



## Application Form

**Sophion Bioscience Early Career Travel Grants** are awarded to students and early career researchers in electrophysiology, drug discovery and related science.

Applicants can be anyone who:

1. Has performed research in the field of ion channels modulation and characterization and would like to present at Sophion Ion Channel Modulation Symposium (Poster/Oral)
2. Has performed research on QPatch or Qube 384 and wish to present data through active participation in any symposium, congress or Sophion User Meeting

The following requirements must be met for your application to be considered for the Sophion Bioscience travel grant:

- Early career includes MSc or PhD students or researchers within five years of BSc/MSc/PhD award
- The grant must cover expenses related to your stay (travel cost, hotel)
- The applicant must be enrolled at a higher education institution
- The applicant must be endorsed by their advisor or a senior researcher at their institution

Full name: .....

Address: .....

Title: .....

City / Postcode: .....

Department / Group: .....

Country: .....

Institution .....

Email address: .....

Which conference would you like to attend?

Preferred presentation format

Application window:

Sophion Ion Channel Modulation Symposium

Oral

February

Other (please specify) .....

Poster

September

Description of research/study area (<300 words):

Presentation title and abstract (<300 words):

Why would you like to go? What is your motivation? (What do you believe you may learn/gain, people to meet, discuss, collaborations, etc.):

Conflict of interest disclosure (if relevant):

\*Endorsement by advisor or senior researcher from institution must be submitted together with this form, and sent to [grant@sophion.com](mailto:grant@sophion.com).