



QPatch Compact Opto 465

Stimulate light-sensitive ion channels

With the **Opto 465 integration kit**, you can take full control of your ion channel experiments—using light to explore cellular excitability, ion channel function, and pharmacology like never before. By pairing optogenetic stimulation with automated patch clamping, you will reveal deeper insights faster. And because QPC Opto 465 is designed specifically for QPatch Compact, you get precise, reproducible stimulation and consistently high-quality data you can trust.

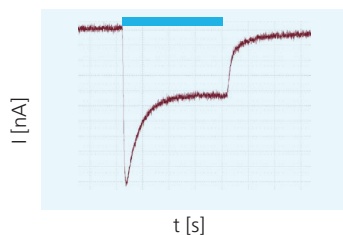
- Perfect for studying light-gated ion channels, light-sensitive chemistry, and light-sensitive internal biological processes
- Well-by-well control for accurate light-to-cell correlation and accurate data analysis
- Adjustable blue-light intensity for precise, reproducible activation
- Guided tracking of stimulation timing to ensure consistency and prevent errors
- Upgradeable wavelengths for future research needs



The custom made **Opto-pen** ensures precise delivery of the light quant without any fake light slipping in. The timing is controlled by the user and is logged by the software in the QPC to make the analysis accurate.

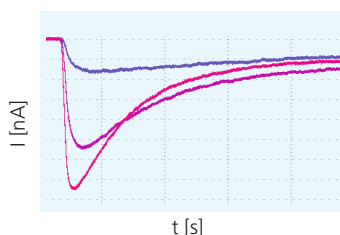
Control both the timing and intensity of light stimulation. Powerful enough to activate intracellular processes.

Optogenetics



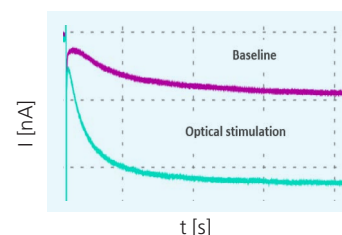
$\lambda = 465 \text{ nm}$
 $t = 200 \text{ ms}$
Channelrhodopsin-2

Optopharmacology



GABAAR
RUBI-GABA

Photoactivation



bPAC/HCN2

Technical specifications

The QPC Opto 465 is compatible with QPatch Compact of any generation and consists of 4 parts

LED for stimulation

Electrical connection

1 meter long pigtailed cable M8 connector
4 positions - male

Dimensions

63 x 59 x 21 mm with baseplate

Weight

100 g

Stimulation activation

Manually, via remote control



465 nm LED for stimulation

CLED_465: Connectorized LED: 465 nm

LED driver

TTL modulation

HI > 2.8 V
LOW < 2.3 V

Analogue modulation

0 - 5 V (400 mA / V)

Output current

40 - 2000 mA or 4 - 200 mA in low power mode
* 0.000 mA when OFF or V-IN < 0.1 V

Rise / Fall time

< 10 μsec

Dimensions

175 x 50 x 105 mm

Weight

580 g

Stimulation activation

Manually, via remote control



LED driver, power source for the LED

LEDD_1: LED driver, 1 channel model
(with adapter for EU, US, UK and AUS)

Opto pen - optic patch cord

Length

50 cm

Material

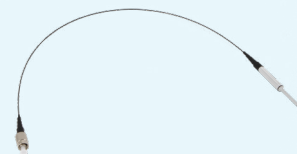
Silica/Silica

Core

400 μm

Jacket

900 μm



Designed for guided insertion
into the QPC manifold
Mono fiberoptic patchcord

Remote control

Stimulation modes

Trigger: Gives one light impulse with
length equal to as long as suppressed
Gate: Is an on/off switch to light stimulation

Logging of stimulation

Timing is guided by QPatch Compact software



External TTL pulser - including USB
power cable and BNC signal cable

GD45922-2