

Ion Channel Modulation Symposium (ICMS) – Japan 2024

Information about “GAKUPOSU” Poster Printing Service

Reception period

May 7th (Tue)JST - May 27th (Mon) JST, 2024

*It will be very crowded near the end of the reception period. We appreciate your cooperation in applying early.

Please note that the deadline cannot be extended.

service

We will print the poster and deliver it to the venue.
(Can be pasted if requested)

Cost and paper type

- | | |
|-------------------|---------------------------|
| ①Semi-gloss paper | 16,500 JPY (tax included) |
| ②Non-woven fabric | 25,300 JPY (tax included) |
| ③LIMEX | 20,900 JPY (tax included) |

Payment method

Credit card payment only

We cannot accept data changes after submission/payment is complete.

Click here to print
your poster

Go to the venue
empty-handed on the
day of the event.



Click here for the
application/submission site ↓

<https://www.kinkosmpm.jp/shopm/icms2024>

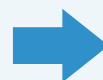
Jump to the dedicated site by clicking the URL
above or the banner on the left.



New user
registration



Login



Submission/
payment

Requests when submitting data

Various manuals and a submission data template here

[Click](#)

①Please submit your application in PDF format.

- When submitting PowerPoint data, characters may be misaligned or errors may occur.

②Printing is available at a size of Width 841mm x Height 1,189mm.

- As a general rule, editing and proofreading will not be performed.
- We will enlarge and print the data you provided to W 841mm x H 1,189mm.
- The long side is finished at 1,189mm in consideration of visibility for the W 850mm x H 1,500mm bulletin board size.
- 5mm margins are automatically adjusted on each of the four sides of the rectangle.

③Please write "Your Name" as the file name.

- We kindly ask for your cooperation in preventing the mixing of customer data and checking the layout location during posting work.
- If the "presentation number" is not listed, you will not be able to post the presentation. Please understand that this is a required item.

■ Contact information regarding poster creation

Kinko's Japan Co., Ltd.

Gakuposu Service Team

E-mail: **icms2024@kinkos.co.jp**

Counter business: Weekdays 10:00-17:00 JST

Email inquiries: Presentation number/name