

The PlateCrane™ from Hudson can be combined with microplate reader systems, washers, and dispensers from Molecular Devices to produce high-capacity workcells.

The PlateCrane allows users to achieve "walkaway automation" of microplate-based assays. Instead of loading and processing microplates one at a time, the user loads stacks of microplates onto the PlateCrane, which then automatically feeds them to the integrated Molecular Devices reader system, dispenser, or washer.

Software interfaces written by Hudson provide communication between the PlateCrane and all attached devices. Automation of large numbers of microplates with these workcells increases productivity in the laboratory. A single process such as plate washing or multiple processes such as washing, dispensing, and reading can be automated. Cell-based or enzyme-based screening assays can be automated in this manner.

PlateCrane Microplate Handling Robot

- Available in 3 models with up to 4 degrees of motion
- Maximum capacity up to 420 microplates
- Works with 96, 384, and 1536-well microplates, deep well plates, and plates with lids
- Up to 8 microplate devices can be configured in a single workcell

PlateCrane Interfaces are available for the following Molecular Devices microplate-based devices.

- SpectraMax Absorbance Microplate Reader Systems
- SpectraMax GEMINI XS Fluorescence Microplate Reader Systems
- CLIPR Chemiluminescence Imaging Microplate Reader System
- FLIPR Fluorometric Imaging Microplate Reader System
- Analyst Multimode Reader
- AQUAmax™ 96-384-1536-well Dispensers
- Embla Plate Washers



Dual device system with SpectraMax reader and Embla washer.



PlateCrane integrated with AQUAmax dispenser.



Close-up of PlateCrane gripper placing a plate in the SpectraMax nest.

Visit www.hudsoncontrol.com for comprehensive information about the Hudson PlateCrane.
 Visit www.moleculardevices.com for comprehensive information about Molecular Devices microplate-based systems.

Hudson Control Group, Inc.
 10 Stern Avenue - Springfield, NJ 07081
 Tel. 973-376-7400 FAX 973-376-8265

Molecular Devices Corporation
 1311 Orleans Drive - Sunnyvale, CA 94089-1136
 Tel. 408-474-1700 FAX 408-747-3601