



WINTER CONFERENCE on Plasma Spectrochemistry

January 9 - 14, 2012 • Tucson, AZ

Booth Exhibition - Agilent booths: 26, 27, 31 and 32

Special Event - Monday 7:00 p.m. (Agilent Booth)

Visit us in our booth and you'll experience the breadth of products in Agilent's Atomic Spectroscopy group including the 7700 ICP-MS, 700 ICP-OES, and all new Agilent 4100 MP-AES. We will also display the latest breakthrough in Atomic Spectrometry technology that we will reveal at our special Monday night event.

Lunch and Learns (registration is required, see below*)

11:30 p.m. – 1 p.m. each day in room **Coronado I**

Monday, January 9

Microwave Plasma – Atomic Emission Spectroscopy, a Revolutionary New Technique and Plasma Source for Atomic Spectroscopy

Tuesday, January 10

Hear the Latest Exciting ICP-MS Developments from Agilent Technologies

Wednesday, January 11

Put the Pedal to the Metal with these Atomic Spectroscopy Tips & Tricks

Customer Appreciation Celebration

Please join us on Tuesday, January 10, from 7 p.m. – 9 p.m. in the Hilton El Conquistador Presidio Foyer for a fun-filled and friendly night. Connect with colleagues and our product experts over refreshments, cocktails, and entertainment.

Informational Posters:

Our informational posters will be on display throughout the event.

For more detailed information about Agilent's posters or technical sessions presented at the 2012 Winter Plasma Conference, please visit our website:

www.agilent.com/chem/wpc

***Registration is required for the Agilent Lunch and Learns and Customer Appreciation Celebration. If you have not yet registered for these events, please stop by the Agilent booth for an admittance ticket.**





Winter Plasma Conference 2012

Agilent Technologies Papers and Posters

Wednesday, January 11, 2012

Oral Presentations

Fundamentals, Instrumentation, and Mechanisms

10:20 W01 Probing an Analytical Microwave Plasma by 2D Imaging and Ion/Atom Thermometry, Richard Morrison, Agilent Technologies Pty Ltd, 679 Springvale Rd, Mulgrave, Melbourne 3170, Australia, richard.morrison@agilent.com; Philip Wilson, Thomas Preuss, Peter Doidge, Mike Hammer

Workshop on New Plasma Instrumentation

3:25 WS102 Evaluation of a Sample Introduction System for ICP-OES That Can Simultaneously Determine Hydride and Non-Hydride Forming Elements with a High Efficiency Nebulizer, Glyn Russell, Agilent Technologies Australia Pty Ltd, 679 Springvale Rd, Mulgrave Victoria 317, Australia, glyn.russell@agilent.com; Dennis Hoobin

4:10 WS105 MP-AES: A Revolution in Atomic Spectroscopy, Michael Hammer, Agilent Technologies Pty Ltd, 679 Springvale Rd, Mulgrave, Melbourne 3170, Australia, mike.hammer@agilent.com

4:25 WS106 The Next Step in Reaction Cell ICP-MS -- Advancing the Technology, Noriyuki Yamada, Agilent Technologies International Japan, Ltd., 9-1, Takakura-cho, Hachioji-shi, Tokyo 192-0033, Japan, oriyuki.yamada@agilent.com; Takeo Kuwabara, Hironori Yamada, Takashi Kondo

4:40 WS107 Eliminating Challenging Interferences on As, Se, and Ti Using ICP-MS Incorporating Novel Next Generation Cell Technology, Amir Liba, Agilent Technologies, 2850 Centerville Road, Wilmington DE 19808, amir_liba@agilent.com

Poster Presentations

Wednesday, January 11, 2012

3:00 WP07 Reaction Cell Frontier: Removing Oxide Polyatomic Ion Interferences Using an Innovative Reaction Cell ICPMS, Katsuo Mizobuchi, Agilent Technologies Japan Ltd., 9-1, Takakura-cho, Hachioji-shi, Tokyo 192-8510, Japan, katsuo_mizobuchi@agilent.com; Naoki Sugiyama

3:00 WP08 New Reaction Cell ICP-MS with Improved Ion Optics Arrangement and Vacuum System, Takeo Kuwabara, Agilent Technologies International Japan, Ltd., Tokyo Analytical Division, 9-1 Takaura-cho, Hachioji-shi, Tokyo 192-0033, Japan, takeo_kuwabara@agilent.com; Noriyuku Yamada, Jun Kitamoto, Hironori Yamada, Takashi Kondo

3:00 WP12 Understanding and Managing the Spectral Background in a N₂ Atmospheric Pressure Microwave Plasma, Richard Morrison, Agilent Technologies Pty Ltd, 679 Springvale Rd, Mulgrave, Melbourne 3170, Australia, richard.morrison@agilent.com; Phil Wilson, Mike Hammer

3:00 WP30 Benefits of the HPLC-ICP-MS Coupling for the Mercury Speciation in Food, Sébastien Sannac, Agilent Technologies, 1, rue Galvani, F-91745 Massy Cedex, France, sebastien_sannac@agilent.com; Yu-Hong Chen, Raimund Wahlen

Thursday, January 12, 2012

3:00 ThP03 Analysis of Wear Metals in Lubricating Oils by Microwave Plasma - Atomic Emission Spectrometry, Craig Taylor, Agilent Technologies Pty Ltd, 679 Springvale Rd, Mulgrave, Melbourne 3170, Australia, craig.taylor@agilent.com

3:00 ThP04 Analysis of Fuel Samples by Microwave Plasma - Atomic Emission Spectrometry, Craig Taylor, Agilent Technologies Pty Ltd, 679 Springvale Rd, Mulgrave, Melbourne 3170, Australia, craig.taylor@agilent.com; Philip Lowenstern, Elizabeth Reisman

3:00 ThP09 Efficient Removal of Polyatomic Ions by ICP-MS Equipped with Novel Reaction Cell: Examples of Highly Purified Chemicals Used for Semiconductor Manufacturing, Junichi Takahashi, Agilent Technologies Inc., 9-1 Takakura-cho, Hachioji-shi, Tokyo 192-0033, Japan, dr-junichi_takahashi@agilent.com; Noriyuki Yamada

3:00 ThP31 Clinical Evaluation of Chromium, Cobalt, and Titanium by ICP-MS for Metal on Metal Implants, Anna Miller, Medtox Scientific, 402 West Country Rd D, New Brighton MN 55112, amir_liba@agilent.com, amiller@medtox.com; Gregory Janis, Amir Liba

3:00 ThP34 Enhancement of Productivity for the Analysis of Food Samples with the 7700x ICP-MS, Sébastien Sannac, Agilent Technologies France, 1, rue Galvani, F-91745 Massy Cedex, France, ebastien_sannac@agilent.com; Jean-Pierre Lener, Jérôme Darrouzes

3:00 ThP35 An Innovative New Plasma Source for Elemental Analysis Using Atomic Spectroscopy, Craig Taylor, Agilent Technologies Pty Ltd, 679 Springvale Rd, Mulgrave, Melbourne 3170, Australia, craig.taylor@agilent.com; Steve Wall, Eric Vanclay, Mike Hammer

3:00 ThP36 Analysis of Food and Beverage Samples by Microwave Plasma - Atomic Emission Spectrometry, Craig Taylor, Agilent Technologies Pty Ltd, 679 Springvale Rd, Mulgrave, Melbourne 3170, Australia, craig.taylor@agilent.com; Tran Nham

3:00 ThP43 Validating ICP-MS Using USP <232>/<233> for Elemental Impurity Analysis in Pharmaceutical Products, Ed McCurdy, Agilent Technologies Ltd, 5500 Lakeside, Cheadle Royal Business Park, Stockport, Cheshire SK8 2GR, United Kingdom, ed_mccurdy@agilent.com; Amir Liba, Samina Hussain

3:00 ThP47 Analysis of Chinese Herbal Medicine Samples by Microwave Plasma - Atomic Emission Spectrometry, Craig Taylor, Agilent Technologies Pty Ltd, 679 Springvale Rd, Mulgrave, Melbourne 3170, Australia, craig.taylor@agilent.com; Chun-hua Wu

Friday, January 13, 2012

Nuclear

3:00 FP09 Reaction Cell Frontier: Analysis of Radionuclides in Environmental Samples Using an Innovative Reaction Cell ICP-MS, Yasuyuki Shikamori, Agilent Technologies International Japan, Letd., 9-1, Takakura-cho, Hachioji-shi, Tokyo 192-0033, Japan, yasuyuki_shikamori@agilent.com; Kazumi Nakano, Naoki Sugiyama, Shinichiro Kakuta

3:00 FP25 Reaction Cell Frontier: Selenium Isotope Dilution Analysis Using Innovative Reaction Cell ICP-MS, Naoki Sugiyama, Agilent Technologies International Japan, 9-1 Takakura-cho, Hachioji-shi, Tokyo 192-0033, Japan, naoki_sugiyama@agilent.com

3:00 FP30 Analysis of Platinum Group Metals in Geological Samples by Microwave Plasma - Atomic Emission Spectrometry, Craig Taylor, Agilent Technologies Pty Ltd, 679 Springvale Rd, Mulgrave, Melbourne 3170, Australia, craig.taylor@agilent.com

3:00 FP38 Chromium Speciation in Drinking Waters Using LC(IC)-ICP-MS, Mina Tanoshima, Agilent Technologies, 9-1, Takakura-cho, Hachioji-shi, Tokyo 192-0033, Japan, mina_tanoshima@agilent.com; Tetsushi Sakai

the **REMARKABLY BETTER**
ATOMIC
SPECTROSCOPY
RUNS revolution
on **AIR**



Introducing the NEW Agilent 4100 MP-AES. Say goodbye to flammable and expensive gases. Say hello to enhanced productivity. The Agilent 4100 Microwave Plasma-Atomic Emission Spectrometer runs entirely on air. It's safer and more cost efficient. It enables superior detection and unattended multi-element analysis. It's the most significant advancement in atomic spectroscopy in decades. In other words:
It's time to run on air.

JOIN THE REVOLUTION. Learn more about the NEW Agilent 4100 MP-AES.
www.agilent.com/chem/RunsOnAir

© Agilent Technologies, Inc. 2012

