



Software validation service

Our on-site SoftMax® Pro GxP Software validation service supports FDA 21 CFR Part 11 guidelines and is conducted by our certified Field Service Engineer (FSE). Each step in the process will be carefully planned and executed.

Key benefits

- Receive a pre-service consultation led by our validation experts
- Get comprehensive software validation testing by our certified FSE
- Receive a complete set of audit-ready documents
- Minimize downtime—leave the validation to us and focus on your research
- Save over a week of your valuable time*

* Estimated time savings

On-site visit

- After completion of all tests, the FSE will provide you with a complete data package:
 - Completed & signed SoftMax Pro GxP validation document (hard & scanned copy)
 - Results cover page with electronic signature
 - Results template with screenshots of each step
 - IQ report for SoftMax Pro GxP Software
 - Copy of the FSE Training Certificate

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	China (Beijing)	+86.10.6410.8669	Japan	+81.120.993.656
United Kingdom	+44.118.944.8000	China (Shanghai)	+86.21.3372.1088	South Korea	+82.2.3471.9531
Europe*	00800.665.32860	Hong Kong	+852.3971.3530		

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