



When Optics Matter.

Reduce impairments and gain deeper insights with advanced spectrophotometric tools and techniques



Register for this free webcast at:

www.spectroscopyonline.com/spec_p/optics

LIVE WEBCAST

Europe:

Wednesday, March 15, 2023
10am BST | 11am CET

North America:

Wednesday, March 15, 2023
1pm PST | 3pm CDT | 4pm EDT

Asia:

Thursday, March 16, 2023
10:30am IST | 1pm SGT | 2pm JST | 4pm AEDT

Presenter



Travis Burt

Global Product Manager –
Cary UV-Vis-NIR Spectrophotometers
Agilent Technologies, Inc.

Moderator



Robert Alaburda

Special Projects Director
Spectroscopy

Event Overview

This webinar will guide you on the best practices, tools, and techniques needed to get the most reliable, verifiable insights into your optical components. Appropriate choice of technique and methodology is critical to getting the information you need to improve design, verify product performance, and accelerate manufacturing.

We will explore the subtleties of reflection and transmission measurements at normal incidence and at oblique angles, emphasizing high absorbance, optical density, linearity, and effects of incident beam collimation. We will discuss diffuse scattering measurements and automation/robotics to broaden quality assurance programs and speed up analysis.

The webinar is relevant to developers of optics, photonics, and electro-optical or semiconductor devices. Improve your understanding of the link between your desired outcomes and the spectrophotometric approaches needed to achieve those insights.

Key Learning Objectives

Attendees will depart with a practical understanding of

- The best spectrophotometric practices to extract deeper insights into the characteristics of optical samples.
- How to overcome the limitations of traditional spectrophotometric tools and verify outcomes.
- Accelerating QA/QC of optics in a production environment.

Who Should Attend

- Commercial or Academic organizations in the fields of; optics, photonics, electro-optical, or semiconductors
- Laboratory Managers
- Scientist/Staff Scientists
- Lab Technician/Technologist
- Analyst
- Project Manager
- SME Directors, Core Lab/ Facility Managers
- Under/Post Graduate Researchers

Sponsored by

Presented by



For questions email
jramesh@mjhllifesciences.com