

Freedom EVOware® Normalization Wizard

Software to help you normalize your DNA samples
to uniform concentration

A wizard to automatically perform the pipetting for quantitation and quality normalization of DNA samples

Tecan offers solutions for high-quality nucleic acid preparation for every throughput demand including PCR setup, reaction cleanup, sample quantitation and normalization. Quantitation and normalization of DNA samples improve the quality of the DNA sequencing data generated from the samples. Freedom EVO® automated workstations are capable of all the pipetting and microplate handling steps needed for sample normalization tasks with the Normalization Wizard when used in con-

cert with Tecan's Freedom EVOware software package. Easily automate sample quantitation and normalization so that samples of varying concentration are diluted to a uniform concentration across the entire microplate. The Quantitation and Normalization Wizards guide the researcher through the normalization process with simple questions and automatically generate a pipetting script.



This Wizard is divided into two components:



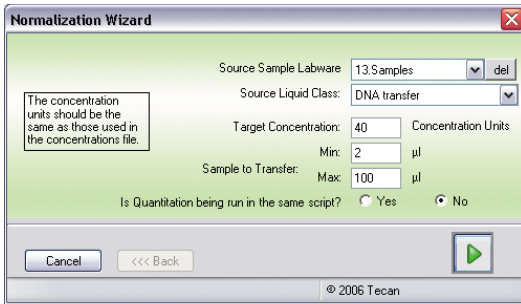
The **Quantitation Wizard** works in conjunction with Tecan microplate readers to quantify samples and the concentration of samples can be determined using Magellan software. The wizard also permits dilution of highly concentrated samples prior to the final measurement.



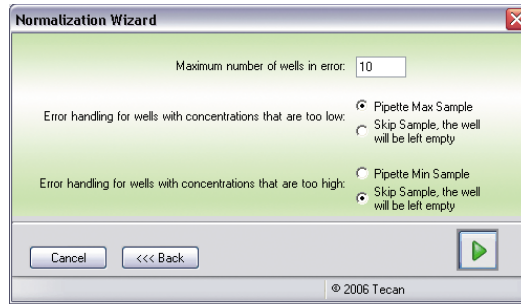
The **Normalization Wizard** allows all the samples to be normalized into a single microplate, with all samples at the same concentration. Data can be imported directly from the Quantitation Wizard during the same script run. Data can also be easily imported from previously generated data files, data files from third party devices such as realtime PCR instrumentation, and data files generated by LIMS.

Product highlights

- Seamless integration of Tecan detection modules and Magellan software package
- The ability to import data from other detection devices and real-time PCR instrumentation
- Simple, wizard-driven interface to combine concentration data, sample data, and pipetting data
- Automated calculations for quantitation and normalization
- Tracking of pipetting information throughout the entire process
- Suitable for normalization applications in the genomic, forensic, protein science, and drug discovery laboratory



Use concentration data from any device.



Built-in error reaction.

Full automation of DNA quantitation and sample normalization

The concentrations of samples are determined using the Quantitation Wizard in conjunction with Magellan software. Concentrations of known samples are used to dynamically generate a worklist for sample normalization with the Normalization Wizard. Data transformations are completed without the need for user interaction during script runtime, enabling a fully automated walkaway system.

Secure data handling and process control

Freedom EVOware is ideal for laboratories working in regulated environments, to provide a full audit trail, electronic records and electronic signatures. The wizard is able to generate printable output files detailing the process data including dispensed volumes and errors. The results are recorded in XML format for easy upload to LIMS systems.

Ideal for forensic applications

Suitable for processing casework samples in the forensic laboratory. Extracted casework or database samples can be processed by the quantitation wizard to prepare RT-PCR quantitation plates. The resulting output files can be imported to normalize sample concentrations prior to STR reaction preparation. Automated solutions result in improved productivity and decreased operator intervention, reducing risks for contamination and for transcription errors.

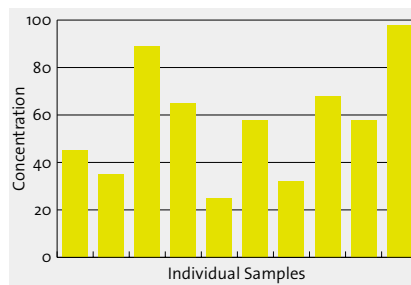
Evolve with Freedom EVOware

Freedom EVOware offers system integrators an extendable library of over 75 drivers for detection, separation, robotics and plate handling and storage devices. System customization is supported through a variety of open interfaces for device drivers, automation, worklisting and messaging, providing flexibility at all levels.

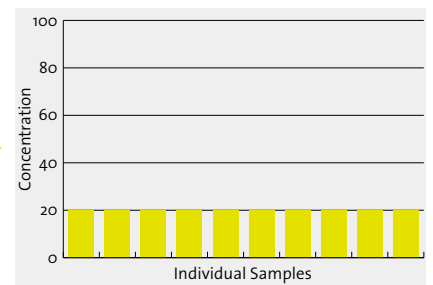
Extraction and Purification



Quantitation



Normalization



via Detection or real-time PCR

Tecan Group Ltd. makes every effort to include accurate and up-to-date information within this publication, however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided in this publication. Changes in this publication can be made at any time without notice. All mentioned trademarks are protected by law. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative. This brochure may contain reference to applications and products which are not available in all markets. Please check with your local sales representative.

© 2006, Tecan Trading AG, Switzerland, all rights reserved. Tecan® is in major countries a registered trademark of Tecan Group Ltd., Männedorf, Switzerland. Freedom EVO® and Freedom EVOware® are registered trademarks of Tecan Group Ltd., Männedorf, Switzerland.

Tecan Asia (Pte) Ltd., T +65 644 41 886, **Tecan Sales Austria GmbH**, T +43 62 46 89 33, **Tecan Benelux B.V.B.A.**, T +32 15 42 13 19, **Tecan Benelux B.V.B.A.**, T +31 18 34 48 17 4, **Tecan Group Ltd., Beijing Rep. Office**, T +86 10 586 95 936, **Tecan Deutschland GmbH**, T +49 79 51 94 170, **Tecan France S.A.S.**, T +33 4 72 76 04 80, **Tecan Italia S.r.l.**, T +39 02 215 21 28, **Tecan Sales International GmbH**, T +43 62 46 89 33, **Tecan Japan Co. Ltd.**, T +81 44 556 7311, **Tecan Nordic AB**, T +46 31 75 44 000, **Tecan Nordic AB, Rep. Office Denmark**, T +45 70 234 450, **Tecan Portugal**, T +351 21 000 82 16, **Tecan Sales Switzerland AG**, T +41 44 922 89 22, **Tecan Iberica Instr. S.L.**, T +34 93 490 01 74, **Tecan UK Ltd.**, T +44 11 89 300 300, **Tecan US Inc.**, T +1 919 361 5200

www.tecan.com

