

## Application form

Do you have an ion channel research project that you cannot start because you lack the resources, time, and funding to move forward?

With the Sophion instrument grant programme, we offer researchers and scientists the unique opportunity to use Sophion's industry-leading automated patch clamp solutions free of charge for one year.

If you are interested in applying for a Sophion instrument grant, please start your registration process by completing the application form below.

You must save the form to your PC and send us your form submission when all required form fields have been completed to [info@sophion.com](mailto:info@sophion.com)

### Basic information:

First name:	Last name:
Job title:	Email:
Company name:	Phone:
Address:	Country / post code:

### Project information:

Primary project title:			
Requested instrument:	QPatch 16	QPatch 48	Qube 384

### Abstract:

**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 are you currently acquiring patch clamp data:**

**How would an automated patch clamp 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.)

**Please list any additional projects, whether your own or lead by another investigator, who would be supported by a grant instrument:**

Additional project title:

Primary principal investigator name:

Primary principal investigator email:

**Abstract:**

**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.

**Detailed project description:**

**Please list any additional projects, whether your own or lead by another investigator, who would be supported by a grant instrument:**

Additional project title:

Primary principal investigator name:

Primary principal investigator email:

**Abstract:**

**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.

**Detailed project description:**



**Please list any additional projects, whether your own or lead by another investigator, who would be supported by a grant instrument:**

Additional project title:

Primary principal investigator name:

Primary principal investigator email:

**Abstract:**

**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.

**Detailed project description:**

**Additional information:**

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