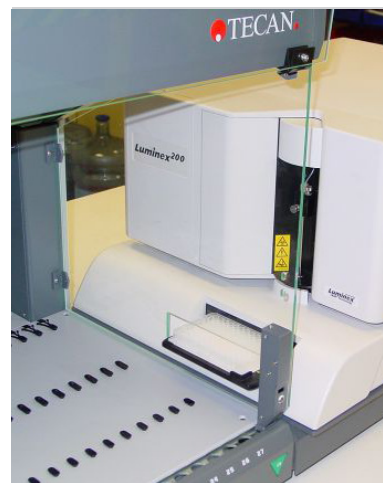


Freedom EVO[®] 75 and Luminex[®] 200 Automation of xMAP[®] biological assays

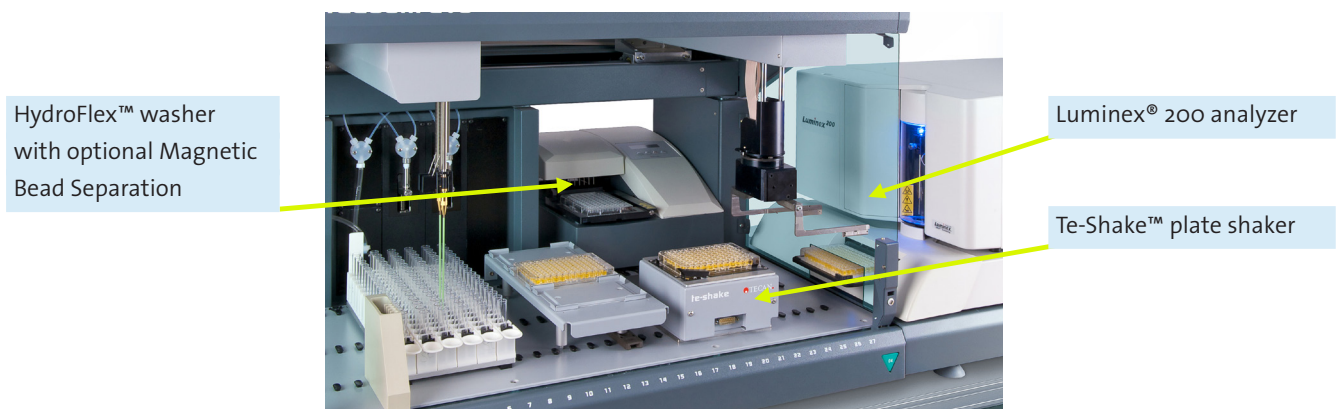
Reference 10003BF100593



The Luminex xMAP multiplexing technology allows researchers to perform numerous tests simultaneously for detection of gene sequences, protein interactions and other biological reactions, on the surface of fluorescent-coded microspheres. The xMAP systems using MagPlex[™] magnetic microsphere to remove unbound material from the assay are especially suited for assay automation.

The Freedom EVO 75 combined with the Luminex 200 analyzer and HydroFlex[™] washer offers a fully integrated platform to carry out the liquid transfers, plate movements, magnetic bead

separation and mixing. The Freedom EVOware[®] software controls all modules and provides dynamic assay scheduling, allowing scalability, from partial plate batch to multi-plate batch. Sample IDs are tracked and the system offers open LIMS integration. The HydroFlex washer ensures efficient plate wash and separation of the MagPlex magnetic beads. This compact and cost-effective system delivers bioassay results with increased reproducibility and significant hands-on-time saving compared to the manual sample preparation.



Suggested configuration

Part Number	Quantity	Product Description
Liquid Handling Hardware and Software		
30025869	1 PC	Instrument base Freedom EVO 75 with scheduling software
30025026	1 PC	Arm liquid handling 2-channel
30025028	1 PC	Robotic arm manipulator
30017037	1 PC	Robotic fingers
10643050	2 PC	Tip, fixed washable
10619532	2 PC	Syringe 1 ml
30025868	1 PC	Wash station for 2 tips
30025024	1 PC	Fast wash pump for tip wash
10613002	1PC	Carrier tubes 13 mm 16 positions (set of 6)
10613006	1 PC	Carrier plate 3-position
10613020	2 PC	Carrier troughs 100 ml 3-positions
Te-Shake Shaker		
10760722	1 PC	Adapter plate Te-Shake front
10760723	1 PC	Shaker orbital base
10760724	1 PC	Holder for microplate
HydroFlex Washer		
30022011	1 PC	Instrument HydroFlex
30022259	1 PC	Housing complete HydroFlex
30022078	1 PC	Arm pivoting
30022335	1 PC	Manifold 8-channel
30023424	1 PC	Option 4 input channels (Option 2 input channels 30023425 may be used instead)
30039258	1 PC	Carrier Smart 2 MBS (Carrier MBS, 30023217, extra strong magnet, may be used instead)
30026398	1 PC	HydroFlex Control Manual
Integration Hardware for HydroFlex and Luminex 200 Analyzer		
30046309	1 PC	Adapter Plate HydroFlex rear
30045976	1 PC	Worktable extension
30045860	1 PC	Adapter plate and right panel with cut-out for Luminex 200

Note 1: Waste and system liquid containers, PC and monitor may also be ordered from Tecan.

Note 2: 96-well flat bottom plates, such as Greiner 655096, should be used for magnetic bead separation on the HydroFlex.

Note 3: The Luminex 200 Analyzer is not supplied by Tecan.

Upgrade possibilities



Heating option for the Te-Shake.
Room Temperature to 80°C ±2°C



Hand-held barcode scanner for identification
of samples and reagents



TouchTools Suite™ including Instant
Pipetting™ and CNS

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

Luminex and xMAP are registered trademarks and MagPlex is a trademark of the Luminex Corporation, Austin, Texas. Tecan, Freedom EVO and Freedom EVOware are registered trademarks and TouchTools Suite, Te-Shake, Instant Pipetting and HydroFlex are trademarks of Tecan Group Ltd., Männedorf, Switzerland.

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