

# SoftMax® Pro 7.6 Software

More control. Clearer records.  
The right fit for your laboratory.



## Introducing GLP Mode

Built for GLP. Not borrowed from GMP.

From early research through regulated workflows, reliable results depend on more than data acquisition alone. Laboratories need the right combination of analysis flexibility, traceability, user control, and documentation—without taking on more compliance infrastructure than their work requires.

SoftMax® Pro 7.6 Software brings plate reader data acquisition, analysis, and workflow controls together in one connected software portfolio. Choose the level of control that fits your laboratory today, with a clearer path forward as your quality requirements evolve.

### New: GLP Mode delivers right-sized compliance

Designed specifically for development and GLP laboratories, GLP Mode provides essential controls for electronic records, traceability, and audit readiness—without imposing the process rigidity of a fully regulated GMP environment.

### Right-sized for the work you are doing now

Get the controls that matter without taking on the administrative burden and process rigidity of a full GMP solution before your laboratory needs it.

**Control access  
with confidence**

**Document review  
in context**

**Trace activity  
more easily**

**Keep data and  
evidence together**

More control now. A clearer path forward.

## Compliance control — without the overkill

A streamlined, GLP-specific administration experience gives your team essential compliance controls without requiring a GxP-oriented administrative framework.



**Audit readiness — without the reconstruction.** Keep audit trails, electronic signatures, review records, and plate reader data together—making it easier to understand what happened, who was involved, and how results were generated.



**Compliance control without overbuilding.** Manage users, permissions, access, and review practices without introducing unnecessary GMP complexity into development-stage work.



**Less risk across users and teams.** Establish more consistent practices across projects, users, and sites while reducing the risk created by disconnected tools and individual workflows.

## One software portfolio. The right level of control.

SoftMax Pro software supports laboratories as their quality requirements evolve:

### Standard Edition

Flexible acquisition and analysis for research laboratories.

**Explore and analyze**

### NEW: GLP Mode

Right-sized controls for development and GLP laboratories.

**Establish control and traceability**

### GxP Mode

Formal, controlled workflows for regulated GMP, QC, and manufacturing environments.

**Support validated, regulated processes**



*Find the right SoftMax Pro solution for your laboratory.  
Explore SoftMax Pro 7.6 Software or talk to a compliance software specialist.*

### 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            | +1.800.266.5338  |

\*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Norway, Portugal, Spain, Sweden, Switzerland and United Kingdom

