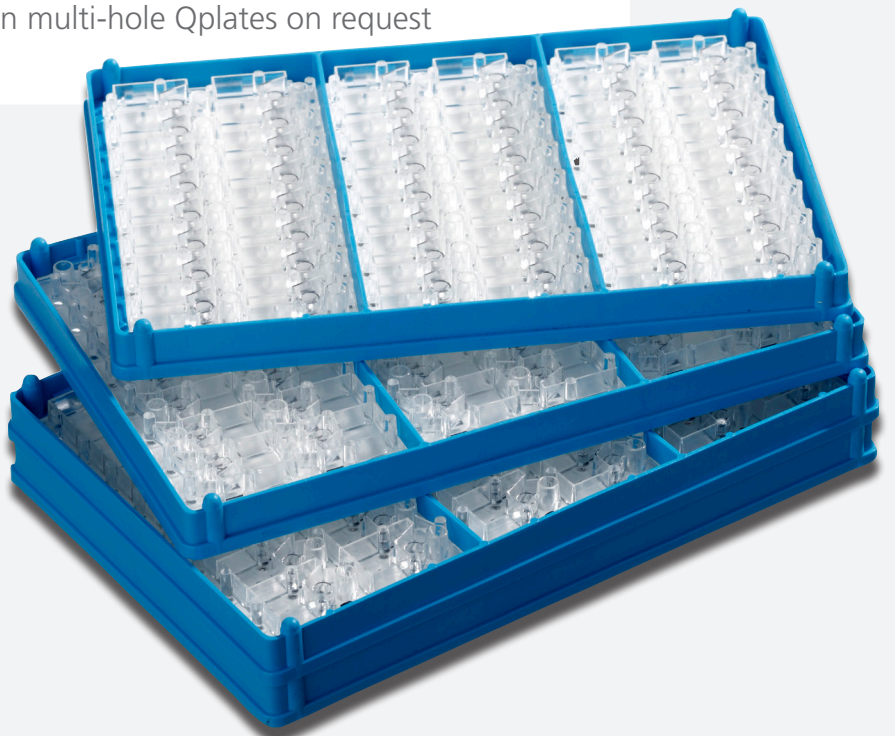


Product Specification

QPlate Measurement Plate

- Genuine giga-ohm seal formation
- Unique flow channel structure, enabling washes and application of multiple concentrations/compounds per well
- Glass-coated flow channels minimize absorption of compounds
- Cell positioning and whole-cell configuration formed by suction
- Both voltage- and ligand-gated applications
- 8, 16 or 48 individually controlled patch clamps measurement sites
- Single- and multihole technology
- Different patch clamp hole-sizes on request
- Different numbers of holes in multi-hole Qplates on request



QPlate - Measurement Plate

The heart of the QPatch technology is the silicon-based patch clamp orifice which replaces conventional glass pipettes.

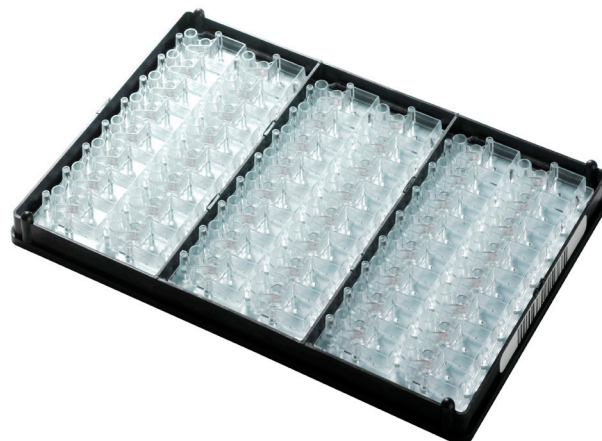
Sophion's proprietary QPlate technology enables the standard formation of genuine classical gigaseals, ensuring uncompromised data quality.

Another unique feature of the QPlate is the compound application principle, which is based on a microfluidic flow channel leading compounds and buffers to the cell. The feature enables the sequential application of a series of compounds or concentrations to the same cell, thus increasing throughput and reducing cost per data point.

Which QPlate to use:

	QPlate 48	QPlate 48X	QPlate 16	QPlate 16X
QPatch 8			•	
QPatch 8X			•	•
QPatch 16			•	
QPatch 16X			•	•
QPatch 48	•	•		
QPatch 48X	•	•		

The unique glass-coated flow channels minimize absorption of compounds. The QPlates exist in both single- and multi-hole versions. Different sizes and numbers of holes (for multi-hole QPlates) are available on request.



QPlate Dimensions:	QPlate 48	QPlate 48X	QPlate 16	QPlate 16X
L x W x H (mm)	128 x 85 x 7	128 x 85 x 7	128 x 85 x 7	128 x 85 x 7
Weight	68 g	68 g	30 g	30 g
Measurement sites	48	48	16	16
Patch clamp holes per site	1	10*	1	10*
Chip material	Si/glass	Si/glass	Si/glass	Si/glass
Chip capacitance	55 ± 5 pF	55 ± 5 pF	55 ± 5 pF	55 ± 5 pF
Orifice resistance	2.0 ± 0.4 MΩ	0.2 ± 0.04 MΩ	2.0 ± 0.4 MΩ	0.2 ± 0.04 MΩ
Max. pipetting vol.	15 µl	15 µl	15 µl	15 µl
Min. pipetting vol.	2 µl	2 µl	2 µl	2 µl
Waste reservoir	250 µl	250 µl	250 µl	250 µl

*QPlates with alternative number of patch clamp holes per measurement site are available on request.

Sophion Bioscience A/S, Baltorpvej 154, 2750 Ballerup, Denmark
Phone: +45 4460 8800 Fax: +45 4460 8899, E-mail: info@sophion.com

sophion.com