

Time & materials planned maintenance services



Lower labor rates. Simpler travel costs. The same trusted Molecular Devices service.

Maximize instrument performance on a limited budget. Powered by PhD level technical support, factory approved repair parts, and OEM certified engineers, Molecular Devices keeps your data precise and your scientific discoveries on track.

Onsite or Depot Preventive Maintenance Services

Protect performance.
Maximize uptime.

- Proactive service that protects your instrument's uptime
- Ensure data integrity through calibrated, optimized optical performance
- Extend instrument life with routine care of critical components
- Improve assay reliability with consistent, reproducible results
- Lower total cost of ownership by avoiding costly repairs and disruptions

**Request a quote for a
PM for your instrument ►**

Onsite Qualification Services

Maximize performance.
Ensure compliance.

- Centralize records with secure, digital documentation for all service activities
- Stay audit-ready with complete validation, maintenance, and repair history
- Meet regulatory standards (FDA 21 CFR and EudraLex Volume 4 guidelines) with certified service engineer execution
- Simplify compliance with standardized IQ/OQ/PM documentation
- Ensure proper installation (IQ) with verified, documented setup
- Validate performance (OQ + PM) with comprehensive testing, calibration, and inspection

**Request a quote for an
IQ/OQ/PMOQ/As Found/post
relocation OQ for your instrument ►**

Validation Plate Recertification Services

Annual recertification ensures plates meet specifications and maintains confidence in instrument performance.

- ISO 17025 accredited (A2LA) for NIST/NMI-traceable accuracy
- Recertified using gold & reference standard readers for full traceability
- Factory-approved repairs with validated parts, when needed
- Fast turnaround ≤5 business days (expedited service available for an additional charge)

**Request a quote for recertification
of your validation plate ►**

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	Taiwan/Hong Kong	+886.2.2656.7585
United Kingdom	+44.118.944.8000	Japan	+81.3.6362.9109
Europe*	00800.665.32860	South Korea	+82.2.3471.9531
China	+86.4008203586	India	+1.800.266.5338

*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Norway, Portugal, Spain, Sweden, Switzerland and United Kingdom

