

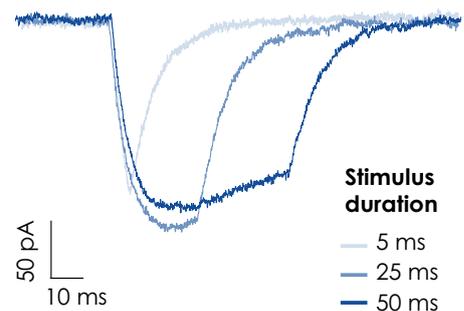
SOL

Optical stimulation for the Port-a-patch

Brighter discoveries start here

- Ideal for light activation of ion channels
- Activation and deactivation of caged or photoswitchable compounds
- Add-on for the Port-a-Patch
- Choose from a range of excitation wavelengths
- Up to 3 independently controlled light sources possible

Light activation of channelrhodopsin 2 (ChR2)



Activation of ChR2 by increasing light pulse length.



Contact us today

SOL for the Port-a-patch

Activate ion channels and compounds using light

Integrate light stimulation with patch clamp recordings

- SOL simply replaces the standard Port-a-Patch lid with an optical lid containing a high-power LED light source.
- Enables simultaneous patch clamp and optogenetic or uncaging experiments.
- Compatible with perfusion and temperature control for fully integrated experiments.
- Synchronizes light activation with voltage or current clamp recordings for precise control.



Achieve precise, programmable light control

- Millisecond-scale timing precision for fast channel or compound activation.
- Control light intensity, timing, and duration directly via software.
- Automate experiments with scriptable light pulses coordinated with recording sequences.
- Swappable LED modules allow flexible wavelength selection for diverse applications.