

## Sharing data between users has never been more secure

- Are **bottlenecks** an issue when multiple users in your lab try to access secure data from a single computer?
- Is **data integrity** a concern but you have limited knowledge or resources to achieve FDA 21 CFR Part 11 or EU Annex 11 compliance?
- Are you spending **valuable time** trying to set up a secure environment instead of doing science?

### If you answered yes, you are not alone.

Many labs grapple with these questions daily. In a quest to become compliant, labs depend on their IT team for setting up a secure networked environment.

Our SoftMax® Pro GxP Software Suite can scale up from a single computer to a global networked environment. Our Professional Services team will partner with your IT team to plan and implement a secure networked SoftMax Pro GxP Software application suite according to your requirements.



### Benefits of a secure networked environment

- Secure, reliable, centralized data repository for easy data sharing—protects against data loss and minimizes failure compared to desktop hardware
- Reduced user management tasks—can control projects, user permissions, and accounts from a single location
- Consolidated, comprehensive system audit trails help prevent audit complications
- Valuable time freed up to focus on research instead of server installation or maintenance
- IT department not required to be a part of the data access process after installation

Installation type	Multi computer
Number of networked computers	Multiple
Number of microplate readers	Multiple
Configuration	SoftMax Pro GxP Admin and database can be installed on different computers
Database	Microsoft SQL Express (SQL Server Standard or Enterprise edition not included)
File sharing	Yes. Files can be shared with any computer connected to the database. The system audit trail remains intact.
User ID and password control	Active Directory Windows Credentials or GxP Admin software
Installer	Script based installation by our Professional Services team
Computer requirements	<p><b>Client computer</b></p> <ul style="list-style-type: none"> <li>• Windows 10 Pro/Enterprise 64-bit</li> <li>• Intel Core i5 or higher with 8 cores</li> <li>• 16 GB RAM (8 GB minimum)</li> <li>• 50 GB of hard drive space</li> <li>• Windows PowerShell 5.1</li> </ul> <p><b>Server computer</b></p> <ul style="list-style-type: none"> <li>• Windows Server 2016 or 2019</li> <li>• Intel Core i5 or higher with 8 cores</li> <li>• 16 GB RAM</li> <li>• 50 GB of hard drive space</li> <li>• Microsoft SQL Express or SQL Server 2016 or 2017</li> <li>• Windows PowerShell 5.1</li> <li>• Configured DNS active for all connected computers</li> <li>• Firewalls and routers configured to allow data transfer using ports 8210, 8211, 1433, and 1434</li> </ul>
Software part number	SMP712GXP-SVR-LIC
Software remote installation service part number	SMPGXP-INSTALLSVR-OS
Customized installation service	SMPGXP-INSTALLADVSVR-OS
Additional client computer installation service	SMPGXP-INSTALLCLT-OS
Additional user license part number	SMP GXP SVR ADD
Software maintenance	<ul style="list-style-type: none"> <li>• SMP CARE-GXP-ESSENTIAL</li> <li>• SMP CARE-GXP-PREFERRED</li> </ul>

The Multi Computer Installation Service includes a consultation call with our Professional Services team to ensure the software and installation meet your IT infrastructure needs. Please begin to discuss the following topics with your IT support team.

- Contact information for IT group if they will be involved with the installation
- Server type: bare metal or virtualized (Hyper-V or VMWare)
- Internet access for remote installation
- Server information
  - SQL database may be colocated on the same server as the GxP Admin software or on separate servers
- SQL database information
- Single site, multisite, or global database access
- Active directory integration
- Installation service package type
  - Standard or advanced

## Contact Us

Phone: +1.800.635.5577  
 Web: [www.moleculardevices.com](http://www.moleculardevices.com)  
 Email: [info@moldev.com](mailto: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, Ireland, Italy, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom