



Hands-Free Software Maintenance

Always have the most up-to-date version of SoftMax® Pro Software with the latest features and benefits. Ensure your lab has the best tools available..

Benefits of the Software Maintenance Plan

- Free feature packed upgrades (6.x to 6.y) for SoftMax Pro Software automatically available to customers
- Significant discount on SoftMax Pro Software major upgrades (6.x to 7.x)
- Additional license seats

New in SoftMax Pro Software 6.2

- Plate Setup Helper to guide your protocol creation
- Print to PDF to enable quick data and report sharing
- Improved well scan feature to display data from each well
- Enhanced curve fit options now with 17 curve fit choices

New in SoftMax Pro Software 6.1

- Redesigned look and feel to have you up and running in minutes
- Clear results display with new 3-D graph feature
- Simplified formula writing with Syntax Helper
- Customized reporting with image import options and live display of mini-graph and tables
- Expanded data analysis tools with plate cloning and additional curve fits
- Data analysis confidence with patent pending curve fit calculation
- Peace of mind with new built-in data recovery feature