

## Preventive Maintenance (PM) Compliance Assurance Digital Installation Qualification (IQ) Operational Qualification (OQ) Services

**PM Service** confirms that your Molecular Devices instrument meets operational specifications.

### Service Description

A factory trained and qualified field service engineer (FSE) performs multi-point inspection procedure on your instrument to verify the instrument meets operational specifications. The rigor of the multi-point inspection procedure allows for proactive identification of replacement parts (as needed) to ensure your instrument is maintained at peak operational performance. At the completion of the PM service, the FSE presents a PM checklist which documents the completion of the multi-point inspection procedure.

**IQOQ and PMOQ Services** certify that your Molecular Devices instrument meets operational specifications.

### Service Description

- The IQ section verifies and documents that all necessary components required for operation are received and properly installed in accordance with Molecular Devices specifications.
- The OQ sections tests the mechanical, electrical, and optical components of each instrument to verify proper operating functionality in accordance with Molecular Devices specifications.
- The IQOQ service is performed for new instrument installations or for instrument relocations.
- The PMOQ service is performed on instruments that have been installed for six months or longer and includes preventive maintenance.

A factory trained and qualified FSE performs operational verification procedure on your instrument, and runs proprietary verification scripts to certify your instrument meets original manufacturer specifications. At the completion of the qualification services, the FSE presents a document which includes specifications testing results and certifies that your Molecular Devices instrument meets original manufacturer operational specifications.

	<b>PM Service</b>	<b>PMOQ/IQOQ Service</b>
<b>Documentation</b>	PM Checklist	Comprehensive digital or printed report, including qualification results and validation certificate to ensure you are audit ready
<b>Execution</b>	Factory trained and qualified FSE	Factory trained and qualified FSE
<b>Fulfillment</b>	FSE confirms that customer's instrument meets operational specifications	FSE confirms that customer's instrument meets operational specifications
<b>Molecular Devices Certification</b>	No	Yes
<b>PM Procedure</b>	Yes	Yes + qualification results and validation certificate
<b>Instrument Operational Procedure</b>	Yes	Yes + qualification results and validation certificate
<b>Proprietary Verification Scripts</b>	Yes	Yes + Performance certification values
<b>Operational Parameter Recording</b>	No	Yes + Performance certification values
<b>FSE Training Record</b>	No	Yes