

Reliable washing of SCHOTT NEXTERION® MTP-96 array plate with Tecan HydroFlex™

SCHOTT Technical Glass Solutions GmbH in Jena, Germany, is a subsidiary of SCHOTT AG, and develops the NEXTERION brand of coated glass products for microarray applications in academic and diagnostic laboratories. The Research and Development group at Microarray Solutions has chosen Tecan's HydroFlex microplate washer for development and quality control of its products.

"At Microarray Solutions, we develop and produce glass surfaces with coatings that have been designed for a range of requirements," explained Dr Rüdiger Dietrich, Director of R&D and Technical Support. "In addition to off-the-shelf products, we also offer volume production for customers who have developed their own working surfaces, developing customized surfaces to specific customer requirements. As well as our standard NEXTERION MPX-16 microscope slide and NEXTERION MTP-96 microtiter plate formats, we can produce coated glass in any format to suit different customer needs."

"With the increasing popularity of the microplate format for microarray applications, we have chosen Tecan's HydroFlex plate washer to run several test applications for automated quality control of NEXTERION MTP-96 format plates. The HydroFlex is used every day, and all the pipetting steps are automated, which



The team at SCHOTT Microarray Solutions

would otherwise be very time-consuming. We are also particularly impressed by its reproducibility; it is essential for us that the process is highly reproducible because, with six people working in the R&D and Technical Support team, several different operators, including newly trained ones, use the instrument."

"The HydroFlex is easy to use, and we have been able to optimize our protocols by adjusting various parameters like liquid volumes and rinsing times. The protocols we develop are shared with the customers, saving them time by helping

them to develop their own wash protocols quickly. We also use a Tecan HS 4800™ Pro hybridization station and a Tecan LS Reloaded™ scanner in our R&D laboratory, and have helped customers with the protocols we have developed with those instruments too."

For more information on Tecan's HydroFlex microplate washer, visit www.tecan.com/hydroflex

For more information on Tecan's HS 4800 Pro hybridization station and LS Reloaded scanner, visit www.tecan.com/hs4800 and www.tecan.com/lr

NEXTERION is a registered trademark of SCHOTT AG, Mainz, Germany.



The NEXTERION microarray slide



The Tecan HydroFlex