

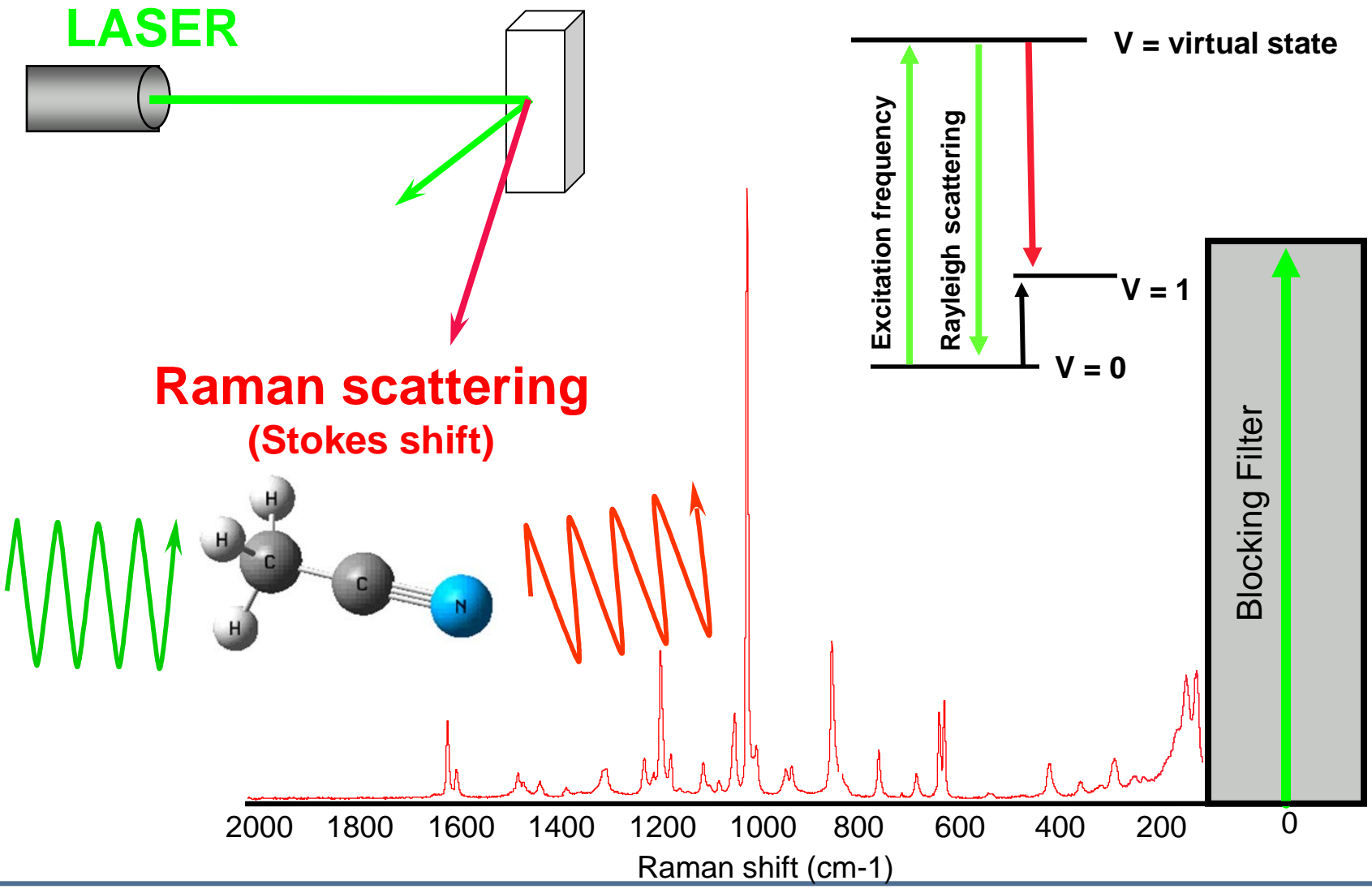
Raman Imaging: A Critical Tool for Realizing Graphene and Graphene Composite Materials

Dr. Mark Wall
Raman Product Specialist
Thermo Fisher Scientific

Overview

- Fundamentals of Raman spectroscopy
- *Modern* instrumentation
- The Raman Characterization of graphene
 - Point measurements
 - Mapping and Imaging
- Raman imaging examples
 - Partial growth graphene
 - Graphene as a protective coating material
- Confocal Raman imaging of graphene
 - Cross sectional and Z-stack measurements

Raman spectroscopy



Modern Raman Instrumentation

- Instruments of the past
 - Large
 - Complex
 - Difficult to maintain
 - Required dedicated operator



- Modern instrumentation
 - Integrated and compact
 - Easy to use
 - Very low maintenance
 - Automated alignment
 - Automated calibration
 - True walk up instruments



Thermo Scientific Family of Dispersive Raman Instruments

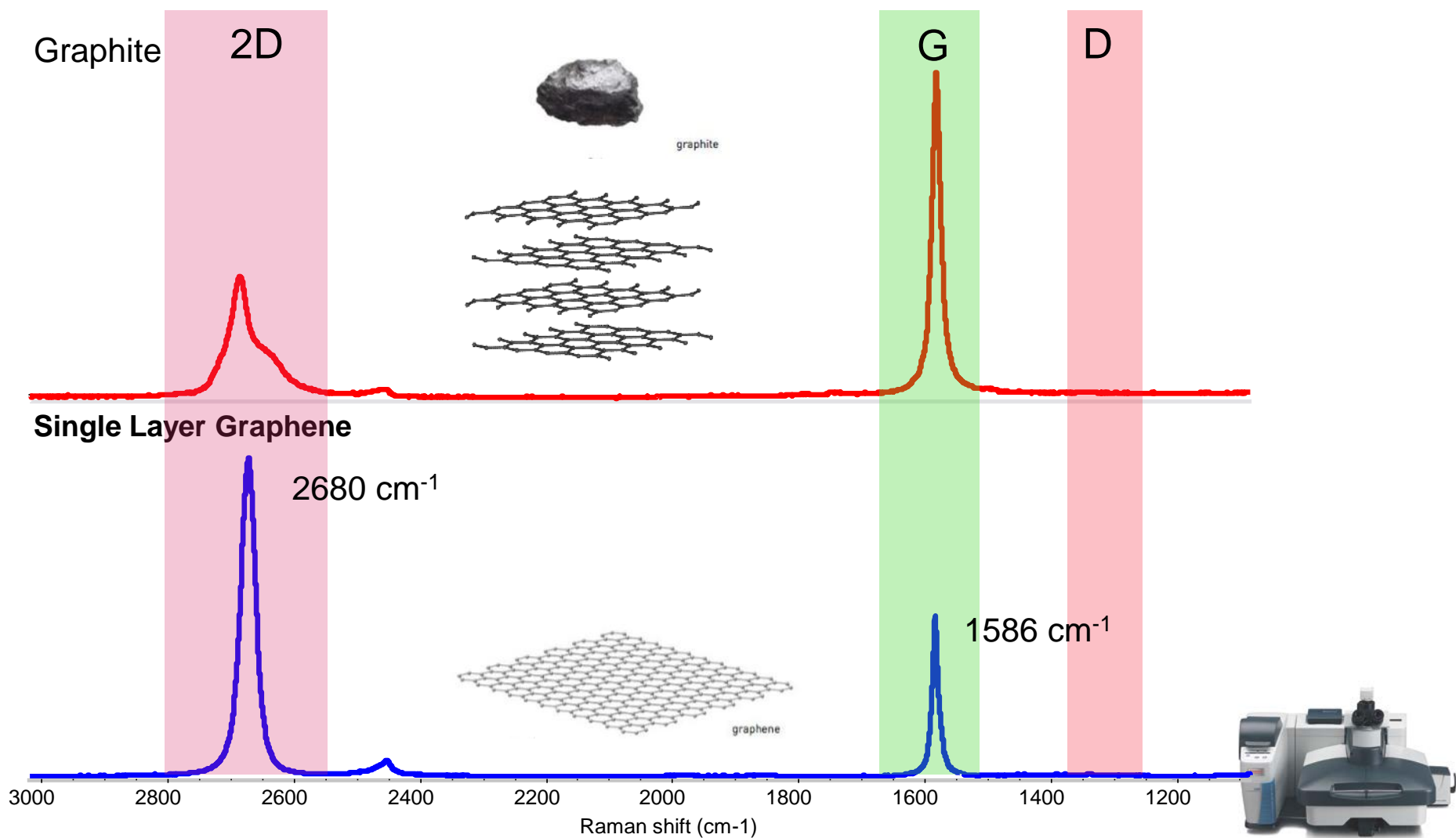
Raman Spectroscopy and Graphene Characterization

- Techniques used for characterization of graphene
 - SEM/TEM
 - Nanometer scale morphological information
 - Sample preparation needed
 - Detailed elemental information lacking in detailed molecular information
 - Requires vacuum
 - Often destructive measurement
 - XPS
 - Nanometer scale surface information
 - Oxidation state and functionalization information
 - Requires vacuum
 - AFM
 - Nanometer scale morphological information
 - Topological, mechanical
 - Lacking in molecular or chemical information

Raman Spectroscopy and Graphene Characterization

- Raman spectroscopy fills the chemical/ molecular “hole”
 - Provides specific and direct chemical information
 - Detailed structural information
 - Sensitive to the local environment
 - Stress and strain
 - No Sample preparation
 - Non destructive
 - No vacuum required
 - Rapid collection times
- Raman microscopy benefits
 - Surface and sub-surface characterization
 - Mapping/imaging capabilities
 - High spatial resolution
 - Sub micron x and y
 - Micron scale in the Z

Graphene: Raman Spectral Features



Raman Spectrum of Graphene – Principle Bands

- G band

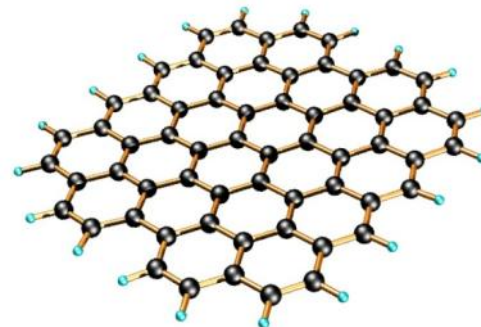
- Peak position and relative intensity is sensitive to
 - Layer thickness
 - Doping
 - Strain

- D band

- Peak intensity is sensitive to
 - Presence of defects or disorder
 - Sampling in proximity of an edge
 - Chemical modification
 - Increase of sp³ hybridized C bonding at the expense sp² hybridized C bonding

- 2D band

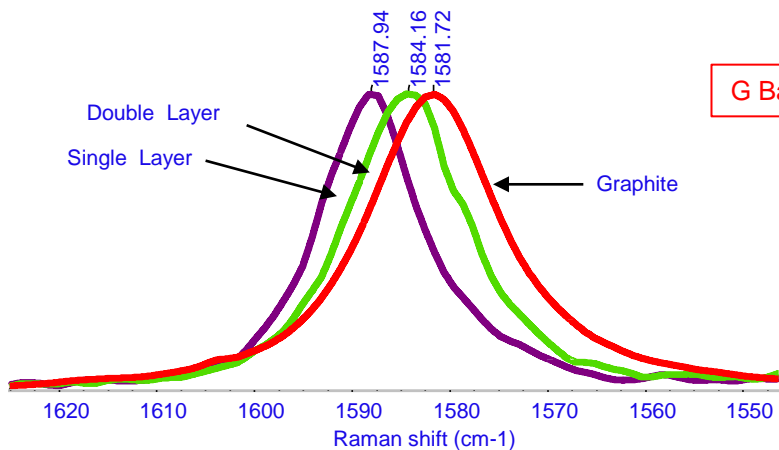
- Peak position, band shape, and intensity sensitive
 - Layer thickness and interlayer orientation
 - Strain



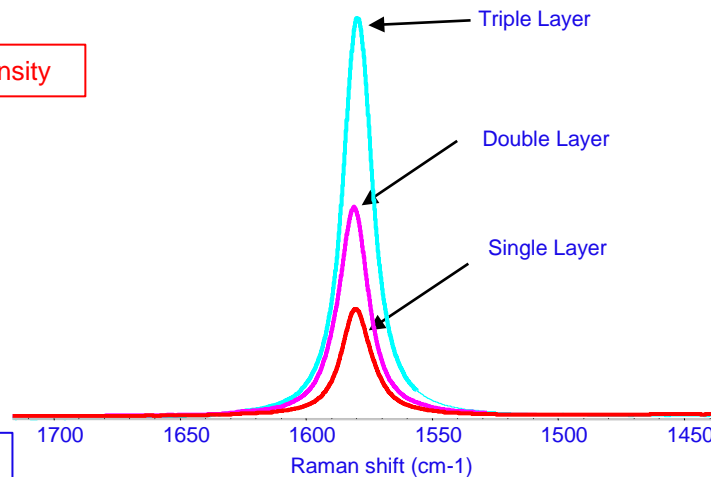
Raman Spectroscopy – Graphene Layer Thickness

$$\omega_G = 1581.6 + 11/(1 + n^{1.6})$$

Linear increase in G band intensity with # of layers present

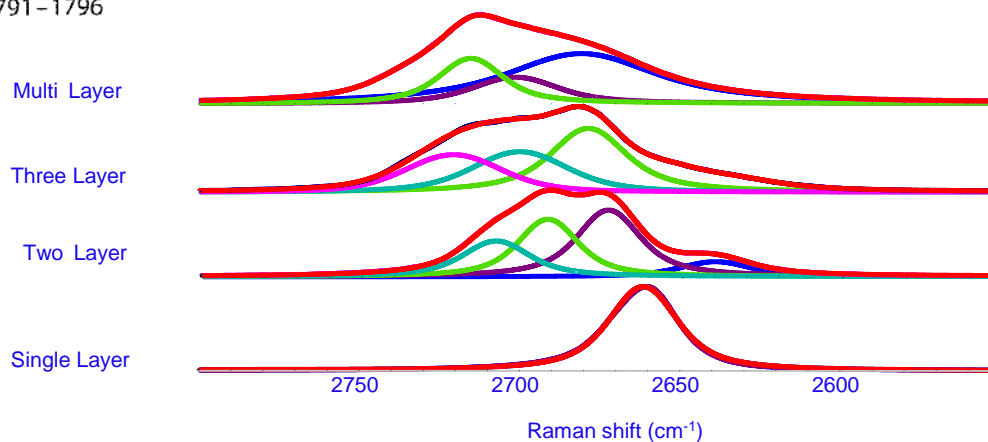


G Band Position and Intensity

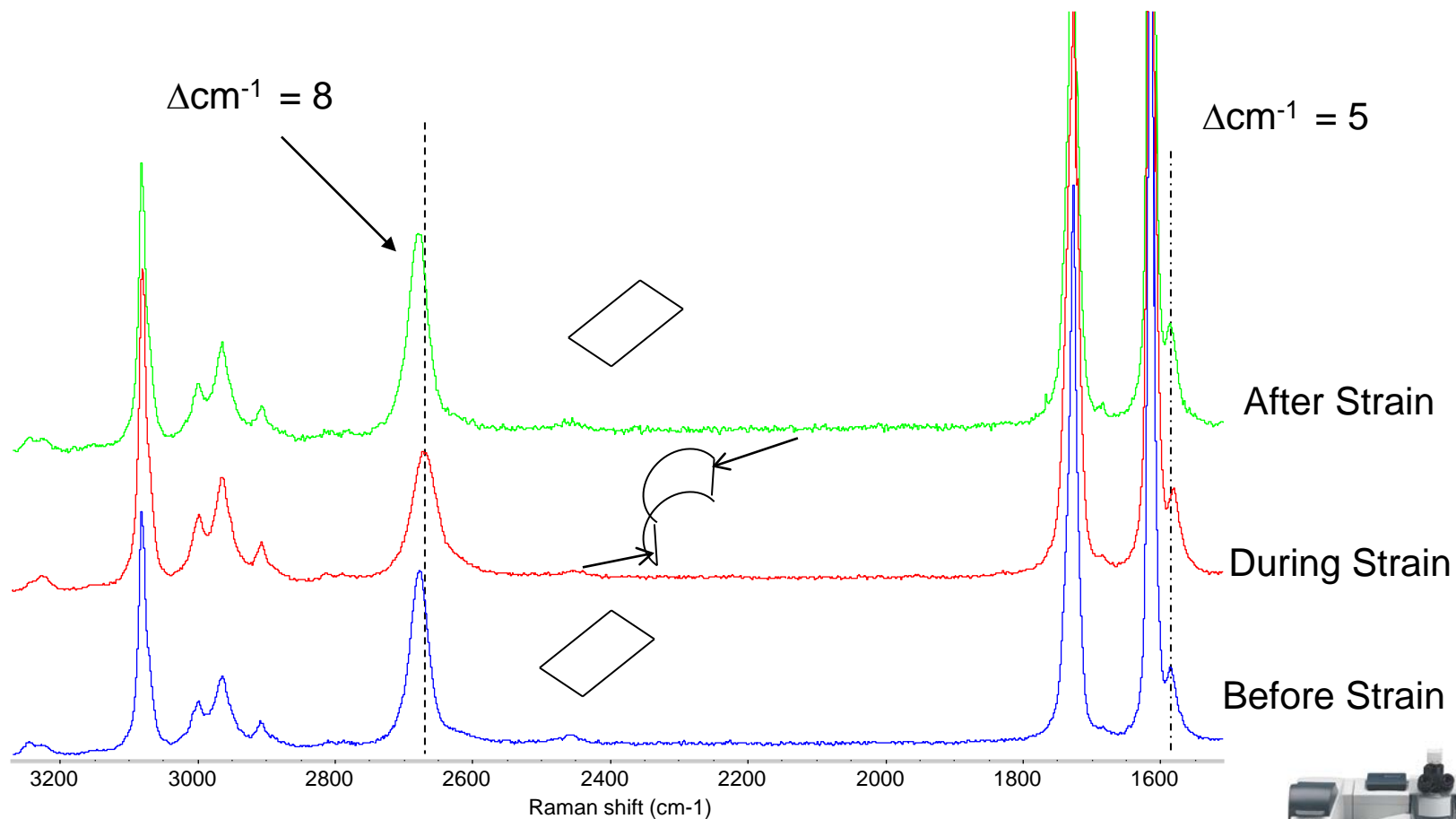


Wang, Hui; Cao, Xuwei; Feng, Min; and Lan, Guoxian
J. Raman Spectrosc. **2009**, *40*, 1791 – 1796

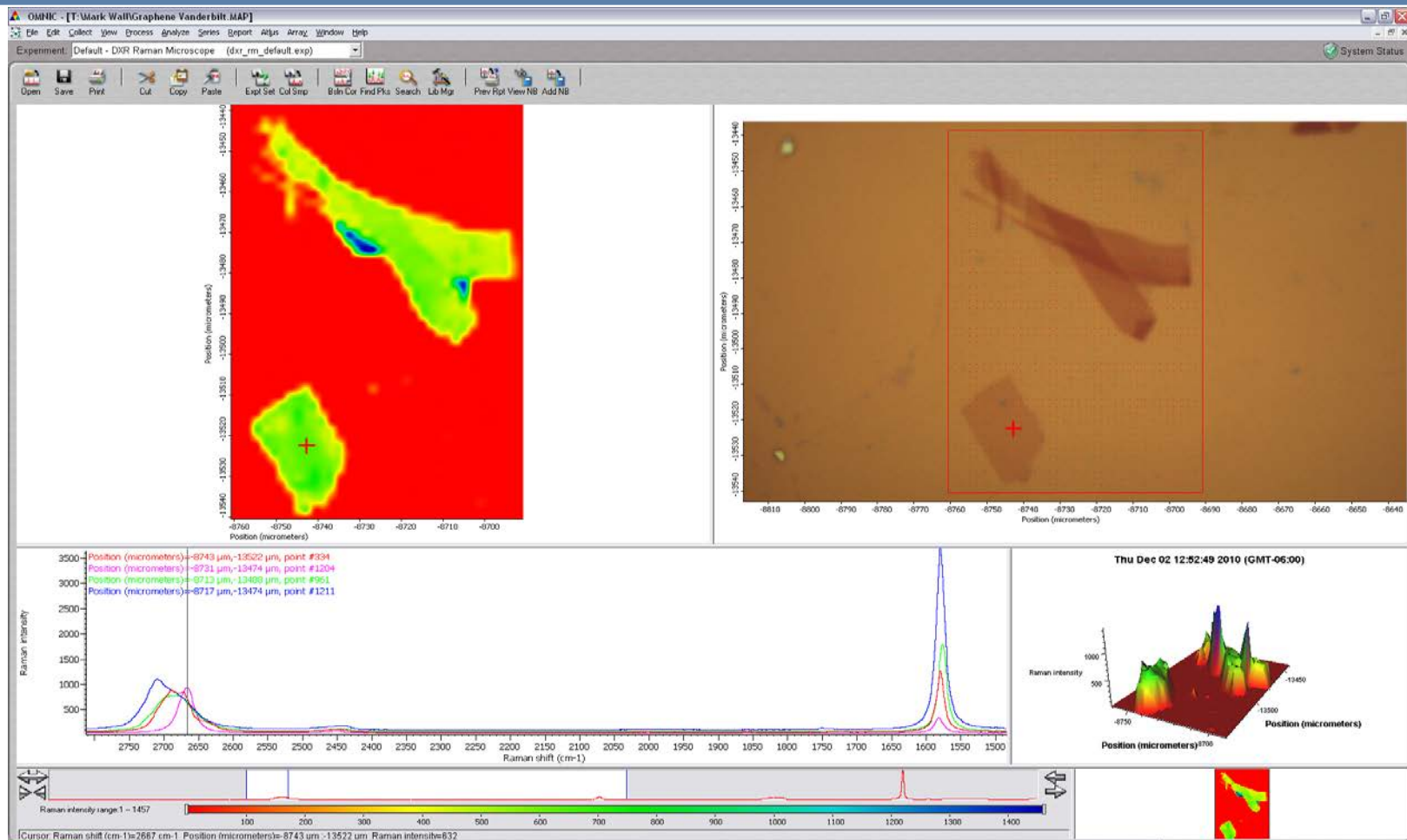
2D Band Shape Analysis



The Effect of Strain on Graphene

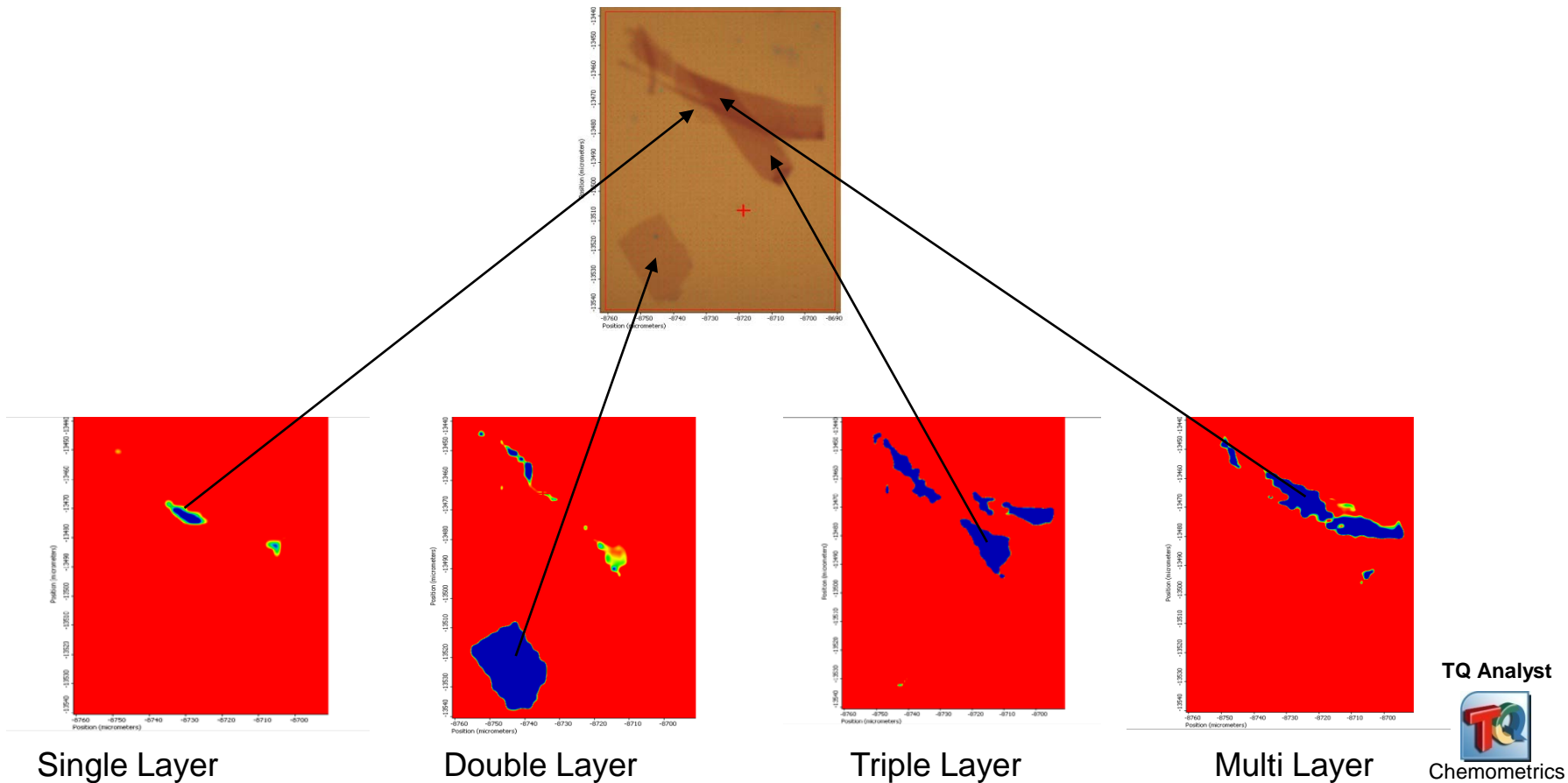


Graphene Characterization – Raman Mapping



Graphene Characterization – Raman Mapping

- Discriminant analysis results based upon the 2D band



The DXRxi Raman Imaging Microscope

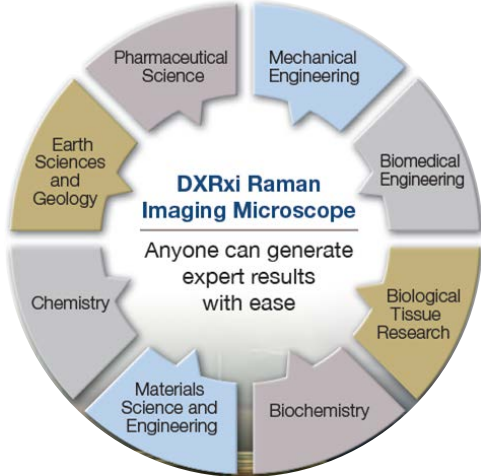
A total imaging system: hardware and software integration combines **powerful performance** with **image-centric** analysis and **ease of use**



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

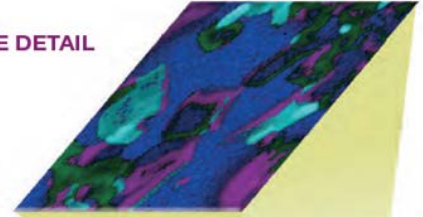
Raman Imaging – A New Tool for Rapid Characterization

The instrument should provide a tool for all users, from basic to advanced

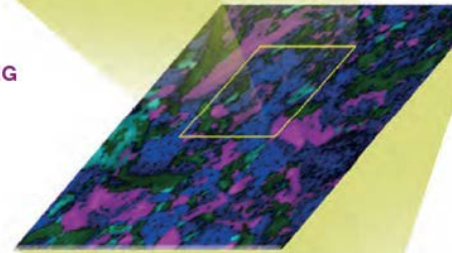


Advanced materials analysis benefits from an image-driven approach

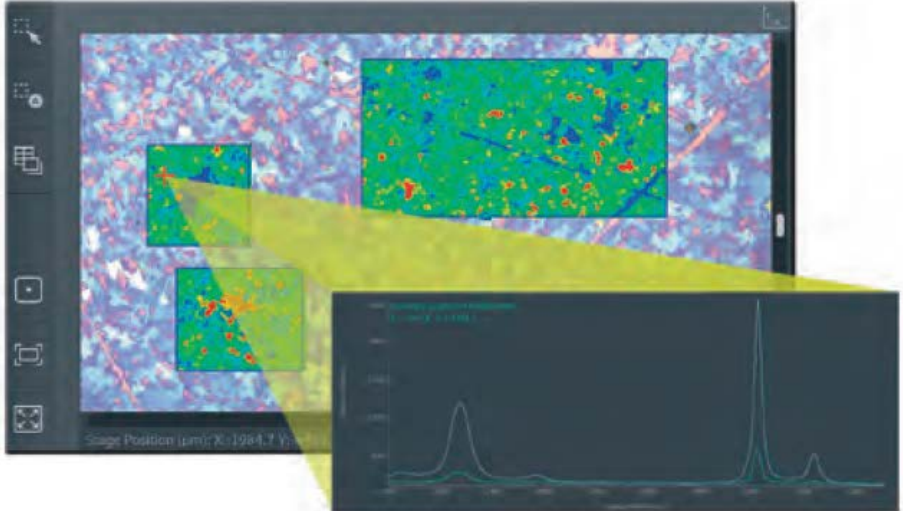
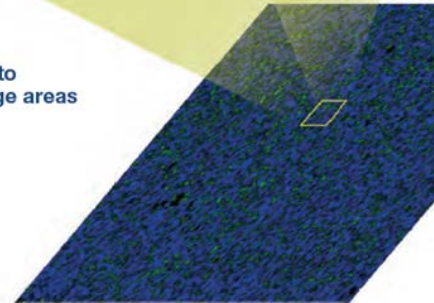
The performance advantage of **FINE DETAIL**



The agility of **RAPID IMAGING**



The **POWER** to look over large areas



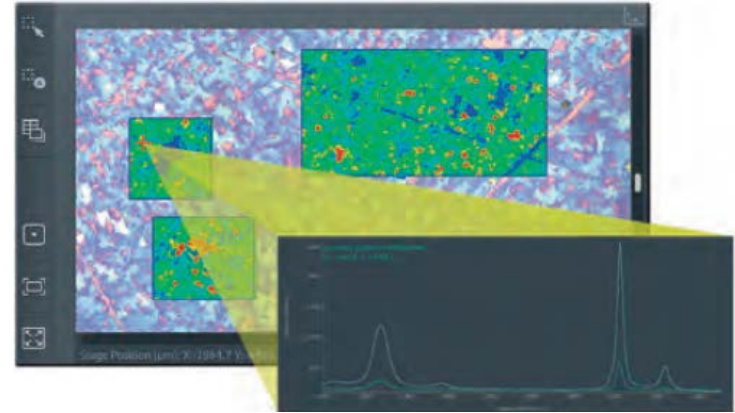
Emphasis on microscopy, powered by spectroscopy keeps the answer in focus

Intelligent Workflow with Excellent Flexibility

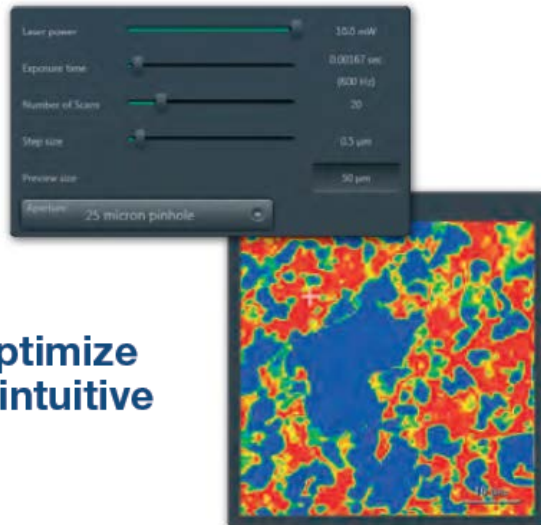
1 Rapid, single-click sample targeting



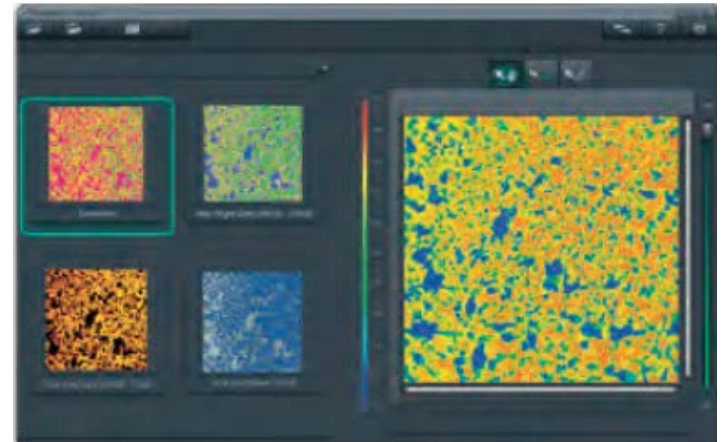
3 Quickly prioritize multiple regions of interest and run



2 Confidently optimize settings with intuitive controls

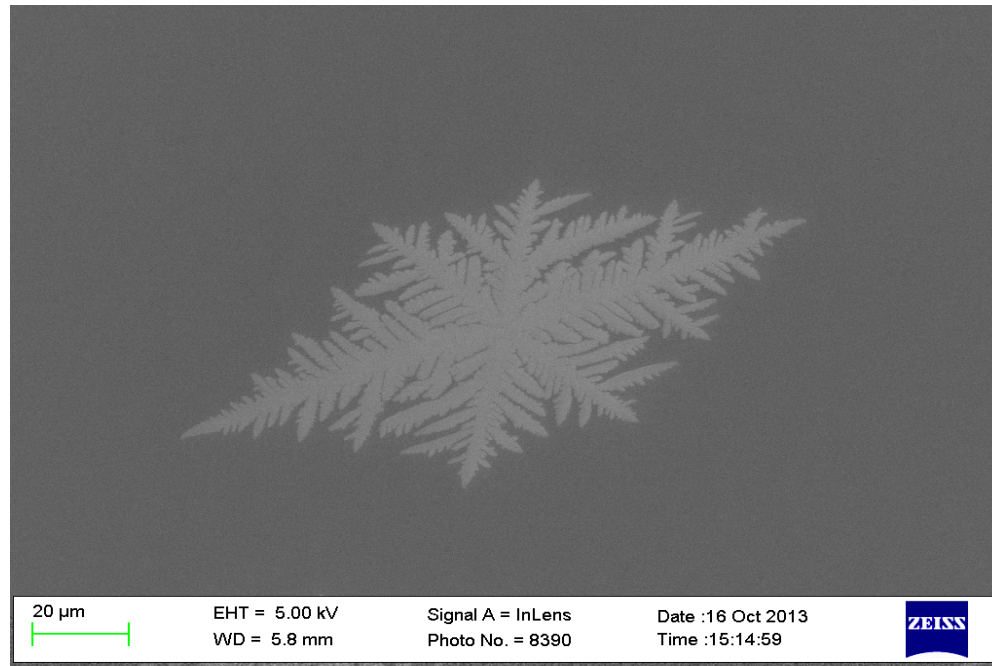


4 Information-rich images reveal a multitude of material characteristics



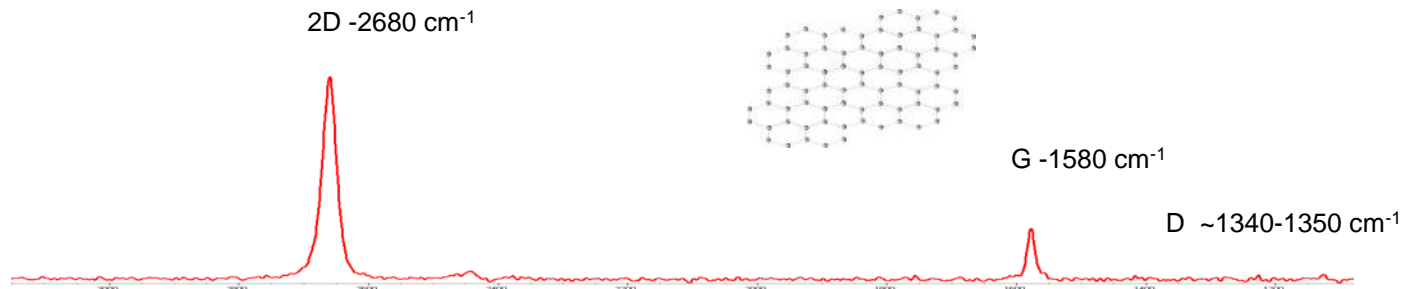
SEM Image of Dendritic Growth Graphene

- SEM illustrates the morphological characteristics that result from particular CVD deposition conditions
- Other techniques are required to obtain deeper insight



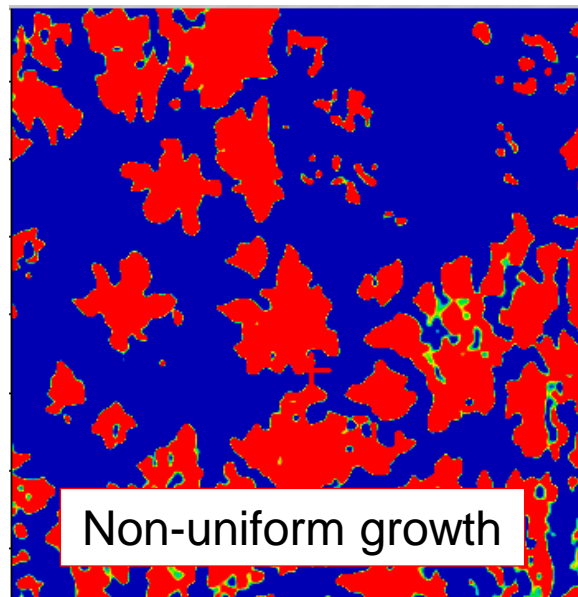
Rapid Raman Imaging of Partial Growth Graphene

- Provides vital information to understand growth dynamics
 - Allows determination of number of layers present
 - Provides information about the density of defects during growth
 - Information can be used to understand critical synthesis parameters and mechanisms that govern nucleation and growth of graphene
- Information from Raman can be used to optimize synthesis graphene
 - Minimize defect densities
 - Drive growth towards single crystal graphene
- Rapid imaging allows for bigger areas to be analyzed far faster than ever



Analysis of Graphene, 532nm laser

Ratio of the 2D to G band of graphene:



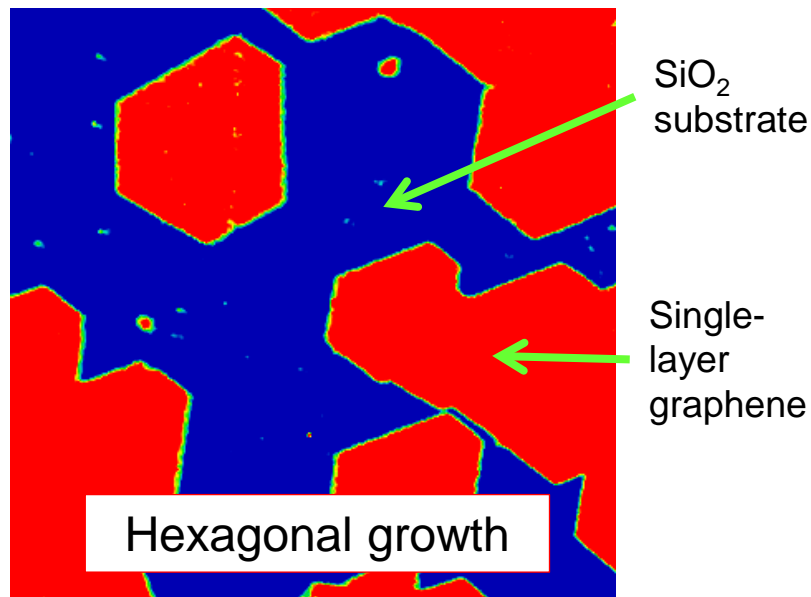
50 x 50 μm surface area

10,000 spectra!!!

Acquisition parameters: 100 Hz (10 ms/spectrum)

~16.7 minute collect time!!

MCR of graphene:



175 x 175 μm surface area

122,000 spectra!!!

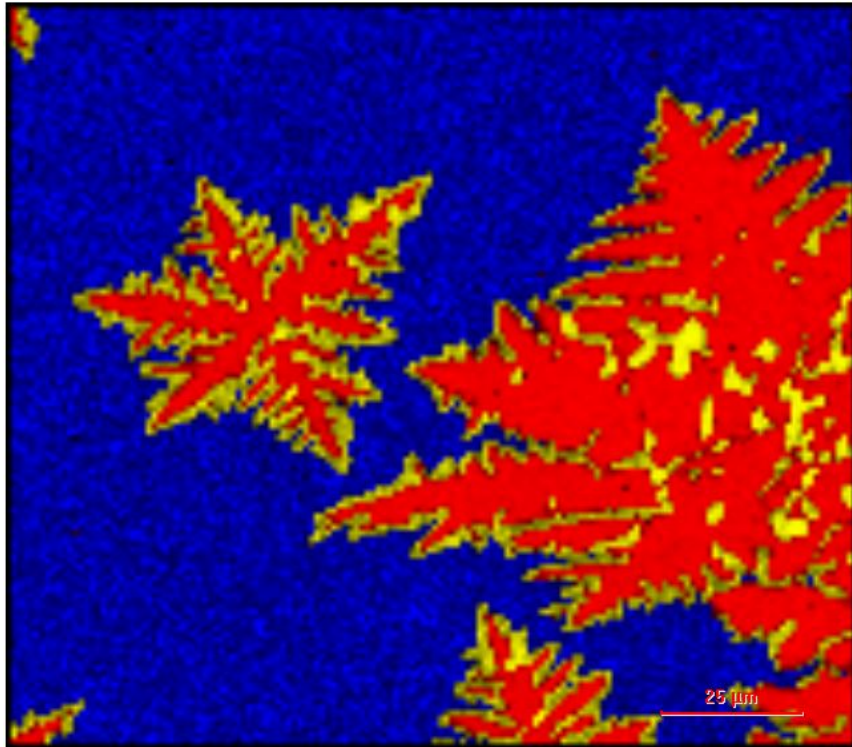
0.5 μm pixel size

Acquisition parameters: 100 Hz, 10 scans

~3.4 hour collect time!!

*Note: This graphene was initially grown on copper and then transferred to a silica substrate because the copper is a significant source of interference

Analysis of Graphene on a Copper Substrate



Graphene grown on copper and analyzed *on the copper* with the 455nm laser

$$I_{scatter} \propto \frac{1}{\lambda_{ex}^4}$$

455nm laser

150 x 150 μm surface area

90,000 spectra!!!

2.0 μm pixel size

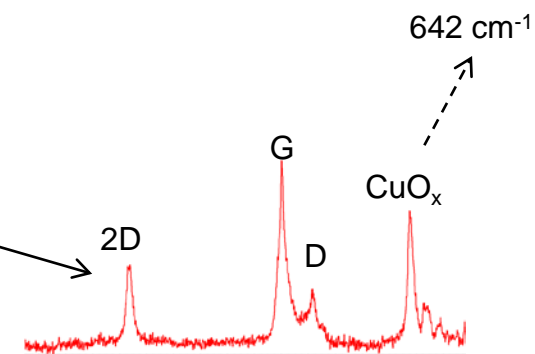
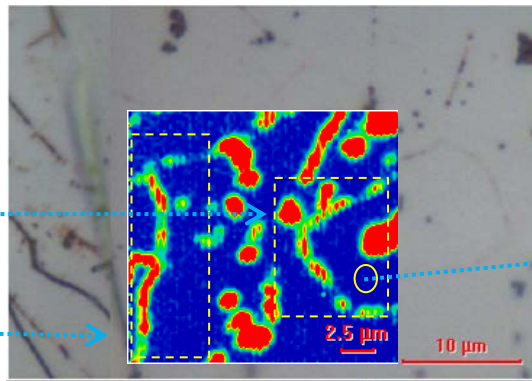
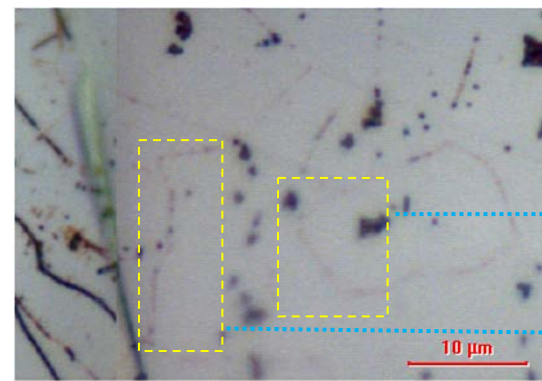
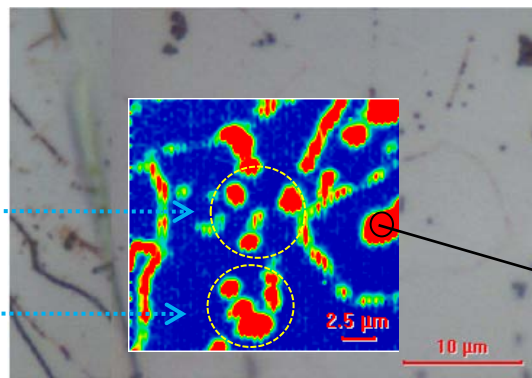
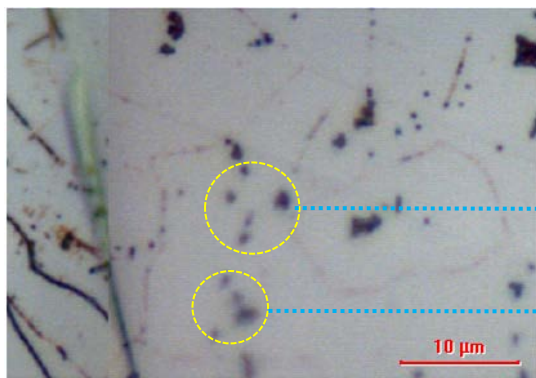
Acquisition parameters: 100 Hz (10 ms/spectrum),
4 scans

■ Copper Substrate ■ Oxidized Copper ■ Single-layer graphene

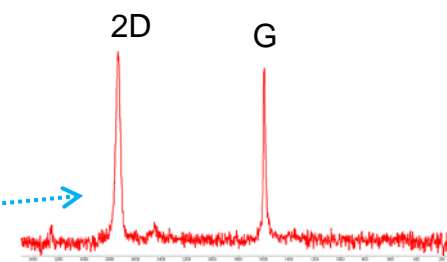


Graphene as a Protective Thin film

Chemical images based on Intensity of 642cm^{-1} band



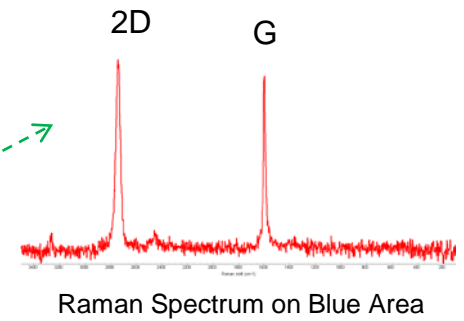
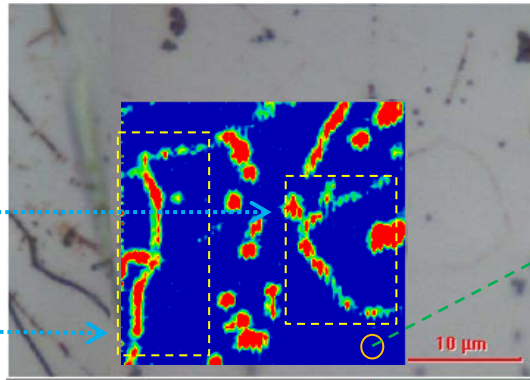
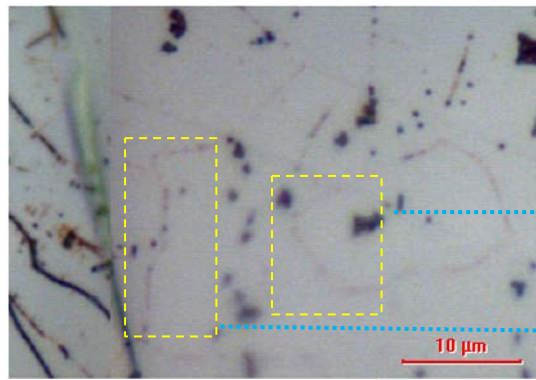
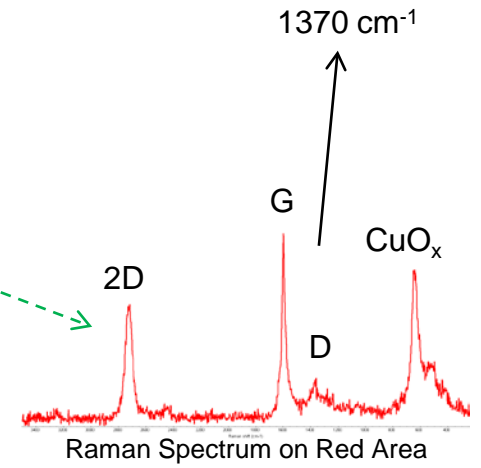
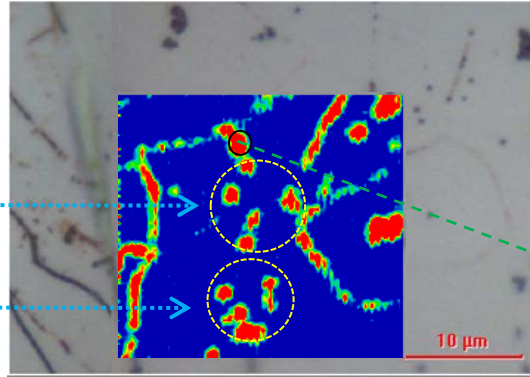
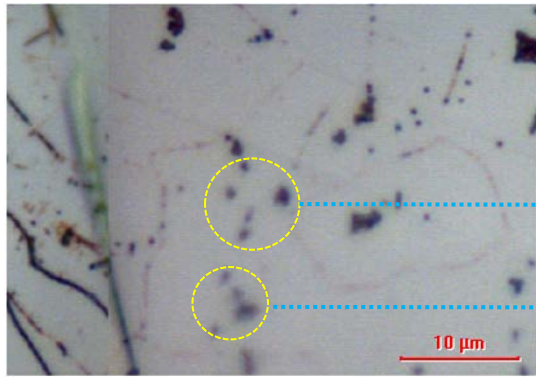
Raman Spectrum Red Area



Raman Spectrum Blue Area



Chemical images based on Intensity of 1370 cm^{-1} band



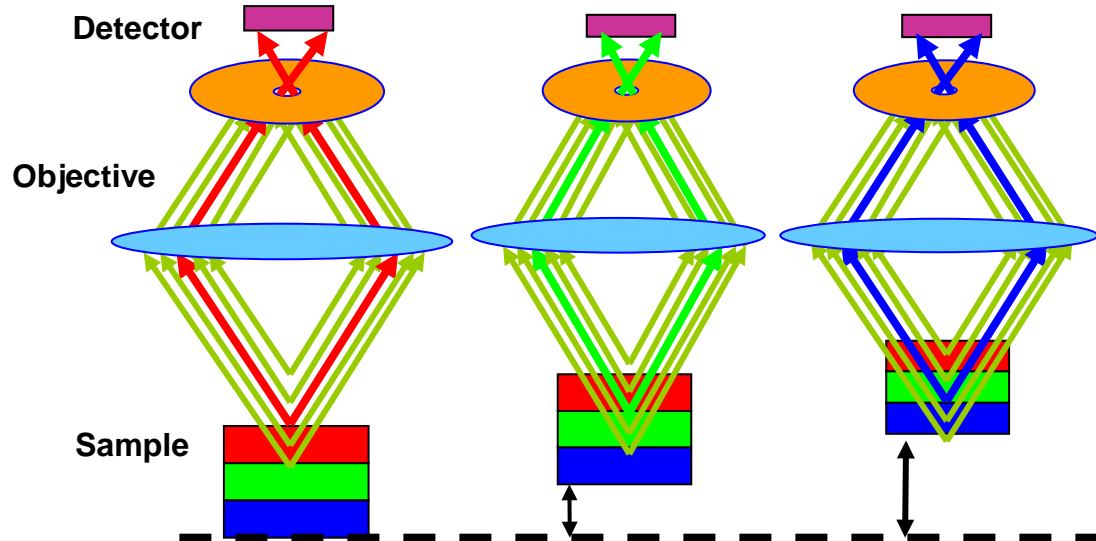
Optical Image Graphene on Cu

Optical Image/ Chemical Image Overlay Graphene on Cu



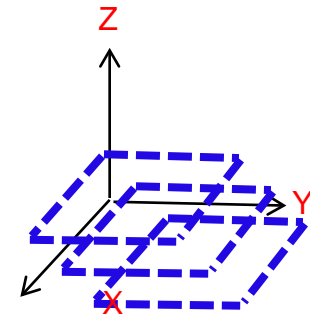
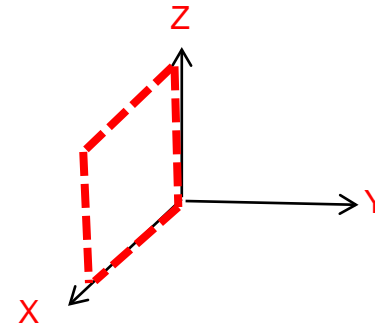
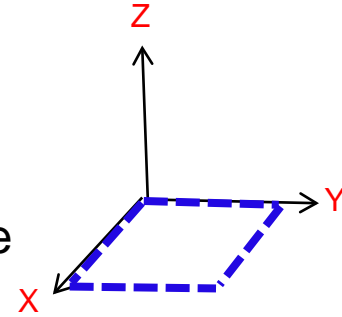
DXRxi Confocal Microscope

- DXRxi is a fully confocal microscope
 - Provides for higher spatial resolution
 - Achieves $<1\ \mu\text{m}$ measured x-y resolution
 - Provides for depth profiling of optically clear samples
 - Achieves $2\ \mu\text{m}$ depth resolution
 - Provides for increased rejection of fluorescence from substrate



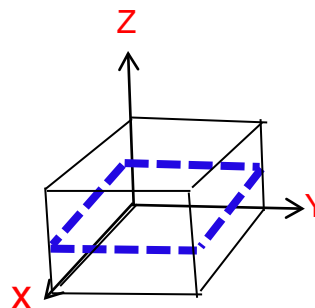
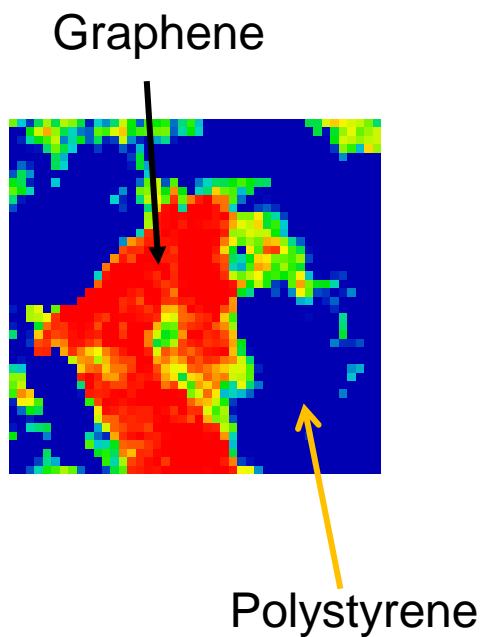
DXRxi Imaging -Sampling

- Raman/Chemical Images – Image Planes
 - XY Image plane
 - Surface image – most common image
 - Subsurface if the z position is below the surface
 - XZ Image plane
 - Confocal
 - Surface and subsurface
 - sample needs to be transparent
 - Also referred to as Cross-section
 - Image stack of XY Image planes
 - Confocal – sample needs to be transparent
 - Z axis image stack
 - Volume rendering capable



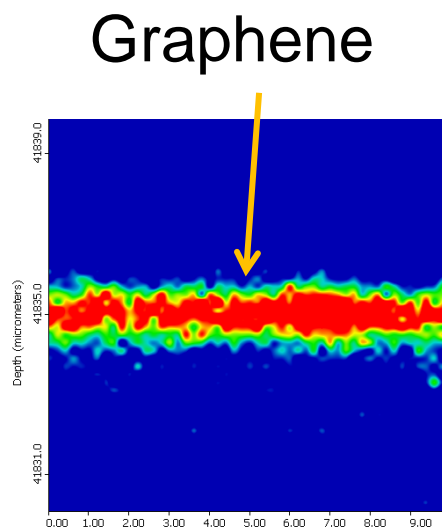
Confocal Raman Imaging Optical Cross Sectioning

Sub-surface xy Raman Image of Multilayer Graphene Flakes in Polystyrene

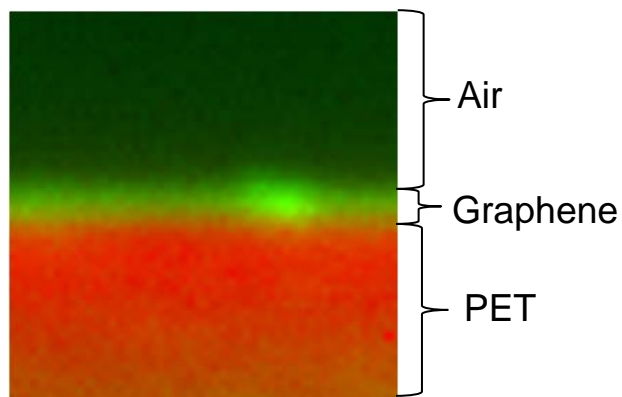


Cross-Section Analysis – Graphene on PET

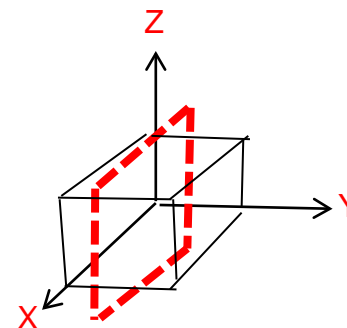
Sub-surface xz Raman Image of Single Layer Graphene on PET



Correlation Profile



MCR Profile

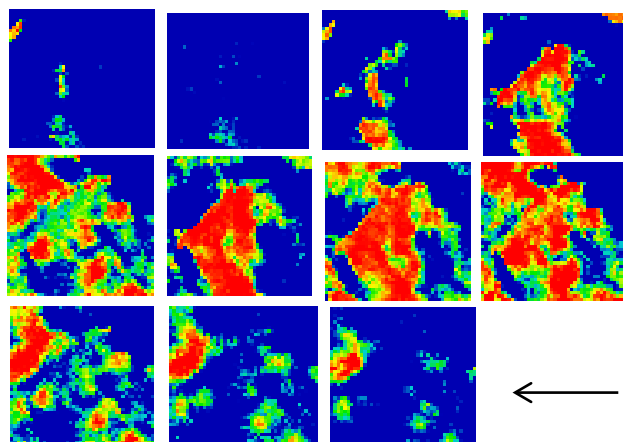


Multilayer Graphene Dispersed in Polystyrene Image Results

- Graphene was dispersed in Polystyrene at 1% w/w
- Confocal Raman images collected on Thermo Scientific DXRxi Raman Imaging Microscope
 - 455 nm excitation 3.7 mW laser power at sample
 - Raman spectra collected at a rate of 35 spectra/sec with 35 co-adds
 - 2 μM x 2 μM image pixel size
 - 11 xy chemical image planes (slices) were collected over an area of 82 μM x 86 μM - 20,000 spectra for the image stack

Top Image →

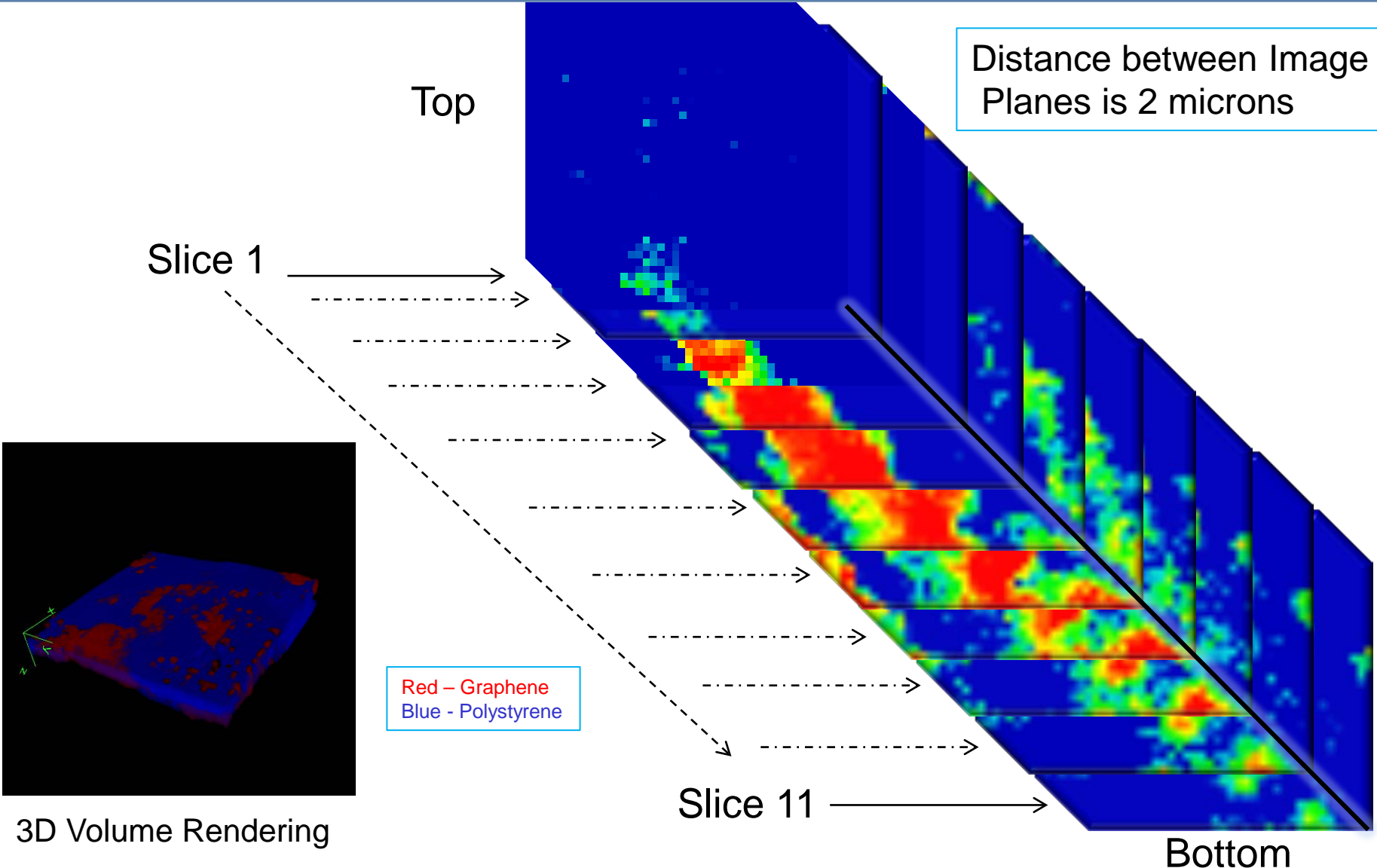
Image Intensity Based on the
2D Band of Graphene at 2711 cm^{-1}



← Bottom Image

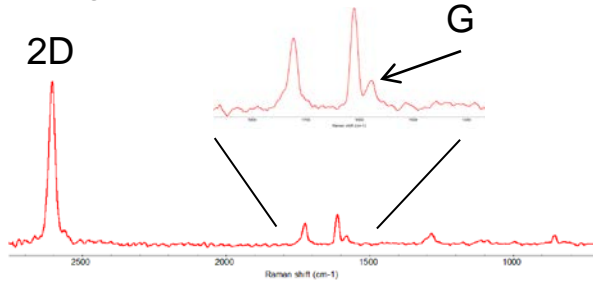


Graphene Dispersed in Polystyrene Z –Stack Slices

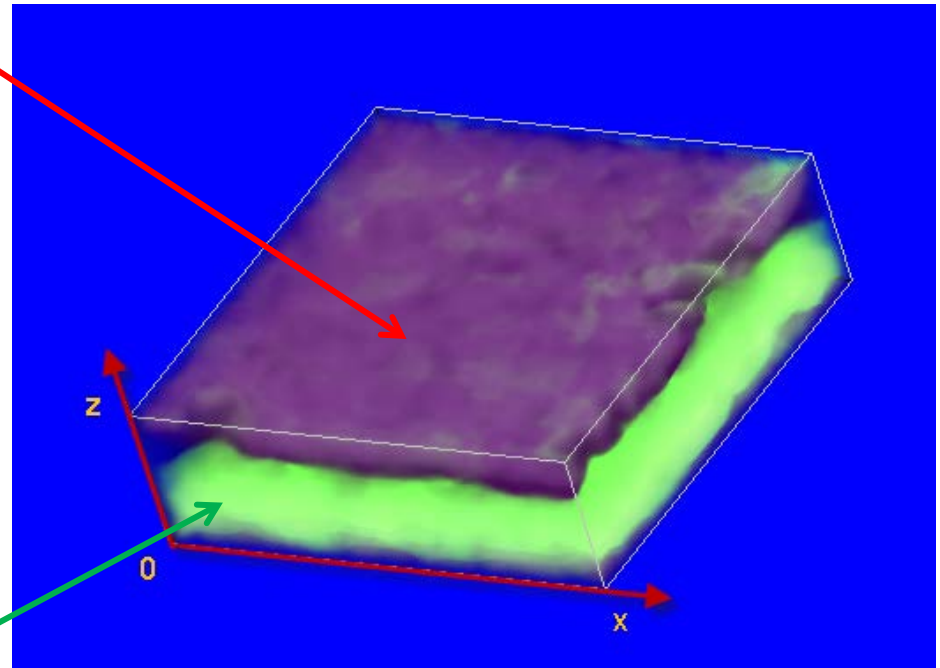
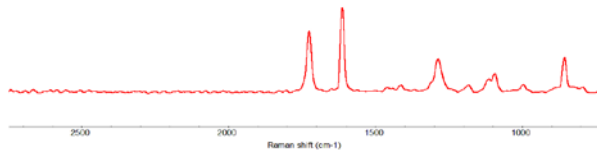


Raman Imaging – Confocal Volume Rendering

Single Layer Graphene



(Poly)ethylene terephthalate



Summary

- *Raman Spectroscopy was presented demonstrating the usefulness of Raman Spectroscopy for characterizing Graphene*
- Modern instrumentation was discussed
- Raman imaging examples were presented
 - Partial growth graphene
 - Graphene as a protective coating material
- Confocal Raman imaging of graphene
 - Cross sectional and Z-stack measurements

Questions?

Thank You for our time and attention!

mark.wall@thermofisher.com