

Join us for a dinner cruise on San Diego harbor



Axion BioSystems and event co-sponsor Nanion Technologies invite you to be our guest:

What A relaxing dinner cruise around San Diego harbor aboard a private yacht

When Monday, November 14th

Where Boarding at the Marriott Marina adjacent to the Convention Center (333 West Harbor Drive)

Time Boarding begins at 5:45pm; yacht sails promptly at 6:00pm

Return to the Marriott Marina at 8:45pm

RSVP dinner-cruise.eventbrite.com
(reservations may include one guest)

Questions? Please contact Catherine Parrish (cparrish@axionbio.com).