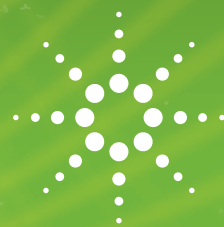


FOOD ANALYSIS

A VALIDATED LC/MS/MS SOLUTION FOR THE ANALYSIS OF PESTICIDES AND OTHER CHEMICALS IN CEREALS



Solutions for Your Analytical Business
Markets and Applications Programs



Agilent Technologies

Authorized Partner Laboratory

Matrices

- Soft Wheat
- Hard Wheat
- Barley
- Oilseed Rape
- Corn
- Soybean
- Peanuts
- Beans
- Sunflower

Target Compounds

- Pesticides
- Growth Regulators

This Solution is dedicated to the analysis of more than 170 pesticides and growth regulators in different cereals.

Pesticides and other chemicals are used widely in agriculture.

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

Agilent offers a turn-key Solution covering:

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



The working range goes from 0.01 mg/kg to 0.10 mg/kg for all analytes. LOQs, LODs, precision and recoveries are in accordance to DG SANCO guidelines.

Sample preparation is done following QuEChERS approach: extraction with organic solvent followed by dispersive SPE with SAMPLIQ material. Sample analysis is performed with an Agilent UHPLC 1260 and an Agilent 6460 triple quadrupole.



Agilent
Guaranteed
Solution

Developed
Installed
Tested
Supported

The Measure of Confidence



Agilent Technologies

© Agilent Technologies, Inc. 2013
Published in USA, October 2013
5991-3461EN