

# Fast structure determination for proteins and small molecules – a MicroED solution


Abhay Kotecha, Bart Buijsse, Michael Janus, Lingbo Yu, Hans Raaijmakers

## ABSTRACT

In this session, you will learn how micro electron diffraction (MicroED) allows fast, high-resolution 3D structure determination of small chemical compounds and biological macromolecules. To efficiently collect diffraction datasets of nano-crystals, a cryo-TEM is equipped with a specially designed diffraction camera (Ceta-D) and a MicroED package. The latter combines the necessary hardware components, as well as optimized optical settings and specialized EPU-D Software for automated data collection. Combined with the intrinsic microscope performance, the data collection is fully automated and can be realized in a matter of minutes.

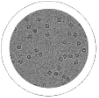
## INTRODUCTION

MicroED is a new technique for structure determination of biological macromolecules and small molecules:



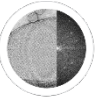
**Cryo-Electron Tomography**

- Whole cells, organelles
- Biomolecules in their native context



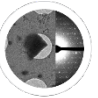
**Single Particle Analysis (SPA)**

- Purified proteins, complexes
- Purified biomolecules, different conformations can be observed



**2D Electron Diffraction**

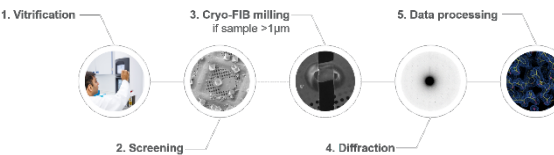
- 2D crystals
- Membrane protein – lipid bilayer interaction can be seen



**MicroED**

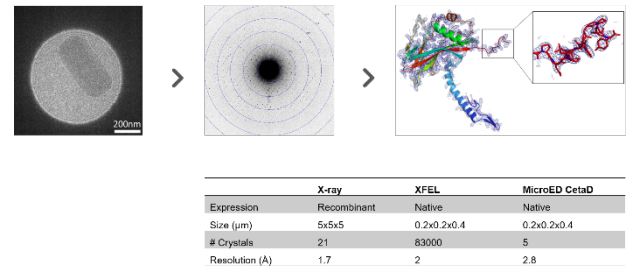
- 3D small crystal
- Highest resolution cryo-EM technique (1-2 Å)

The MicroED workflow consists of four or five steps depending on the size of the crystals under investigation:



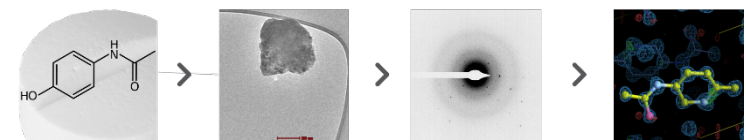
## EXAMPLES

Figure 1. MicroED of nanocrystals - Granulovirus



Granulin, is a small protein that forms the coat of granulovirus. This protein is of size 29.4 kDa (245 amino acids) and forms nanocrystals (200-400nm in size) containing on average 9000 unit cells, with cubic symmetry and a unit cell of 103-Å. This corresponds with a volume of the diffracting body of less than 0.02 µm<sup>3</sup>. This is an order of magnitude smaller than what was possible so far using protein nanocrystals in electron diffraction and has only been achieved using XFEL or large recombinant crystals by X-rays.

Figure 2. MicroED of small molecules and natural compounds



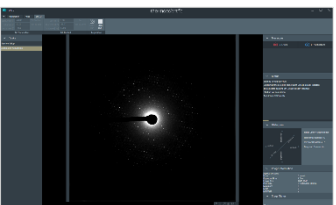
MicroED of small molecules is fast, orthogonal to NMR and mass spectrometry, and sample prep is very easy (~10<sup>-12</sup>g of sample, even mixtures). Shown here, the structure of paracetamol was obtained from a ground-up tablet. The contained crystals, diffracted to 0.8 Å and were collected in less than 2 min. The structure was obtained by direct phasing.

## EQUIPMENT

Table 1. Existing microscopes can be upgraded easily

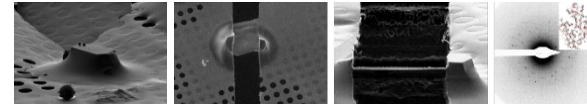
	Glacios/Talos	Krios
<b>1 Ceta-D camera</b>		
• Optimized for diffraction application: increased accuracy and sensitivity	✓	✓
• Compatible with SPA screening requirements	✓	✓
• Compatible with bottom-mount filter (refractable)	✓	✓
<b>2 MicroED package</b>		
Data acquisition SW	✓	✓
Modified beam stop, optimized for MicroED application	✓	✓
Optimized C2/SA aperture set	✓	✓
MicroED lens series	✓	✓
90° rotation projection system	✓	✓

Figure 3. EPU-D software for MicroED



Dedicated MicroED software with a focus on automation and through-put with the same ease-of-use EPU is known for.

Figure 4. Cryo-FIB milling of microcrystals



Intermediate-sized crystals (1-50 µm) can be thinned to make them suitable (150-600 nm) for MicroED while preserving the internal order of the crystal lattice. Lysozyme, used as a test system, resolves to 1.9 Å in cryo-FIB milled crystals.

## CONCLUSIONS

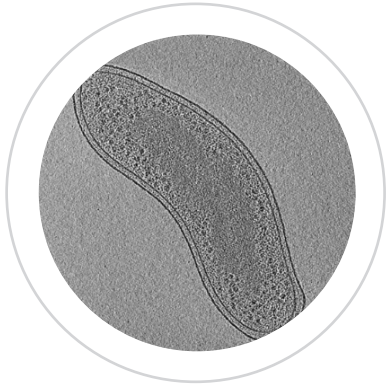
- Fast atomic-resolution 3D structural information.** Diffraction data from nanocrystals in minutes.
- Instant productivity.** Nanocrystals as small as 100 nm can be readily analyzed, removing the burden of growing large crystals (as with X-ray crystallography). Also reduces the amount of sample material required. Mixtures of different polymorphs and compounds can be analyzed.
- Complete turnkey solution.** Including hardware, software and support from one single vendor. Acquired data can be readily processed using established reconstruction packages for X-ray crystallography.
- 2-in-1 solution.** MicroED and single particle analysis (SPA) can be performed on the same cryo-electron microscope. This solution is compatible with new microscopes but is also retrofittable on existing units.

## FURTHER INFORMATION

Please scan this QR-code to get to the website containing the MicroED introduction video as well as a download link for the PDF describing the MicroED package and the Ceta-D camera.

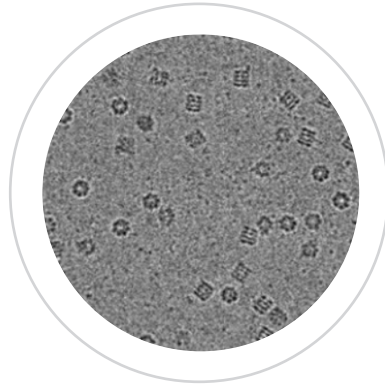


# MicroED is a New Technique for Structure Determination of Biological Macromolecules



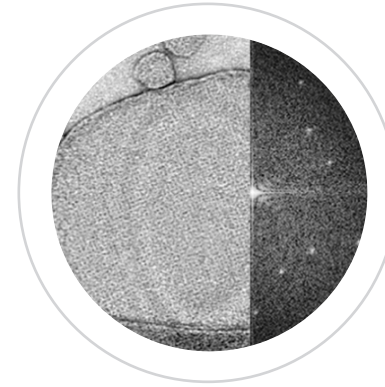
## Cryo-Electron Tomography

- Whole cells, organelles
- Biomolecules in their native context



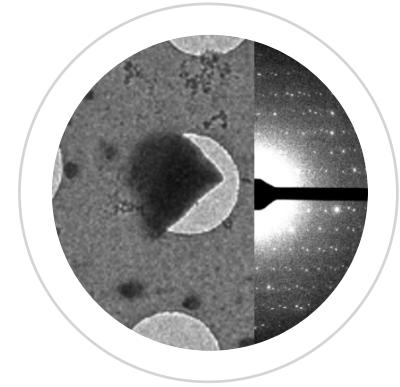
## Single Particle Analysis (SPA)

- Purified proteins, complexes
- Purified biomolecules, different conformations can be observed



## 2D Electron Diffraction

- 2D crystals
- Membrane protein – lipid bilayer interaction can be seen

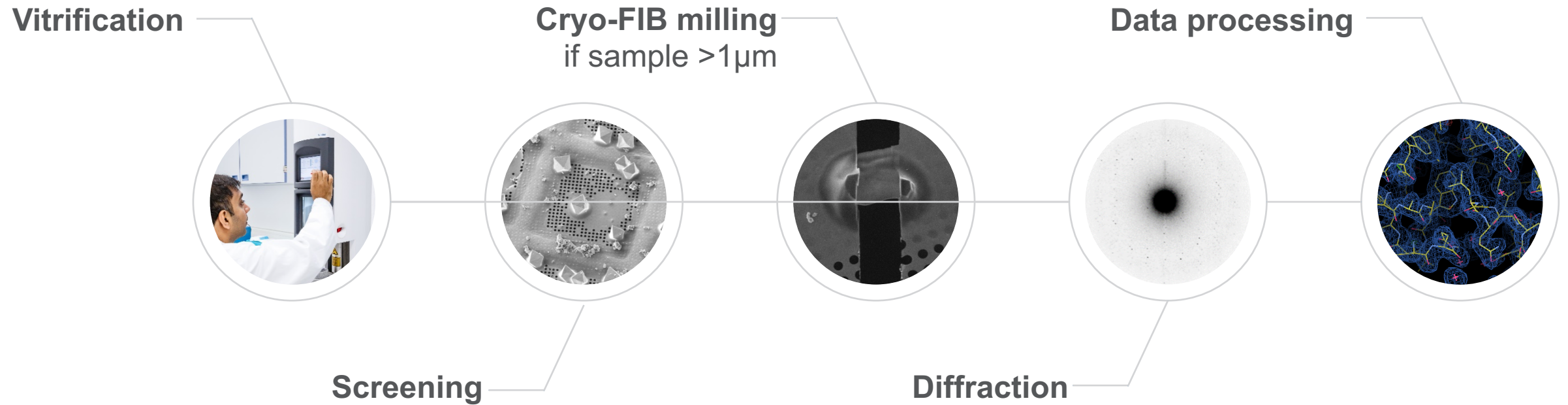


## MicroED

- 3D small crystal
- Highest resolution cryo-EM technique (1-2 Å)



# The MicroED Workflow

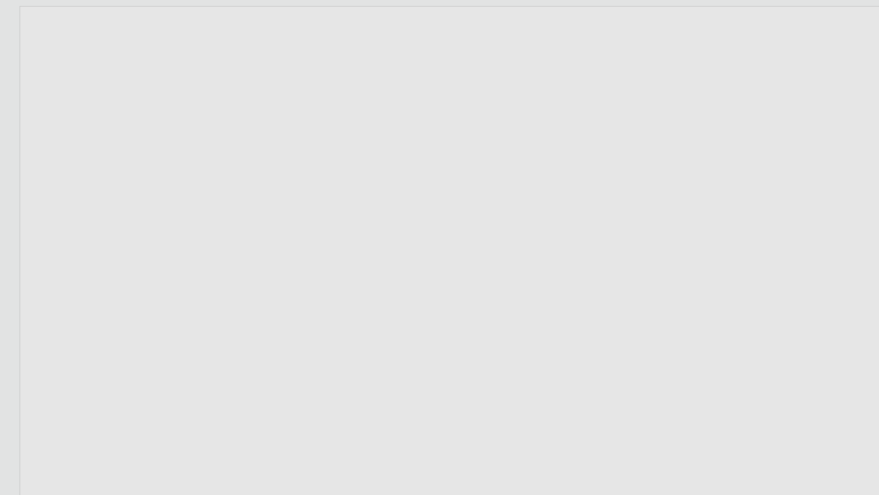
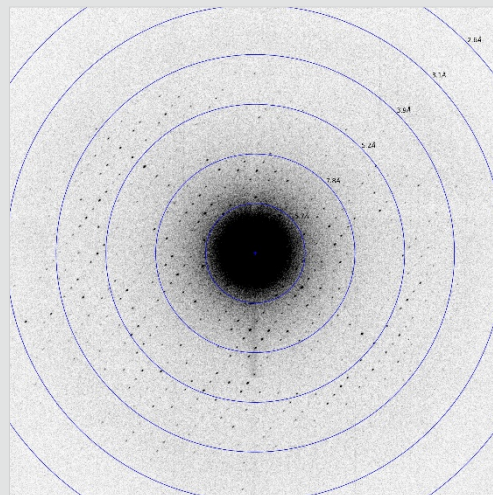
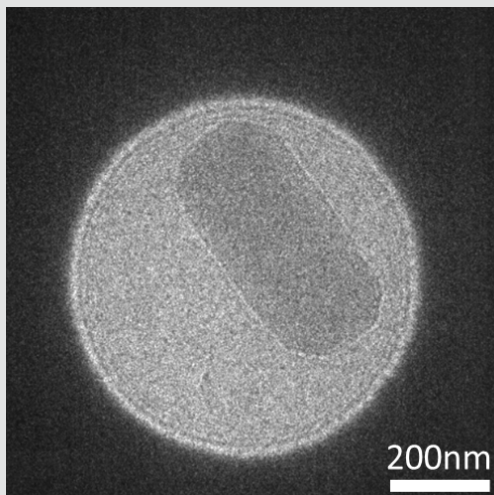




## MicroED offers unique advantages

- Provides **high resolution**
- **Fast** method
- **Small crystals**
- Needs **few crystals**
- **Mixtures** allowed

# Resolving Nano-crystals of Granulovirus

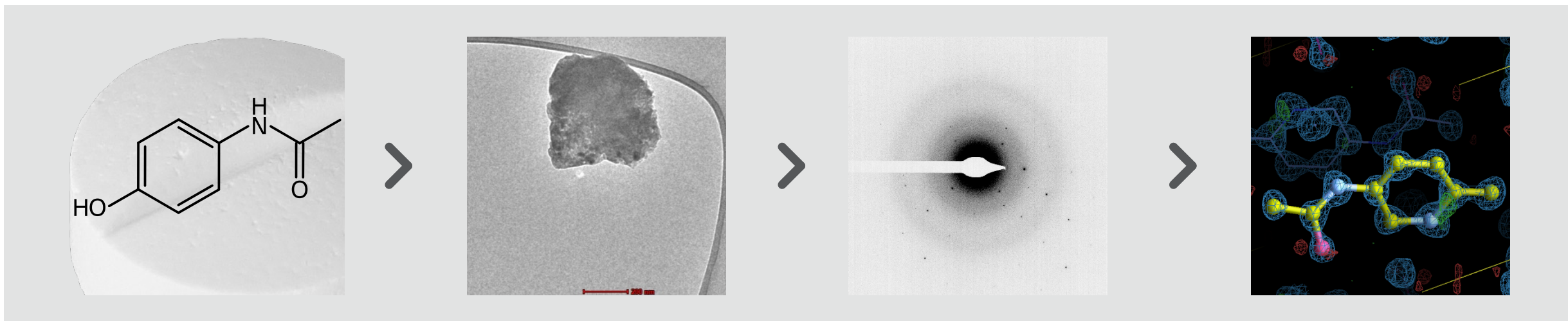


## Granulin

Protective virus coat  
**29 kDa** protein  
9,000 unit cells/virion  
**0.5 hr data collection** on  
5 crystals

	X-ray	XFEL	MicroED CetaD
Expression	Recombinant	Native	Native
Size (μm)	5x5x5	0.2x0.2x0.4	0.2x0.2x0.4
# Crystals	21	83000	5
Resolution (Å)	1.7	2	2.8

# Structure Determination of Drugs and Natural Compounds Using MicroED



## Sample

Grinded paracetamol

## Prep.

Lacey carbon grid,  
cryo temperature

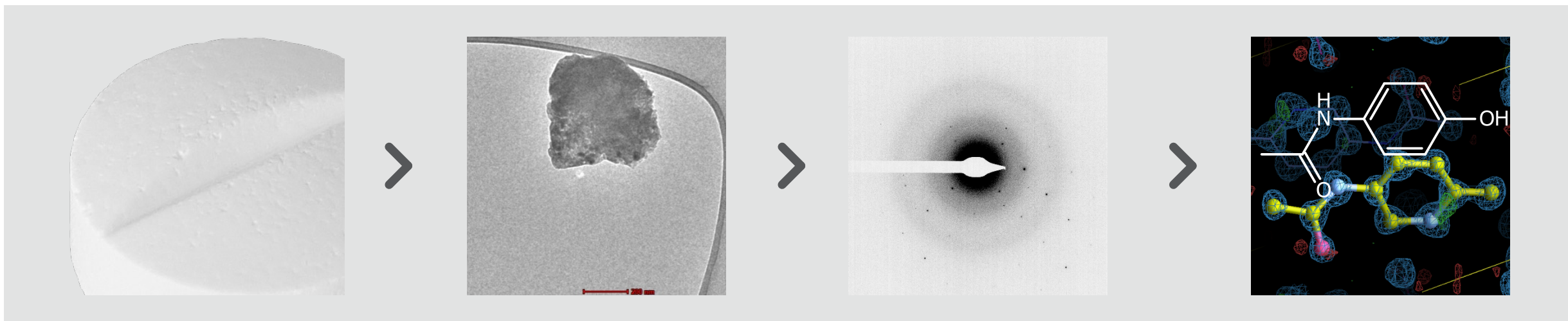
## Acquisition

160 x 1.0° x 1sec  
0.88Å

## Analysis

Dials → shelxt  
70% complete

# Structure Determination of Drugs and Natural Compounds Using MicroED



## Sample

Grinded paracetamol

## Prep.

Lacey carbon grid,  
cryo temperature

## Acquisition

160 x 1.0° x 1sec  
0.88Å

## Analysis

Dials → shelxt  
70% complete



## Small molecule structure determination

- Orthogonal to NMR, mass spec etc.
- Easy sample prep
- $\sim 10^{-12}$ g of sample
- Data collection <3 min
- Direct phasing

# MicroED is Enabled by the Ceta-D Camera and the MicroED Package

Existing microscopes can be upgraded easily

## ① Ceta-D camera

- Optimized for diffraction application: increased accuracy and sensitivity
- Compatible with SPA screening requirements
- Compatible with bottom-mount filter (retractable)

## ② MicroED package

EPU-D: Data acquisition SW  
Modified beam stop, optimized for MicroED application  
Optimized C2/SA aperture set  
MicroED lens series  
90° rotation projection system

**Glacios/Talos**

**Krios**

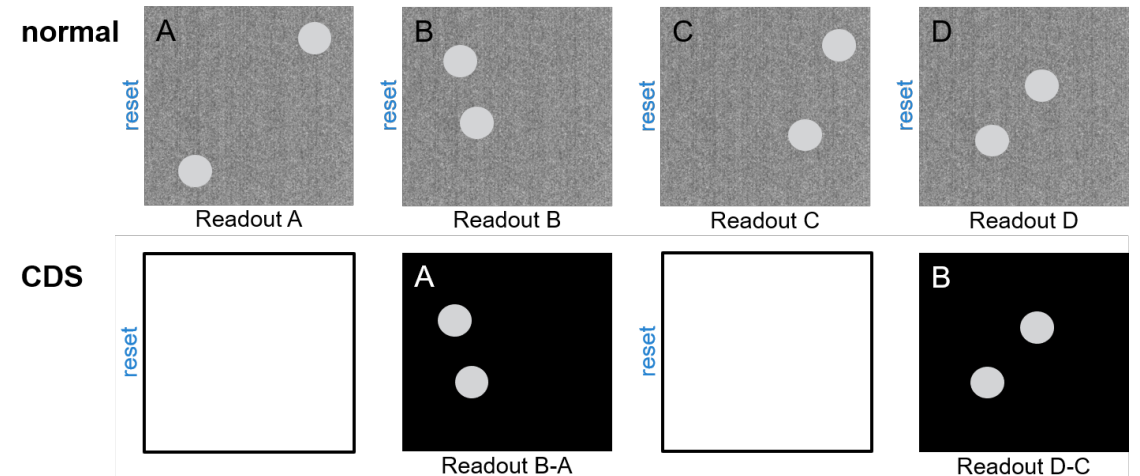
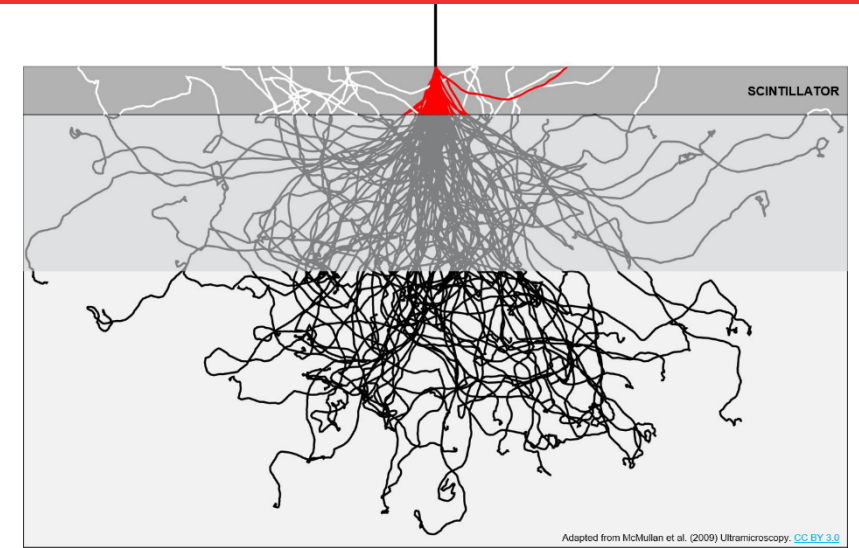


# Ceta-D Camera

Scintillator-based detection with a CMOS sensor **designed for MicroED**

**The Ceta-D offers improved signal-to-noise and enables continuous tilt diffraction**

- **4x higher sensitivity compared to Ceta:** Optimized scintillator thickness for enhanced signal-to-noise ratio and improved DQE at low frequency.
- Movie mode is enabled for **continuous-tilt diffraction**.
- **Correlated Double Sampling (CDS)** in movie mode **decreases noise 2.5x**.

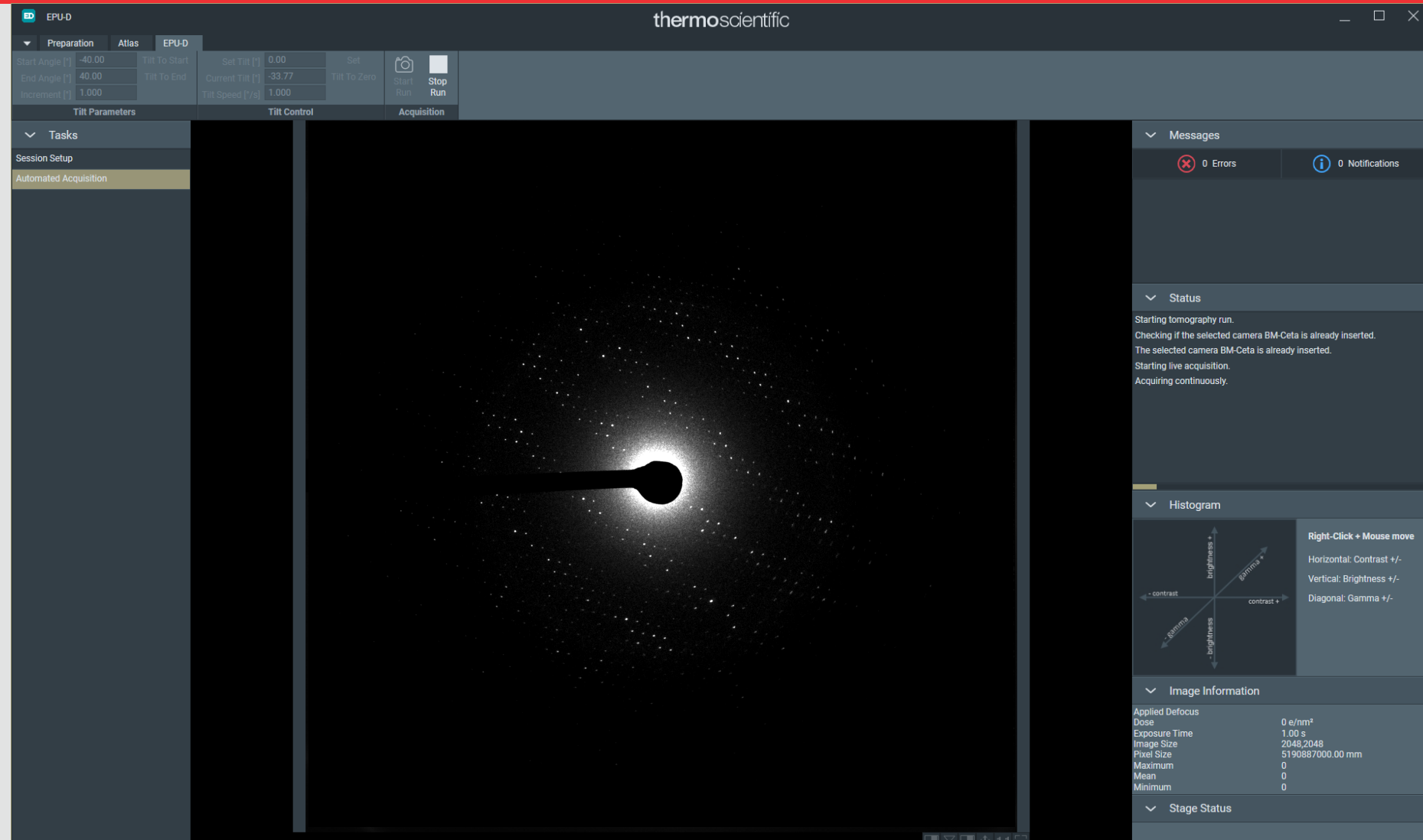


# EPU-D Acquisition Software

Dedicated MicroED software with a focus on **automation, usability and throughput**

**EPU-D facilitates MicroED with the same ease-of-use EPU is known for**


- Intuitive workflow **GUI** for navigation, set-up and acquisition.
- **Fast and easy setup** of MicroED experiments
- **Continuous tilt series** acquisition.




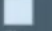
Preparation

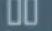
Atlas

EPU-D









Start Position

Close to centre

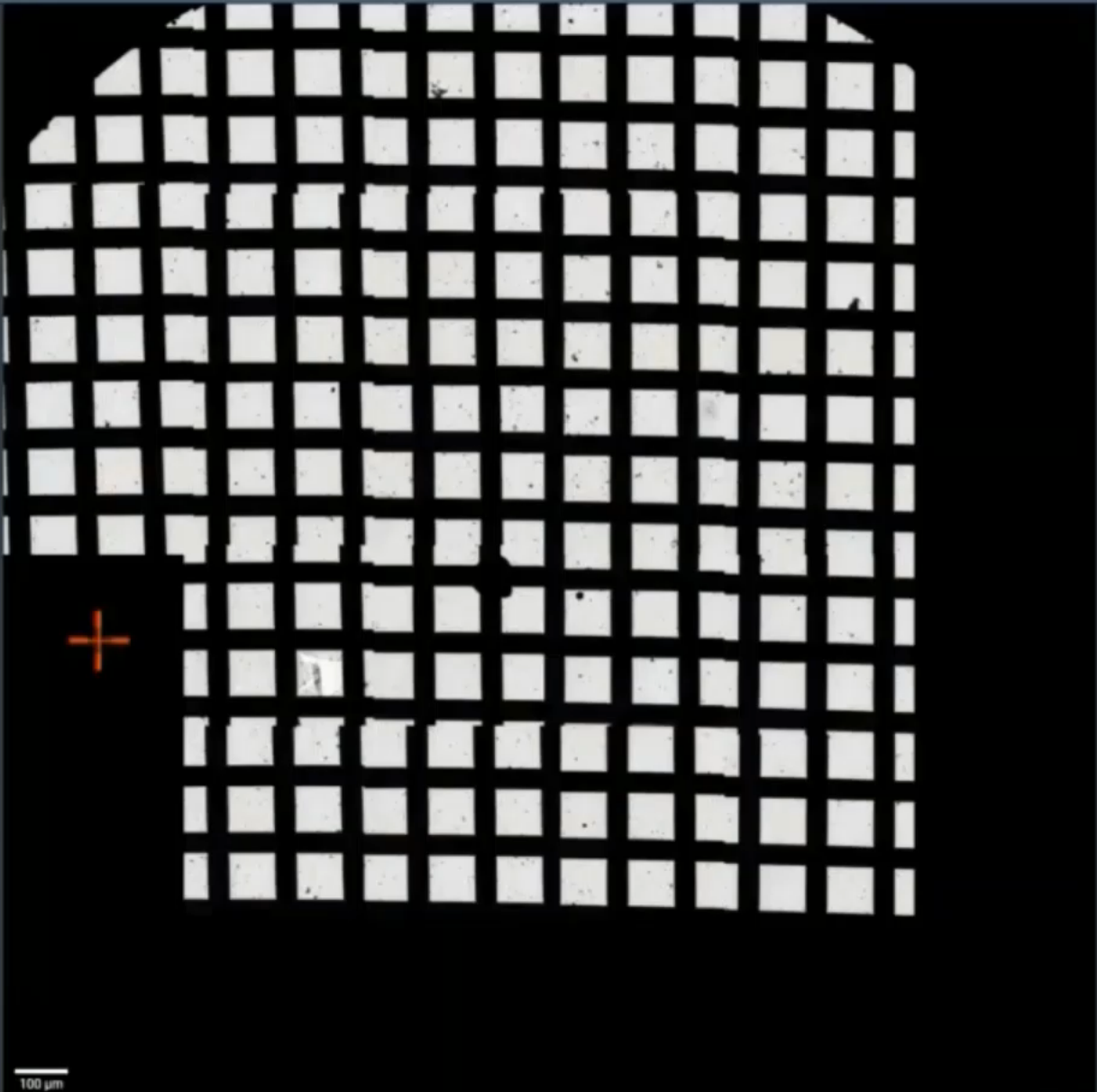
Acquisition

Atlas Settings

Tasks

Session Setup

Atlas Acquisition



Messages

 0 Errors

 0 Notifications

Status

- Postprocessing tile
- Moving stage to next position
- Set the defocus
- Acquiring tile
- Reset the defocus
- Postprocessing tile
- Moving stage to next position
- Set the defocus
- Acquiring tile
- Reset the defocus
- Postprocessing tile
- Moving stage to next position

Histogram

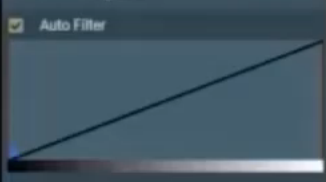


Image Information

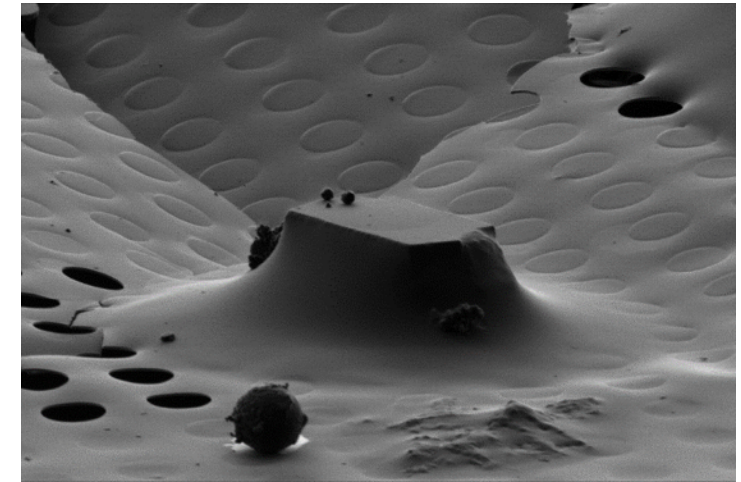
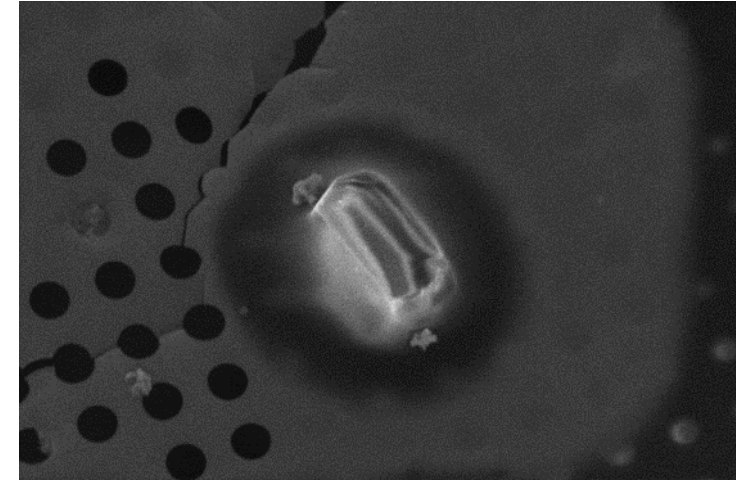
Applied Defocus	-500.00 μm
Dose	0
Exposure Time	2.00 s
Image Size	4096,4096
Pixel Size	505.00 nm
Maximum	562
Mean	6.31
Minimum	0

# Aquilos Cryo-FIB

Cryo-FIB milling makes micro-crystals usable for MicroED

**Intermediate-sized crystals can be thinned to make them suitable for MicroED while preserving the internal order of the crystal lattice.**

- Crystal thickness for MicroED is limited to ~700nm.

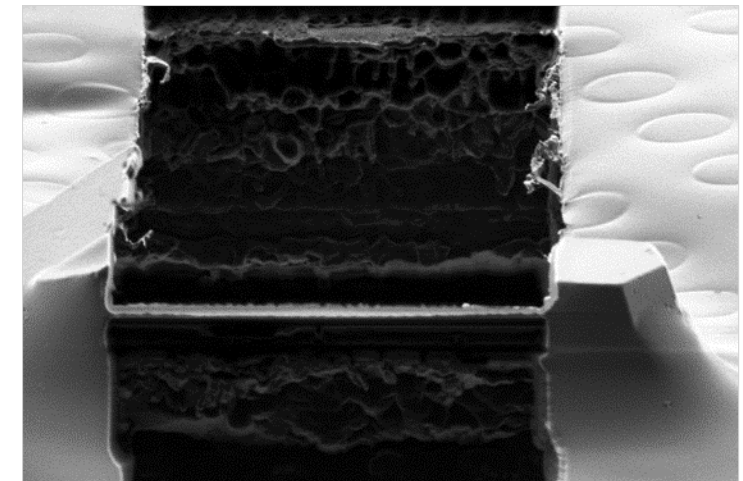
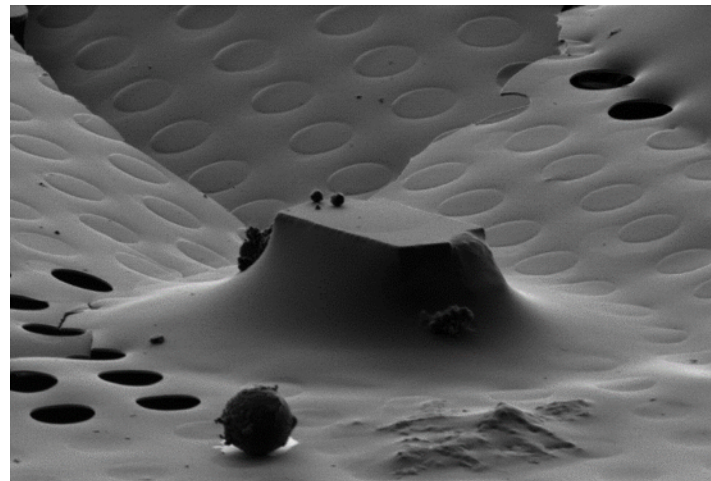
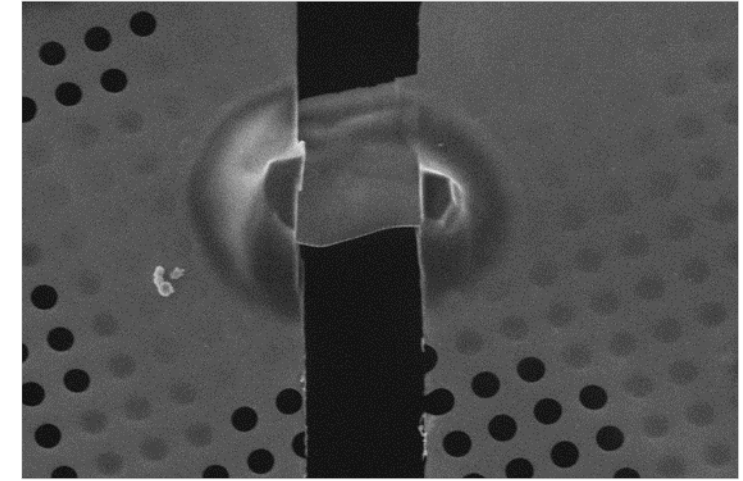
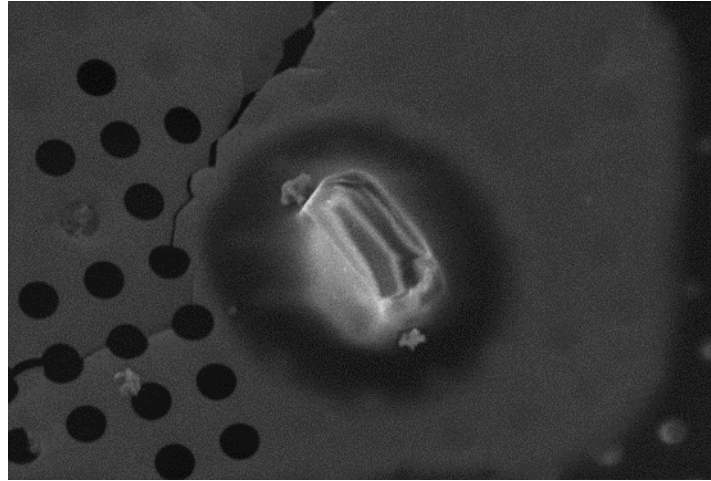


# Aquilos Cryo-FIB

Cryo-FIB milling makes micro-crystals usable for MicroED

**Intermediate-sized crystals can be thinned to make them suitable for MicroED while preserving the internal order of the crystal lattice.**

- Crystal thickness for MicroED is limited to ~700nm.
- Crystals ranging from 1 $\mu$ m to 50 $\mu$ m thickness can be thinned 150-600nm using cryo-FIB milling.

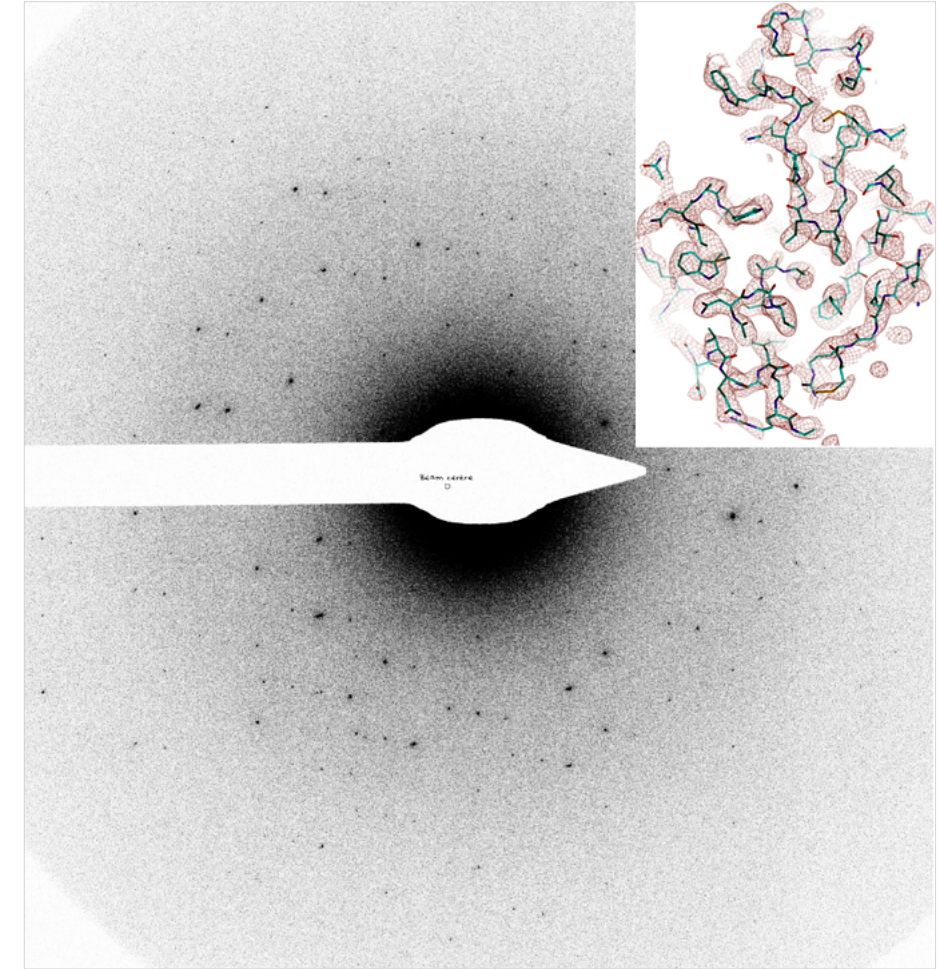


# Aquilos Cryo-FIB

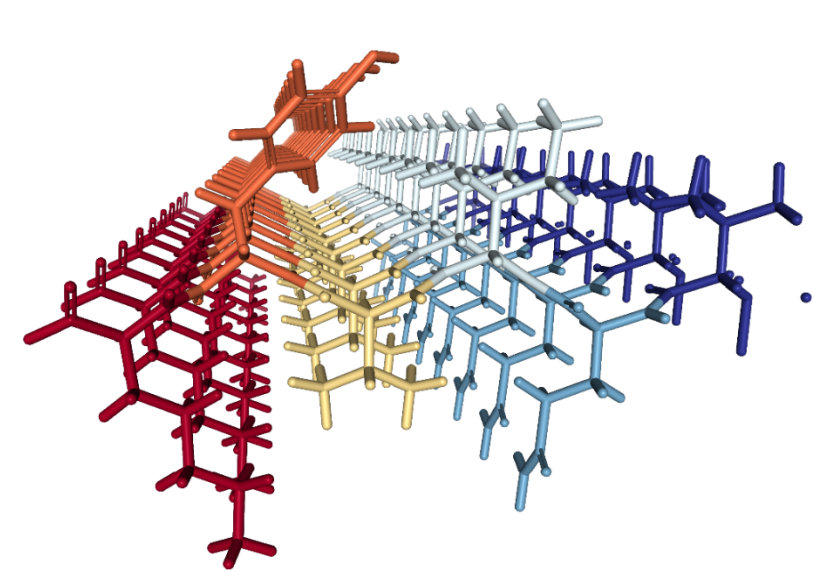
Cryo-FIB milling makes micro-crystals usable for MicroED

**Intermediate-sized crystals can be thinned to make them suitable for MicroED while preserving the internal order of the crystal lattice.**

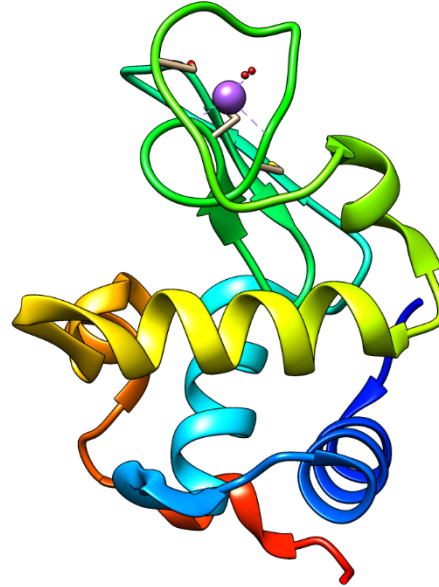
- Crystal thickness for MicroED is limited to ~700nm.
- Crystals ranging from 1 $\mu$ m to 50 $\mu$ m thickness can be thinned 150-600nm using cryo-FIB milling.
- Lysozyme, used as a test system, resolves to 1.9 Å in cryo-FIB milled crystals.



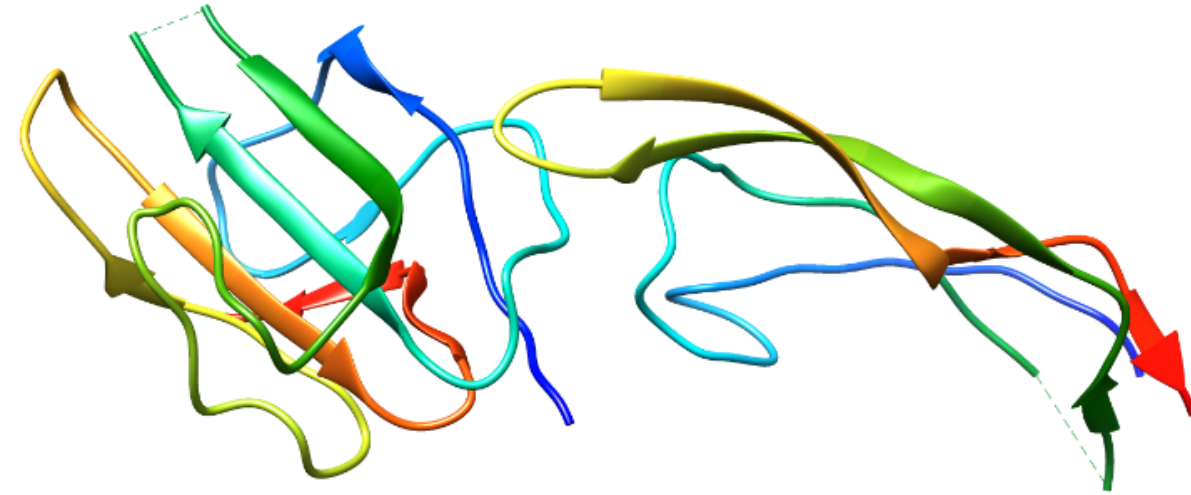
# Summary



Tau peptide



Lysozyme



TGF- $\beta$ m:T $\beta$ RII

Images of PDB 5K7N; 5K7O; 5TY4; 5K7P; 5K7Q;  
5K7R; 5K7S; 5K7T created with Chimera



MicroED can make use of  
crystals too small for XRD



High resolution, high throughput  
technique



Components can be retrofitted  
(depending on configuration)



Cryo-FIB opens a path for  
intermediate-sized microcrystals



**Thank you!**