

# AGILENT WEBINAR:

## Ultra-fast Analysis of Small Molecule Analytes in Urine and Blood-based Matrices Using an SPE/MS/MS System

The Measure of Confidence



February 16, 2012

Agilent is hosting a webinar that will demonstrate the ability of an ultra-fast SPE/MS/MS system to analyze small molecule analytes in human urine or blood-based matrices with much faster sample cycle times and similar analytical results compared to LC/MS/MS assays.

<b>Presenter:</b>	Dr. Vaughn Miller RapidFire Applications Manager, Agilent Technologies, Inc.
<b>Date:</b>	February 16, 2012
<b>Time:</b>	8:00 am (PST)/11:00 am (EST)
<b>Location:</b>	Your Desktop
<b>Register:</b>	<a href="http://www.agilent.com/chem/SPMSMS_webinar">www.agilent.com/chem/SPMSMS_webinar</a>

### Join our Webinar, where we will:

- Analyze clinical research and forensic toxicology samples 10X faster than traditional methods
- Investigate a variety of matrices in a high-throughput manner
- Streamline workflows with RapidFire to achieve results that correlate with LC/MS/MS

We look forward to your participation!

Register today at:

[www.agilent.com/chem/SPMSMS\\_webinar](http://www.agilent.com/chem/SPMSMS_webinar)

### Analytes Assessed

- benzodiazepines
- marijuana metabolite
- cocaine metabolite
- tacrolimus
- levetiracetam
- hydroxymidazolam

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. 2011

C-00005816



Agilent Technologies