



PANalytical
get insight



Cryostream Plus Compact from Oxford Cryosystems enables XRD experiments in transmission geometry at temperatures from 227 °C (500 K) down to -193 °C (80 K).

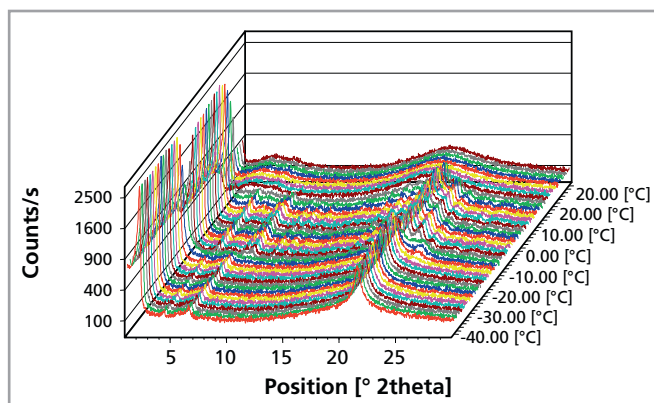
Non-ambient attachment for XRD

Cryostream Plus Compact

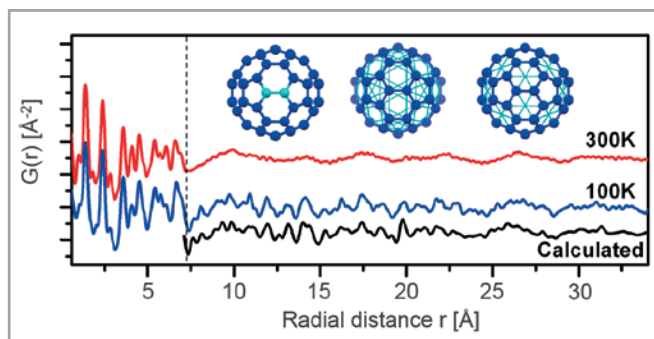
Benefits

- Unique device for low- and medium-temperature experiments in capillary geometry:
 - minimizes preferred orientation
 - excellent low-angle data
 - allows study of samples sensitive to air or humidity
- Wide temperature range
- Laminar gas flow significantly reducing ice formation
- Fast cool down to 100 K in just 20 min
- Stable temperature regime in excess of 0.1 K
- Low and constant N₂ consumption
- Optional automatic refill unit available, allowing experiments of infinite duration

Application examples



In situ monitoring of phase transformation of LOL (a fat) recorded 'live' at a temperature ramp of 2 deg/min

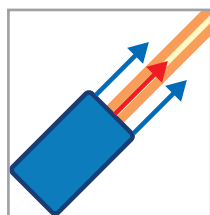


Pair distribution function of C₆₀ (fullerene) as a function of temperature

Cryostream Plus Compact

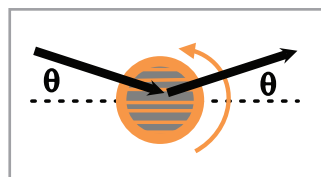


Features



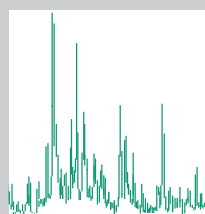
Gas stream cooling/heating

80 K – 500 K
Cool-down time to 100 K: 20 min

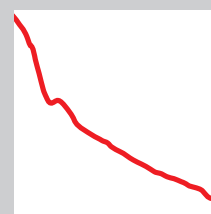


Capillary spinner in transmission geometry and glass capillary

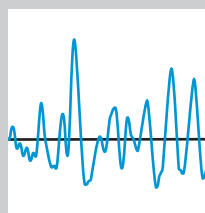
Applications



Powder XRD



Small-angle X-ray scattering*



Pair distribution function*

* Application developed by PANalytical

Conclusion

The Cryostream Plus Compact is an ideal attachment for *in situ* XRD applications at variable temperatures requiring capillary transmission geometry. It is suitable for XRD studies of both organic and inorganic powders.