



BECOME A SPONSOR

We invite you to become part of the Ion Channel Modulation Symposium (ICMS®), Japan hosted by Sophion Bioscience. The ICMS will take place Thursday the 30th and Friday the 31st of May 2024 and the event will be held at Tokyo University, Japan. The ICMS brings together a broad range of people from academia, CROs and the pharmaceutical industry to discuss the science of ion channel modulation. Over 20 speakers from labs around the world will present at the symposium to around 130 attendees.

Join us in advancing the science of Ion Channel Modulation

Why sponsor ICMS in Tokyo?

Exposure:
Showcase your brand to a global audience of ion channels experts and decision makers.

Custom Packages:
Tailored sponsorship options to meet your objectives and requirements.

Innovation:
Stay ahead with insights into the latest research and technology.

Networking:
Connect with industry leaders and scientists for potential collaborations.

Sponsorship packages	Platinum ¥700,000/€5,000	Gold ¥420,000/€3,000	Silver ¥70,000/€500
Company logo on the ICMS 2024 Japan website and in printed program	●	●	●
Table top exhibition space in exhibition area	●	●	
Roll up panel/zap stand (one per stand) to be used in the exhibition area	●	●	
One poster presentation	●		
Full symposium attendance for two persons	●		
Full symposium attendance for one person		●	
Invitation to the drinks reception	●	●	
Option to have literature available to attendees on shared table top in exhibition area			●

Please fill in the below form and send it by mail to Erina Katayama eka@sophion.com. Once the application has been approved we will inform you by email. We will need a purchase order from you in order for us to issue an invoice. Your sponsorship is not secured until we have received the your payment. We reserve the right to reject any sponsorship applicant.

Company	
Address 1	
City and Post code	
Contact person	
Mail address contact person	
Phone no.	
Type of sponsorship	Platinum: <input type="radio"/> Gold: <input type="radio"/> Silver: <input type="radio"/>
PO number:	
Signature & date:	