

AIR ANALYSIS EDUCATIONAL WEBINARS

ACCELERATE RESULTS IN EVERY ENVIRONMENT

Regulatory agencies globally are increasing requirements and rigor for the monitoring of air to protect human health and the environment. Ever-changing regulations, emerging contaminants, method variation, and difficult matrices means your laboratory testing capabilities must be agile.

Indoor and outdoor air monitoring requires accurate, reliable analysis. The Air Analysis on-demand webinars are designed to help you overcome these common challenges and improve efficiencies in your lab.

Who Would Benefit from Watching these Air Analysis Webinars?

- Laboratory Analysts
- Air Quality Scientists and Chemists
- Laboratory Managers and Directors

Advantages of Passive Air Monitoring



This webinar examines the methodology and benefits of passive sampling for monitoring volatile organic compounds in air. EPA Method 325 will be used as an example to demonstrate how passive monitoring is an accurate, reliable and

easy way to collect air samples over both short and long-term periods, allowing efficient, cost effective characterization of air quality.

<https://www.perkinelmer.com/passiveairmonitoringwebinar>

Air Sampling Techniques Examined



This webinar highlights different techniques for air compound monitoring and the applications in which they are utilized. You will learn about optimum analytical parameters and how to manage the vast data typically collected for ozone precursor monitoring.

<https://www.perkinelmer.com/airsamplingtechniqueswebinar>

Learn How to Analyze Both Volatile and Semi-volatile Compounds in a Single Air Analysis



In this webinar, best practices on how to streamline collection and analysis of volatile and semi-volatile organic compounds in air will be covered. The focus of the discussion is on two main applications, Soil Gas and Manufacturer Gas Plant Sites. Learn how to

analyze these compounds in one method recovering compounds through the boiling point of nC44 while retaining the most volatile compounds required by regulations.

<https://www.perkinelmer.com/simplifiedVOCandSVOCairanalysis>

Lab Quality Analysis in Continuous Ambient Air Monitoring



This webinar examines quality control practices and optimized processes necessary for VOC data collected, from automated gas chromatography systems (AutoGCs) at Photochemical Assessment Monitoring Sites (PAMS), that will be utilized in ambient air modeling. Learn how Orsat has

helped to establish needed automation to maintain high quality hourly data that has been involved in the PAMS monitoring program for the Texas Commission on Environmental Quality (TCEQ) for over 20 years.

<https://www.perkinelmer.com/continuousambientairmonitoringwebinar>

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com


PerkinElmer
For the Better

For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2020, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.