

Lab shakers

Best practices for using adhesive mats on orbital shakers

Thermo Scientific™ adhesive mats are reliable and easy to use when properly handled on any of our orbital shakers.

When working with our adhesive mats, the top three best practices are:

1. Choose the appropriate adhesive mat to fit your vessel type and run speed.
 - Adhesive mats can be used to secure a wide range of tube racks, Erlenmeyer flasks and other flask types, multiwell plates, wash basins, and more to an orbital shaker platform. The maximum load and speed are dependent on the type of adhesive mat chosen, the base area-to-height ratio of the desired labware, and the material that comprises the labware.
2. Clean and dry the adhesive mat at least once each week.
 - At least once each week or before attaching the adhesive mat to the shaker platform, clean the mat and the shaker platform thoroughly to help ensure they are free of oil, fingerprints, dust, or dirt. Use mild dishwashing soap without skincare additives, then rinse with 70% ethanol to remove any residues left by the cleaner. After the mat is completely dry, press firmly to attach to the shaker platform.



3. Ensure that the vessels are properly attached to the adhesive mat.
 - Attach each vessel to the mat by pressing firmly and then apply a gentle clockwise twist. This twisting motion secures the vessel and locks it into place. Ensure that the entire bottom surface of the vessel is firmly attached. To remove, gently tip the vessel to one side and then pull to that same side.

Ordering information

Description	Maximum rpm*	Cat. No.
High Adhesion Sticky Mat, 200 mm x 200 mm	400 rpm	75004126
High Adhesion Sticky Mat, 280 mm x 170 mm	400 rpm	75004127
Low Adhesion Sticky Mat, 200 mm x 200 mm	200 rpm	75004111
Low Adhesion Sticky Mat, 280 mm x 170 mm	200 rpm	75004117
Sticky Mat, 140 mm x 140 mm	225 rpm	88881126
Sticky Mat, 230 mm x 230 mm	250 rpm	300349
Adhesive Mat Rolling Applicator	N/A	75004124

* See manual for suitability of your labware material at maximum shaking speed.

Tips and tricks for adhesive mats

Our adhesive mats are a practical alternative to using and storing stainless steel shaker clamps. To fully realize the adaptability of our shaker mats, check out the tips and tricks listed below:

- Use the mats at temperatures between 0°C and 40°C; higher temperatures will not allow proper adhesion
- The mats can be cut to size and fitted to any shaker platform, or you can use multiple mats to cover the entire surface area of larger platforms
- Do not exceed the maximum load size and speed—the Thermo Scientific™ Low Adhesion Sticky Mat is suitable for all vessels up to 200 rpm, while the Thermo Scientific™ High Adhesion Sticky Mat is suitable for some vessels up to 400 rpm
- Our adhesive mats are color coordinated by adhesion level for easy recognition; each part number is a different color to ensure quick identification in the laboratory
- Adhesive mats can be applied by hand or using the Thermo Scientific™ Adhesive Mat Rolling Applicator—the applicator is an optional tool to help press the adhesive mat to the platform
- Make sure the mat and platform are completely dry—they are slippery when wet; regular cleaning is critical to proper use because any dust, dirt, oil, or fingerprints will block adhesion and put your vessels at risk

Summary

When adhesive mats are properly cleaned, maintained, and operated within published guidelines, they provide a flexible, convenient, and easy-to-use alternative for attaching a variety of vessels to any orbital shaker.



Thermo Scientific™ sticky mats are flexible, convenient, and easy to use for attaching a wide variety of vessels to an orbital shaker.

Learn more at thermofisher.com/shakers

thermo scientific