains unclear, there is strong evidence that dietary factors play a role in the onset of this disease. Previous studies have indicated that cinnamon may mimic insulin effects and thus may improve glucose uptake.<sup>1-2</sup> Considering the complex composition of cinnamon, it is important to be able to isolate and characterize the possible active components in it. A range of reports have been devoted to the studies of this topic. However, most work has been mainly focused on the studies of monomer and dimers. The study of trimer or even larger compounds using on-line LC/MS<sup>n</sup> approach has not been reported.

The fast cycle time and excellent MS<sup>n</sup> sensitivity of the LCQ Fleet ion trap mass spectrometer provide enhanced analytical throughput for complex natural product analysis. In this report, profiling and structural elucidation of polyphenol polymers, the possible active compounds in cinnamon, are carried out using an LC/MS<sup>n</sup> methodology using the LCQ Fleet ion trap mass spectrometer. The application of pulsed-Q dissociation (PQD<sup>TM</sup>) and high resolution isolation (HRI) facilitates the analysis of this complex sample.

## Methods

An aliquot (0.5 g) of cinnamon was ground and extracted in 8 mL 0.1 N acetic acid using a sonic bath for 3 minutes. The supernatant liquid was cooled to room temperature and diluted in mobile phase. The final solution was filtered prior to HPLC separation. LC separation was performed using a C18 column and detected at 280 nm wavelength. Sample fractions CB4 with insulin enhancing activity were collected, and the powders were dissolved in water:MeOH: ACN = 25:50:25 solvent for LC-MS<sup>n</sup> analysis using the LCQ Fleet ion trap mass spectrometer.

LC instrument: Surveyor Plus LC system

Column: Hypersil GOLD (50 × 2.1 mm, 3 μm particle size, Thermo Fisher Scientific)

Mobil phase: A: Water with 0.1% formic acid

B: Acetonitrile with 0.1% formic acid

Flow rate: 400 μL/min

t (min)	A%	B%
0.00	95	5
0.20	95	5
3.20	65	36
3.30	65	35
3.31	95	5
4.00	95	5

**Mass Spectrometer:** The LCQ Fleet ion trap mass spectrometer was operated in positive electrospray mode. The electrospray voltage was 5 kV. The capillary temperature was 200 °C and the sheath gas flow was 30 units. An isolation width of 2 Da was used with a 30 msec activation time for MS<sup>n</sup> experiments. All scan events were acquired with a 200 ms maximum ionization time.



Figure 1: Picture of the LCQ Fleet Ion Trap Mass Spectrometer with Surveyor Plus LC System

## Key Words

- LCQ Fleet<sup>TM</sup>
- Surveyor Plus<sup>TM</sup>
- Hypersil GOLD<sup>TM</sup> Column
- MS<sup>n</sup>
- Mass Frontier<sup>TM</sup>
- Natural Product Analysis

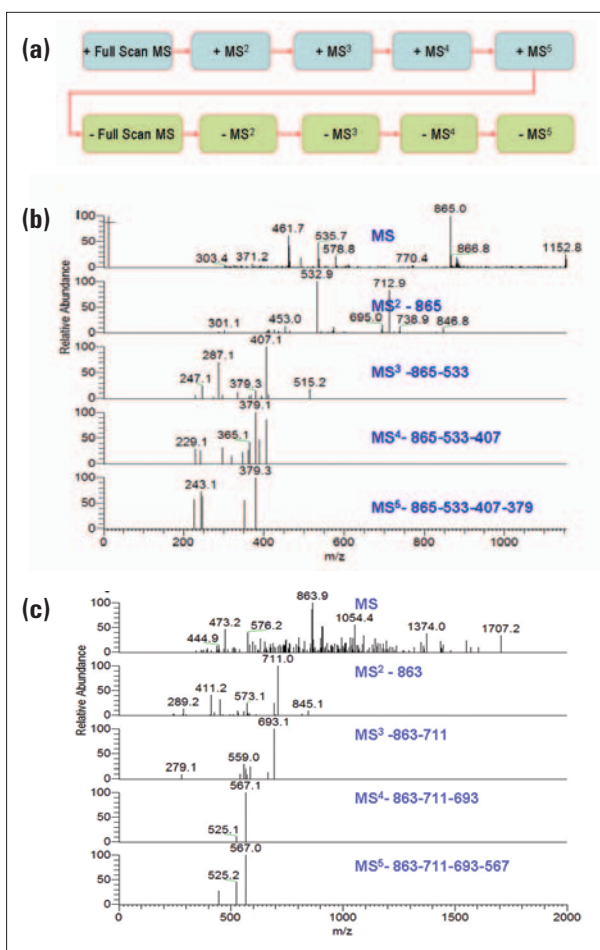


Figure 2: LC-MS<sup>n</sup> Analysis of doubly linked A-type Procyanidin from Cinnamon. (a) Ten-scan-event MS<sup>n</sup> method setup: MS<sup>n</sup> was performed on the most abundant ions from the previous level of analysis; (b) MS to up to MS<sup>5</sup> data for the target compound with  $m/z$  865 in positive polarity; (c) MS to up to MS<sup>5</sup> data for the target compound with  $m/z$  863 in negative polarity. (The data was collected from fraction CB4)

## Results

It has been reported that a group of polyphenolic polymers found in cinnamon may function as antioxidants to potentiate insulin action, and therefore, may be beneficial in the control of glucose intolerance and diabetes. Two major types of polymeric procyanidins have been observed including A-type and B-type linkages. Among them, A-type has two interflavan linkages (IFL) while B-type has a single IFL, which results in a 2 amu unit difference. Studies have been performed on both

types of moneric, dimeric and trimeric procyanidins. However, detailed structural elucidation results for higher orders of polymerization are still rare, particularly using an on-line LC/MS<sup>n</sup> approach.<sup>1,3</sup> In this work, a range of procyanidins from cinnamon extract fractions have been identified and characterized within a rapid ten-scan-event LC-MS<sup>n</sup> analysis (Figure 2a). The fast positive/negative switching the capability of the LCQ Fleet mass spectrometer ensured the comprehensive information collection from the samples. Figure 2b and Figure 2c demonstrated the MS<sup>n</sup> (up to MS<sup>5</sup>) data of one trimer procyanidin in both positive (Figure 2b) and negative (Figure 2c) polarities from fraction CB4. The ability to perform MS<sup>n</sup> analysis on precursor ions (isolated using HRI) using the ion trap enables the structural elucidation of these compounds with less interference from the matrices. The presence of  $m/z$  289 in negative MS/MS spectrum is considered to be specific for a C-O-C IFL in the polymers.<sup>4,5</sup>

Mass Frontier software contains a number of tools for structural identification. Its fragment prediction module is very useful for accurate structural characterization from MS<sup>n</sup> data. As an example, the positive MS/MS spectrum of type A procyanidin trimer was processed with Mass Frontier, and all the major fragments were assigned with a predicted sub-structure (Figure 3). This feature can be further coupled with the database searching of Mass Frontier for compound identification.

Mass Frontier also predicted that the ion at  $m/z$  287 could be generated from singly linked B-type procyanidin oligomers, which is consistent with the observations in this work. The fragmentation of the ion at  $m/z$  867 resembles that of  $m/z$  865 with  $m/z$  287 as a common fragment ion (data not shown). Therefore, the ion at  $m/z$  287 may not be a good indicator ion for the presence of A-type procyanidin oligomers, as noted in a previous report.<sup>1</sup>

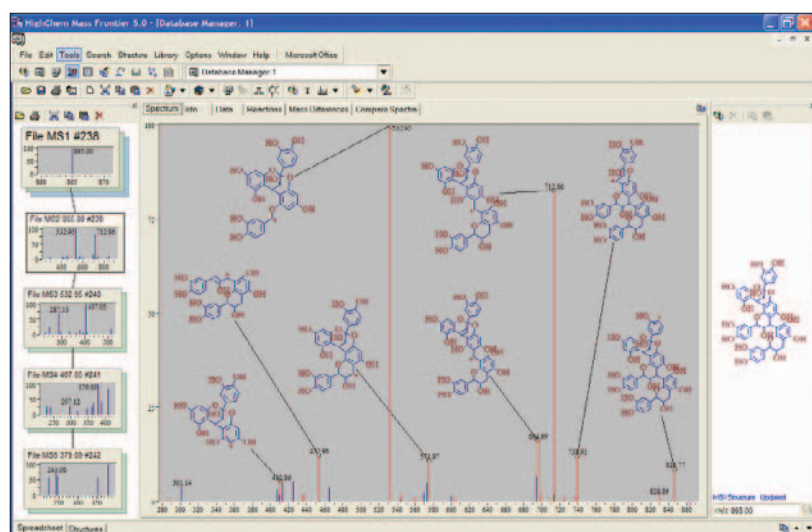


Figure 3: Fragmentation prediction for the identification of A-type procyanidin trimer using Mass Frontier

Dimers and tetramers have been found, in addition to the trimers, in this fraction with up to MS<sup>5</sup> data for identification purpose in both polarities. With the traditional CID method, the *m/z* 289 from MS/MS spectrum of *m/z* 1151 (the tetramer) in negative polarity is not observable in an ion trap mass spectrometer due to the low mass cutoff (Figure 4a). In order to observe this indicator ion, an MS<sup>3</sup> analysis is usually required (see insert in Figure 4a), which may be impacted by both cycle time and signal strength. With the new dissociation method, Pulsed Q Dissociation (PQD), it is possible to access directly the low mass product ions in the ion trap without the need for the further step of MS<sup>n</sup> analysis. As demonstrated in Figure 4b, the ion at *m/z* 289 was observed directly from the MS/MS spectrum of the ion at *m/z* 1151.

## Conclusions

A rapid and rugged LC-MS<sup>n</sup> method has been developed to characterize the various procyanidin oligomers in the Chinese cinnamon extract. Both A-type and B-type polymers were found. The fast cycle time and excellent MS<sup>n</sup> capability of the LCQ Fleet enabled the confident identification of the components of interest. Comprehensive information was obtained in both positive and negative polarities. Use of PQD provided diagnostic low mass product ions without the need for higher order MS<sup>n</sup> experiments.

## References

- Anderson, R.A.; Broadhurst, C.L.; Polansky, M.M.; Schmidt, W.F.; Khan, A.; Flanagan, V.P.; Schoene, N.W.; Graves, D. *J. Agric. Food Chem.* 2004, 52, 65.
- Mang, B.; Wolters, M.; Schmitt, B.; Kelb, K.; Lichtinghagen, R.; Stichtenoth, D.O.; Hahn, A. *Euro. J. Clin. Invest.* 2006, 36, 340-344.
- Lampke, S.M.; Pavlovich, J.; Polansky, M.; Anderson, R.; Graves, D. Poster on Experimental Biology, 2006.
- Sun, W.; Miller, J.M. *J. Mass Spectrom.* 2003, 38, 438.
- Guyot, S.; Vercauteren, J.; Cheynier, V. *Phytochemistry*, 1996, 42, 5.

## Acknowledgements

We would like to thank Dr. James Pavlovich from the University of California, Santa Barbara for helpful technique discussions.

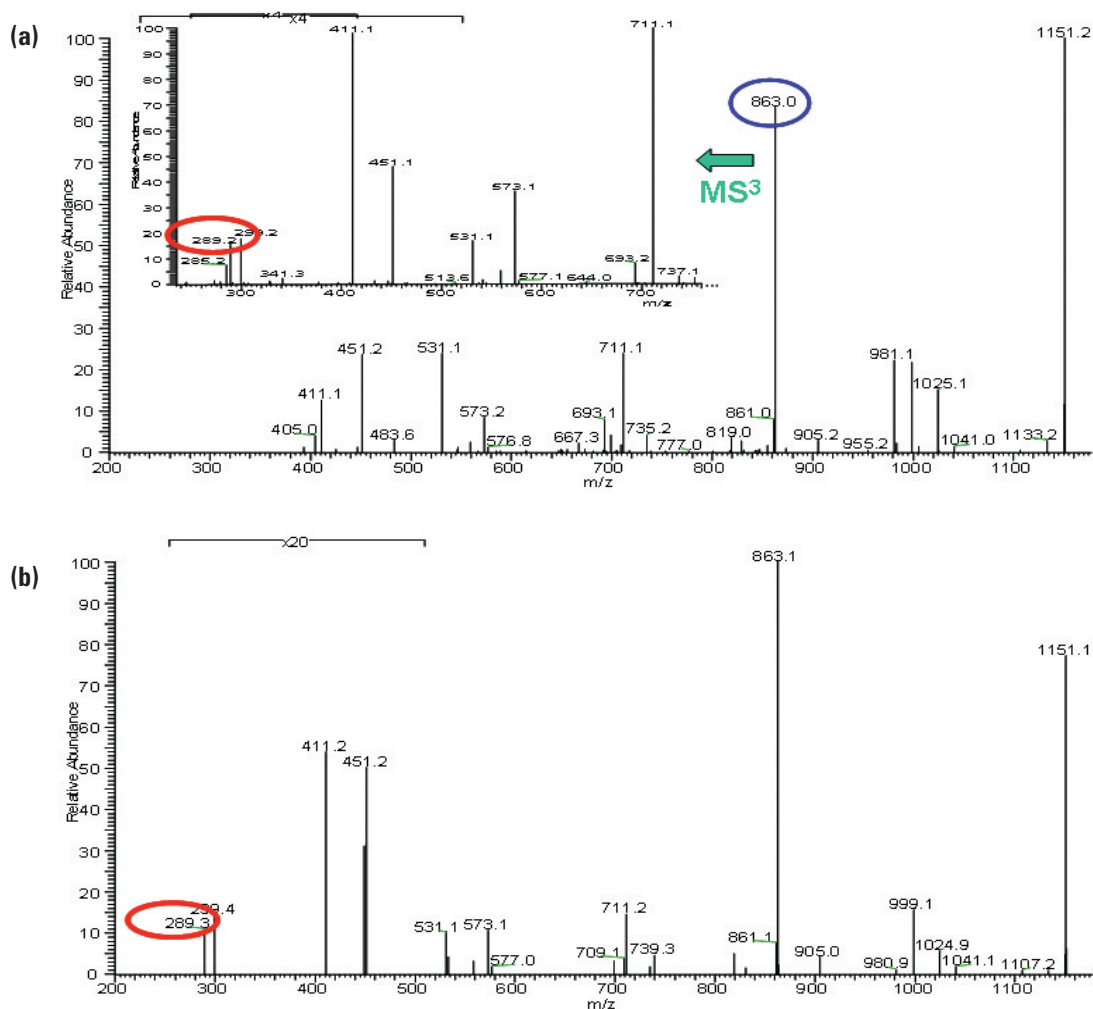


Figure 4: Negative product ion spectra of the tetramer ion (*m/z* 1151) (a) CID spectrum, the insert is the MS<sup>3</sup> data of *m/z* 863; (b) PQD spectrum.

## Laboratory Solutions Backed by Worldwide Service and Support

Tap our expertise throughout the life of your instrument. Thermo Scientific Services extends its support throughout our worldwide network of highly trained and certified engineers who are experts in laboratory technologies and applications. Put our team of experts to work for you in a range of disciplines – from system installation, training and technical support, to complete asset management and regulatory compliance consulting. Improve your productivity and lower the cost of instrument ownership through our product support services. Maximize uptime while eliminating the uncontrollable cost of unplanned maintenance and repairs. When it's time to enhance your system, we also offer certified parts and a range of accessories and consumables suited to your application.

To learn more about our products and comprehensive service offerings, visit us at [www.thermo.com](http://www.thermo.com).

In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

**Australia**  
+61 2 8844 9500  
**Austria**  
+43 1 333 50340  
**Belgium**  
+32 2 482 30 30  
**Canada**  
+1 800 532 4752  
**China**  
+86 10 5850 3588  
**Denmark**  
+45 70 23 62 60  
**France**  
+33 1 60 92 48 00  
**Germany**  
+49 6103 408 1014  
**India**  
+91 22 6742 9434  
**Italy**  
+39 02 950 591  
**Japan**  
+81 45 453 9100  
**Latin America**  
+1 608 276 5659  
**Netherlands**  
+31 76 587 98 88  
**South Africa**  
+27 11 570 1840  
**Spain**  
+34 91 657 4930  
**Sweden/Norway/  
Finland**  
+46 8 556 468 00  
**Switzerland**  
+41 61 48784 00  
**UK**  
+44 1442 233555  
**USA**  
+1 800 532 4752

[www.thermo.com](http://www.thermo.com)

### Legal Notices

©2007 Thermo Fisher Scientific Inc. All rights reserved. Mass Frontier is a trademark of HighChem, Ltd. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific Inc. products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.



Thermo Fisher Scientific,  
San Jose, CA USA is ISO Certified.

View additional Thermo Scientific LC/MS application notes at: [www.thermo.com/appnotes](http://www.thermo.com/appnotes)

AN62438\_E 08/07S