



Example integration

# plate::handler Connect

Automate. Integrate. Scale.

Unlock the full potential of your AssayMate™ liquid handling workstation, increase productivity, and reduce costs. plate::handler™ Connect converts your AssayMate liquid handling workstation into a powerful, modular, integrated workflow automation solution. Broaden your application range by combining our trusted Revvity automation solutions in one compact, cost-effective package.

## Benefits

- Affordable, future-ready
- Adds sample tracking, off-deck labware storage, random-access plate incubation and integration capabilities (reader, washer, etc.) to the AssayMate liquid handler
- Powerful, dynamic scheduling software supporting real-time decision making
- Sample tracking through built-in barcode reader

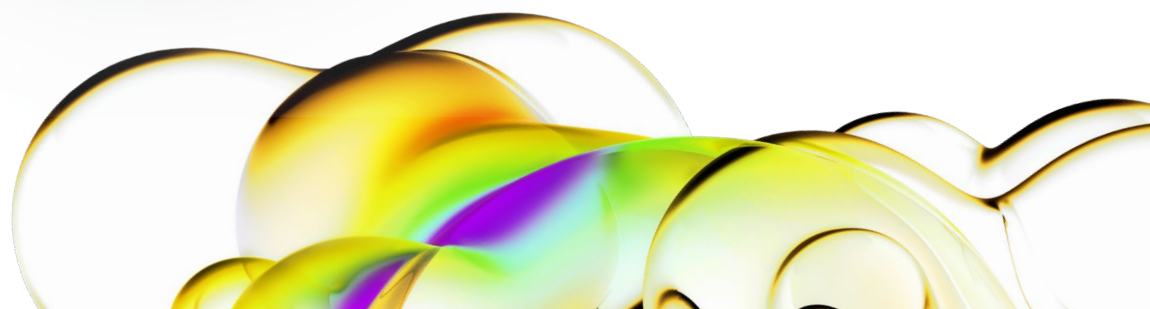
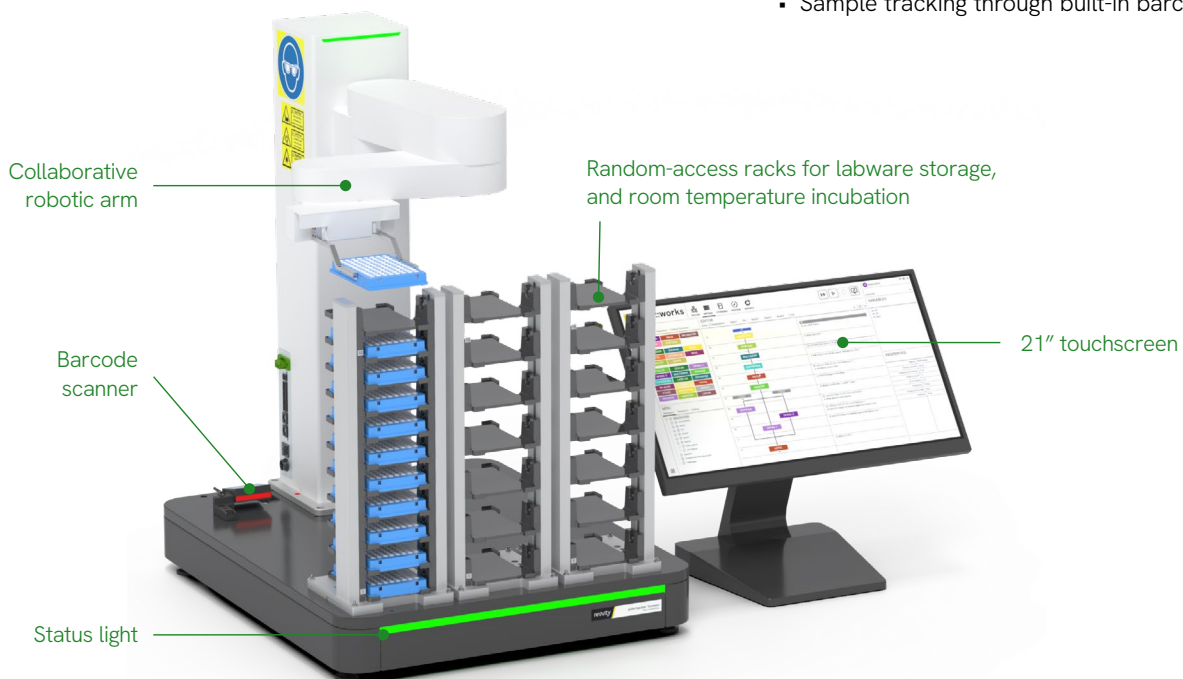
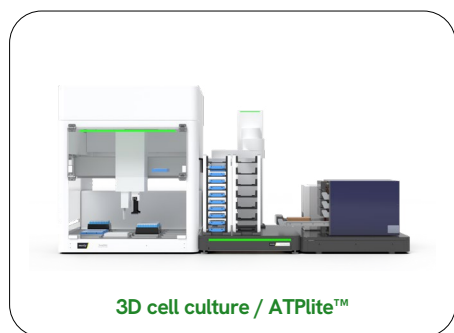


plate::handler Connect specifications	
Plate Handling	Collaborative, four-axis SCARA robot supporting side-by-side human-robot cooperation, built-in servo gripper enabling microplate handling in both landscape and portrait orientation
Labware storage, incubation (RT)	Up to 3x random-access plate racks, capacity depending on labware dimensions. Rack design supporting easy labware loading
Barcode reading	SICK CLV601, for reliable barcode reading in laboratory automation applications
Scheduling software	plate::works™ software V6.35, job::manager process planning add-on
Control PC/monitor	HP® EliteDesk 8 Mini PC (5i, 16GB RAM, 512GB) or similar PC build into robot base, 21" touchscreen monitor, optional hardware switch/firewall for secure image/data upload
Electrics	1-phase 230V/50 Hz (110V/60Hz)
Dimensions (robot only) (W x D x H)	530 mm x 650 mm x 780 mm

### plate::handler Connect with the AssayMate liquid handler



### Other example configurations

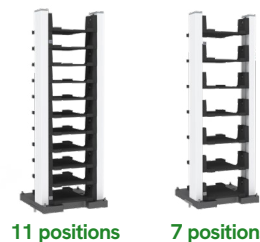


Part number	Description
HH20260001	plate::handler Connect, robotic plate mover with barcode reader, control PC, 21" touchscreen monitor and plate::works scheduling and control software.  Racks for labware storage and room temperature incubation to be quoted separately.



### Labware storage (add up to 3x to quote, same type or mixture)

HH20260002	11-position random access rack for SW plates
HH20260004	7-position random access rack for DW plates
HH20260003	6-position random access rack for tip boxes
HH20260005	Random access rack with custom spacing



### Baseplate extensions

HH20260006	Baseplate extension S, including mounting bracket
HH20260007	Baseplate extension M, including mounting bracket



### Instrument integration

Please submit IS request form to obtain a quote

[Request a quote](#)

### Installation

HH20110069	Installation on site (by local Revvity Service Team, 1 day including travel)
------------	--

### Application development

2770001	Level 3 support application development, 1 day
---------	--



[Learn more](#)

Supporting labs do more with less.

revvity