

FOOD ANALYSIS

A VALIDATED LC/MS/MS SOLUTION FOR THE ANALYSIS OF MYCOTOXINS IN CEREALS



Solutions for Your Analytical Business
Markets and Applications Programs



Matrices

- Corn
- Soft Wheat
- Hard Wheat

Target Compounds

- Aflatoxins
- Fumonisin
- Ochratoxin
- Deoxynivalenol
- Zearalenone

This Solution is dedicated to the analysis of Aflatoxins, Fumonisin, Ochratoxin, Deoxynivalenol and Zearalenone in corn, soft wheat and hard wheat

Moulds can develop on food commodities of plant and animal origin. These moulds can, in suitable environmental conditions, produce chemical toxic compounds known as mycotoxins.

European Community regulations determine the maximum allowed amounts of mycotoxins in different matrices for its member states.

Agilent offers a turn-key Solution covering:

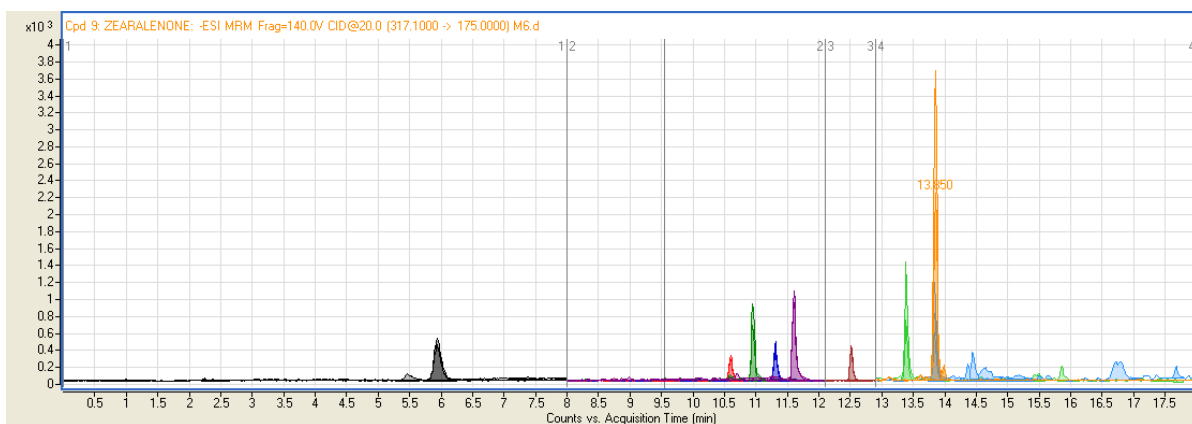
- Sample preparation
- Sample analysis
- Data analysis
- Full Agilent support



Recoveries, precision, LODs and LOQs are in accordance with COMMISSION REGULATIONS (EC) No 401/2006, No 1881/2006, No 105/2010 and No 165/2010.

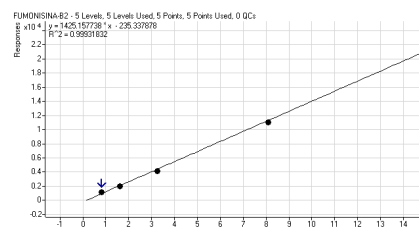
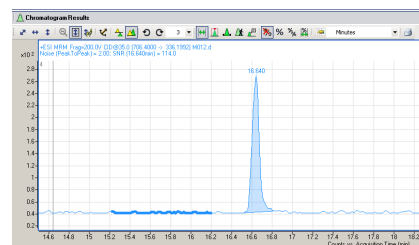
Sample preparation consists of one single extraction with organic solvent and water for all matrices. Sample analysis is performed with an Agilent UHPLC 1260 and an Agilent 6460 triple quadrupole. All mycotoxins are determined in a single analytical run.





The Agilent Mycotoxins Solution for cereals comprises the following:

- Complete hardware set-up (Agilent 1260 UHPLC and Agilent 6460 triple quadrupole) with dedicated analytical column;
- Standard operating procedure (SOP) with detailed descriptions of the analysis and validation procedures;
- Method of analysis (a DVD/CD containing sample preparation and sample analysis methods, recommended consumables and materials);
- One week training at Laboratorio Químico-Microbiológico (Murcia, Spain) on sample prep, validation procedures and analytical methods;
- On-site Training and Support (at customer site) covering Qualitative and Quantitative analysis;



STANDARD OPERATING PROCEDURE

Rev. 00

DATE: 2011/02/01

REVIEWED: 2011/02/01

APPROVED: 2011/02/01

MADE BY: [Name]

REVIEWED BY: [Name]

APPROVED BY: [Name]

DATE: 2011/02/01

DATE: 2011/02/01

DATE: 2011/02/01



maps_agilent@agilent.com
www.agilent.com/chem/solutions_app

Agilent Products are for Research Use Only.
 Not for use in diagnostic procedures.
 Information, descriptions and specifications in this
 publication are subject to change without notice.

© Agilent Technologies, Inc. 2013
 Published in USA, October 2013
 5991-3458EN

The Measure of Confidence



Agilent Technologies