

# Analysis of Patulin in Fruit Juices and Extracts Using Liquid Chromatography Triple Quadrupole Mass Spectrometry

Claudia P. B. Martins, Cristina C. Jacob, Michael Volný, Thermo Fisher Scientific, 355 River Oaks Parkway, San Jose, CA, 95134, USA

## ABSTRACT

**Purpose:** Patulin is a mycotoxin produced by different types of fungi as their secondary metabolite. The regulatory authorities have imposed restrictions on maximum patulin levels in fruit products, which creates a need for sensitive analytical methods.

**Methods:** Two different LC-MS/MS methods were tested using Thermo Scientific™ TSQ Fortis™ Triple Quadrupole Mass Spectrometer and Thermo Scientific™ Vanquish™ UHPLC system. SPE was used to extract patulin from fruit juices.

**Results:** SPE LC-MS/MS method for quantification of patulin was developed and tested

## INTRODUCTION

Patulin, 4-hydroxy-4H-furo[3,2-c]pyran-2(6H)-one (CAS#149-29-1), is a polyketide produced as a mycotoxin by several fungi, namely *Aspergillus* and *Penicillium*. Fruits that were damaged or improperly stored are susceptible to the growth of patulin-producing molds. If these fruits are used to make further consumer products, patulin can be present as a toxic contaminant. Because patulin is a heat-stable lactone that resists thermal denaturation, the normal pasteurization treatment is not sufficient to decompose it. The regulators in different jurisdictions around the world have imposed restrictions on maximum patulin levels in different products, especially in apple juice. These levels differ depending on the product and the country, but predominantly were set in the concentration range between 5-100ppb (50ug/kg in the US and EU).

## MATERIALS AND METHODS

### Chemicals and Sample Preparation

Patulin was obtained from Sigma Aldrich.; all solvents and reagents were obtained from Fisher Scientific. For method development the samples were prepared as neat in 0.1% acetic acid in the concentration range 0.5-500ppb. To prepare matrix samples, apple juice was purchased in a local market and spiked with patulin.

### Mass Spectrometry

The TSQ Fortis Mass Spectrometer was used for all the examples described in this work.

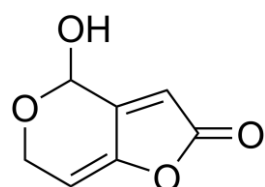
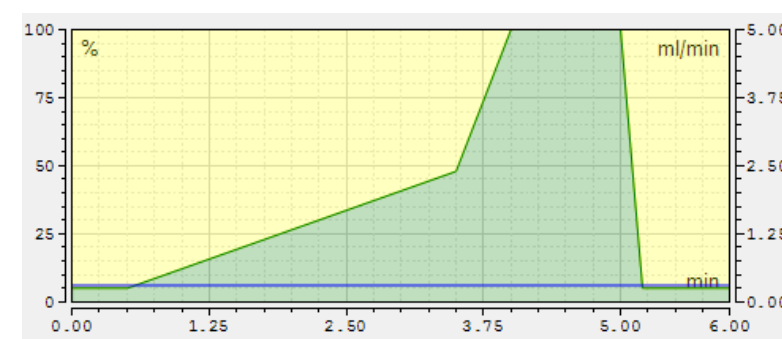


Table 1. Patulin SRM parameters

Precursor m/z	Product m/z	CE (V)	Tube lens (V)	
153.0	108.9	8	76	Quan
153.0	81.2	11	76	Confirm

### LC methodology



[A] water  
[B] acetonitrile  
Injection volume: 10uL  
Temperature: 30°C  
Column: HSS T3 1.8µm; 2.1x100mm



[A] 0.2mM NH<sub>4</sub>F in water  
[B] acetonitrile  
Injection volume: 10uL  
Temperature: 30°C  
Column: HSS T3 1.8µm; 2.1x100mm

### Solid-Phase Extraction

#### Format

**Sample pre-treatment**  
Condition  
Equilibration  
Sample load  
Wash 1  
Wash 2  
Elution  
**Post elution and reconstitution**

Oasis Max 6cc  
Spike apple juice at 1; 2.5; 5; 10; 50; 100 and 200 ppb (2 mL)  
6 mL of methanol  
6 mL water  
load 2 mL of spiked apple juice  
3 mL 5mM ammonium acetate  
3 mL water  
Eluate patulin with 4 mL Methanol  
The extracts were evaporated to dryness under N<sub>2</sub> stream and reconstituted in 1 mL 0.1% acetic acid

### Data Analysis

Thermo Scientific™ TraceFinder™ and Thermo Scientific™ FreeStyle™ software were used for data processing.

Table 2. Results overview for LC-MS/MS method #1 (left) and #2 (right); based on 3 replicate injections

Sample	Average diff. from theoretical value [%]	RSD [%]	Sample	Average diff. from theoretical value [%]	RSD [%]
0.5 ppb	6.1	11.93	0.5 ppb	9.3	14.8
1 ppb	13.8	3.65	1 ppb	6.0	3.9
2 ppb	6.5	7.29	2 ppb	3.1	1.7
5 ppb	13.8	11.8	5 ppb	3.7	4.8
10 ppb	6.5	2.7	10 ppb	3.4	2.3
20 ppb	9.7	1.1	20 ppb	3.3	1.2
50 ppb	5.9	4.0	50 ppb	3.5	3.9
100 ppb	0.8	1.5	100 ppb	4.8	1.1
200 ppb	2.6	0.7	200 ppb	1.5	2.0
500 ppb	4.0	0.9	500 ppb	1.4	2.0

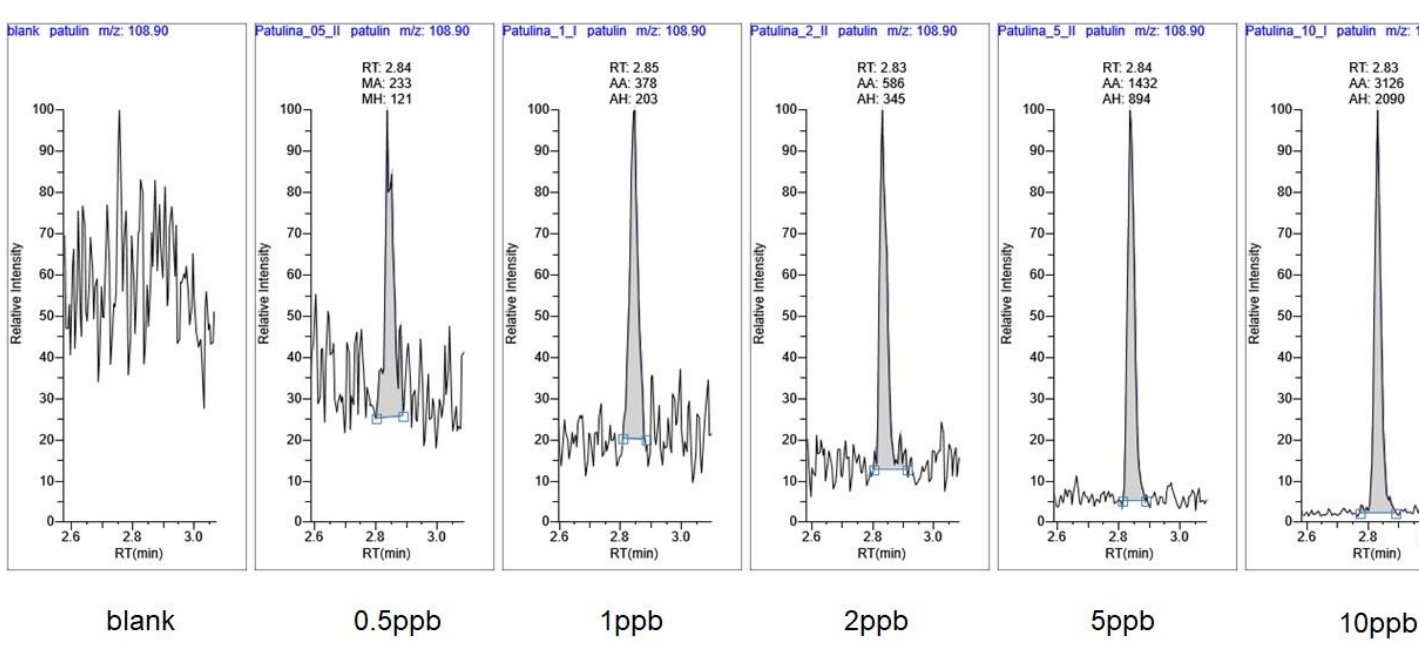


Figure 5 Extracted Ion chromatograms for lowest calibrators in neat solvent and solvent blank; LCMS method #1

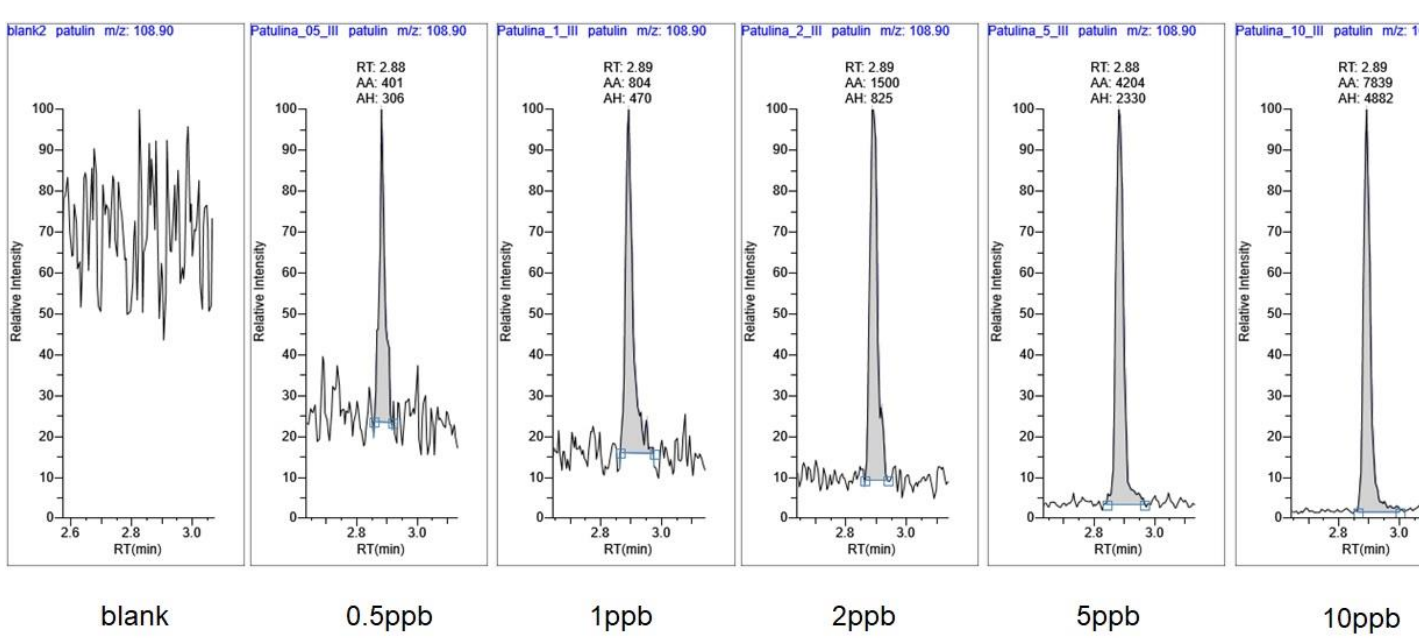


Figure 6 Extracted Ion chromatograms for lowest calibrators in neat solvent and solvent blank; LCMS method #2

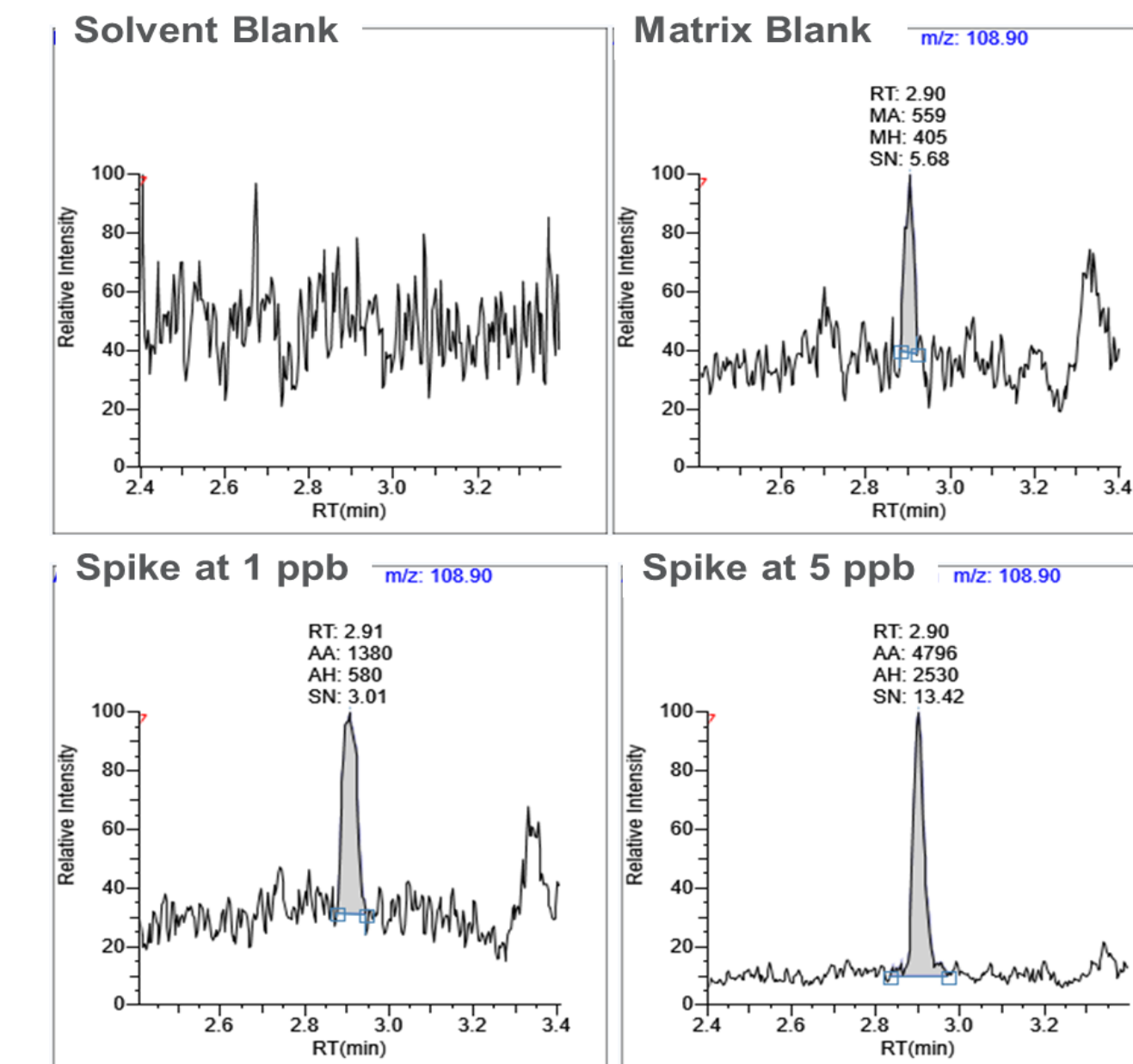


Figure 7 Patulin Extracted Ion Chromatograms in apple juice; LC-MS method #2 was used for apple juice matrix samples

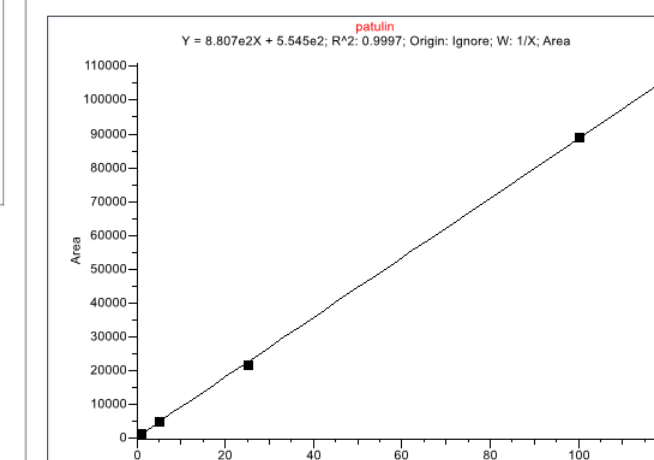


Figure 8 Patulin calibration curve in apple juice matrix.

Table 3. Overview of Patulin calibrators in apple juice

Calibrators in apple juice	Calculated Conc. (ppb)	CV %	Ion ratio
1 ppb	1.005	0.53%	36.38
5 ppb	5.11	2.17%	40.54
25 ppb	24.14	-3.45%	37.34
100 ppb	100.748	0.75%	37.49

Table 4. Sample preparation accuracy and precision \*

Apple juice Samples N=4	Average concentration /ppb	Precision %	Accuracy %
Spiked at 1 ppb	1.01	10%	101%
Spiked at 5 ppb	4.90	9%	98%
Spiked at 25 ppb	28.37	11%	113%
Spiked at 100 ppb	97.02	12%	97%

\*Isotopically labeled internal standard is commercially available for patulin, but was not used in this work

## CONCLUSIONS

- Method for LC-MS/MS quantitation of Patulin in negative ion mode was developed using TSQ Fortis mass spectrometer and Vanquish-Flex UHPLC system.
- Two different LC methods were evaluated, one with addition of NH<sub>4</sub>F as modifier
- TSQ Fortis showed linearity with R<sup>2</sup> >0.999 in the 0.5-500 ppb concentration range in both neat and apple juice matrix
- Method #2 demonstrated better sensitivity, but required longer column equilibration than simpler method #1, which was based on the Chinese regulation GB5009.185-2016 (ref 6)
- Calibrators were analyzed in triplicates, all RSD were <15%
- Both methods showed very good precision and accuracy even without using any internal standard
- The SPE method provided a recovery of 80-90% with a signal suppression of ca.14%, affording lower limit of quantitation (LLOQ) of 1 ppb for patulin in apple juice (mean accuracy 101%; CV=10%) using LC-MS method #2 (with NH<sub>4</sub>F in the mobile phase)



## REFERENCES

- Patulin in Apple Juice, Apple Juice Concentrates and Apple Juice Products, Compliance Policy Guidance, FDA, September 2011
- FDA Memorandum, "Hazards or Patulin in Apple Juice." May 5, 1994.
- Brain, P. W. (1956) Production of patulin in apple fruits by *Penicillium expansum*, Nature (London) 178: 263.
- Becci, P. J., Hess, F. G., Johnson, W. D., Gallo, M. A., Babish, J. G., Dailey, R. E., & Parent, R. A. (1981) Long-term carcinogenicity and toxicity studies of patulin in the rat. J. Appl Toxicol 1(5):256-261.
- GB 2761-2011 Food Safety National Standard for Maximum Levels of Mycotoxins in Foods, National Health and Family Planning of People's Republic of China (NFHPC), China, 2011
- GB5009.185 – 2016 National food safety standard Determination of patulin in in apple and hawthorn products, National Health and Family Planning of People's Republic of China (NFHPC), China 2016

## TRADEMARKS/LICENSING

© 2019 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.

PO65492-EN0419S

**ThermoFisher**  
SCIENTIFIC

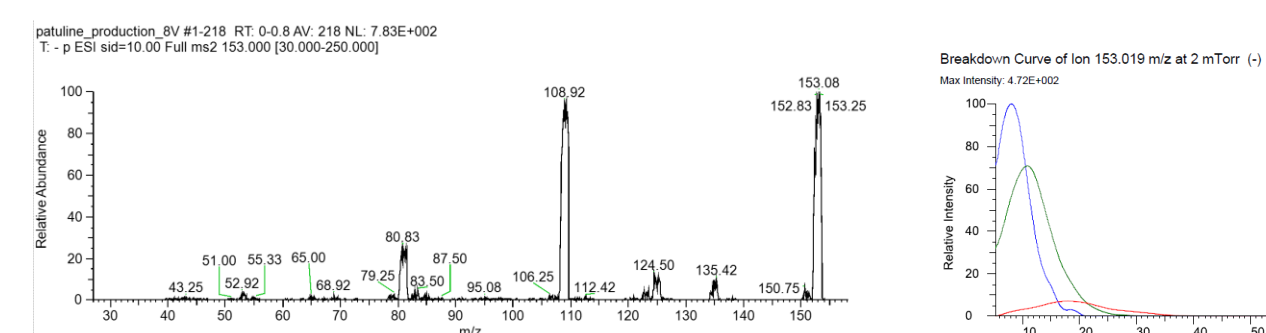


Figure 3 Product spectrum of patulin at 10V CID

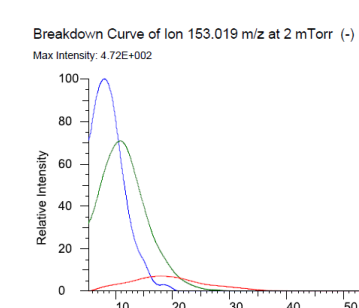


Figure 4 Patulin CID breakdown curves