

Automated Plate Feeding for Operetta CLS™ High-Content Analysis System & Opera Phenix® Plus High Content Screening System

To automate Operetta® CLS™ High-Content Analysis System & Opera Phenix® Plus High Content Screening System, PerkinElmer can offer a two turn-key plate::handler™ FLEX plate feeding solutions designed to provide walk-away automation for **Fixed Cell Screening** and **Live Cell Screening**. With plate::handler™ FLEX plate feeding solution using the same robotic arm, PC and control software as our larger explorer™ G3 workstation ensures full upgradeability incl. the possibility to reuse existing components towards the build of a larger automation platform.

Robotic Arm	Collaborative, four-axis SCARA robot with built-in safety features enabling side-by-side human-robot cooperation; Control PC with 2x SDD in RAID 1 configuration running Windows 10 Enterprise LTSB
Storage Capacity	Fixed Cell Screening Package: 42x 96/384w (random access) Live Cell Screening Package: 44x 96/384 in Liconic® STX44 IC SA robotic incubator (32-50°C, 0-10% CO ₂ , max. 95%RH), 42x 96/384w in random access racks
Control Software	plate::works™ software 6.25 or better
Application Templates (incl. in delivery)	<ol style="list-style-type: none"> 1. Reading a variable number of plates using a fixed experiment (selected from drop-down menu) 2. Reading a variable number of plates with two subsequent experiments (PreciScan™ plug in – where supported by imager) 3. Reading a variable number of plates using plate specific experiments; with information on which experiment to be used for which plate to be taken from an Excel worklist
Options & Accessories	Larger racks, different size incubator job::manager workflow managing software (multiuser)

Operetta® CLS™
High-Content Analysis System
or Opera Phenix® Plus
High Content Screening System
on heavy duty lab table

plate::handler™ FLEX
Plate Feeding Solution

plate::handler™ Flex 750
collaborative robot arm

3x 14-position random
access plate racks &
barcode scanner

Workstation Extension

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.

FIX CELL IMAGING

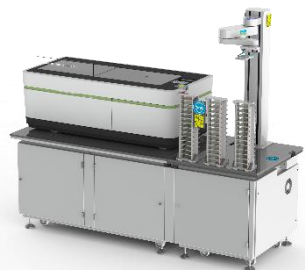


Benchtop Fix Cell Imaging for Operetta® CLS™ High-Content Analysis System

HH2020 0010

Benchtop Fixed Cell Imaging Package for Operetta® CLS™ High-Content Analysis System

- plate::handler™ Flex collaborative robot
- up to 42x96/384w plates stored in random access plate::racks
- barcode reader



Fix Cell Imaging for Operetta® CLS™ High-Content Analysis System*) or Opera Phenix® Plus High Content Screening System*)

HH2020 0011

Standalone Fixed Cell Imaging Package for Operetta® CLS™ High-Content Analysis System or Opera Phenix® Plus High Content Screening System

- plate::handler™ Flex collaborative robot
- up to 42x96/384w plates stored in random access plate::racks
- barcode reader

with onboard liquid handling**)



Fix Cell Imaging for Opera Phenix® Plus High Content Screening System*) with onboard liquid handling option**)

HH2020 0013

Standalone Fixed Cell Imaging Package for Opera Phenix Plus® Plus High Content Screening System with onboard liquid handling option

- plate::handler™ Flex robot
- up to 63x96/384w plates stored in random access plate::racks
- barcode reader
- Safety shielding

LIVE CELL IMAGING



Live Cell Imaging for Operetta® CLS™ High-Content Analysis System*) or Opera Phenix® Plus High Content Screening System*)

HH2020 0012

Standalone Live Cell Imaging Package for Operetta® CLS™ High-Content Analysis System or Opera Phenix® Plus High Content Screening System

- plate::handler™ Flex robot
- Liconic® STX44 IC, 44-position plate incubator (32-50°C, high RH, 0-10% CO₂)
- up to 42x96/384w plates stored in random access plate::racks
- barcode reader

with onboard liquid handling**)



Live Cell Imaging for Opera Phenix® Plus High Content Screening System*) with onboard liquid handling option**)

HH2020 0014

Standalone Live Cell Imaging Package for Opera Phenix Plus® Plus High Content Screening System with onboard liquid handling option

- plate::handler™ Flex robot
- Liconic® STX44 IC, 44-position plate incubator (32-50°C, high RH, 0-10% CO₂)
- up to 63x96/384w plates stored in random access plate::racks
- barcode reader
- Safety shielding

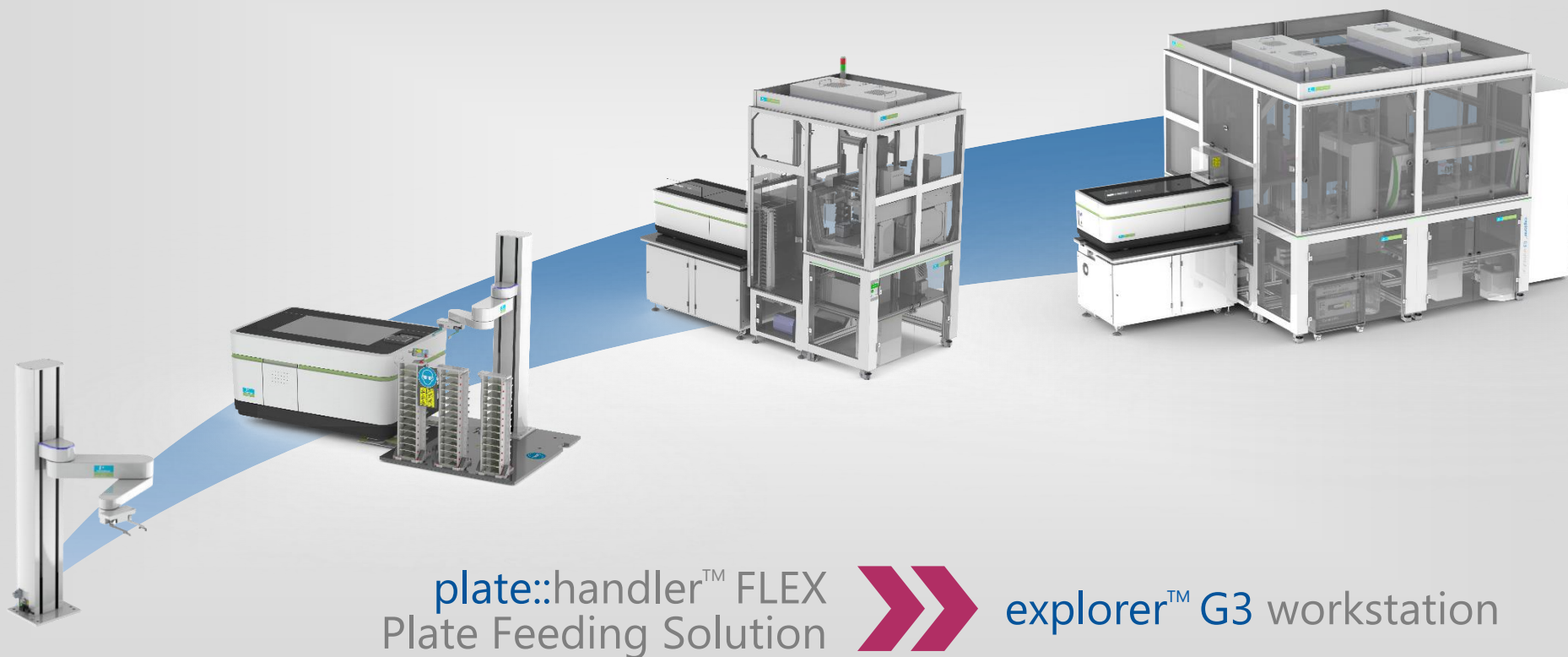
*) Imager on heavy-duty labcart, **) Standard pHF Package to supporting loading Opera Phenix® Imager with imaging plate only. Please contact PerkinElmer for configurations supporting storage and automated loading of compound plates and tip boxes

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.

LINEAR UPGRADE PATH TO *explorer™ G3* workstation
FULL WORKFLOW AUTOMATION SOLUTIONS



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.