

Detailed project description:

Background and scientific relevance:

(describe the scientific background of your project and explain why it is relevant to your laboratory).

Feel free to link to relevant publications.

How would the QPatch Compact system help you achieve your research goals:

What technologies do you currently have access to that will aid your use of an automated patch clamp system:
(i.e. alternative techniques or platforms, cell lines or expression systems, etc.)

Organizational information:

Project staff: Please provide information on the researchers, who will be working on the project and their qualifications.
(describe the ability of your group or department to support the project and instrumentation in the short and long term).

Collaborators: Please list any collaborators involved in the project or who would be interested to utilize our technology if you are awarded the QPatch Compact instrument grant.

Organization funding: Please describe the current funding sources used by your department for ion channel research and the approximate amount of funding.

Additional information:

Feel free to add any additional information you think is relevant to us, here: