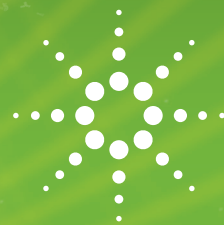


FOOD ANALYSIS

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



Solutions for Your Analytical Business
Markets and Applications Programs



Matrices

- Soft Wheat
- Hard Wheat

Target Compounds

- Chlormequat
- Mepiquat

This Solution is dedicated to the analysis of Chlormequat and Mepiquat in different cereals.

In agriculture, the quaternary ammonium herbicides Chlormequat and Mepiquat are used as plant growth regulators either individually, as mixtures or with other pesticides.

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

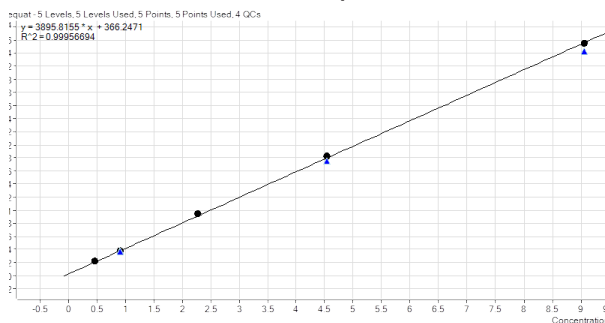


The working range goes from 0.01 mg/kg to 0.10 mg/kg for all analytes. LOQs, LODs, precision and recoveries are calculated in accordance to DG SANCO guidelines and used whenever possible the MRLs established in the Regulation (EC) No 396/2005 of the European Parliament and the Council on maximum residue levels (MRLs) in products of plant and animal origin.

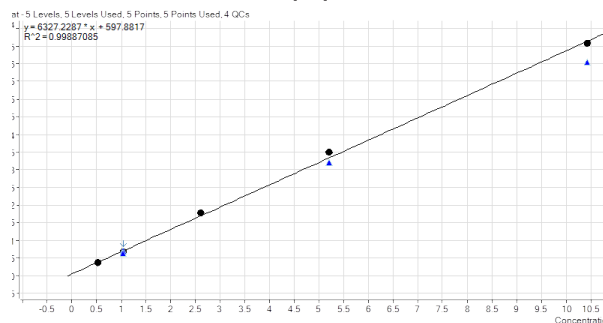
Sample preparation is done with a single solvent extraction. Sample analysis is performed with an Agilent UHPLC 1260 and an Agilent 6460 triple quadrupole.



Chlormequat



Mepiquat

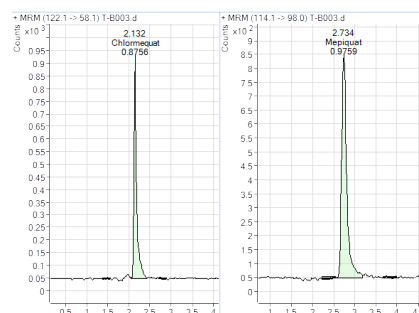


Straight calibration from 0,5 ppb to 10 ppb

The Agilent Pesticides 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;

Chlormequat and Mepiquat at 1 ppb



STANDARD OPERATING PROCEDURE

Rev. 00

SMALL CHANGES CONTROL

Date	Page(s)

MADE BY: [Signature]
 REVIEWED: [Signature]
 APPROVED: [Signature]

Date: 01/10/2011 Date: 01/10/2011 Date: 01/10/2011

STANDARD OPERATING PROCEDURE

1. PURPOSE

2. SCOPE

3. REFERENCES

4. GENERAL

5. ANALYTICAL PROCEDURE

6. VALIDATION

7. RESULTS

8. APPENDICES

9. REFERENCES

10. APPENDICES

11. APPENDICES

12. APPENDICES

13. APPENDICES

14. APPENDICES

15. APPENDICES

16. APPENDICES

17. APPENDICES

18. APPENDICES

19. APPENDICES

20. APPENDICES

21. APPENDICES

22. APPENDICES

23. APPENDICES

24. APPENDICES

25. APPENDICES

26. APPENDICES

27. APPENDICES

28. APPENDICES

29. APPENDICES

30. APPENDICES

31. APPENDICES

32. APPENDICES

33. APPENDICES

34. APPENDICES

35. APPENDICES

36. APPENDICES

37. APPENDICES

38. APPENDICES

39. APPENDICES

40. APPENDICES

41. APPENDICES

42. APPENDICES

43. APPENDICES

44. APPENDICES

45. APPENDICES

46. APPENDICES

47. APPENDICES

48. APPENDICES

49. APPENDICES

50. APPENDICES

51. APPENDICES

52. APPENDICES

53. APPENDICES

54. APPENDICES

55. APPENDICES

56. APPENDICES

57. APPENDICES

58. APPENDICES

59. APPENDICES

60. APPENDICES

61. APPENDICES

62. APPENDICES

63. APPENDICES

64. APPENDICES

65. APPENDICES

66. APPENDICES

67. APPENDICES

68. APPENDICES

69. APPENDICES

70. APPENDICES

71. APPENDICES

72. APPENDICES

73. APPENDICES

74. APPENDICES

75. APPENDICES

76. APPENDICES

77. APPENDICES

78. APPENDICES

79. APPENDICES

80. APPENDICES

81. APPENDICES

82. APPENDICES

83. APPENDICES

84. APPENDICES

85. APPENDICES

86. APPENDICES

87. APPENDICES

88. APPENDICES

89. APPENDICES

90. APPENDICES

91. APPENDICES

92. APPENDICES

93. APPENDICES

94. APPENDICES

95. APPENDICES

96. APPENDICES

97. APPENDICES

98. APPENDICES

99. APPENDICES

100. APPENDICES



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

The Measure of Confidence



Agilent Technologies

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
 5991-3459EN