



Technical Specifications:

QPatch - automated patch clamp system

Performance/features	QPatch	
Hardware modules (standard)	Automatic cell preparation QPlate and compound plate stacking	
Hardware modules (optional – all can be retrofitted)	Temperature control; heating / cooling at recording site	
Unattended operation	>4 hours, up to 15 QPlates	
Parallel test capacity	16	48
Number of amplifiers	16	48
Number of pipettes	4 or 8	8
Target throughput per month	<10,000	<30,000
Pipette types	TiO ₂ coated on both inner and outer surfaces	
Stimulation/clamp modes V, L, P, and C are individual for each site	Voltage-gated Ligand-gated Pressure Current clamp (optional) Temperature (optional)	
V _{xx} adaptive voltage-clamp	√	
I _{adapt} adaptive current-clamp (requires current clamp)	√	
Recording configuration	Whole-cell / Perforated patch	
Cell types applicable	Cell lines Stem cells Primary cells iPSC	
QPlate compatibility	Single-hole Multi-hole Custom-size QPlates	
Maintenance of electrodes / User maintenance of instrument	None / None	
No. of electrodes per recording site	Individual pairs for all sites	
Electrode stability	Drift in voltage offset (V _{off} ~ 0.005 mV/min)	
Seal quality	True GΩ-seal in physiological solutions	
Need of seal fortifying agents	None	
Test compound consumption, per well	2-15 μL User-configurable and can be broken into multi-spit	
Total possible liquid addition, per well	250 μL	
Liquid added accounting	Automatic	
Liquid journey to cell	Individual microfluidic channels	
Liquid exchange rate	τ < 40 ms	
Compound plate formats	MTP-96 (SBS standard)	
Fast R _{series} compensation, τ < 400 μs	√ (optional)	
R _{series} compensation, τ > 400 μs	√	
C _{total} , C _{fast} & C _{slow} , compensation	√	
Analyzer software	√ (unlimited licenses)	
Bandwidth	20 kHz	

RMS noise	< 14 pA @ BW=20 kHz < 4 pA @ BW= 5 kHz <1.6 pA @ BW= 1 kHz
Sampling rate, maximum	500 kHz, digitally down sampled to 50 kHz, 16 bit
Filter options	Butterworth; 2, 4 & 8 th order Bessel; 2, 4 & 8 th order
Current range	±25 nA, ±50 nA, ±100 nA Largest range is automatically forced with multi-hole QPlates
Resolution	16 bit
Data storage and security	External database, automatic backup, full log of activity, user-hierarchy
External database operating system and version	Windows 10 Pro Oracle 19c
Internal PC operating system	Windows 10 LTSC

Dimensions & Requirements	QPatch 16	QPatch 48
Width	104 cm	
Depth	78 -121 cm (open)	
Height	174 - 199 cm (open)	
Weight	350 kg (772 lbs)	
Point pressure	10 kg/cm ³	
Foot print	0.55 m ²	
Power supply	AC 100-240 V 50-60 Hz Max 6 A	
Pressure	4 - 8 Bar	
Vacuum	0.7 - 0.9 Bar (1 m ³ /hour)	



Sophion Bioscience A/S
 info@sophion.com
 sophion.com