

**Freedom EVO – HSM workstation includes:**

- 1 Compact Instrument:
  - Freedom EVO 100 on support table
  - 120 cm (w) x 110 cm (d) x 190 cm (h)
- 2 Promega HSM 2.0 (heater, shaker, magnet) integration includes worktable cut-out and support frame
- 3 Liquid system for fast bulk reagent dispensing, including reagent rack, six-way valve and 5 ml syringes
- 4 Aspiration system for fast removal of supernatant, including vacuum pump, eight-channel manifold, electronic valves for each channel, separate tubing and separate active wash station
- 5 Racks for 16 mm or 13 mm Vacutainers® and 2 ml Sarstedt, or similar sized, tubes
- 6 Carriers for troughs, microplates and disposable tips
- 7 Computer with flatscreen monitor, touchscreen interface and wireless hand-held barcode scanner
- 8 Software: Freedom EVOware® Standard & TouchTools Suite™



**Additional options:**

- Dust cover (shown in picture)
- Primary tube barcode reading (not shown)
- Base plate (shown in white on bottom)

**Items to be purchased from Promega:**

XAT1005 Installation package, consisting of:

- Promega HSM 2.0
- Integration and programming
- Biological sample demonstration
- Installation/training

A2751 ReliaPrep™ Large Volume HT gDNA Isolation System Reagents\*

\*For research use only

**Service and Support**

Tecan and Promega offer joint support for the Freedom EVO - HSM workstation through one point of contact. As the chemistry provider, Promega Corporation will provide the first point of contact for support of the Freedom EVO - HSM workstation with the ability to call on resources from Promega's automation team as well as Tecan instrument support for off-site and on-site trouble shooting and maintenance.

**Training**

We offer fast, convenient, effective training and certification for everyone who uses Tecan products. Freedom EVO - HSM workstation user training, including maintenance and routine use of the system, is provided as part of the installation.

Australia +61 3 9647 4100 Austria +43 62 46 89 33 Belgium +32 15 42 13 19 China +86 21 220 63 206 Denmark +45 70 23 44 50 France +33 4 72 76 04 80  
Germany +49 79 51 94 170 Italy +39 02 92 44 790 Japan +81 44 556 73 11 Netherlands +31 18 34 48 17 4 Singapore +65 644 41 886 Spain +34 93 490 01 74  
Sweden +46 8 750 39 40 Switzerland +41 44 922 89 22 UK +44 118 9300 300 USA +1 919 361 5200 Other countries +43 62 46 89 33

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. In general, the trademarks and designs referenced herein are trademarks, or registered trademarks, of Tecan Group Ltd., Männedorf, Switzerland. A complete list may be found at <http://www.tecan.com/trademarks>. Product names and company names that are not contained in the list but are noted herein may be the trademarks of their respective owners. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative.

Tecan is in major countries a registered trademark of Tecan Group Ltd., Männedorf, Switzerland.

© 2016 Tecan Trading AG, Switzerland, all rights reserved.

For research use only. Not for use in diagnostic procedures.

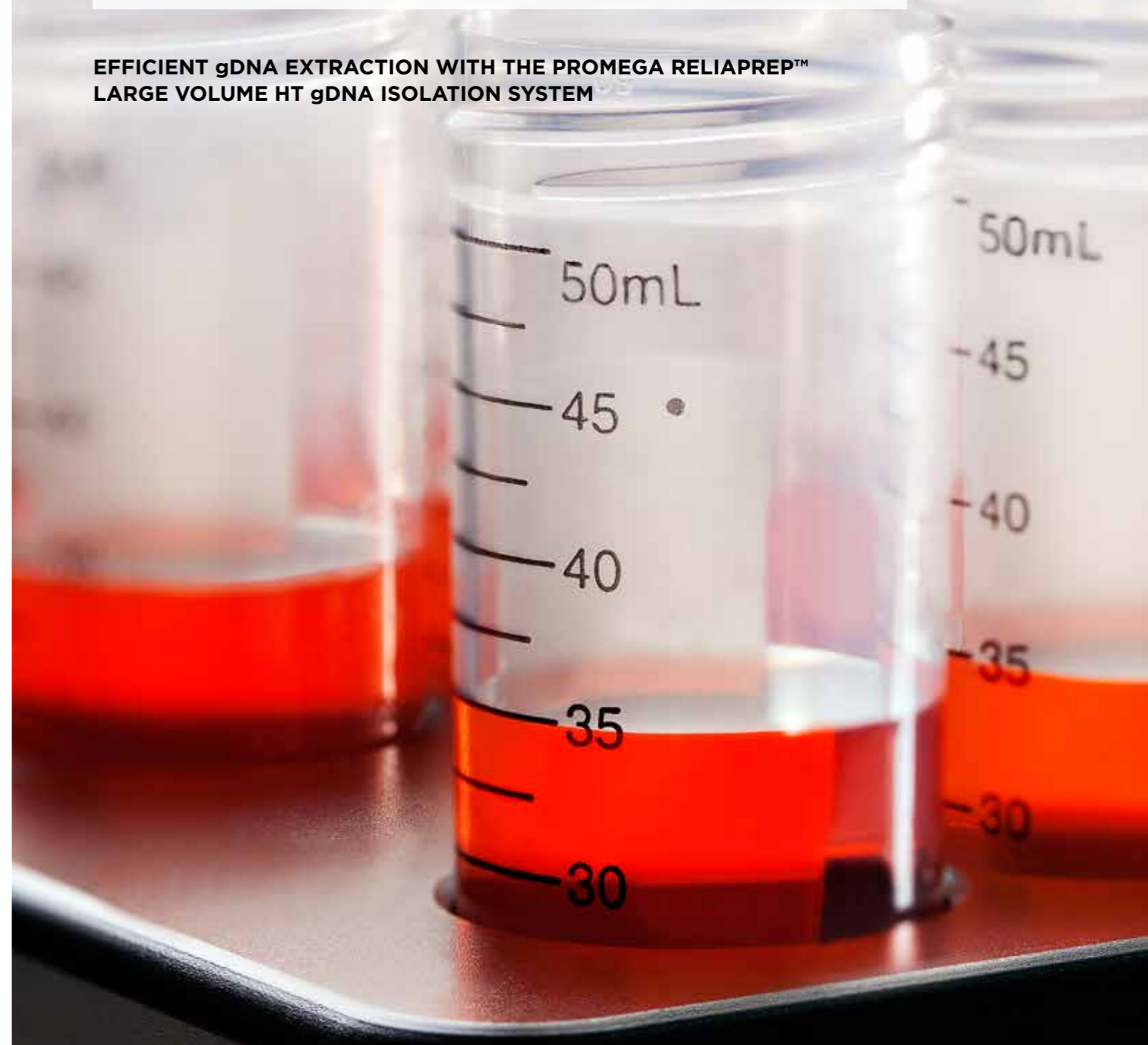
[www.tecan.com](http://www.tecan.com)



# Freedom EVO®

# HSM workstation.

**EFFICIENT gDNA EXTRACTION WITH THE PROMEGA RELIAPREP™  
LARGE VOLUME HT gDNA ISOLATION SYSTEM**



397618 V2.0 03-2016 30111830

## Efficient gDNA extraction with the Promega ReliaPrep™ Large Volume HT gDNA Isolation System – tailored to your DNA biobanking needs

Tecan and Promega are partnering to provide a uniquely efficient and reliable system for isolating genomic DNA from large volumes (1-10 ml) of blood, typically drawn into a standard 10 ml Vacutainer tube. The combination of the Freedom EVO, the ReliaPrep™ HSM and the ReliaPrep™ Large Volume HT gDNA Kit allows automated recovery of pure gDNA from up to 32 blood samples (each up to 10ml) in four hours.

This workstation resolves many of the challenges encountered when processing large volume samples. For instance, there is no risk of sample pellet loss during decanting or manual resuspension during precipitation-based methods. As the entire extraction process is performed in one 50 ml conical tube, there is minimal plastic waste and all hazardous waste is collected in one bottle, eliminating dangerous handling of open containers.

### PROMEGA RELIAPREP™ LARGE VOLUME HT gDNA ISOLATION SYSTEM

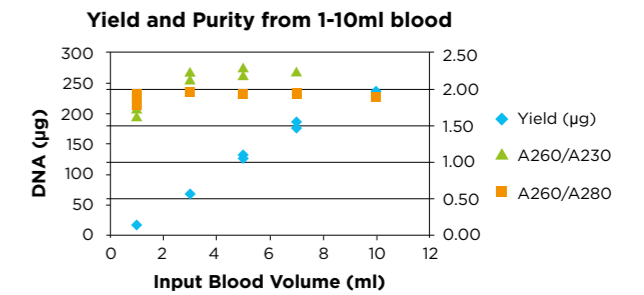


#### Intelligent one tube extraction

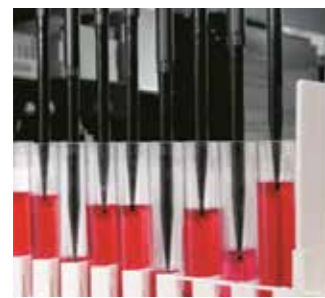
- The ReliaPrep™ HSM instrument, with its shaking, heating and magnetizing functions, enables gDNA extraction in one 50 ml conical tube
- Minimal plastic waste; one 50 ml tube, five tips per sample and one plate per run
- No need to move reagent filled tubes – reduces the risk of errors and cross-contamination

#### Highly pure gDNA for all your downstream needs:

- 1-10 ml input samples
- Purity specifications:  
 $A_{260/280} > 1.7$ ,  $A_{260/230}$  1.8-2.2
- High molecular weight DNA, ideally suited for applications requiring longer fragments, such as NGS
- Similar yields from fresh and compromised samples



### FREEDOM EVO – HSM WORKSTATION

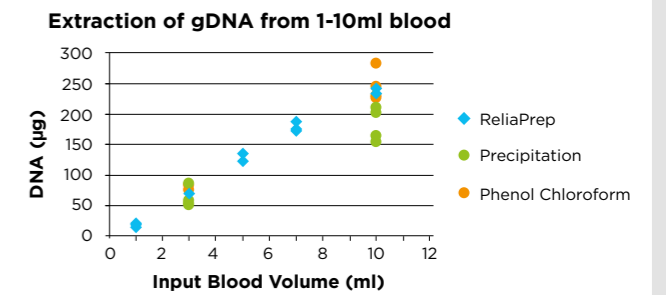


#### Optimized for each individual sample

- Optional sample transfer from primary tubes, or manual scanning of barcodes and transfer to 50 ml tubes
- Detection of sample volume and automatic reagent volume scaling for each sample conserves reagent and allows maximum yield and high purity from each sample, irrespective of the input sample volume
- DNA is eluted in water ready to use or store

#### Scalable chemistry:

- Any sample volume from 1 to 10 ml of blood can be processed in a single run using the same method
- No reagent is wasted as the exact volume, needed for each sample, is used

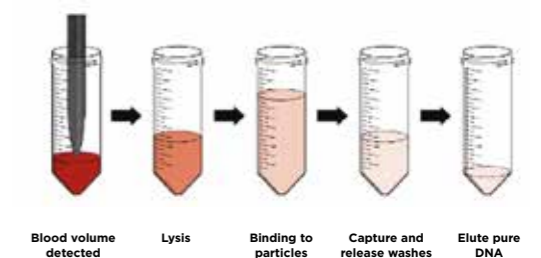


#### Designed for rapid large volume processing

- Bulk reagent dispensing, for fast reagent addition
- Eight-channel vacuum manifold for highly efficient large volume waste removal after each magnetic separation step
- Compact footprint 120 cm (w) x 110 cm (d) x 190 cm (h)

#### Optimized throughput:

- 4.5 - 6.5 hours for 32 x 10 ml whole blood samples, depending on workflow options.
- Shorter run times for smaller sample volumes
- Option to include primary sample barcode scanning and sample transfer from primary tubes to 50 ml tubes in the HSM instrument
- Minimal space requirements



### TOUCHTOOLS SUITE INTERFACE



#### Interactive touchscreen for intuitive set-up

- Pre-installed protocols
- Simple guided set-up
  - Selection of sample input and output format
  - Selection of sample number (1-32)
  - Selection of elution volumes (0.5 ml to 1.50 ml)
  - Sample and reagent barcode options

#### Simplified set-up:

- Pre-installed protocols allow you to get your first samples processed within days
- Easy guided set-up minimizes the training needs for your laboratory staff
- Customizable report generated to document each purification run

