

nanji[on

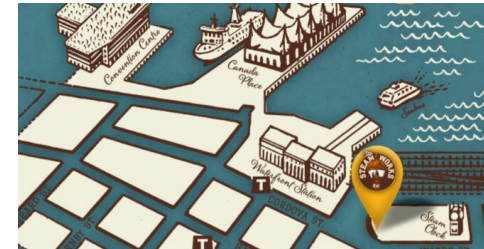
Meet Nanion at the SPS 2016: Nanion's Networking Reception:

Global VIP meet & greet – CiPA highlight dinner

Directly after our hosted company session at the SPS we move on to our after dark event!

Enjoy the great networking opportunity with Nanion's scientists and friends.

Monday 19th of September, 19:30 pm,
Steamworks Brew Pub, Wine room,
375 Water Street, Vancouver
GPS: 49°17'06.6"N+123°06'38.9"W
(600 m distance from Convention Center)



Meet Dr. Sonja Stoelzle-Feix, Director, Scientific Affairs at Nanion and in the Leadership team of CiPA High throughput screening (HTS) team

Atsushi Ohtsuki, Customer Support Manager Nanion, Japan

Corina Bot, Senior Scientist at Nanion Inc., USA



Don't miss:

Nanion's Hosted Company Session
Nanion's Poster on cardiac safety
Nanion's Networking Reception