

CUSTOM, MODULAR AUTOMATION SOLUTIONS

Automation will be key to increase lab efficiency and to drive scientific advancement. **explorer™ G3 workstation** comprises a hardware and software platform which can be freely configured to create modular, bespoke automation solutions which provide turnkey automation solutions for virtually all scientific workflows.

WHY AUTOMATE YOUR SCIENCE?



IMPROVE EFFICIENCY



INCREASE PRODUCTIVITY



REDUCE HANDS-ON TIME



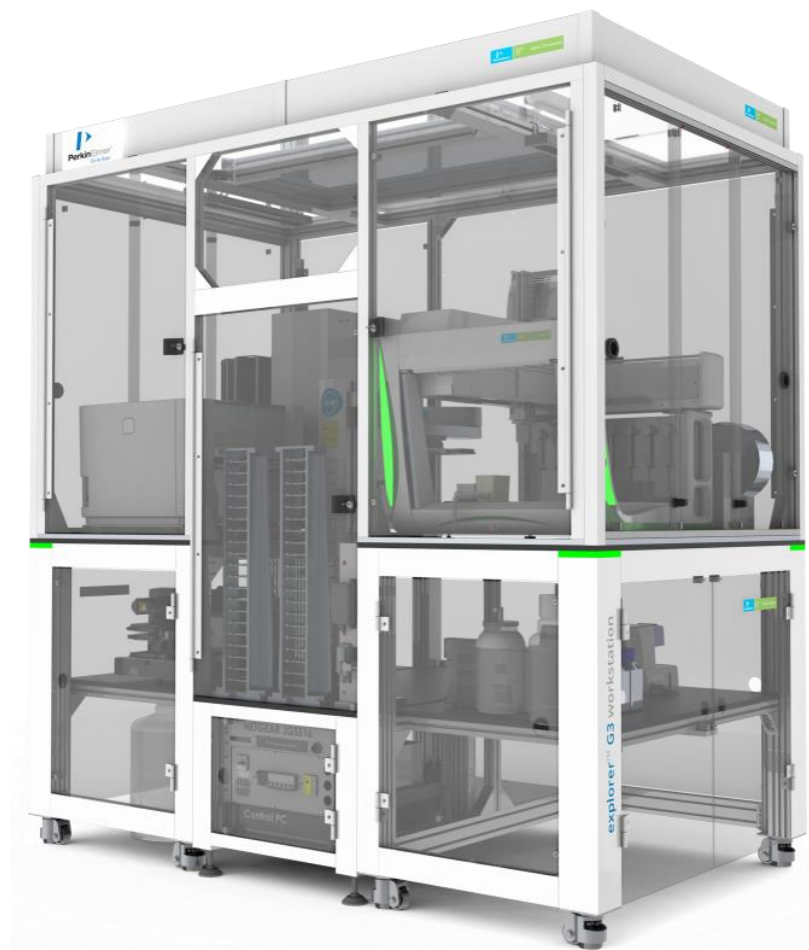
OPERATORS AND ROBOTS
SHARING SAME WORKSPACE



STANDARDIZATION, FAIR DATA



MODULARITY & SCALABILITY



LEARN MORE ABOUT INTEGRATED LABORATORY AUTOMATION SOLUTIONS

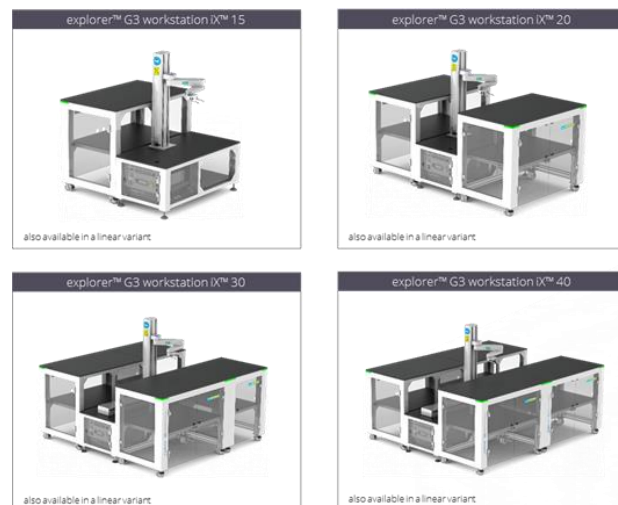
For research use only. Not for use in diagnostic procedures.

© 2021 PerkinElmer. All Rights Reserved. Schnackenburgallee 114 | 22525 Hamburg | perkinelmer-appliedgenomics.com/home/products/integrated-lab-automation/
All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

explorer™ G3 workstation

PLATFORM

Center piece of explorer™ G3 workstation platform are a range of standardized Instrument and Robot Tables which can be assembled to create modular, easy-to-upgrade workstations of different shapes and sizes.

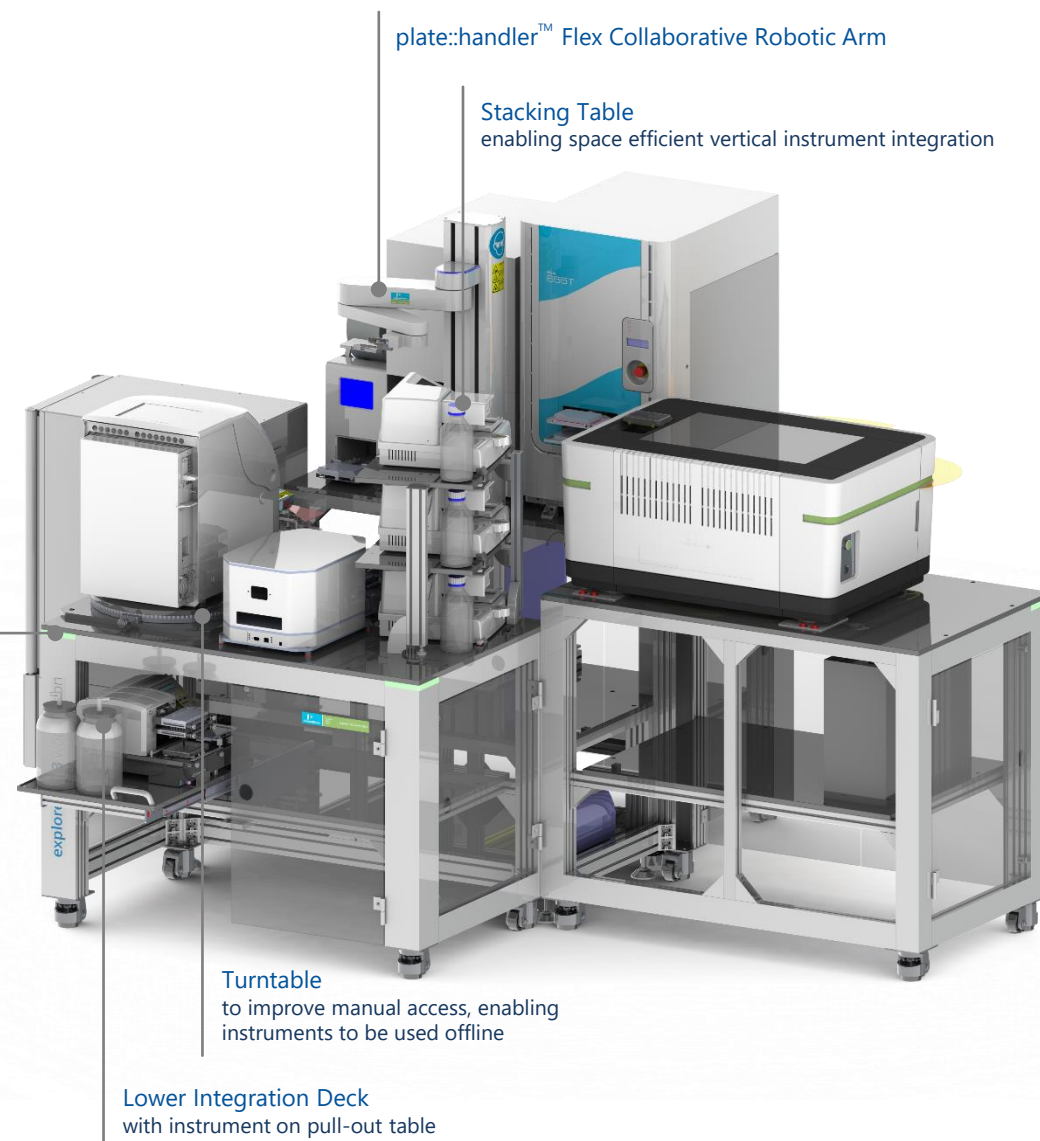


SPACE SAVING VERTICAL INTEGRATION CONCEPT




For research use only. Not for use in diagnostic procedures.

© 2021 PerkinElmer. All Rights Reserved. Schnackenburgallee 114 | 22525 Hamburg | perkinelmer-appliedgenomics.com/home/products/integrated-lab-automation/
All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.



LEARN MORE ABOUT INTEGRATED LABORATORY AUTOMATION SOLUTIONS



The image displays three PerkinElmer Flex robots, labeled FLEX 400, FLEX 750, and FLEX 1160, arranged from left to right in increasing order of height. Each robot is a white, four-axis SCARA robot with a built-in servo gripper. The FLEX 400 is the shortest, FLEX 750 is medium height, and FLEX 1160 is the tallest. All three robots have a similar design with a vertical column and a horizontal arm.

- + Collaborative, four-axis SCARA robot with built-in safety features enabling side-by-side human-robot cooperation; no safety shielding needed
- + Full metal casing, space saving design with motion controllers build into the structure of the robot
- + Fast, fluid and quite movements
- + Build-in Servo gripper enabling robot to grip plates on either long or short side
- + Hand guided teaching
- + Robot available in 3 different heights (400, 750 and 1160mm), two different arm length with option to increase lateral reach by placing robot on a linear track

LEARN MORE ABOUT INTEGRATED LABORATORY AUTOMATION SOLUTIONS

For research use only. Not for use in diagnostic procedures.

© 2021 PerkinElmer. All Rights Reserved. Schnackenburgallee 114 | 22525 Hamburg | perkinelmer-appliedgenomics.com/home/products/integrated-lab-automation/
All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

plate::works™ SCHEDULING SOFTWARE

BUILDING ON 20 YEARS OF EXPERIENCE

Since its launch in 1997, plate::works™ scheduling and control software has been used to automate a great variety of customer workflows, with numerous installations spread over all five continents.

FLEXIBLE

plate::works™ software empowers operators to create their own methods with the plate::works™ event-based scheduling model to support even the most challenging workflows. Flexible control elements to set-up sample and/or plate specific processes, with plate::works™ software to support multiple workflows being processed in parallel.

EASY-TO-USE

Software guiding operators through the steps to set-up an automated process. All labware movements handled automatically by the scheduler with robot to use transport speeds and gripping positions based on information stored in a central labware database.

RELIABLE

Software to make every conceivable attempt to recover from an error situation, making a failed run an exceptional event. Advanced error handling routines guiding operators through the steps and options to get the system quickly back into operation.

> 150 INSTRUMENT DRIVERS AVAILABLE

Extensive library of readily available instrument drivers. Drivers for PerkinElmer Instruments to be provided free of charge.



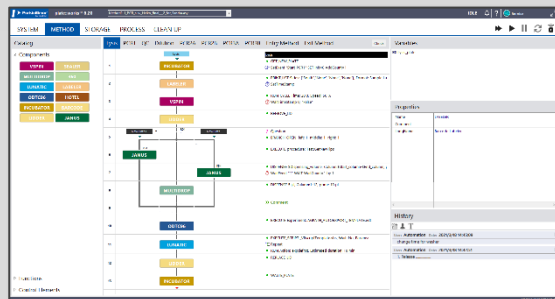
LEARN MORE ABOUT INTEGRATED LABORATORY AUTOMATION SOLUTIONS

For research use only. Not for use in diagnostic procedures.

© 2021 PerkinElmer. All Rights Reserved. Schnackenburgallee 114 | 22525 Hamburg | perkinelmer-appliedgenomics.com/home/products/integrated-lab-automation/
All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

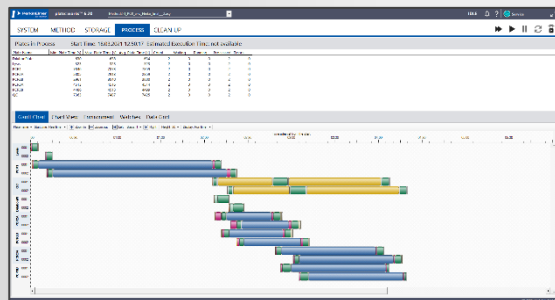
plate::works™ SCHEDULING SOFTWARE

METHODS EDITOR



STORAGE EDITOR

GANTT CHART



KEY FEATURES

Event Driven Scheduling

plate::works™ software to combine advantages of dynamic and static scheduling enabling operators to control and fine-tune dynamic scheduling by adding constraints and/or controlling elements

Real Time Decision Making & Re-Scheduling

Scheduler to support on-the-fly re-scheduling allowing critical parameters to be updated at any time during a run and plate processing to respond to external data or events (e.g. results, conditions, LIMS, scripts, ...)

Parallel Methods

Scheduler to support multiple independent methods being executed in parallel

Continuous / On-Demand Processing

Scheduler to support continuous plate processing allowing new plates and labware to be added to an already running process as well as on-demand plate processing with system to process plates when they become available

Pooling

Up to 3 identical instruments to be treated as one logical instrument (for easier programming and added redundancy)

Simulations

To quickly optimize workflows (test different process variants and conditions) and to check for correct execution prior to committing time and reagents

21CFR11 Support

plate::works™ software to support setting-up regulated processes by providing user rights management and by logging changes being made to methods

Worklist Support

Plate/sample specific parameters or conditions (e.g. incubation times, dispense volumes, ...) can be read from worklists. Support for cherry picking, normalization and other tasks relaying on external information

Scripting Support

Enabling operators to add own functionality to scheduling process

Offline Use

plate::works™ software to support operators taking critical detection instruments off-line and to use manually up till the point where instrument is been needed to support automated process.

Add-ons / Options

job::manager™ Process Planning and Workflow Scheduling software
wellmap::creator™ Transfer Map Editor

LEARN MORE ABOUT INTEGRATED LABORATORY AUTOMATION SOLUTIONS

For research use only. Not for use in diagnostic procedures.

© 2021 PerkinElmer. All Rights Reserved. Schnackenburgallee 114 : 22525 Hamburg : perkinelmer-appliedgenomics.com/home/products/integrated-lab-automation/
All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

CUSTOM, MODULAR, SCALABLE WORKFLOW AUTOMATION SOLUTIONS

**DESIGNED TO
RUN 24/7**

**COMPACT, SPACE
SAVING DESIGN**

**REAL-TIME
DECISION MAKING**

**> 150 INSTRUMENT
DRIVERS AVAILABLE**

**> 20 YEARS EXPERTISE IN BOTH
INSTRUMENTS & INTEGRATIONS**



Please contact us to discuss your automation needs and how PerkinElmer can help you automating your science

LEARN MORE ABOUT INTEGRATED LABORATORY AUTOMATION SOLUTIONS

For research use only. Not for use in diagnostic procedures.

© 2021 PerkinElmer. All Rights Reserved. Schnackenburgallee 114 | 22525 Hamburg | perkinelmer-appliedgenomics.com/home/products/integrated-lab-automation/
All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.