



May 4, 2021

Subject: Notification of Global Discontinuation of SpectraMax 340PC384

Dear Valued Customer,

You are receiving this email because our records indicate that you have a SpectraMax® 340PC384 Microplate Reader. After careful consideration, we have decided to cease manufacturing this reader. This letter serves as an Advanced Notice of product discontinuation.

Last day to purchase a SpectraMax 340PC384 Microplate Reader: December 31, 2021

Service and Support: We will continue to offer repair and maintenance services for the next 5 years.

Last day to purchase a service plan: December 31, 2025.

All service plan coverage will end by December 31, 2026, after which our ability to perform repairs and provide support will be on a “reasonable effort” basis.

The discontinuation of the SpectraMax 340PC384 reader is in keeping with our practice of continuous improvement and providing our users with new products that have improved design and functionality. [SpectraMax ABS and ABS Plus](#) Absorbance Microplate Readers replace the SpectraMax 340PC384 reader.

If you have any questions, please contact Technical Support:
North America: 1.800.635.5577 or email support@moldev.com.
EU: 00800.665.32860 or email techsupport.eu@moldev.com.
For all other regions, [click here](#).

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

Thanks!

Laxmi Chokhani

Sustainment Product Manager