



January 2022

Subject: Discontinuation of VersaMax Microplate Reader

Dear Valued Customer,

Thank you for being a valued customer of Molecular Devices. We are focused on continuously providing new innovative products with improved security, design, and functionality in support of your scientific research. After careful consideration, we have decided to cease manufacturing of our VersaMax™ Microplate Reader, which was launched over 20 years ago, to align with this strategy. This letter serves as an Advanced Notice of Discontinuation to help you plan accordingly.

Important dates for support and purchase:

- Last day to purchase a VersaMax Microplate Reader is December 30, 2022. After this date, it will no longer be available for purchase.
- Last day to purchase a service plan for VersaMax Microplate Reader is December 31, 2027. All service plan coverage will end by December 31, 2028. After this date, all service and support will be on a reasonable effort basis.

At this time, we are happy to help you upgrade to the SpectraMax® ABS or ABS Plus Microplate Readers which contain new features, extended dynamic range, and compatibility with SpectraTest® ABS2 Validation Plate. SpectraMax ABS and ABS Plus readers are compatible with SoftMax® Pro 7.1.2 Software – which is available in standard and GxP options – contain new features, offers improved security, and supports current Windows operating systems. SoftMax® Pro 7.1.2 GxP Software also addresses the updated guidance released by the FDA in 2016 regarding electronic signature and record-keeping requirements in [21 CFR Part 11](#).

Additional resources for SpectraMax ABS and ABS Plus readers are available [here](#).

For more information, please contact our team:

North America: 1.800.635.5577 or email support@moldev.com

Europe: 00800.665.32860 or email infoboxeu@moldev.com

For all other regions, [click here](#).

As always, we remain committed to providing you with the highest quality of products and services.

Please let us know if we can answer any questions.

Sincerely,

Laxmi Chokhani

Product Manager