

Tecan supports ELRIG sponsorship of 'TechnoBotts'

Tecan continues to encourage the growth of innovative technology through a donation to help European Laboratory Robotics Interest Group (ELRIG), at the ELRIG Drug Discovery conference, gold-sponsored by Tecan, in Liverpool, UK in September 2009. Tecan contributed to the sponsorship of a promising group of young automation specialists named TechnoBotts, a UK-based team of teenage robotics developers. With funding from Tecan and other vendor companies, a check for £5,000 was donated to support the team in its venture to enter RoboCup 2010, an international robotics competition in Singapore in June 2010.

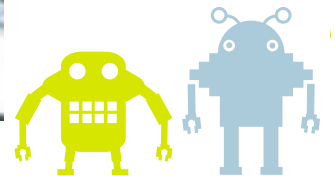
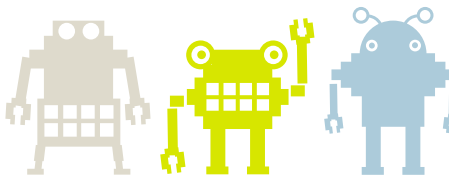
As team TechnoBotts³, Andrew Wightwick (14), Oliver Wightwick (12), Eric Wieser (15) and Alex Gilday (14) enjoyed great success in 2008 – 2009, becoming the UK and Ireland FIRST LEGO® League (FLL) champions. They also represented the UK in the FLL World Festival in Atlanta, Georgia in April 2009,

coming 4th in Robot Performance and winning the first place trophy for Innovative Design. For 2009 – 2010, Andrew, Oliver and Eric have re-formed as TechnoBotts⁴ and are looking for further success in RoboCup Junior, in both the Rescue and Soccer categories.

LEGO is a registered trademark of the LEGO Group.



(l to r) Kevin Moore (Tecan), the TechnoBotts team, Adrian Kinkaid (ELRIG), Robin Grimwood (FluidX)



Tecan to supply fully automated solution for Hologic Cervista® HPV tests

Tecan has made a global OEM agreement with Hologic, Inc., a leading developer, manufacturer and supplier of premium diagnostics and surgical products for the healthcare needs of women, to supply a fully automated solution for Hologic's Cervista molecular diagnostics human papillomavirus (HPV) tests. The system is based on Tecan's Freedom EVO® liquid handling platform with an integrated Tecan detection device. Prototypes of the automated solution have been tested successfully with select Hologic customers. Validation for regulatory approval of the automated system has started before the end of 2009, and Tecan expects to start shipping the Cervista system in the US in

2010. Tecan will also supply Hologic with plastic consumables as an integral part of the validated system.

HPV is the most common sexually transmitted disease in the US and is recognized as the cause of most cervical cancers. To help prevent the onset of disease, the American College of Obstetrics and Gynecology (ACOG) suggests routine Pap and HPV testing for women over the age of 30 to identify individuals most likely to develop cervical cancer.

Hologic and Cervista are trademarks and/or registered trademarks of Hologic, Inc. and/or its subsidiaries in the US and/or other countries.

