



Molecular Devices recommends annual recertification of your validation plates in order to ensure that they meet specifications and to give you confidence in the performance of your instrument. Only Molecular Devices has the necessary knowledge and equipment to recertify SpectraTest® validation plates.

Why recertify your validation plate with Molecular Devices?

- Molecular Devices is **ISO 17025** Certified by American Association of Laboratory Accreditation (**A2LA**), which requires rigorous outside auditing on a regular basis so that you can be assured of the accuracy of your certificate with respect to **NIST** traceable standards.
- Gold Standard and Reference Standard Molecular Devices plate readers are used for recertification so that your certificate values are traceable back to factory standards.
- If repair is needed, only new factory-approved parts that meet our internal specifications for these plates are used.
- Rapid turnaround time. Typically 5 business days or less. Expedited service is available as needed.

Benefits

- Ensure data accuracy of your plate reader
- Detect instrument problems quickly and easily when they happen so that product quality is not compromised
- Minimize downtime by quickly being able to distinguish between instrument failure and assay failure
- Maintain conformance to your validation protocols

You can trust Molecular Devices to provide the highest quality service at the best value.



Call or e-mail us to schedule your plate recertification.

USA and Canada+1.800.635.5577 ext. 1815

Email: servicelogistics@moldev.com

United Kingdom+44.118.944.8000

We developed your validation plates. We are your best choice for recertification service.