

SyncroPatch 384 NPC-384 Chips

Consumables tailor-made for your experiments

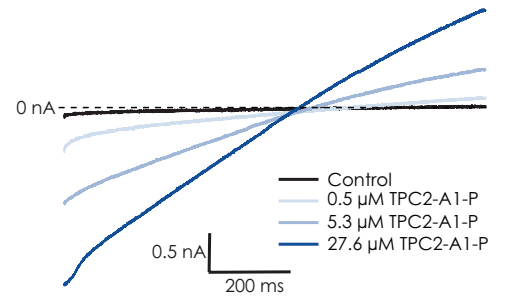
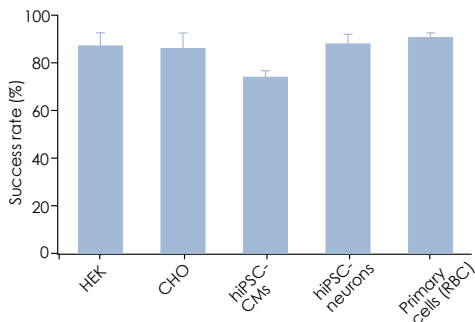
- 1 Borosilicate glass substrate
- 3 Large range of applications
- 5 Low adsorption
- 2 Made in-house
- 4 Customizable
- 6 Cost efficient

Chip type	Number of holes	Standard cell lines	iPSCs and primary cells	Mechano-stimulation	Organelles	Small cells (e.g. RBC, T-Cells)	Fast ligand	Fluoride-free
NPC-384T S	1x, 4x, 8x	●	●	●		●	●	
NPC-384T L	1x	●	●	●			●	
NPC-384T S F/M-STIM	1x, 4x, 8x	●	●	●		●	●	
NPC-384T L F/M STIM	1x	●	●	●			●	
NPC-384T microS	1x, 4x				●	●		
NPC-384T nanoS	1x, 4x				●	●		
NPC-384 FF high*	1x, 4x					●		●
NPC-384 FF medium*	1x, 4x							●
NPC-384 FF low*	1x, 4x							●

* Available as kit with solutions and chips

NPC-384 chips

For all your patch clamp needs



Lysosome-TPC2

Standard experiments

Excellent success rates for cell lines, iPSCs and primary cells.

Recording organelles

Lysosomes, mitochondria or other intracellular organelles.

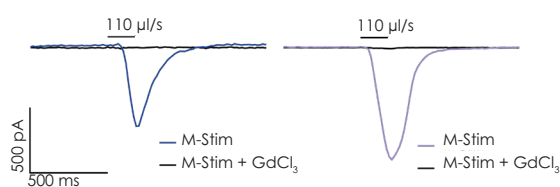


Mechano-activated ion channels

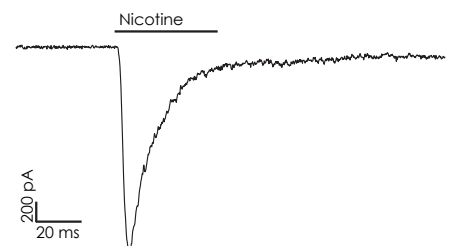
Mechanical activation of piezo (or other) channels.

Fast ligand-gated ion channels

Reliably record fast desensitizing ligand-gated ion channels.



HEK-Piezo1



HEK-nAChR α 7



Ready to start your SyncroPatch 384 experiment?

Scan the QR code & order your chips!