

## Discontinuation Notice

### Service plans for SpectraMax i3 and i3-R Multi-Mode Microplate Readers

Dear Valued Customer,

You are receiving this email because our records indicate that you have a SpectraMax i3 or i3-R Multi-Mode Microplate Reader. After careful consideration, we have decided to cease service plans coverage for our SpectraMax® i3 and i3-R Multi-Mode Microplate Readers. This letter serves as an Advanced Notice of Service Plans Discontinuation.

#### **Last day to purchase a service plan: December 31, 2021**

All service plans coverage will end on December 31, 2022, after which our ability to perform repairs and provide support will be on a “reasonable effort” basis.

Please note: Discontinuation of these service plans is specific to the SpectraMax i3 readers only. It does not in any way affect service for the SpectraMax i3x readers.

The discontinuance of the SpectraMax i3 readers is in keeping with our practice of continuous improvement and providing our users with new products that have improved design and functionality. As such, our **SpectraMax® i3x Multi-Mode Microplate Reader** replaces the SpectraMax i3 reader.

If you have any questions, please contact Technical Support:

- North America: 1.800.635.5577 or email [support@moldev.com](mailto:support@moldev.com).
- EU: 00800.665.32860 or email [techsupport.eu@moldev.com](mailto: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.

Thank you,

**Laxmi Chokhani**

Sustainment Product Manager



[Privacy Policy](#) | [Terms & Conditions](#) | [Trademarks & Logos](#)

For Research Use Only. Not for use in diagnostic procedures. The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners.

©2021 Molecular Devices, LLC 3860 N First Street, San Jose, CA 95134 USA

[View email in browser](#), [manage your subscriptions](#), or [unsubscribe](#) from future newsletters.

Please be advised that this email may contain confidential information. If you are not the intended recipient, please notify us by email by replying to the sender and delete this message. The sender disclaims that the content of this email constitutes an offer to enter into, or the acceptance of, any agreement; provided that the foregoing does not invalidate the binding effect of any digital or other electronic reproduction of a manual signature that is included in any attachment.