



Hotel Nikko Tokyo in Odaiba, overlooking Tokyo Bay. The name 'Odaiba' originated from the late Edo period (1603-1867) and derives from the government-built artillery platforms (daiba) to protect Edo, the old Japanese capital that is now Tokyo, from attack by foreign ships.

Photo provided courtesy of Hotel Nikko Tokyo

Asia Pacific Sales Meeting in Japan

Tecan's Asia Pacific 2008 Sales Meeting was held from February 18 to 19, at the Hotel Nikko (Japan Airlines) Tokyo in Odaiba, located in the Tokyo Bay area. The annual sales meeting, co-hosted by Tecan Japan and Tecan Asia Pacific, brought together Tecan's regional offices and sales agencies, and was attended by participants from Australia, China, India, Korea, Malaysia, Singapore, Taiwan, Thailand and Japan.

The meeting was opened by Carl Severinghaus, Senior Vice President, followed by a review of 2007 and future plans for 2008 in Asia Pacific and Japan and summaries by managers of each business unit. There were announcements and information about new products from head office representatives and, with a presentation from each regional office and sales agency, it was an excellent opportunity to exchange information among the Asian countries. Five section meetings were held for liquid handling, detection, REMP sample management, components, and service and consumables, during which the opportunity for hands-on training for new products was enthusiastically received.

On the evening of the first day, attendees enjoyed an evening boat cruise and Japanese cuisine such as tempura,

sashimi and chanko-nabe (one-pot stew), gazing upon the night view of the seaside city sub-center from a yakata-bune – traditional Japanese cruising boat – floating in Tokyo Bay. Further entertainment during the feast came from performances by dancers, called 'furusode-san' ('furusode' are long sleeved kimono, traditional Japanese formal attire), and the evening was perfect for members of the team to catch up with each other.



A section meeting in progress



Attendees at the Asia Pacific 2008 Sales Meeting



'Furusode-san' entertained on the first evening