

HydroFlex™ – ELISA configuration

The reliable microplate washer for ELISA assays & cells



High wash efficiency

Building on a decade of experience in microplate washing, the HydroFlex configured for ELISA-washing & cell washing, sets a new standard in performance and reliability.

Its advanced design allows you to combine full control over critical wash parameters with advanced options for online process control. For a minimum residual volume of only 2µl per well, the HydroFlex provides a sophisticated crosswise aspiration function, using multiple aspiration points per well.

To achieve high wash-efficiency without compromising throughput, overflow washing replaces tedious, sequential aspiration- and dispense-steps with a single procedure.

High reliability

For uncompromised performance and high long-term stability, the HydroFlex uses advanced pump technology.

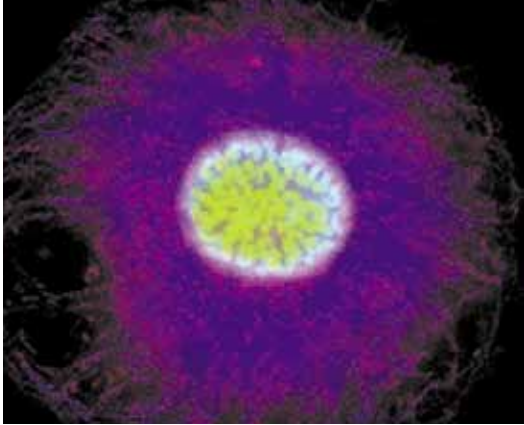
The HydroFlex provides a unique proof of quality via an MTTF-record (mean time to first failure record) of over 40,000 plates, typically corresponding to several years of operation.

Regulatory compliance

The HydroFlex configured for ELISA washing, has been designed to meet the European directive 98/79/EC for in-vitro diagnostic products. Additionally, the HydroControl™ software fulfils the requirements for electronic records and electronic signatures, and has been designed to meet the FDA regulation 21 CFR part 11.

Product highlights:

- Low residual volume of only 2µl/well
- Multiple aspiration points per well
- High wash efficiency via overflow procedure
- Drop-wise dispensing for gentle cell washing
- Automated buffer switching
- Proven reliability via MTTF-benchmark value
- Advanced process control options
- IVDD 98/79/EC compliant for ELISA washing
- Ready for robotic integration



Cell washing

For gentle washing of adherent cells for research use, the HydroFlex provides an extra-gentle drop-wise dispense mode.

Additionally, a set of tunable wash parameters including aspiration speed and aspiration height is available, to minimizing cell detachment even with weakly adherent cell lines.

Available options

Process control option:

This advanced function provides online process control of wash steps in overflow position for process safety and traceability.

Liquid Level control option:

This process-control function allows online monitoring of filling volumes in the entire bottle set of the HydroFlex, including bottles for wash buffer(s) and waste.

Specifications

Parameters	Characteristics
General	
Display unit	Liquid Crystal Display with two rows of sixteen digits
Keyboard	Four key membrane keyboard
Number of dispensing channels	1-, 2- or 4-channels
Manifold types	8- or 16-way for parallel processing of 8- or 16-wells, 8- or 16-way Process control manifold
Variable	
Volume of solution dispensed	50 – 3000 µl in 50 µl steps for washing 50 – 400 µl in 50 µl steps for dispensing
Dispensing accuracy	<= 2 % (300µl dispense volume of wash-buffer, dispense rate: 3, 8-way manifold, Greiner 96-well flat bottom plate).
Dispensing uniformity	CV: <= 4% over the plate (300µl dispense volume of wash-buffer, dispense rate: 3, 8-way manifold, Greiner 96-well flat bottom plate).
Residual volume	Crosswise aspiration: <= 2 µl per well for 96-well flat bottom Greiner plate (asp. rate 3, asp. time 4 sec., head speed 10mm/sec, 8-way manifold, wash buffer in all wells)
Interface	USB
Shaking	Linear shaking; high-, medium- and low-setting
Physical	
Outside dimensions for plate washing	Width: 27.5 cm, Depth: 36.6 cm, Height: 18.1 cm
Weight	6.6 kg

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.

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

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

HydroFlex and HydroControl are trademarks of Tecan Group Ltd. Männedorf, Switzerland.

Austria +43 62 46 89 33 **Belgium** +32 15 42 13 19 **China** +86 10 5869 5936 **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 48 174 **Portugal** +351 21 000 82 16 **Singapore** +65 644 41 886 **Spain** +34 93 490 01 74 **Sweden** +46 31 75 44 000
Switzerland +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **ROW** +41 44 922 8125

www.tecan.com/HydroFlex

395 957 V1.0, 04-2009

