

Creative customer competition

Earlier this year we launched a competition inviting you to show us how much you love your Tecan and win a party for your team. We have received some great, creative photos showing just how important Tecan systems are to your teams, a selection of which are illustrated. We hope you enjoy working with your instruments as much as we love making them for you.

Burak Tiryakioglu,
EUROMMUN, Turkey



Kristen Tamburro,
UNC, USA

We only have one winner & prize... but Agata Favon-Gorecka and the Institute of Pharmacology, Polish Academy of Sciences, Poland was rated very highly



WE OUR TECAN



WINNER
Gema Santamara,
Pharma Mar, Spain





Natalya, FCI Gianniah, Russia



Andrew Flatley, Helmholtz Zentrum München, Germany

Bree Rumballe, Institute for Molecular Bioscience, Australia



Rosemary Nelson, Stryker Biotech, USA



Stefan Alatalo, Octapharma, Sweden



Also highly commended was Yener Selcut, MD GmbH, Germany



Jeanette Wu, Novartis Institute for Tropical Diseases, Singapore



Thomas Edmonds, Institut De Recherche Servier, France