

Analysis of Vomitoxin (DON) and T-2 Toxin by LC/MS

Chromatography and Mass Spectrometry Application Note

John H. Moncur, Thermo Electron Corporation

Introduction

Grain and grain products may be susceptible to a variety of fungal infections. *Fusarium* species fungi are known to produce a number of mycotoxins including Vomitoxin (DON) and T-2 Toxin. These mycotoxins may evolve pre-harvest in the field or post harvest during processing, storage and transportation. The presence of these mycotoxins can lead to human and animal toxicology and in the severest cases, death. The US FDA (United States Food and Drug Administration) has issued 'advisory' limits for Vomitoxin in grain and grain products as follows:

- 1 ppm—Finished wheat products for human consumption.
- 5 ppm—Grain and grain by-products destined for swine and other non ruminant species.
- 10 ppm—Grain and grain by-products for ruminating beef and feedlot cattle older than 4 months.

Additional restrictions on percentage make up of the diet are also advised.

Fusarium infections of grain products affect not only the vigor of the crop itself leading to lower yields, but also the quality of the crop as a feedstuff. The severity of effect of the mycotoxin contamination is to a large extent dependent on the age and species of the particular livestock. In particular, pigs are susceptible to even low levels of Vomitoxin in their feed.

Goal

This report describes a simple LC/MS method for the determination of fungal mycotoxins. It will allow analysts to develop and optimize sample preparation, extraction and clean-up techniques for these important contaminants. The flexibility of simultaneous SIM and scanning acquisition will facilitate the development of quantitative and qualitative methodologies.

Key Words

- Fungal Infections
- Pseudomolecular Ion Species
- LC/MS
- Finnigan™ Surveyor™ MSQ

Chromatographic Conditions

LC Column: 150×2.1 mm
Packing: 5 µm C18
Flow Rate: 200 µL/min
Mobile Phase: Water/Methanol
Gradient: 80/20 to 20/80 in 10 min,
held 5 min.
Injection Volume: 10 µL
Column Temp: 40 °C

Mass Spectrometer Conditions

Instrument: Surveyor MSQ
Ionization: Electrospray
Polarity: Positive
Needle Voltage: 3 keV
Probe Temp: 350 °C
Cone Voltage: 90 eV
Scan Range: 100 – 650 Da
Scan Rate: 0.8 seconds per scan
Data Acquisition: Centroid
Selected Ions: m/z 319 (M+23) and 489 (M+23)
Dwell Time: 100 ms

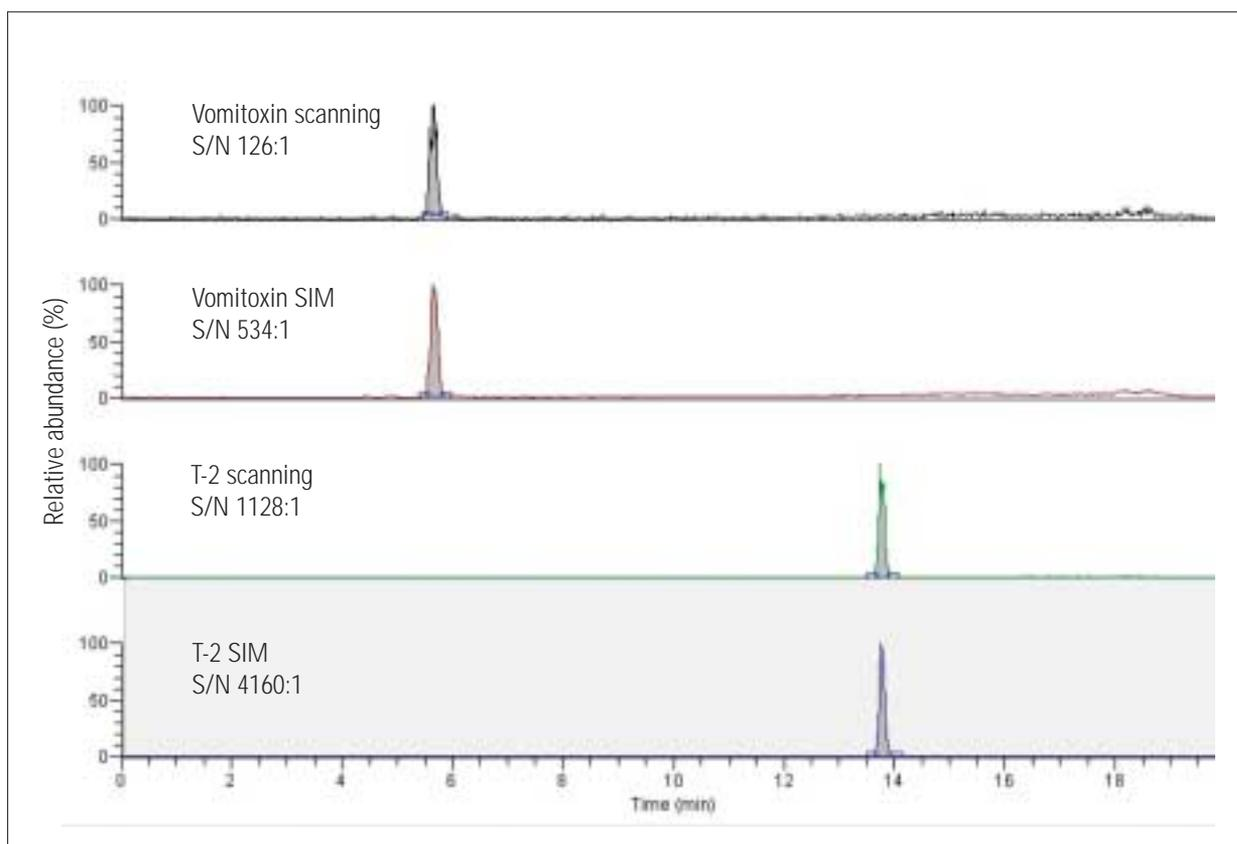


Figure 1. SIM and Scanning Mass Chromatograms can be run simultaneously.

Results and Discussion

Figure 1 shows the unique capability of the Surveyor MSQ to collect selected ion monitoring (SIM) and scanning data from the same run. The improvement in signal to noise from the scanning (upper) to the selected ion (lower) acquisitions is shown on the pseudomolecular ion traces for both Vomitoxin (DON) and T-2 Toxin respectively. This allows low level quantitative data and qualitative survey data to be obtained from the same injection.

The full-scan mass spectral data obtained from this run is shown in Figure 2. Pseudomolecular ion species, presumably sodium ion adducts, were observed at $[M+23]$ for both the Vomitoxin (upper) and the T-2 Toxin (lower).

The calibration data from 5-500 $\text{pg}/\mu\text{L}$ for Vomitoxin is shown in Figure 3. The selected ion monitoring chromatogram for the 5 $\text{pg}/\mu\text{L}$ standard shows a signal to noise ratio of 28 to 1 and the correlation co-efficient (R^2) was 0.9994. For T-2 Toxin a similar calibration line was observed (not shown) giving a signal to noise of 7 to 1 for the 5 $\text{pg}/\mu\text{L}$ standard level and a correlation coefficient (R^2) of 0.9982.

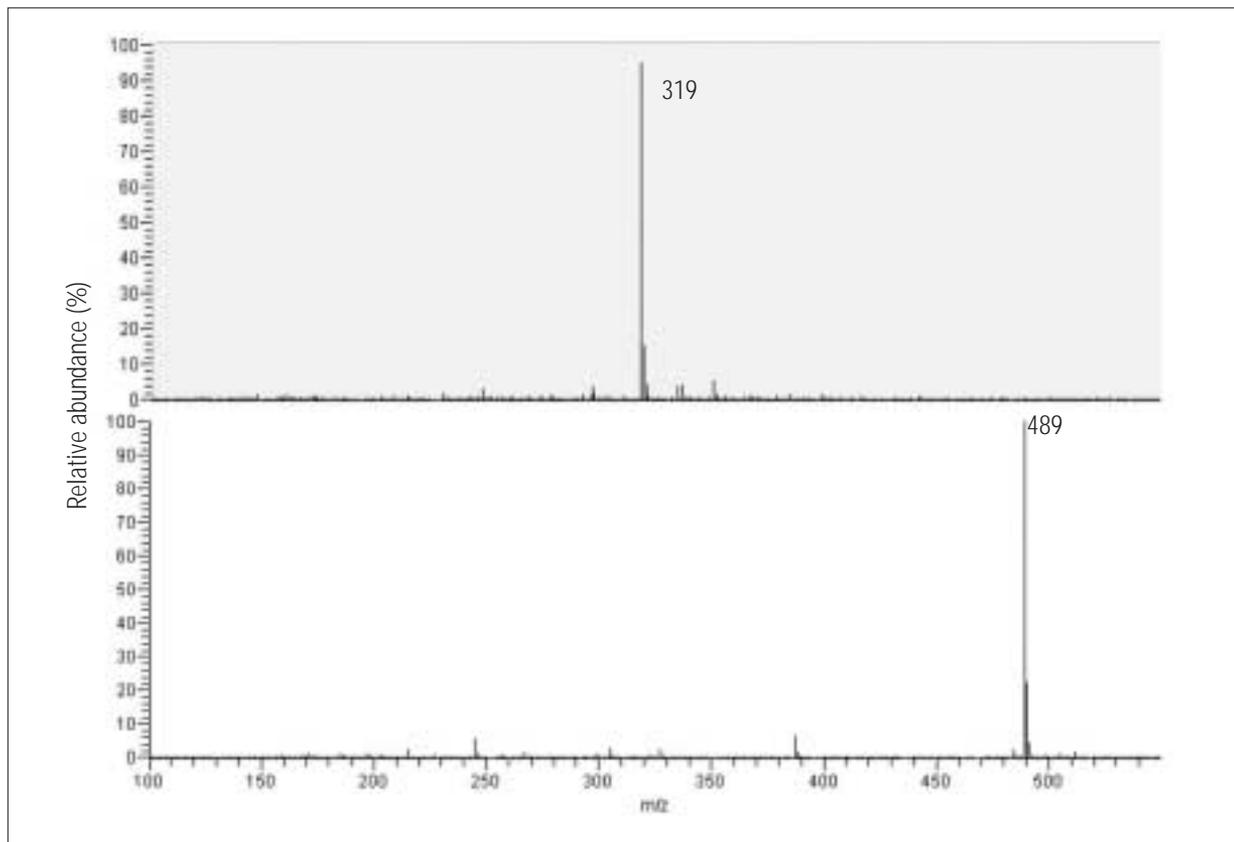


Figure 2. Vomitoxin (upper) and T-2 Toxin (lower) mass spectra

Conclusions

A simple LC/MS method for the separation and analysis of these important fungal metabolites is shown. The data show excellent linearity and sensitivity for these compounds that would readily allow determination on contamination to below ppm levels in grain samples when combined with a suitable extraction technique. The unique capability of the Surveyor MSQ to acquire both SIM and scanning data from the same sample injection is demonstrated

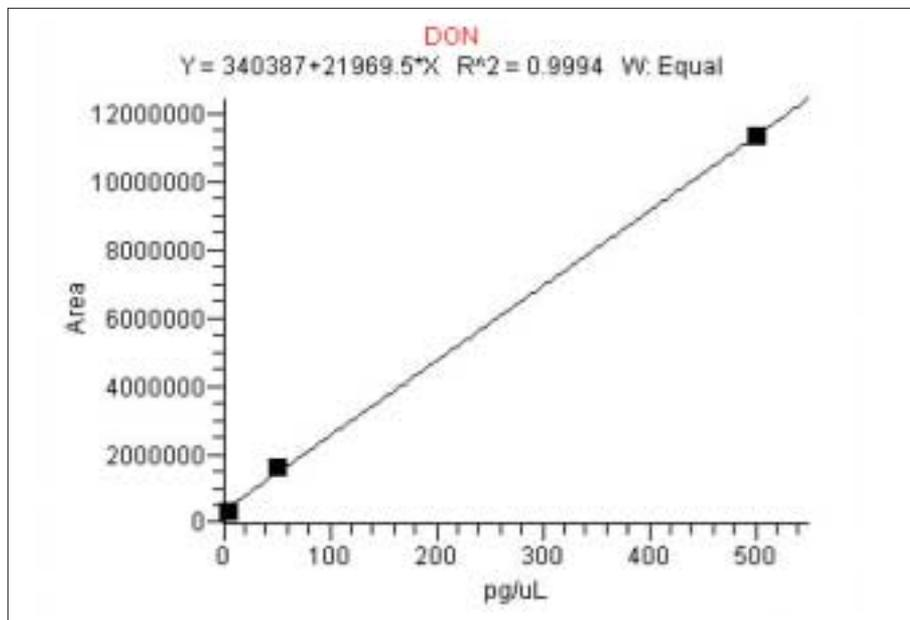


Figure 3. Vomitoxin calibration data

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

Australia
+61 2 9898 1244

Austria
+43 1 333 50340

Belgium
+32 2 482 30 30

Canada
+1 800 532 4752

China
+86 10 5850 3588

France
+33 1 60 92 48 00

Germany
+49 6103 4080

Italy
+39 02 950 591

Japan
+81 45 453 9100

Latin America
+1 512 251 1503

Netherlands
+31 76 587 98 88

Nordic
+46 8 556 468 00

South Africa
+27 11 570 1840

Spain
+34 91 657 4930

Switzerland
+41 61 48784 00

UK
+44 1442 233555

USA
+1 800 532 4752

www.thermo.com



Thermo Finnigan LLC,
San Jose, CA USA
is ISO Certified.

AN61568_E 03/04S