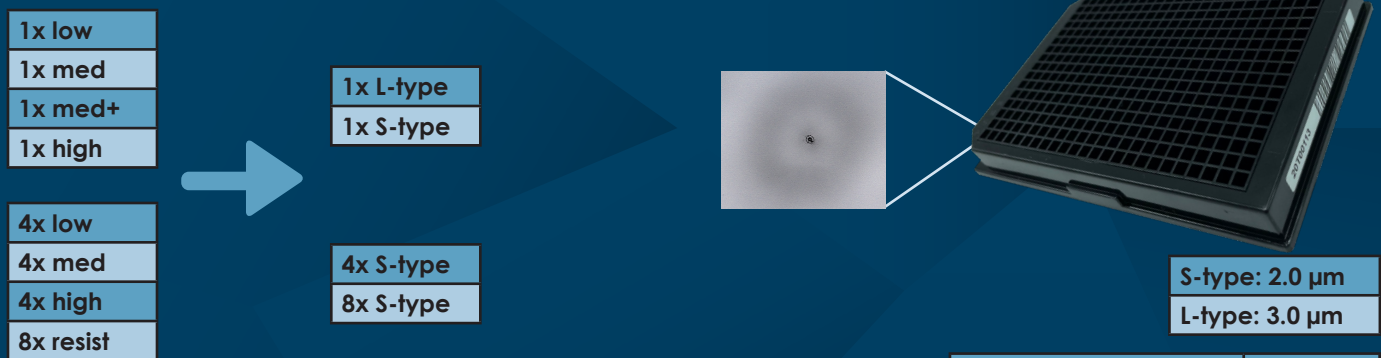


## New Consumables: NPC-384T

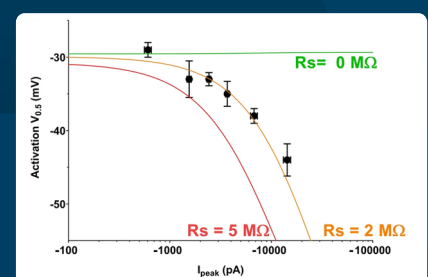
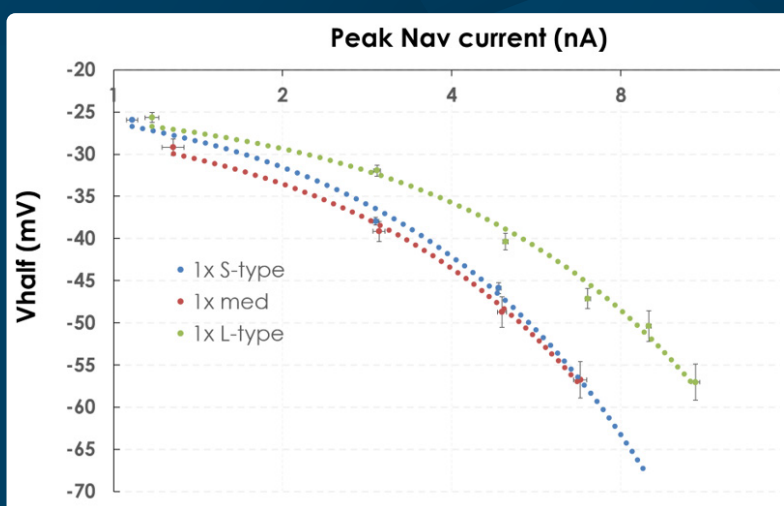
- NPC-384T chips are made from a thinner version of the same borosilicate substrate used for other NPC chips.
- Similar or higher success rates obtained by internal and external evaluations when using NPC-384T compared to current standard chips. Automated recognition by software and only minor adjustments in the experimental settings required.
- The thin glass reduces the access resistance and improves voltage control, representing the future version of consumable for the SyncroPatch 384.

### Standard Chips Types / NPC-384T Equivalents



Chiptype	Order #
NPC-384T 1x S-Type	22 2101
NPC-384T 1x L-Type	22 2102
NPC-384T 4x S-Type	22 2401
NPC-384T 8x S-Type	22 2801

### Voltage Clamp Properties



Montnach J et al. Sci Rep. 2021 Feb 8;11(1):3282.  
doi: 10.1038/s41598-021-82077-8

- 1x S-Type - and 1x med standard chips show similar voltage clamp properties
- 1x L-Type chips allow for less Vhalf deviation over a broader range of current amplitudes