



Molecular Devices, now a part of MDS Analytical Technologies has discontinued the following products:

- Skan Washer 300/400 Microplate Washer
- SkanStacker
- EMBLA 96/384 Well and Cell Washers
- Cell Harvesters
- AquaMax DW4 Dispenser/Washer\*

Customers with active service contracts will continue to receive support until July 31, 2013. Post 2013, MDS Analytical Technologies may attempt to support discontinued products but do not guarantee such services due to obsolescence of parts and technologies.

We recommend our customers to migrate to the new generation [AquaMax 2000 and 4000 series Microplate Washers](#) with the following performance enhancements:

- Modular design for 96 and 384 capability
- 4-Fluid inlets for multiple assay or multi user labs
- One Touch cleaning and priming utilities enhancing robustness and reliability
- Integrated stacker for up to 50 microplates

Information concerning the replacement products can be found at:

<http://www.moleculardevices.com/pages/instruments/aquamax.html>

We want to thank you for your business and look forward to supporting your future research needs. Please contact [support@moldev.com](mailto:support@moldev.com) for further assistance.

\*Customers with active service contracts will continue to receive support until May 31, 2014.