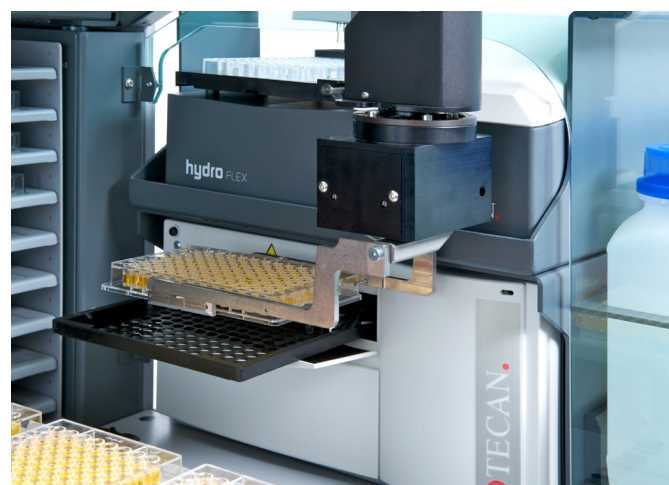


# The Freedom EVO® 75 ELISA configuration

## Fully automated ELISA assays

Reference 10003BF100596

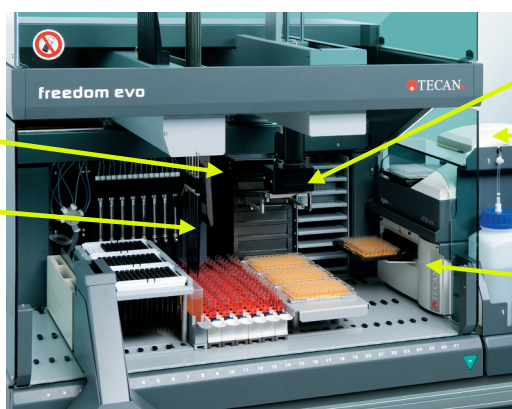


The Freedom EVO 75 workstation configured with the 8 Plus 1 Access™ Arm and Robotic Manipulator Arm economically enables the full automation of ELISA assays. Time consuming reagent addition steps are reduced by more than 80% compared to those using a single tip instrument. The system can process multiple assays per plate and up to three plates in a batch. The EVO 75 provides the flexibility of an open system; as a result you are not locked into a single reagent provider. The integrated HydroFlex™ plate washer and Sunrise™ plate reader help ensure the consistent and reliable microplate washing and reading steps with no user intervention required.

You will be up and running quickly with the easy user interface. All modules are controlled by the Freedom EVOware® Plus software, enabling hands-free automation from distribution of the samples to detection of the assay signal. The dynamic scheduling capability of the software ensures optimal usage of the robotic resources and exact incubation times. 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 to the acquisition of the detection data. A sample oriented wizard allows easy definition of tests, plate layouts and dilution schemes. Assays can be run in batch mode or with a job-list that specifies the test and dilution for each sample.

Heated incubator  
available with or without  
shaking

8 Plus 1 Access liquid handling arm



Robotic plate manipulator arm

HydroFlex plate washer

Sunrise plate reader

## Suggested configuration

Part Number	Quantity	Product Description
30025019	1 PC	Instrument Freedom EVO 75 base unit
30031708	1 PC	Liquid Handling arm 8 Plus 1 Access Arm
30025028	1 PC	Robotic Plate Manipulator arm
30017037	1 PC	Set of eccentric fingers for Robotic Manipulator
30025024	1 PC	Fast Prime pump
20734805	8 PC	Syringe 1000 µl
10650037	1 PC	Disposable tip waste, trough carrier 3-position and wash station
10613022	1 PC	Carrier for disposable tips, 3-position
10613020	1 PC	Carrier for troughs 100 ml, 3-position
various	–	Disposable tips
10613048	1 PC	Troughs 100 ml (108 pieces) used with 10613020
various	1 PC	Carrier for tubes, 16-position
10613006	2 PC	Carrier for microplates, 3-position (or use flat carrier 10612624 instead, to access lower slots of the Incubator)
various	1 PC	Monitored Incubator Option, holds 6 plates
10612674	1 PC	Hotel, holds 9 plates
30039408	1 PC	Kit reader, washer, worktable extension and scheduling software

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

## Other possibilities



2-channel liquid handling arm with fixed washable tips



Hand-held barcode scanner for positive identification sample tubes



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, 8 Plus 1 Access, Freedom EVOware, TouchTools Suite, Instant Pipetting, HydroFlex and Sunrise 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](http://www.tecan.com), p. 2/2