

Magellan™ offers new features for ELISA

m magellan
data analysis software

PLAY&WIN

To be in with a chance of winning Magellan v6.6 software, pre-installed on a netbook, go to www.tecan.com/magellan-package-deal

As from September 15

Tecan has set a new standard for ELISA assays with its Magellan v6.6 data analysis software. This latest version of Magellan supports eight languages to offer even more user-friendly operation, including Chinese, English, French, German, Italian, Portuguese, Russian and Spanish.

Magellan v6.6 is available as part of several packages designed to provide customers with attractive solutions for ELISA applications, even on low cost netbooks.

To find out more about these special offers, visit www.tecan.com/magellan-package-deal



Welcome
Magellan speaks
your language!

Tecan's roadshow is coming to the UK!

Tecan will be touring the UK from the end of August until the end of October, visiting 13 universities, research institutes and science parks to demonstrate the capabilities of the Freedom EVO® 100 automated liquid handling workstation and the Infinite® 200 series of modular, multimode plate readers.

Everyone is welcome to visit us at any of the venues and see our instruments in action. Take this opportunity to see Tecan systems in action and speak to our representatives to learn how automation can help you to achieve reliable results and make tedious manual handling a thing of the past.

For more information, please visit www.tecan.com/ukroadshow

