



Jan 14, 2022

Subject: Update regarding 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 Our older products often rely on outdated and unavailable third party parts, posing a challenge to providing exceptional support to our customers.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. Advanced Notice of this Product Discontinuation was sent to you in December 2020. Below is an important update in support of the transition to SpectraTest ABS2 Validation Plate.

Key dates to remember:

Dates	Actions
6/1/2022	All services, including PM, PMOQ, and IQOQ will be performed using SpectraTest ABS2 Validation Plate.
6/1/2022 to 6/1/2023	We recommend users evaluate data comparisons between ABS1 and ABS2 validation plates. To enable this, we will offer temporary complimentary service of validating instruments with ABS1 validation plates during PM, IQOQ, and PMOQ.
6/2/2023 onwards	Customers can still request services with ABS1 plate for an additional charge.
12/29/2023	Last day to purchase SpectraTest ABS1 validation plates.
Until 12/29/2028	Continued service and support for ABS1 plates including recertification and repair. After this date, all service and support will be on a reasonable effort basis.

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 [here](#).
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 of products and services.

Thanks!

Laxmi Chokhani

Sustainment Product Manager