

Time & materials repairs



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

We don't just repair instruments—we protect your science. 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.

What does this mean for you:



Lower, market aligned labor rates make urgent service needs easier to justify and approve internally.



Faster decisions and reduced downtime when service is needed.



Simpler, predictable travel fees with flat, zone based pricing—not influenced by traffic or distance, no surprises.



Clearer, more transparent quotes that are easy for you to review and approve.



The same trusted service quality, delivered by OEM certified engineers using proven diagnostics and factory approved parts.



Modernized pricing aligned to customer expectations, without reducing scope, expertise, or service standards.

Request a quote for your instrument repair.

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

