



Application Form

Research Grants for MSc or PhD students & early career researchers within five years of BSc/MSc/PhD award in electrophysiology, drug discovery & related science.

The grant can be used in connection with an MSc, PhD or post doc or as Technology Access Program for existing research projects and usually includes a one to 10 weeks stay at our R&D laboratories in either Tokyo, Boston or Copenhagen with accommodation and lab expenses included.

An application can be made by students or early career researchers performing research in the field of ion channels that would benefit from access to high performance, high throughput automated patch clamp instrumentation.

Full name:

Address:

Title:

City / Postcode:

Department / Group:

Country:

Institution / Company:

Email address:

Where would you like to go?

☐ Sophion Copenhagen (headquarters)

☐ Sophion Boston

☐ Sophion Tokyo

When would you like to go?

☐ Quarter ☐ Year

Estimated duration:

Weeks:

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

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

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.