

HUMAN HEALTH

ENVIRONMENTAL HEALTH

# AUTOMATE YOUR OPERETTA FOR GREATER PRODUCTIVITY

## Automated Workstations for Operetta High Content Imaging System

Simply choose the automation application solution that meets your lab's needs and increase the productivity of your Operetta® High Content Imaging System by up to five times, saving time and money.

As this is a complete solution from PerkinElmer, you do not have to be concerned about dealing with different vendors, and you can rest assured that our global service and applications teams will be on-hand to support you.

### Live Cell Imaging

Improve cell health and viability over prolonged periods with the addition of a dedicated incubator:

- Reduce well-to-well cell variance with a temperature, humidity and CO<sub>2</sub> controlled incubator
- Improve productivity by up to five times with automated plate loading
- Ensure high data integrity with barcode tracking throughout the whole process.

### Staining and Fixation

Improve your lab's reproducibility and productivity by automating cell staining:

- Run staining and Operetta imaging processes in parallel
- Fix your cells and remove formaldehyde aerosols automatically
- Move to higher densities, such as 384- or 1526-well plates, and save reagent costs
- Reduce staining variability
- Improve consistency through automated procedures, leading to improved results

## Screening Process Solutions

Automate the whole siRNA or compound screening process for maximum productivity:

- Reduce the risk of contamination through the use of HEPA filtered air
- Increase accuracy and reduce variability of compound addition through the use of automation
- Increase your flexibility by combining live and fix cell imaging, staining and fixation as well as siRNA and compound screening processes in one system
- Vastly improved screening time lines through an integrated, automated solution



Operetta workstation for automating the iRNA screening process

Application System	Fixed Cell Imaging	Live Cell Imaging	Cell Signal Induction	Fixed Cell Staining	Cell Fixation	Live Cell Staining	Antibody Addition	Compound Addition	3D Micro Tissue Handling	HEPA Filtered Air	Operetta TCO
Plate Handling (PH)	•										
Live Cell Imaging (LI)		•									•
Life and Fix Cell Imaging (LIF)	•	•									•
Life Cell Imaging and Signal Induction (LISi)		•	•								•
Staining and Imaging (SI)	•			•	•*						
Staining, Imaging and Cell Fixation (SIF)	•	•	•	•	•	•	•**				•
Antibody Addition and Staining (AS)	•		•	•	•*		•				•
Antibody Addition, Staining Compound and 3D tissue Screening (ASCD)	•	•	•	•	•*	•	•	•	•	•	•

\* Might require additional fume removal to prevent formaldehyde exposure

\*\* User has to accept higher dead volumes due to use of dispenser

Can't find the automation workstation to suit your application and workflow here?  
We also offer tailored automation configurations to cover every lab's unique requirements.

Please contact us for more information and expert advice to help you choose the right automated workstation for your needs.

PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright © 2013, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.