

Cleaning Instructions for MicroMounts™, MicroLoops™, MicroMeshes™, MicroGrippers™ and MicroTools™

The best and easiest way to clean MiTeGen's tools is to soak them in a bath of a detergent-containing solution. When mounted in MiTeGen's Base Holders (Cat.# GB-BH-x), large quantities of mounts can be cleaned at one time.

To remove protein residues, we recommend an enzyme-containing detergent such as Alconox's Tergazyme or Decon's Contrex EZ; otherwise, a standard laboratory glassware/plasticware cleaner like Alconox is adequate. We recommend an initial 10 to 15 minute soak, followed by a quick rinse in a bath of water to remove the detergent. The optimal length of the soak needed will be determined by what residual materials are left on the mount or tool (protein, etc.) and how long it has been since they were last used. The sooner you clean them after use the easier it will be.

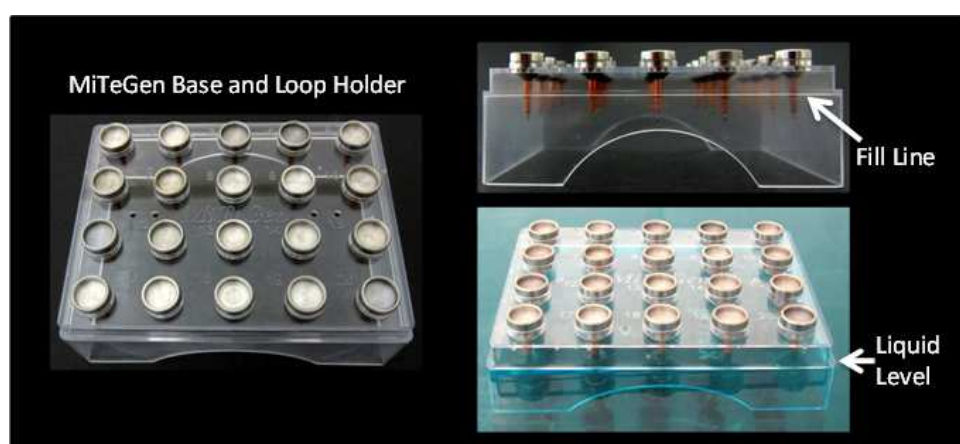


Figure 1: Base Holders and fill line for soaking MiTeGen's mounts and tools e.g. in detergent containing solution

Step by Step Instructions for Cleaning:

1. Load your loop-containing goniometer bases into the Base Holder.
2. Place the Base Holder into a plastic or glass tub or dish.
3. Mix your detergent-containing solution according to the instructions on the detergent box.
4. Then pour the detergent-containing solution into the tub until the level rises to the fill line on the Base Holder. This will ensure that the loop tips are soaked, while keeping the stainless steel part of the goniometer bases (caps) out of the solution and preventing corrosion.
5. Incubate for 10 - 15 minutes.
6. Replace the detergent-containing solution in the tub by ultrapure water to rinse the loops. This step should be repeated 1 - 2 times until the detergent is completely removed.
7. Let the loops air-dry at a clean place for about 15 - 30 minutes.

In the rare case when soaking doesn't remove deposits, you can use a Medium (M) or Fine (F) paper wick. For easiest handling, cut back from the widest end using a razor blade so that the diameter fits into a 0.7 mm mechanical pencil. Then dip the other end in water, a detergent solution or in isopropanol, and gently stroke the polymer from base to tip.



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