



IT BENEFITS

SoftMax Pro 7.1 GxP Software

Address your compliance needs

SoftMax® Pro 7.1 GxP Software is the latest, most secure software to achieve full FDA 21 CFR Part 11 compliance with streamlined workflows to ensure data integrity. Every step is optimized to simplify analysis and reporting to support our microplate readers. Our expert team will partner with you to establish environments to conduct GMP/GLP (GxP) compliant work.

Solutions for your IT compliance issues



Microsoft SQL Express database

- Enterprise-level document sharing with complete control over file access permissions
- Windows folder permission system is no longer needed
- Can be easily upgraded later to a regular SQL database
- Reduced support from IT departments



Windows Active Directory

- Direct connection to your company's Windows Active Directory
- More control over password requirements and reduced IT support needed to reset passwords
- Includes a classic built-in user admin system for setting password strength, auto lock, and password aging
- Can simultaneously use both Windows Active Directory system and the classic built-in GxP Admin system for external auditors or users
- For current SoftMax Pro GxP customers, our GxP Admin EDB Converter tool quickly and easily imports .edb user account files from GxP Admin Software version 2.x



Projects feature

- Can quickly create various project teams and increase usability
- Project team members only have access to projects they are assigned to, eliminating the need for IT to create Project folders with custom permissions

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.