

## MetaMorph® 7.0 4D Viewer Video Card Requirements

### OVERVIEW

This document describes the video card requirements for the optional 4D Viewer dropin in MetaMorph 7.0 and above. The 4D Viewer enables you to view multi dimensional data sets, stacks, and sequential images. With it you can simultaneously view multiple z sections, wavelengths, time points, and positions from a single viewing window.

To take full advantage of the 4D Viewer's advanced 3D rendering and stack playback capabilities, the computer running MetaMorph must use a modern video card with sufficient memory. The 4D Viewer was optimized to work with following video card:

ATI Radeon® 9600 Pro AGP card with 256 Megabytes (MB) DDR SDRAM

Other video cards with similar or higher specifications should provide equal performance as the Radeon 9600.

### NOTES:

- The 4D Viewer will not work with a video card that has less than 128 MB of memory.
- Molecular Devices Corporation (MDC) recommends installing the latest version of your video card's driver from the manufacturer's web site before using the 4D viewer.
- If you are using an ATI Radeon 9600 series video card, you must have ATI Catalyst™ Control Center driver software version 5.10 or greater installed.
- Your video card must support the OpenGL API. Most modern video cards include support for OpenGL® with the card's drivers.
- If you are experiencing performance issues when viewing a 3D volume, try lowering the detail level from the 4D Viewer window.

**ARTICLE #**  
T20052

**PRODUCTS**  
MetaMorph® 7.0 and  
above with optional 4D  
viewer dropin

**CREATED**  
23-March-2006

**LAST UPDATED**  
27-March-2006