

Automated Software Qualification Tool

For Functional Testing of SoftMax Pro GxP Software Suite

Our newest service innovation for software functional testing features an automated testing tool that streamlines the manual entries needed to configure software settings and step-by-step execution.

Key features

- Reduces validation time
- Automates the manual steps at all test cases
- Generates a system audit trail report which documents step-by-step validation activity with accurate time/date stamps
- Generates a Validation Report with select screenshots
- Follows a non-scripted testing approach to Computer Software Assurance (CSA)

Reduced validation time

Functional specification testing (negative & positive tests) of the user administration of permissions, roles, and projects between the GxP Admin Portal software and their expected behavior in SoftMax Pro GxP is a manual and repetitive process. Physical transcription of test results and generation of screenshot evidence can take up to four days to complete the entire Validation Test Suite. With the automation of the manual steps to execute these functional test cases, the full end-to-end session using the automated testing tool is reduced to less than two hours.

Intended purpose

A qualified Field Service Engineer (FSE) will execute and monitor the progress of the automated software qualification (ASQ) tool and will provide the auto-generated system audit trail report and the Validation Report complete with screenshots and pass/fail criteria.

Service offerings

Part Number: **SMPGXP-OQ-VALSVC-OS**

- Operational Qualification [OQ]: Automated functional testing of SoftMax Pro GxP Software Suite: GxP Admin Portal and SoftMax Pro GxP

Part Number: **SMPGXP-PQ-VALSVC-OS**

- Performance Qualification [PQ]: Manual execution of performance tests of data analysis features: curve fitting graphing interpolations, and connectivity to the GxP Admin server.
- Available option for qualifying additional clients connected to a validated single server (multi PC) environment.

Software validation services are available for SoftMax Pro GxP Software version 7.2 and above. Based on your needs, services can be purchased separately or bundled together.

Contact Us

Phone: [+1.800.635.5577](tel:+18006355577)
 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	+91.73.8661.1198

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

