

Oxford's PheniX is a closed-cycle helium cryostat designed for powder X-ray diffraction at the temperature range from 25 °C (298 K) down to - 261 °C (12 K)

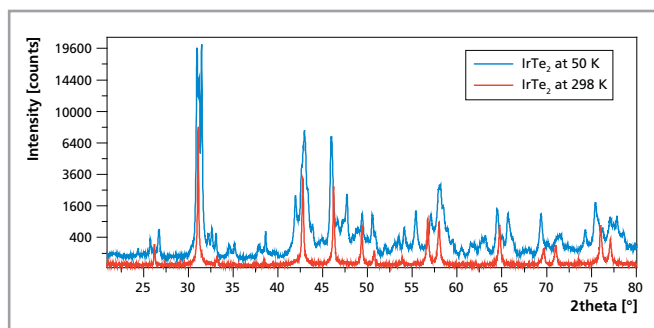
## Non-ambient attachment for XRD

# PheniX – low-temperature cryostat

### Benefits

- Extremely low temperatures
- High temperature uniformity and stability across the sample (0.1 K at 12 K)
- Fast cool-down and warm-up of the sample
- Easy to use and operate due to the compact design of the closed-cycle cryostat
- Virtually no helium consumption
- Easy sample preparation and mounting

### Application example



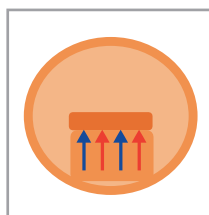
*The structure of the low temperature phase of IrTe<sub>2</sub> was solved using combined single crystal neutron diffraction and powder X-ray diffraction data.*

H. Cao et al. Phys. Rev. B 88, 115122 (2013) Courtesy of Huibo Cao, Oak Ridge National Laboratory

# PheniX

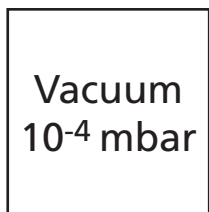
## low-temperature cryostat

### Features

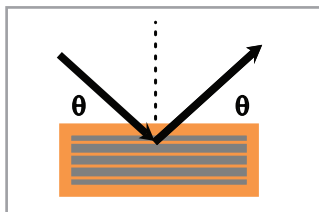


Heating/cooling (plate)

12 K – 290 K (vacuum)  
Cool-down time to 100 K: 40 min  
Cool-down time to 12 K: 60 min  
Warm-up time from 12 K: 45 min



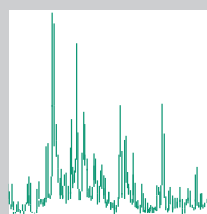
Atmospheres



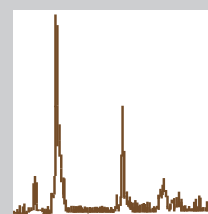
Flat plate reflection geometry.  
Sample holders made of Cr-Cu alloy.



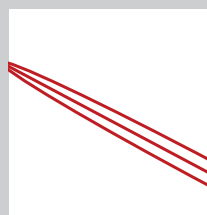
### Applications



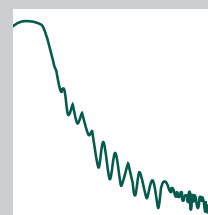
Powder XRD



Basic grazing incidence XRD\*



Basic stress\*



Basic reflectivity\*

\* Limited sample alignment options (no tilt and rotation axis)

### Conclusion

The PheniX cryostat is an ideal choice for in situ powder XRD studies of organic and inorganic materials at extremely low temperatures, down to 12 K (-261°C).