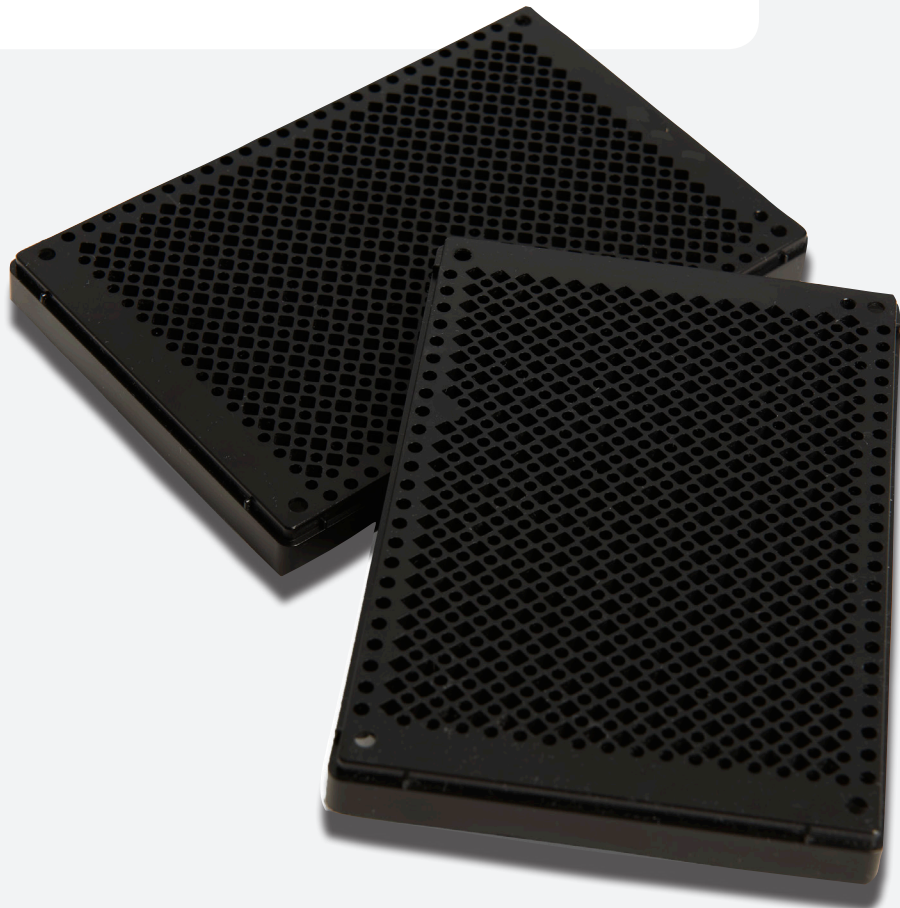


Product Specification

QChip 384 Measurement Plate

- 384 individual patch clamp sites with integrated flow channels
- Maintenance free integrated individual electrodes
- Both voltage- and ligand-gated applications
- Single-, 10-hole and variable number of holes are standard
- Custom number of holes on request
- Custom size of holes on request



QChip 384 Measurement Plate

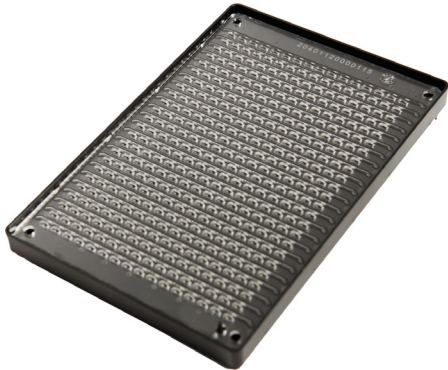
One of the core technologies in Qube is the QChip 384. QChip 384 integrates 384 individual patch clamp into one plate.

QChip 384 has built-in microfluidic flow channels that ensure fast and complete exchange of liquid. These channels make it possible to measure reliably on ligand- as well as voltage-gated ion channels while also adding sequentially an unlimited series of compounds to the same site.

Features & outcome

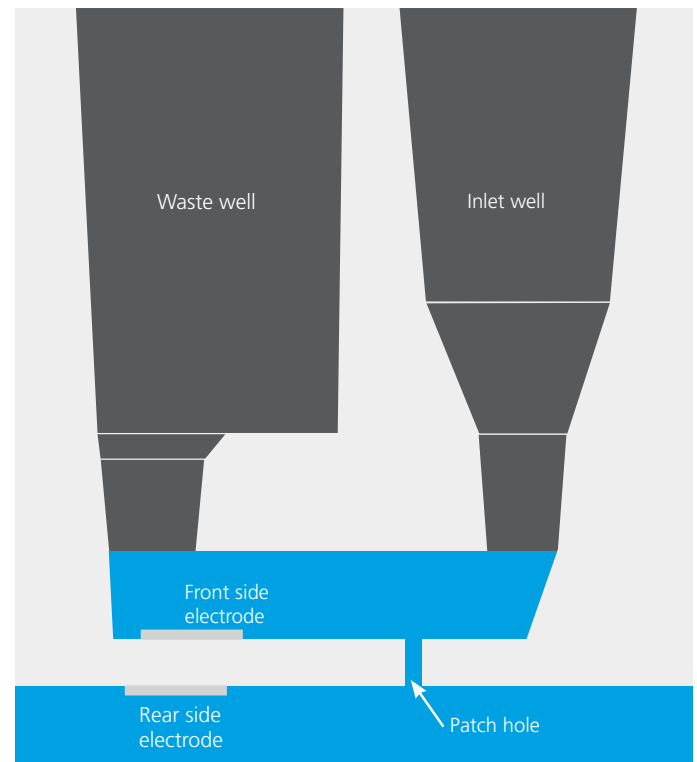
Features	Outcome
Integrated individual flow channels	Precise, fast and complete liquid exchange allowing for infinite liquid additions
Integrated individual electrodes	Stable recording over long periods of time without the need of maintenance
Barcode	Makes Qube recognize type of QChip e.g. for employing Rs-compensation, current clamp* or not and for ensuring full traceability of results

*Optional feature for Qube



QChip 384 has conventional Ag/AgCl-electrode pairs for each site allowing for stable recordings over a long period of time. There is no need for electrode maintenance and the QChip is always ready to use

QChip 384 exists in versions with both a single-hole (QChip 384), a 10-hole (QChip 384X), a variable number of holes (QChip 384D) and custom number of holes.



Dimensions	QChip 384	QChip 384X	QChip 384D	QChip 384 Custom
L x W x H (mm)	128 x 86 x 11	128 x 86 x 11	128 x 86 x 11	128 x 86 x 11
Weight	74 g	74 g	74 g	74g
Measurement sites	384	384	384	384
Patch clamp holes per site	1	10	1, 2, 6, 10, 16 & 36	Optional (=nn)
Site resistance	2.0 ± 0.4 MΩ*	0.2 ± 0.04 MΩ*	[0.06 - 2] MΩ*	Depending on nn
Volume surrounding cell	2 µl	2 µl	2 µl	2 µl
Waste reservoir	26 µl	26 µl	26 µl	26 µl
Possible liquid additions	∞	∞	∞	∞

* Depends on solutions

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

sophion.com