

Features

- **Easy manipulation and mounting** of crystals from 700 μm to less than 5 μm .
- **Rigid yet thin construction**: minimal bending in the cryostream and in viscous solutions.
- **Convenient scoop-like action** for retrieving crystals and transferring them between solutions, reducing chance of crystal loss or damage.
- Gold hue provides **good mount-crystal contrast** during alignment.
- **Low background scatter** – much less than 20 micron nylon – increasing achievable resolution with small crystals.
- **Ideal for use at microfocus beamlines.**
- **Completely non-magnetic** → no accidental pick-up by magnetic tools and manipulators.
- **Allows more complete optimization of cryoprotection and flash cooling.**
- **Low thermal mass** → rapid cooling.
- **Crystal aperture can be matched to crystal size and shape**, minimizing surrounding liquid.
- **Easy removal of excess liquid.**
- **Less liquid** → less background scattering, easier cryoprotection, better resolution.
- Designed specifically for **easy automated crystal alignment** in the X-ray beam. Crystals can be aligned without ever optically observing them.
- **Fully compatible with existing crystal mounting hardware.**
- **Custom designs** available on request.