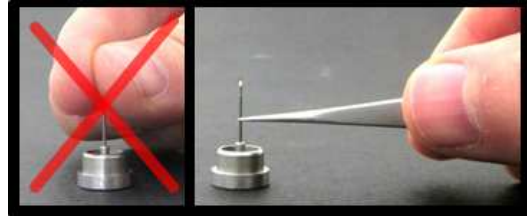


Handling Do's and Don'ts

The microfabricated tips of MiTeGen's MicroMounts, MicroLoops and other tools are made from the most durable unoriented polymer available, for ultra-low background X-ray scatter. Compared with an oriented fiber like nylon, unoriented polymers can be more easily torn or cut if used improperly. Follow the instructions below to get the maximum life out of our products.

1. **Never** touch the tips with your hands. Always use our heavy duty tweezers along the pin shaft, not the tip.



2. **Never** push the beveled steel pin into the tip against a hard surface.

3. **DO:** Clean the tips by soaking in detergent while held in a MiTeGen Base Holder (Cat.# GB-BH).

DO NOT: Sonicate for extended times. Sonication for a few minutes or so is ok, but sonication for long periods can damage some tips. Using an enzyme-containing detergent, sonication should be unnecessary.



4. **DO:** Remove stubborn residues using MiTeGen's Medium (M) or Fine (F) Paper Wick.

DO NOT: Use cotton or foam-tipped swabs. Swabs can catch and tear the polymer.

5. To dry the tips after cleaning, let them air dry. You can briefly dip in isopropanol for faster drying.

Avoid blow drying. If the nozzle is placed too close to the tip, the resulting turbulent drag force can damage the tip. Keep the nozzle at least 20 cm away from the tip.



Please contact xtals@jenabioscience.com with comments or suggestions.