



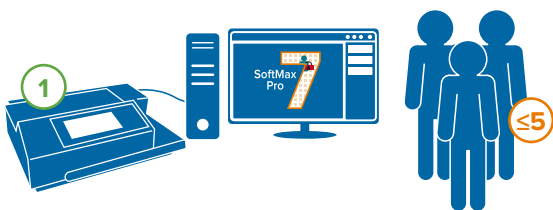
PURCHASE GUIDE

SoftMax Pro GxP Software Suite

The SoftMax® Pro GxP Software Suite can scale from a single computer to a networked environment of multiple computers. Consult with your Molecular Devices representative and your network administrators to plan and scale your implementation according to your requirements.

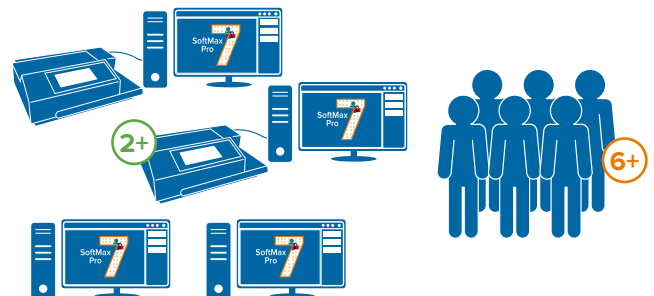
Single Computer Setup

- Best for small labs with all users sharing **one** microplate reader
- All software components are installed on a single computer, which is used by all users for data acquisition, analysis, and user administration



Multi Computer Setup

- Best for labs with all users sharing **multiple** microplate readers through a server on a corporate network
- Data acquisition, analysis, and user administration may occur on multiple computers, which all connect to a database hosted on a central server



SoftMax Pro GxP Software Suite components

- **GxP Admin Software**—Database and security backend application that you install on a dedicated, centrally located server with a static IP address. This server must be accessible to all computers that run the GxP Admin Portal Software and the SoftMax Pro GxP Software.
- **GxP Admin Portal Software**—The user interface that you use to interact with the GxP Admin Software. Install the GxP Admin Portal Software on a computer or computers that have access to the dedicated server on which you install the GxP Admin Software.
- **SoftMax Pro GxP Software**—Controls the microplate reader for data acquisition and statistical data analysis. These computers must have access to the dedicated server on which you install the GxP Admin Software.

Installation type	Single computer	Multi computer
Concurrent users	One	Multiple
Number of readers	One	Multiple
Configuration	SoftMax Pro GxP, Admin and database installed on the same computer	SoftMax Pro GxP, Admin and database installed on different computers
Database	Microsoft SQL Express with 10 GB limit	Microsoft SQL (not included)
File sharing	No	Yes. Files can be shared with any computer connected to the database. The system audit trail remains intact.
Database backup	Tool supplied	Not included
User ID and password control	GxP Admin software	Active Directory Windows Credentials or GxP Admin software
Installer	Simple wizard	Script based installation
Computer requirements	<ul style="list-style-type: none"> Windows 10/11 Enterprise or Windows 10/11 Pro (all 64-bit) Intel Core i5 or higher with 8 cores (minimum 4 cores to install) 16 GB RAM (8 GB minimum) 50 GB of available space, or more Windows PowerShell 5.1 Microsoft SQL Express will be installed by the wizard (10 GB storage limit) 	<p>Client computer</p> <ul style="list-style-type: none"> Windows 10/11 Enterprise or Windows 10/11/Enterprise Windows Server 2016 or 2019 (all 64-bit) Intel Core i5 or higher with 8 cores 16 GB RAM (8 GB minimum) 50 GB of available space, or more Windows PowerShell 5.1 <p>Server computer</p> <ul style="list-style-type: none"> Windows Server 2016 or 2019 (all 64-bit) Intel Core i5 or higher with 8 cores (minimum 4 cores to install) 16 GB RAM 50 GB of hard drive space Microsoft SQL Express or SQL Server 2016 or 2019 Windows PowerShell 5.1 must be 64-bit. Must allow script execution with the execution policy set to at least Remote Signed Configured DNS must be active for all connected computers
GxP Software	Latest version of SoftMax Pro 7 GxP Software Suite includes: 1 software activation for each user license, GxP Admin Software, self-guided documentation for software IQ/OQ validation package, user license certificate, compliance certificate, Windows 11 compatible	
Software installation service	Standard server installation (Applies to initial purchase only)	
Customized installation service	Custom server installation for multi computer setup only	
Additional user license	Additional user license and software activation purchases	
Software maintenance	Annual maintenance and technical support plan for SoftMax Pro 7.1.1 GxP Software or higher (Required for first year of new license purchase)	
Software validation service	Onsite software validation for SoftMax Pro GxP 7.1.1 or higher	

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
- Internet access required for remote installation by Teams, Webex, or Zoom
- 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
 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