



ICMS '25 SAN FRANCISCO
ION CHANNEL MODULATION SYMPOSIUM

SEPTEMBER 23-24, 2025
UNIVERSITY OF CALIFORNIA,
SAN FRANCISCO, CA, USA

BECOME A SPONSOR

Join us in advancing the science of Ion Channel Modulation

We are happy to invite you to become part of the **Ion Channel Modulation Symposium (ICMS®)** in the USA, hosted by Sophion Bioscience. The ICMS will take place on **23 -24 September 2025** at University of California, San Francisco. The ICMS brings together a broad range of people from academia, CRO's and the pharmaceutical industry to discuss the science of ion channel modulation. More than 20 speakers from labs around the world will present at the symposium to around 200 attendees.

Why sponsor ICMS in San Francisco, USA?

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 € 5,000 \$ 5,400 ¥ 820,000	Gold € 3,000 \$ 3,200 ¥ 500,000	Silver € 500 \$ 550 ¥ 82,000
Company logo on the ICMS 2025 USA website and in printed program	•	•	•
Table top exhibition space in exhibition area	•	•	
Roll up panel (one per booth) 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			•

Other opportunities	Half page (A5 landscape)	Full page (A4 portrait)
Advertisement in printed programme	€ 500 / \$ 550 / ¥ 82,000	€ 1,000 / \$ 1,050 / ¥ 160,000

Please fill in the below form and send it by mail to Duncan Jarman dja@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 zip code	
Contact person	
Mail address contact person	
Phone no.	
Type of sponsorship	Platinum: <input type="radio"/> Gold: <input type="radio"/> Silver: <input type="radio"/> Half-page ad: <input type="radio"/> Full-page ad: <input type="radio"/>
PO number:	
Signature & date:	