## Visualizing, Characterizing, and Analyzing Pharmaceutical Constituents with Raman Imaging

Robert Heintz, Ph.D. Senior Applications Specialist <u>Robert.heintz@thermofisher.com</u>

The world leader in serving science

#### Agenda

• Brief overview of Raman spectroscopy & Raman imaging

- Introducing the Thermo Scientific<sup>™</sup> DXR<sup>™</sup>xi Raman imaging microscope
- Raman imaging for pharmaceutical products
  - Examples
    - Pharmaceutical Tablet homogeneity and content uniformity
    - Low Dose Tablet distribution of polymorphs
    - Hot Melt Extruder Products component characterization



## What is Raman Spectroscopy?

- Complementary technique to infrared (IR) spectroscopy
- Uses light to probe covalent chemical bonds by looking at vibrations
- Provides detailed molecular information: sensitive to even slight changes in bond angle or strength
- Useful for identifying unknown solids and liquids, including both inorganic and organic materials
- Can also detect sensitive changes in structure, morphology, and even temperature!



#### Raman Spectroscopy – The Raman Effect





## What Can Raman Imaging Do?

- Extends the advantages of Raman analysis across the sample
- Rapid collection of vast amounts of spectroscopic data
- Provides visual images depicting differences in molecular structure and chemical environment
- Raman images provide views of the samples that are not always apparent in the visual images





#### **Application Areas for Raman Imaging**

#### **Pharmaceuticals**





#### **Other Application Areas**

Polymers and Packaging Semiconductors and Thin Films Carbon Nanomaterials Geology / Mineralogy Life Sciences



#### Introducing the Thermo Scientific DXRxi Raman Imaging Microscope

A total imaging system: hardware and software integration combines **powerful performance** with **image-centric** analysis and **ease of use** 



A completely **new approach** to Raman imaging!



## Intelligent Workflow with Excellent Flexibility





2 Confidently optimize settings with intuitive controls

#### Quickly prioritize multiple regions of interest and run









#### No Raman Expertise Required to get the Best Results

Visual controls and instantaneous, continuous visual feedback

- NO lengthy trial and error
- NO guesswork
- You can see when parameters are optimal
- Focus quickly on the problem, not the technique

Laser power	5.9 mW			
Exposure time	0.02000 sec (50 Hz)			
Number of Scans	25			
Step size	0.5 µm			
Aperture 50 micron slit	J			
	0	<u>_</u>		
	3200 3000 2800 2600	2400 2200 2000 1800 1600 140 Raman shift (cm-1)	0 1200 1000 800 600	400 200
				>>



#### **Raman Image Preview**

- Just like using an visual image to inspect the sample now it is possible to use a Raman image preview of the sample.
- Don't waste time guessing at your region of interest
- Don't wait until an image is collected to learn if parameters were ideal
- Rapidly see and identify constituents and domains without the wait





#### Get There Faster By Getting Just What You Need

- Optimize *image* collection, not individual spectra
  - Quickly and visually balance image collection time with necessary detail level
  - Remove unanticipated results somewhere in an image
  - Stop any time if results are good enough rather than wait for multiple scans of each point in entire image to finish, one at a time



20 micron image pixels 1900 spectra, 2 scans 1 minute



5 micron image pixels 30,000 spectra , 2 scans 4 minutes



1 micron image pixels 191,000 spectra , 2 scans 25 minutes



#### Image-centric vs. spectral-centric at same "spectral speed"

100 x 100 microns 1 micron spacing (10,000 spectra); 500 Hz data collection; 10 "co-adds"

**DXRxi**, rastering entire images to desired quality level, like other microscopes



Single scan of entire image with MCR, 10,000 spectra

1 minute



3 scans of entire image with MCR 200 seconds





**Other Raman** Imaging Systems, building images one spectrum at a time



1000 co-added spectra, no useful image information



3000 co-added



10 scans of entire

image with MCR

10,000 spectra, no useful spectra image information collected

MCR Processed Map



#### Built-In Expertise: Profiles and MCR (Multivariate Curve Resolution)

- Standard profiles (correlation, chemigram, peak area, peak height, peak ratios, peak shift) applied immediately via graphical interface
- Component analysis calculated in real-time

One-click application of image profiles is a unique concept – adds value at every step of the workflow







## Data Processing Concurrent With Data Collection

- Raman images with component identification are created in real time
  - Without configuring a spectroscopic method
  - Without prior operator knowledge of what's in the sample
  - Without waiting for an entire image data set to be collected
- Instant and obvious interpretation even if you don't know what you're looking for





#### Integration and Cross-Compatibility

- Access to raw data
  - All data (chemical image, video image, spectra) can be quickly exported using a full array of formats
  - HDF5 provides open-source solution for compatibility with third party packages
  - Send data to OMNIC and Specta with a single click!













## **Microscopy Options**

- Supports a wide variety of sample measurement options, including:
  - Single and dual microscope slide holders
  - Heating and cooling stages even during imaging!
  - Rotating stage insert
  - Industry standard wafer holders and SEM accessories
  - 'Breadboard'-style holder for custom configurations
  - Holder for the Thermo Scientific K-Alpha XPS!
- Integrated Olympus research grade optics for peak performance and stability:
  - High NA and long working distance objectives
  - Optional brightfield and darkfield optics
  - DIC and visual polarizers for more challenging samples
  - Available with transmission illumination











## **Pharmaceutical Formulations**

- Typically complex multi-component mixtures
- Need to identify and verify components
  - Known components
  - Impurities
  - Identify changes in components during processing
- Distribution of components
  - Homogeneity
  - Particle size
  - Content uniformity









#### **Tablet Imaging Example**



Video Mosaic Image (10X objective, 100X total magnification)

# Migraine Relief Tablet

11 mm diameter, 676 mg

APIs Acetaminophen Aspirin Caffeine

250 mg (37%) 250 mg (37%) 65 mg (9.6 %)

#### Inactive

corn starch, microcrystalline cellulose, sodium lauryl sulfate, sodium starch, glycolate, crospovidone, polyethylene glycol, polyvinyl alcohol, povidone, stearic acid, talc, titanium dioxide



## Imaging the Whole Tablet

Raman MCR Image



Area Imaged - 11 x 11 mm<sup>2</sup> 10X objective Image Pixel Size - 25 μm 226,000 spectra Exposure Time 1.8 ms (550 spectra per s) 532 nm laser,

#### 8 minute collect time!!



## Higher Resolution Image – Whole Tablet



Area Imaged - 11 x 11 mm<sup>2</sup> 10X objective Image Pixel Size - 5 μm 5.4 million spectra Exposure Time 1.8 ms (550 spectra per s) 532 nm laser 36 GB file Size only computer limited (128 GB RAM) 6 hour collection time (3 hr estimated)

#### Image Analysis % Area of Particles

Component	Calculated % (Surface Area)	Reported %
Aspirin	38.6	37
Acetaminophen	35.4	37
Caffeine	7.7	9.6



#### Tablets Components – From Multivariate Curve Resolution (MCR)



**Titanium Dioxide** 



Acetaminophen



#### Caffeine



## Small Area – Longer Exposure Time





1.6 x 1.7 mm<sup>2</sup>
50X Objective
5 micron image pixel size
116000 spectra
Exposure time 5 ms (200 spectra per s)
532 nm laser
5 averaged scans
55 minutes

#### <u>500 μm</u>

Blue – Caffeine, Green -Acetaminophen, Yellow – Aspirin, and Red – starch



#### Smaller Area, Higher Resolution, Longer Exposure Time



# 225 x 250 μm<sup>2</sup> 100X Objective 0.5 micron image pixel size 229000 spectra Exposure time 10 ms (100 spectra per s) 532 nm laser 5 averaged scans 3 hr collect

100 μm

Blue – Aspirin, Green – Acetaminophen, Yellow – Caffiene, Red – starch, Fuchsia – microcrystalline cellulose, Orange – sodium lauryl sulfate.



## Summary of Tablet Imaging

- Possible to Image an entire 11 mm diameter tablet in 8 minutes
- Higher resolution images on whole tablets are possible but may not be necessary and there are other alternatives to imaging the whole tablet (select regions, multiple regions)
- Raman imaging can give spatial distribution of components including particle size estimates and relative percentages based on areas occupied by different components
- APIs tend to be strong Raman scatterers. Weaker excipients may require longer exposure times and possibly better spatial resolution to differentiate them.



#### Low Dose Tablet Example



Video Mosaic Image (10X objective, 100X total magnification)

#### **Tibolone Tablet**

3% Tibolone

A synthetic steroid used in hormone replacement therapy

6 mm diameter

Polymorphs (monoclinic, triclinic)





#### Raman MCR Image of Low Dose Tablet

#### Raman MCR Image



5.7 x 5.7 mm<sup>2</sup> area
10X objective
25 μm image pixel size
52000 spectra
Exposure Time 20 ms (50 spectra per s)
532 nm laser
10 averaged scans
3 hr collect

Tibolone was not readily differentiated using MCR



Starch 📕 Lactose 📕 Fluorescent Compound

#### A Peak Height Profile Readily Shows the Location of the Tibolone





#### Smaller Area, Higher Resolution, More Scans



1.1 x 1.6 mm<sup>2</sup> area
10X objective
5 μm image pixel size
75000 spectra
Exposure Time 20 ms (50 spectra / s)
532 nm laser
25 averaged scans
10 hr collect

Tibolone is one of the components defined by the MCR analysis



## Distribution of Polymorphs of Tibolone





## Summary of Low Dose Tablet Imaging

- Lower dose tablets are commonly encounter pharmaceutical products
- Other profile options (peak height etc.) might be better choices than MCR for displaying the spatial distribution of low concentration components.
- Multiple options are available in the OMNICxi software for generating different types of Raman images based on different aspects of the spectra.
- Whole tablet imaging is possible
- Evaluation of the presence and distribution of polymorphs is possible.



## Raman Imaging of Hot Melt Extruder (HME) Samples

Hot melt extrusion is a novel way of formulating solid dosage pharmaceutical products (tablets, granules, pellets, and transdermal films)

Has been used extensively for a long time in the plastics industry

API and other components are combined with a pharmaceutically approved thermoplastic polymer (usually at higher temperatures).

Screw threads control the mixing and transport properties at various stages.

Final form depends on the die and post extruder processing.





## **HME General Processing Advantages**

- Continuous process inline monitoring and control
- Establish stable solid solutions
  - Increase the availability of poorly soluble ingredients
- Flexibility to easily produce different dosage products
- Availability of time release forms
- Taste masking
- Special dosage form designs (films, rods, etc...)
  - Die change provides different shapes for special applications
- Reduce the consumption of solvents
  - Compare to wet granulation process



## Raman Imaging of HME Products

- Hot melt extrusion produces new forms using new processes
- Processing can effect the components
- Monitoring components
  - Changes in molecular structure
- Spatial distribution and particle size of components
- Identification of unknowns (impurities and defects)



## Example of Raman Imaging of HME Products\*



Video Mosaic Image (10X objective, 100X total magnification) Darkfield Illumination HPMCAS (hydroxypropyl methyl cellulose acetate succinate or hypromellose acetate succinate) polymer carrier

Ibuprofen (25-33%) and ibuprofen (25-33%) + D-mannitol (7-15%).

Cross-sections mounted in epoxy for analysis

To evaluate the spatial distribution of components and to look for any unforeseen changes caused by processing conditions

\* HME Samples Provided by Dr. Adrian Kelly, School of Engineering, Design and Technology, Bradford University, UK



## Raman Image of HME Product



Raman MCR Image purple – HPMCAS, green –ibuprofen, yellow – epoxy, red – cyanoacrylate, blue – inorganic impurity



#### Smaller Area, Higher Resolution, Raman Imaging



(50X objective - brightfield illumination)

Raman MCR Image blue – HPMCAS, green –ibuprofen , yellow – ibuprofen

50X objective, 780 nm laser, 24 mW, 687 x 423  $\mu$ m area, 3.0  $\mu$ m image pixel size, 32289 spectra, 0.0100 s exposure time, 100 scans



## Subtle Differences in Ibuprofen Spectra





## Differences in the Raman Spectra of Ibuprofen

- Ibuprofen is a mixture of stereoisomers
  - Not distinguishable with this type of experiment (ROA required)
- The active form is S (+) ibuprofen
- Different polymorphs of ibuprofen have been reported
  - Phase I (thermodynamically stable) & Phase II (metastable)
- Degree of crystallinity effects the Raman spectra
  - Crystallinity versus amorphous
- Co-crystallization of ibuprofen with other components can alter the Raman spectrum
- Ibuprofen association with other carriers (polyvinyl pyrrolidone (PVP) can cause slight differences in Raman spectra
- The observed differences in the ibuprofen spectra in these products does not match with any of these effects
  - However it does illustrate how sensitive Raman imaging can be



#### Raman Imaging – From Whole Samples to Small Particles



#### Millimeters to Microns





#### Hot Melt Extruder Product – Ibuprofen & D-Mannitol



6.65 x 4.65 mm<sup>2</sup> area
10X objective
25 μm image pixel size
49476 spectra
Exposure Time 10 ms (100 spectra /s)
780 nm laser
100 averaged scans

Raman MCR Image blue – HPMCAS, green –ibuprofen, yellow – epoxy, orange – cyanoacrylate, fuchsia – D-mannitol



#### Summary of Raman Imaging of Hot Melt Extruder Products

- Raman imaging allows evaluation of the spatial distribution of components in hot melt extruder products. This is not generally available with the typical inline monitoring of the products.
- Raman can be used to monitor any changes in molecular structure and chemical environment including molecular associations that might be induced during the HME processing.
- There are many options for Raman imaging from imaging whole samples down to small particles (millimeters to microns). This is important for these types of samples where there can be a significant range in particle sizes.



## Let the Power of Raman Imaging Work for You

- Raman imaging is clearly a very useful analytical tool for evaluating pharmaceutical products
- Raman imaging extends the power of Raman spectroscopy across greater areas and further expands the utility of Raman spectroscopy
- Raman imaging gives you the ability to identify materials and to assess subtle differences in molecular structure and chemical environment
- The DXRxi Raman imaging microscope will help you:
  - Get results and solve problems quickly with exceptional performance
  - Allow more people to solve problems without the need for a central expert
  - Walk up and use the system within minutes, anytime, without a significant learning curve





#### Thank you!

- Thank you for attending!
- Visit <u>thermoscientific.com/DXRxi</u> for more information on the Thermo Scientific DXRxi Raman imaging microscope

# Accelerate your work Visualize your answers



