



QPrep cell preparation unit

Enhance your ion channel research with simple and consistent cell preparation

Reduce variability in suspension-based e-phys assays and gain reliable, consistent data with our new cell preparation device. The compact, easy-to-use design streamlines your workflow and saves space. The combined cell preparation and cell hotel deliver efficient, scalable results on QPatch Compact or any e-phys device dependent on homogeneous single cell preparations.



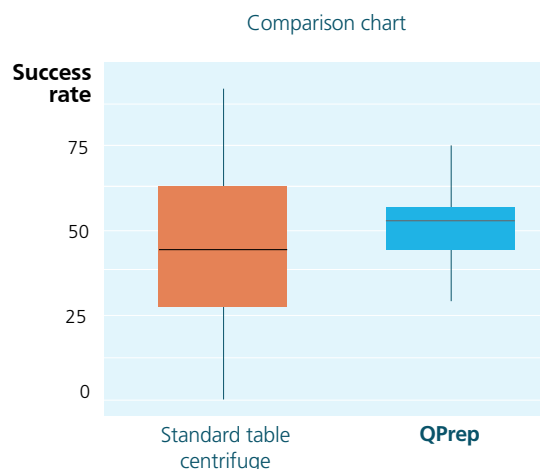
The centrifuge insert holds 1.5 mL lidless tubes (max spinning volume 1.0 mL) and operates at speeds up to 200 G. There's no maximum time limit, allowing flexible, ad hoc usage. The built-in magnetic stirrer ensures efficient cell stirring during processes.

- Cells kept gently in suspension with onboard QStirrer
- Consistent cell handling through precise pipetting guidance
- Unique 90° swing-out centrifuge ensures homogeneous cell separation
- Preset settings for reproducibility, good for GLP environment
- Real-time status is shown using indicator lights
- Clean and easy to operate
- Built with QPatch Compact workflows in mind

Sophion QPrep performance comparison

Proven superior, consistent cell handling

Built on 20 years of experience from QPatch® automatic cell preparation, the QPrep delivers more reproducible cell preparation than using a standard table centrifuge where success rates varies.



Technical specs

Width x Depth x Height (mm)	160 x 323 x 152
Touch screen size (mm)	95 x 65
Weight	5.95 kg / 13.1 lbs
Footprint	517 cm ²
Power supply	100-240 V, 50-60 Hz, Max 0.48 A
Centrifuge positions	1
Centrifuge insert volume, lidless tubes	1.5 mL (max spinning volume 1.0 mL)
Centrifuge speed	Up to 200 G
Centrifuge time	Ad hoc, no max limit, steps in seconds
Cell stirrer	Magnetic stirrer built-in
Magnet stirring bar	30 x 3 mm
Cell stirrer cup volume	60 mL
Connecting to PC/QPC	Prepared for software upgrades
Programs	Ad hoc. Possibility to save up to 5 presets
Tools	4 mm Allen key
Pipette tips to be used	ARTTM Wide bore filtered tips, 1,000 µL, cat.no 2079G
Light-signaled status:	
White:	Idle (breathing) or after spin cycle (constant)
Green:	Completed cycle (constant) waiting for user acknowledgement
Blue:	+/- Accelerating (two rolling) or steady speed (one rolling)
White/Red:	Emergency brake engaged or other fault (rolling)
Light intensity	Adjustable
Safety	Build in emergency brake, triggered by opening lid
Certificates	CE



GD43556-6