



Dec 9, 2020

Subject: Discontinuation of SpectraTest ABS1 Validation Plate

Dear Valued Customer,

We are continually innovating new products and streamlining our existing product portfolio. Our older products often rely on outdated and unavailable third party parts, posing a challenge to providing exceptional support to our customers. After careful consideration, we have decided to cease the manufacturing of SpectraTest® ABS1 Validation Plates. This letter serves as an Advanced Notice of Product Discontinuation.

Last day to purchase a SpectraTest ABS1 Validation Plate : December 29, 2023

Continued Service and Support: We will continue to recertify, repair, and support the SpectraTest ABS1 Validation Plates until **December 29, 2028**

Obsolescence: The ability to perform recertifications, repairs, and provide technical support will be on a “best effort” basis after **December 29, 2028**

SpectraTest ABS2 Validation Plate (P/N 0200-6191) is our next generation, automated, absorbance validation tool, replacing the ABS1 plate. It is compatible with SpectraMax® iD3, iD5, i3x, i3, M2, M2e, M3, M4, M5, M5e, Plus 384, ABS, ABS Plus and FlexStation® 3 microplate readers. For additional information on SpectraTest ABS2 Validation Plate, visit our website at <https://www.moleculardevices.com/products/microplate-readers/gxp-compliance-solutions/spectratest-validation-plates>. If you have any questions, please contact Technical Support at 1800-635-5577 or email support@moldev.com

Please contact your Molecular Devices sales representative for quotes on SpectraTest ABS1 or ABS2 Validation Plates.

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

Thanks!

Laxmi Chokhani

Sustainment Product Manager