



QC 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 QC compliance issues



System audit trail

- The system audit trail tracks and records all actions for easy identification
- Electronic signatures and date/time stamps provide paperless documentation
- Filtering information simplifies quality control and handling during audits making data export to Excel unnecessary



Microsoft SQL Express database

- Centralized Microsoft SQL database provides complete control over file administration and file access permissions
- Can be upgraded to a regular SQL database



Document workflow and status system

- The paperless, document status system maintains data integrity with control over document workflows
- Project teams can track documents while they are moving through development, review, release, and usage in a controlled environment



Improved auto-save functionality

- Documents can be defined at the design stage
- Documents are automatically saved before and after a read to prevent the loss of data

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.