

EVO Logic™ simplifies Genesis™ trade-in

The Genesis workstation has been superseded by our more powerful Freedom EVO® platform and, from January 2012, we will no longer be providing support for the Genesis. Last year, many customers took advantage of our special offer to make the transition to Freedom EVO workstations, and our network of service and application specialists has been working closely with these customers to ensure this change has been as smooth as possible.

To further simplify switching to the Freedom EVO platform without disrupting existing laboratory processes, Tecan has developed EVO Logic. Designed to run alongside the new platform's Freedom EVOware® software, this open and flexible

solution is intended for laboratories currently using the Logic software for the Genesis platform, and helps to ensure full compatibility with existing working practices, protocols and LIMS interfaces. EVO Logic's sample-driven architecture makes it easy to transfer existing protocols and begin using your Freedom EVO workstation, without the need for extensive process validation or staff training.

To find out more on Tecan's EVO Logic software, visit

www.tecan.com/evologic

