

Quick Start Guide

AAW™ Flex Gripper

99900244



AAW™ Flex Gripper

The AAW™ Flex Gripper is an instrument that can pick up labware and move it around the deck automatically. Adding a gripper to the AAW™ Automated Assay Workstation (AAW™ Workstation) helps improve automated workflows in your lab.

User Guide

Read the AAW™ Workstation user guide before attempting to operate. It can be downloaded from the product page at SigmaAldrich.com. The user guide contains detailed information about this product, modules, and accessories.

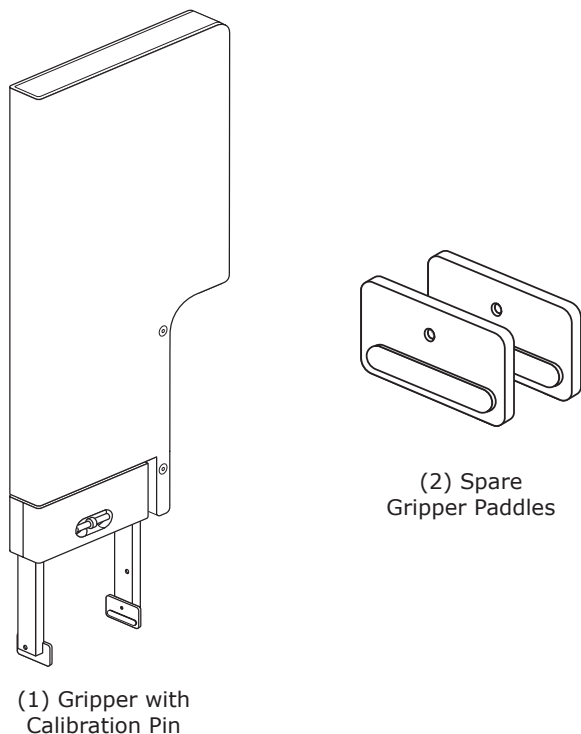
Labware Compatibility

The gripper is compatible with labware of the following types:

- Certain full-skirted 96-well plates, including Opentrons Tough PCR plates
- Certain deep well plates
- All Flex tip racks

The AAW™ Flex Gripper may work with other ANSI/SLAS compliant labware, but this is not recommended.

Box Contents



Installation

For assistance with installation, contact SigmaAldrich.com/TechService.

Attach

Instructions on the touchscreen, or in the Opentrons App, will guide you through the attachment process. On the touchscreen, tap Instruments and then tap Extension Mount. To attach your gripper:

1. Clear the deck area so it is free from obstructions. The gantry will move during the installation process.
2. Connect the gripper to the right side of the gantry arm by aligning the mounting holes on the back of the gripper with the pins on the side of the gantry arm.
3. Secure the gripper by tightening the two mounting screws with a 2.5 mm hex screwdriver. These are captive screws recessed into holes located on the front of the gripper.

Calibrate

The gripper requires calibration before use. Instructions on the touchscreen, or in the Opentrons App, will guide you through the calibration process.

To calibrate your gripper:

1. Clear the deck area so it is free from obstructions. The gantry and pipette will move during calibration.
2. Attach the calibration pin to the front jaw of the gripper and follow the instructions on the touchscreen. Repeat the calibration process for the rear jaw.
3. Remove the calibration pin and store it in the magnetic storage area when calibration is complete. Your gripper is ready for use.

Clean

Clean the outside of the gripper with a soft cloth and isopropyl alcohol or ethyl alcohol. Do not use acetone or other aggressive solvents. For more cleaning and decontamination information, see the AAW™ Workstation User Guide.

Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

Technical Assistance

Visit the tech service page on our web site at SigmaAldrich.com/TechService.

Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at SigmaAldrich.com/Terms.

Contact Information

For the location of the office nearest you, go to SigmaAldrich.com/Offices.

Merck, Millipore, AAW, and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.
© 2025 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.

