

Gambarou Nippon! Let's do our best together Japan!

In the aftermath of the deadly earthquake, tsunami and radiation threat in March, there was widespread fear that this disaster would also impact on Japan's national economy, with the country being left behind in the global market. For the first couple of months after the disaster, most of the tradeshows, festivals and concerts were cancelled, with International BioEXPO – the annual life sciences trade show held in Tokyo in June – among those threatened. Making a success of this show had special significance for Tecan Japan, helping to create a positive impression and further encouragement for the Country's industry, and we are pleased to report that the show was highly successful, attracting similar visitor numbers to last year.

"The devastation caused to the earthquake and tsunami-hit area is huge, and far beyond any description," commented Miyako Sakai, Marketing Communications at Tecan Japan. "With more than 20,000 dead and missing, many lost their loved ones, houses and jobs. This tragedy has reminded the Japanese

people of their fortitude in the face of adversity and of the strong ties we have with the rest of the world; we appreciate very much the support and prayers we have received from across the globe. Life goes on and we are hanging in there with our favorite slogan: 'Gambarou Nippon!' Let's do our best together Japan!"



Tecan exhibits at International BioEXPO 2011

Tecan and its employees worldwide have contributed a total of CHF 50,000 to the relief effort through Swiss Solidarity, a foundation that raises funds for humanitarian projects and works closely with the SRG SSR (Swiss Broadcasting Corporation) and relief organizations around the world.

がんばろう 日本



Tecan-sponsored graduate team triumphs in Californian sun



TMP project team

Tecan is pleased to announce that a Company-sponsored team has won the Keck Graduate Institute (KGI) Team Masters Project (TMP) Collaborative Excellence Award, part of the Institute's Master of Bioscience (MBS) degree program focusing on business and

management skills within the life science sector. The TMP is the capstone of the MBS program, and is designed to encourage innovative, application-focused thinking as multidisciplinary teams of postgraduate students work with industry figures to solve real problems using novel and existing technologies.

Each year, leading companies from throughout the life sciences industry are invited to sponsor individual teams, helping to investigate how cutting-edge technologies could be applied to novel or previously unexplored market sectors. This year's Collaborative Excellence Award was won by students Jessica Chang, Neelima Kumar,

Alice Lai, Jessica Lin, Sergio Sanchez and Jeffrey Weiss, who worked with Marc Feiglin and Tony Mamone from Tecan to explore new areas for the application of Tecan's technology portfolio. Craig Adams, KGI Faculty Supervisor for the team, commented: "This was an exceptional team; they worked closely together while maintaining their analytical objectivity and adding their individual perspective into each market analysis. Finally, Marc and Tony facilitated such a great project that we are thrilled Tecan has chosen to sponsor another TMP during the 2011-2012 academic year."