



Extended Uniformity of Dosage Unit Testing of Pharmaceutical Tablets, Using Transmission Raman Spectroscopy

LIVE WEBCAST

Airing 1 - Europe:

Tuesday, March 28, 2023

1pm BST | 2pm CET

Airing 2 - North America:

Tuesday, March 28, 2023

11am PDT | 1pm CDT | 2pm EDT

Airing 3 - Asia:

Wednesday, March 29, 2023

10:30am IST | 1pm SGT | 2pm JST | 4pm AEDT

Presenter



Anders Sparén

Associate Principal Scientist
AstraZeneca Gothenburg, Sweden

Moderator



Megan Manzano

Senior Editor of Special Projects
Spectroscopy



Register for this free webcast at:

www.spectroscopyonline.com/spec_p/dosage-unit

Event Overview

In this webinar, Anders Sparén, Associate Principal Scientist at AstraZeneca, will present work utilizing Transmission Raman spectroscopy (TRS).

The increased sampling volume minimizes sub-sampling and makes the technique ideal for the assessment of assay, uniformity of dosage unit, and polymorphs in tablets and capsules. Due to its non-destructiveness and high sample throughput, transmission Raman spectroscopy can be applied to extended uniformity of dosage unit testing and enhanced control strategies aiming at real-time release for batch and continuous operations.

Key Learning Objectives

- Learn how AstraZeneca utilizes transmission Raman analytical workflows in PAT
- Learn different approaches how to interpret and evaluate the spectral data
- Find out more about the features and benefits of transmission Raman towards an enhanced control strategy
- Find out how rapid, nondestructive technology can benefit your lab

Who Should Attend

- Pharmaceutical scientists specializing in solid oral dose formulations and manufacturing
- Lab managers with interest in tools to improve analytical workflows and operational efficiencies
- Spectroscopists and process analytical technology (PAT) scientists interested in vibrational spectroscopy

Sponsored by



Presented by

Spectroscopy

Pharmaceutical
Technology

For questions email
JdelaBandera@mjhlifesciences.com