



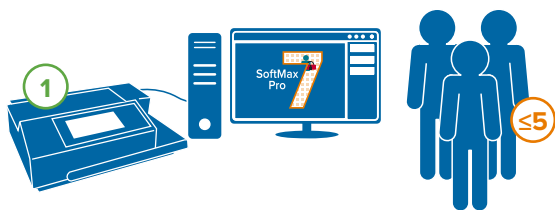
PURCHASE GUIDE

SoftMax Pro 7.1.1 GxP Compliance Software Suite

The SoftMax® Pro 7.1.1 GxP Compliance Software Suite can scale from a single computer to a multi-computer networked environment. Consult with your Molecular Devices representative and your network administrators to plan and scale your implementation according to your requirements.

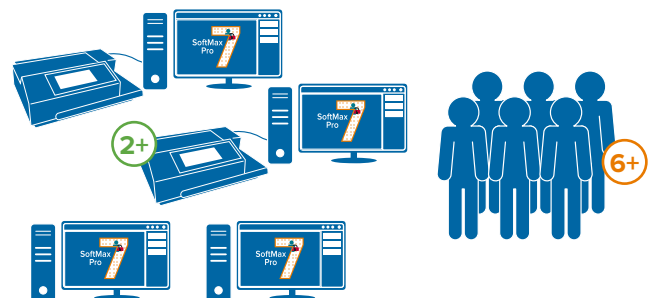
Single Computer Setup

- Best for small labs with **five** or fewer users on **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 **six** or more users on **multiple** microplate readers
- Data acquisition, analysis, and user administration may occur on multiple computers, which all connect to a database hosted on a central server



SoftMax Pro 7.1.1 GxP Compliance 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 Pro/Enterprise 64-bit Intel Core i5 or higher with 8 cores 16 GB RAM (8 GB minimum) 50 GB of hard drive space Windows PowerShell 5.1 Microsoft SQL Express 	<p>Client computer</p> <ul style="list-style-type: none"> Windows 10 Pro/Enterprise 64-bit Intel Core i5 or higher with 8 cores 16 GB RAM (8 GB minimum) 50 GB of hard drive space Windows PowerShell 5.1 <p>Server computer</p> <ul style="list-style-type: none"> Windows Server 2016 or 2019 Intel Core i5 or higher with 8 cores 16 GB RAM 50 GB of hard drive space Microsoft SQL Express or SQL Server 2016 or 2017 Windows PowerShell 5.1 Configured DNS active for all connected computers Firewalls and routers configured to allow data transfer using ports 8210, 8211, 1433, and 1434
Software part number	SMP7X GXP SINGLE COMP	SMP7X GXP SERVER
Software remote installation service part number	SMPGXP-INSTALL1COMP-OS	SMPGXP-INSTALLSVR-OS
Customized installation service	None	SMPGXP-INSTALLADSVR-OS
Additional user license part number	SMP GXP ADD	SMP GXP SVR ADD

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 by 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
 - Self-install requires knowledge of PowerShell script-based installation

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 (Osaka)	+81.6.7174.8331
United Kingdom	+44.118.944.8000	China (Shanghai)	+86.21.3372.1088	Japan (Tokyo)	+81.3.6362.5260
Europe*	00800.665.32860	Hong Kong	+852.3971.3530	South Korea	+82.2.3471.9531

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