

The Freedom EVO® 75
Nucleic Acid Extraction configuration
Magnetic bead separation with a flexible
input for 1 to 24 samples

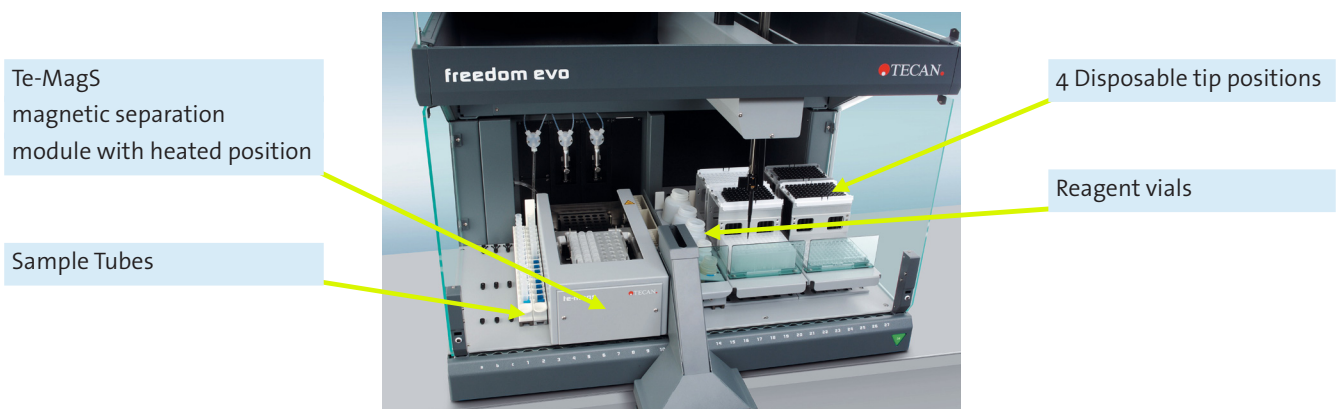
Reference 10003BF100598



The Freedom EVO® 75 workstation compact pipetting platform can be configured for extracting genomic DNA, viral DNA or RNA from 1 to 24 samples.

The instrument is equipped with a 2-channel liquid handling arm and the Te-MagS™ magnetic bead separation module; a barcode scanner can also be integrated with the system for sample identification. The automated steps of the extraction system include sample and reagent pipetting, heating, incubation, magnetic capture, washing and elution.

All modules are controlled by Freedom EVOware® software, enabling hands-free automation. In addition, the Freedom EVOware® Sample Tracking option offers the ability to track the pipetted volumes of each sample or plate throughout the entire process.



Suggested configuration

Part Number	Quantity	Product Description
30025019	1 PC	Instrument Freedom EVO 75 base with Freedom EVOware, Sample Tracking embedded and Sample Oriented EVOware
30025021	1 PC	Safety panel, right side
30025027	1 PC	Panel Black (required if no fast priming pump is configured)
30025026	1 PC	Liquid handling arm 2-channel
10619532	2 PC	Syringe 1000 µl
10612502	2 PC	Disposable tip option
30025029	1 PC	Lower disposable tip eject option
10650037	1 PC	Disposable tip waste and wash station
30028830	1 PC	Tube Carriers set: - one carrier for tubes 13 mm, 16 Pos. - one carrier for tubes 13mm, 14 Pos & 2 controls
10650010	2 PC	Carrier for microplates and disposable tip boxes
10290317	2 PC	Tip box for 1000 µl disposable tips
10290318	2 PC	Tip box for 200 µl disposable tips
Various	1 PC	Carrier for reagent bottles
30028829	1 PC	Output block for 1.5 ml tubes, 24 positions
30018042	1 PC	Instrument Te-MagS
30028419	1 PC	Te-MagS Plate carrier, 24 position
30028417	1 PC	Te-MagS Heating block, 24 position
10760701	1 PC	Te-MagS Magnet block, 48 position

Note: Waste and system liquid containers, PC and monitor may also be ordered from Tecan

Upgrade possibilities



Hand-held barcode scanner for positive identification of sample tubes



Membrane pump for fast priming of the liquid system



TouchTools Suite™ for instrument operation through a touchscreen

Scientific instrumentation; not for use in human, clinical or diagnostic procedure.

Tecan is in major countries a registered trademark and Freedom EVO, Freedom EVOware, TouchTools Suite, Instant Pipetting, Te-MagS are trademarks of Tecan Group Ltd., Männedorf, Switzerland.

© 2009, Tecan Trading AG, Switzerland, all rights reserved.

Please contact your local dealer for more information. www.tecan.com, p. 2/2