



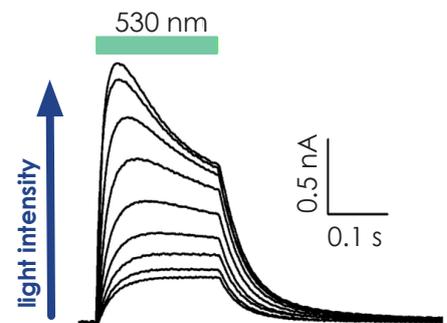
SOL

Optical stimulation for the SyncroPatch 384

Light up your experiments

- Ideal for light activation of ion channels
- Activation and deactivation of caged or photoswitchable compounds
- Add-on for the SyncroPatch 384
- Activate with light in 384 wells simultaneously
- Dual wavelength excitation possible
- Choose from more than 10 different wavelengths

Light activation of channelrhodopsin



Activation of ChR by increasing light intensity, adapted from [Govorunova et al, 2022](#).



Contact us today

SOL for the SyncroPatch 384

Activate ion channels and compounds using light

Choose from more than 10 different wavelengths

- Couple voltage protocols and light activation.
- Each Illuminator can be equipped with two independently controlled wavelengths.
- Dual wavelengths can be used simultaneously or separately.
- Choose and combine different wavelengths to find the right setup for your experiments:

Stimulate 384 wells simultaneously

- Base designed to fit on top of the SyncroPatch 384 measurehead.
- Use one to four illuminators depending on your assay format and throughput needs.
- Temperature-stable operation ensures reliable optical power during recordings.
- Easy to remove for experiments not requiring light stimulation.

Light type	Dominant or Peak wavelength (nm)	
	Minimum	Maximum
 UV*	380	390
 Visible	-	-
 Royal Blue	440	460
 Blue	460	480
 Cyan	490	510
 Green	520	540
 Amber	588	592
 Red-Orange	610	620
 Red	620	645
 Deep Red	650	670

*Other wavelengths for UV LEDs available. Custom wavelengths also available. Please ask us for details.

