



Streamline your compliance journey in GMP/GLP labs



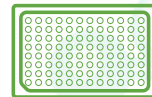
READERS & WASHERS

Best-in-class microplate readers and washers for all your assay needs.



SOFTMAX PRO GxP SOFTWARE

Our most secure software to achieve full FDA 21 CFR Part 11 compliance.



VALIDATION PLATES

Validation plates test the performance of your microplate reader using traceable materials for reliable results.



IQ/OQ & PM/OQ SERVICES

Comprehensive OEM operational qualification and repair coverage for microplate readers and washers.



INSTALLATION SERVICES

Software installation services verify and document that required components are installed to operational specifications.



VALIDATION SERVICE

On-site software validation service is conducted by our certified Field Service Engineer.



RECERTIFICATION SERVICES

Validation plates are cleaned, calibrated, and recertified according to ISO 17025, and returned to you with a new certificate of calibration.

Contact Us

Phone: +1.800.635.5577 Web: www.moleculardevices.com Email: info@moldev.com

Check our website for a current listing of worldwide contact information.