



MetaMorph NX

Makes Imaging Easier

MetaMorph[®] NX Microscopy Automation and Image Analysis Software

Hundreds of Hardware Drivers

- Supports the broadest range of microscopy components to enable your unique application
- Continuous driver development targets state-of-the-art peripheral hardware and camera drivers to keep you ahead of the curve
- In-house driver and component testing gives you confidence that your system will work every time

Multi-Threaded Application

- Acquisition and analysis can now be accomplished simultaneously
- Asynchronous control of peripheral hardware shortens time-lapse intervals to improve temporal resolution
- Device streaming captures images at the maximum hardware rate for high-speed device synchronization during acquisition