



Why recertify your validation plate?

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

- Molecular Devices is ISO 17025 accredited* 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 and NMI traceable standards.
- Gold Standard and Reference Standard Molecular Devices microplate readers are used for recertification so that your certificate values are traceable.
- If validation plate repair is needed, only factory-approved parts that meet our internal specifications for these plates are used.
- Rapid turnaround time of five business days or less. Expedited service is available as needed.

*[View our ISO 17025 certificate of accreditation.](#)

Benefits

- Ensure data accuracy of your reader
- Monitor reader performance on demand to ensure data quality is not compromised between PM/OQ maintenance services
- Minimize downtime by quickly being able to distinguish between reader and assay issues
- Maintain conformance to your validation protocols

We developed your validation plate—you can trust us to provide you with high quality service at the best value.

Email us to schedule your plate recertification.

North America: Customer.Relations@moldev.com

China: Support.China@moldev.com

Europe: Service.EU@moldev.com

Rest of World: [Please contact your local sales representative.](#)

Contact Us

Phone: +1.800.635.5577

Web: www.moleculardevices.com

Email: info@moldev.com

Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada +1.800.635.5577

United Kingdom +44.118.944.8000

Europe* 00800.665.32860

China (Beijing) +86.10.6410.8669

China (Shanghai) +86.21.3372.1088

Hong Kong +852.3971.3530

Japan +81.3.6362.9109

South Korea +82.2.3471.9531

*Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden and Switzerland