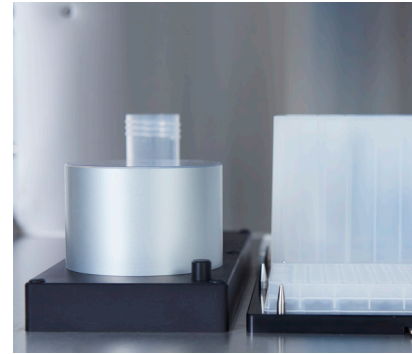
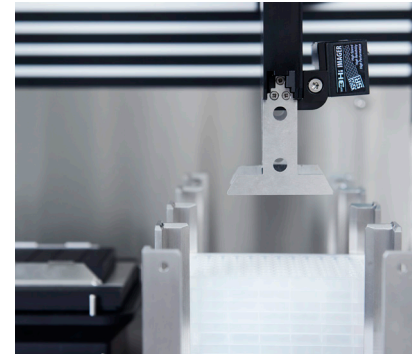


QPatch has been the benchmark automated patch clamp solution for cardiac safety and advanced electrophysiology since 2005. QPatch II builds on the legacy of its predecessor but is completely redeveloped to fulfil the requirements of the efficient ion channel laboratory of tomorrow.

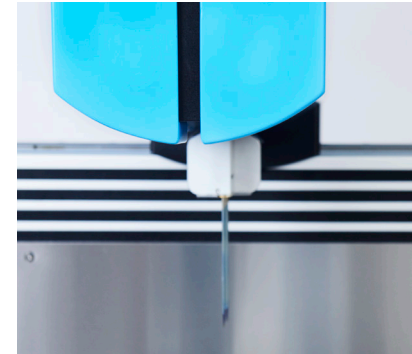
- Faster, easier to use and with fewer cells required providing same proven high-quality data
- The new intuitive user interface reduces human errors and enable easier staff rotation
- 100% correlation to QPatch. No re-validation of assays required



The compact cell hotel and cell preparation station ensures standardization. The compact design allows multiple cell hotel possibilities.



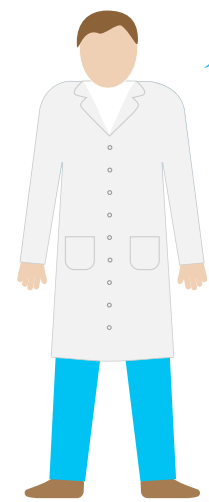
Stacking of plates on the work plane and a barcode scanner on the gripper arm provides faster and more robust assays.



Individual pipetting and pressure system is still the hallmark of the QPatch II operation, ensuring that all wells are handled individually.

We asked ion channel experts worldwide:

What are the main challenges in your lab today and in the future?



Pressure to increase laboratory efficiency

Need robust full automation with unattended use

Difficult to find qualified staff

Fast and robust

Greater flexibility with the intelligent assay scheduler
Barcode reader on the gripper arm saves time
New QPlate locking system provides greater QPlate position accuracy
Embedded QPlate electrodes ensure no electrode drift or maintenance

CiPA and cardiac safety

Pre-installed CiPA protocols
Prepared for multiple cell hotels
Glass coated microchannels reduce compound adsorption
Temperature controlled measurement site

Optimized cell handling

Patented automated cell preparation unit reduces user variation
Decreased cell consumption
Cell lines, stem cells and primary cells
Automated cell preparation ensures consistency

All cells treated individually

Individual pipetting and pressure system
Cherry picking of compounds
Adaptive whole-cell protocols
Adaptive voltage protocols
Individual estimated $V_{1/2}$

Service and support

Unrivalled application support
Trained and certified staff
Continuous support and training

True walkaway operation

6 hours or 10 plate unattended runs
QPlate stacking on the work plane, greater simplicity and robustness
Saves time for data analysis or assay development

Ease-of-use equals efficiency

Learn how to operate QPatch II in 10 minutes
Parallel analysis of data while running
Enables staff rotation
Free up time for assay development and analysis

New core engine with more power

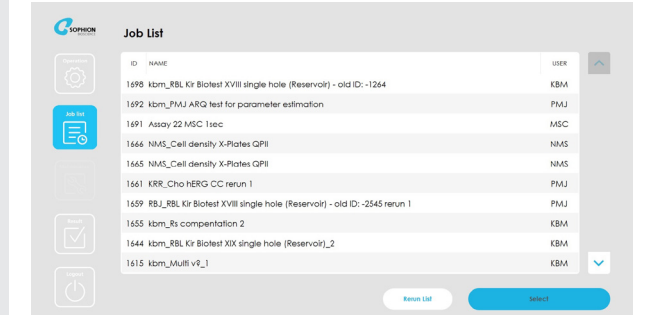
More powerful data engine
Larger internal memory
Easy access to USB ports
16 or 48 channels

Health and safety

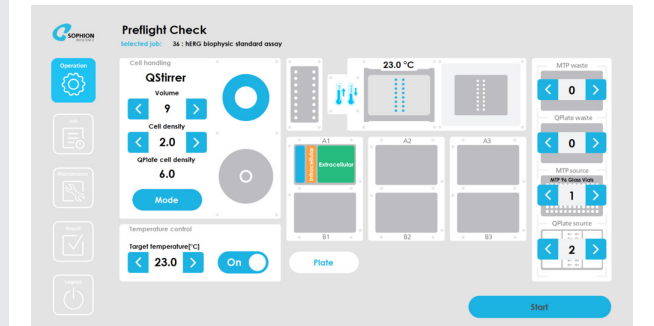
Adjustable height of work plane and touchscreen
Touch screen on arm or on table
Full visibility of work plane
Low waste volume



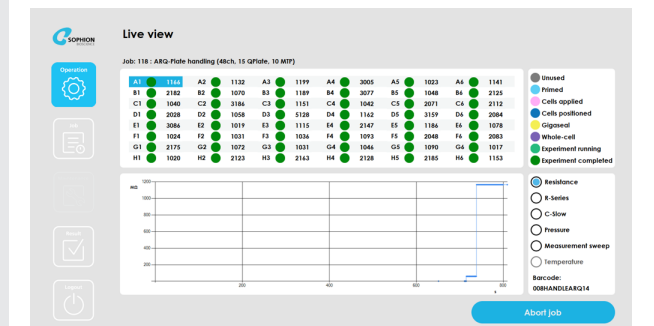
It takes less than 10 minutes to learn the intuitive user interface and how to operate QPatch II



Reduce human errors using the simple job list overview and re-run possibility



The work plane instructions are easy to follow when setting up your next job



During run informative live view data is displayed along with useful assay information

Easy to use

Simple and intuitive user interface reduces human errors
Learn to operate in less than 10 minutes
Pre-defined assays and re-runs directly from the touch screen

Efficient

Ease of use enables staff rotation
Overnight unattended operation
Automated data analysis and reporting

Versatile

Proven on all commonly used ion channels
Use for cardiac safety, profiling or small/medium library screens
Giga-seal quality without seal enhancer

Standardized

High reproducibility and repeatability
Temperature controlled testing environment
QPatch has been the benchmark for cardiac safety since 2005

Compliant

Electronic data records compliant with 21 CFR Part 11
Designed for Good Laboratory Practice (GLP)
Windows 10 compliant

QPatch II

AUTOMATED PATCH CLAMP SYSTEM

Cardiac safety

Compound characterization

Medium throughput screening

Advanced electrophysiology

The benchmark solution for cardiac safety and advanced electrophysiology since 2005

