

Guidelines for abstracts submission

You are invited to submit an abstract on original work for presentation as a poster. All abstracts submitted will be reviewed by the committee and you will be notified end of September 2024 if your abstract has been accepted.

You do not need to be registered for the conference before abstract submission, but once your abstract has been accepted, the presenting author must register and be present at the meeting. Fill in the form, save it and send it to [Sophion by email](#).

Poster title:

- Keep it brief - max. 40 words

Authors and affiliations:

- List authors and co-authors
- Full name
- Comma separated names,
- No academic degrees

Presenter name & affiliation:

Abstract:

- 500 word limit excl. title, authors etc.
- Do not include figures, graphics or references

Best Poster Award



Each year the ICMS® Scientific Board acknowledges the most outstanding poster presented at a specific Ion Channel Modulation Symposium in either Europe, United States or Japan.

The evaluation criteria for identifying the best poster include:

1. Background provides appropriate perspective/context for the subject
2. Objectives/research questions are clearly stated
3. Research design/methods/modeling is appropriate and transparent
4. Data sources and/or sampling procedures are clear and appropriate
5. Research objectives are met/addressed
6. Factual information is kept separate from interpretations or implications
7. Abstract is presented in an unbiased manner
8. Clarity of presentation

PLEASE NOTE: The winner **MUST** be present in the auditorium when the award winner is announced.