

# And we have a winner!



*HydroFlex and Sunrise in the different colors available*

Discussion continues to rage over whether orange, blue, red or bright green is the best color for our new HydroFlex™ 3-in-1 platform and the popular Sunrise™ microplate reader. Many of you have joined in with your thoughts and for a chance to win one of these instruments in one of the available colors, or an Apple iPod nano 8GB.

Mrs Annamari Heiskanen of Glykos Finland Ltd in Helsinki, Finland, is the first winner of the iPod nano from the October draw. Annamari voted for red as her favorite colored Sunrise reader and won herself a shiny iPod in magenta.

Cell biologist Annamari Heiskanen works as a project manager at the biotech company Glykos Finland Oy in Helsinki. Glykos Finland Oy specializes in glycobiology and carbohydrate chemistry, offering product development and services to multinational customers in the field of functional foods and pharmaceuticals. The company's innovative development of bioactive carbohydrate molecules has led to a

variety of applications, such as biomarker screening from cell and cancer glycomes, as well as a product for treating gastric ulcers and development of functional food components. "The company is growing rapidly so we need new, reliable workhorse instruments for our laboratories," said Annamari. "It is a bonus if they also look inspiring!"

There is still time to join the debate. Visit us at [www.tecan.com/colors](http://www.tecan.com/colors) for your chance to win one of these prizes. A draw will be held at the end of every month until March 2008 for the iPod nano, and a separate draw in January 2008 for the HydroFlex or Sunrise.

*Mrs Annamari Heiskanen of Glykos Finland, receives her new iPod nano from Hanna Laiho at PLD Finland Oy*

