



Example integration

# explorer G3 SmartBench workstation

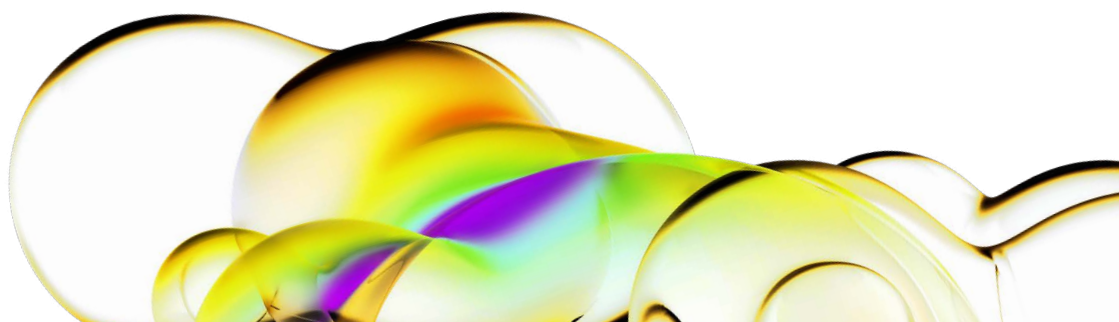
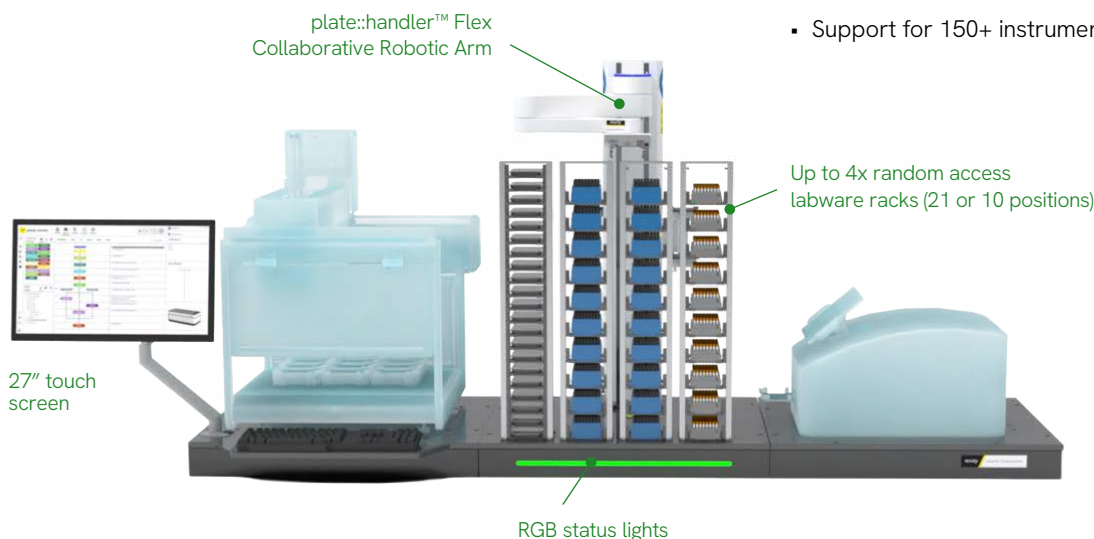
## Integrated Lab Automation

**Modular, scalable, all-in-one solutions to automate your workflows**

Elevate your productivity with **explorer™ G3 SmartBench**, our cost-effective solution to seamlessly integrate instruments with advanced robotics to create modular workflow solutions.

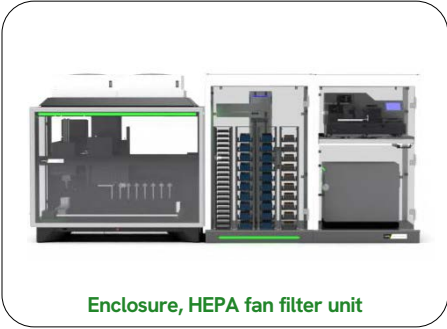
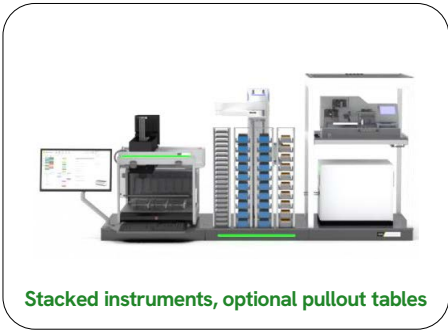
### Benefits

- Affordable, future-proof
- Built-in support for Liquid Handling instruments: 2x active de-lidding positions, passive positions for plate alignment and orientation change
- Powerful dynamic scheduling, software supporting real-time decision making
- Sample tracking through built-in barcode reader
- Support for 150+ instruments



explorer G3 SmartBench specifications	
Platform	explorer G3 SmartBench workstation
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. Maximum reach: 731 mm; maximum z-travel: 750 mm; repeatability: 0.2 mm
Barcode reading	SICK CLV601, for reliable barcode reading in laboratory automation applications
Lidhandling	2x lidhandling stations mounted to robotic arm, option to increase to 4x
Scheduling software	plate::works software V6.35, job::manager process planning add-on
Control PC / network	19" rack mountable Industrial PC running Windows® 11 Enterprise LTSC version, Hardware Switch/Firewall for secure image/data upload
Electrics	1-phase 230V/50 Hz (110V/60Hz) with central mains switch
Dimensions	2400 mm x 800 mm, height to depend on configuration

Configurations and options





[Learn more](#)

Supporting labs do more with less.

