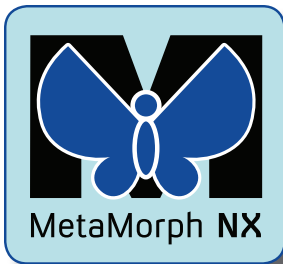


Export

Rotate

Multiple Channels

Filmstrip



4D Viewer for MetaMorph[®] NX Microscopy Automation and Image Analysis Software

Experience It

- View your data in 3-D without leaving the MetaMorph NX Software workspace—the 4D Viewer is seamlessly integrated into the user interface.
- Quickly switch between datasets and time points in 3-D using the built-in filmstrip.

Explore It

- Readily observe the spatial relationships between objects in 3-D by displaying multiple fluorescent channels at the same time.
- Investigate your data from all angles by spinning, rotating, and zooming into your images.

- Peer into the various structures and layers of your images by interactively slicing through your dataset.

Share It

- Effortlessly create a presentation with your data by exporting your 3-D display directly to Microsoft PowerPoint[®].
- Quickly prepare data for publication by saving your images with publication-specific DPI settings.
- Easily share your data with others by creating a movie of your dynamic 3-D renderings.