



# MetaMatters

## ***You wanted it, you got it!***

### *MetaMorph® Software*

*...now compatible with*

*Microsoft Windows® 7!*

- Click **Here** to download a FREE copy of Windows 7 compatible MetaMorph Software
- **OR** Contact your local MetaMorph distributor to obtain the NEW MetaMorph Software version 7.7:

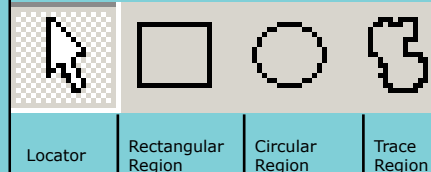
[www.moleculardevices.com/pages/distributors.html](http://www.moleculardevices.com/pages/distributors.html)

- Download the Windows XP compatible MetaMorph Software version 7.7 here:  
[www.moleculardevices.com/pages/MM-new/meta\\_smp.html](http://www.moleculardevices.com/pages/MM-new/meta_smp.html)
- MetaMorph Software version 7.7 is the **ONLY** version compatible with the new Windows 7
- 32-bit and 64-bit Windows 7 supported
  - If interested in the 64-bit version, ask your local sales representative whether your hardware and additional modules are supported in that environment.

#### **Inside this issue:**

MetaMorph Software News	1, 4
MetaMorph Software Incentives	2
FOCUS: Graphing Astrocytic Calcium Intensity with MetaMorph Software	2, 3
Upcoming Events	4

#### MetaTool Tips: Region Shortcuts



## Amazing Software Incentives!

---

Submit a 500 - 800 word MetaMorph® Software method and receive a FREE software upgrade or application module.

---

Submit a one paragraph MetaMorph® Software tip and receive a FREE 12 month software maintenance agreement.

The method description and tip paragraph will be published in *MetaMatters*.

Do you have an amazing image that is a "Work of Art"?

If yes, share your artistry with microscopists world wide and receive  
FREE MetaMorph Software!

In exchange for giving Molecular Devices the right to use your fully credited image in advertisements and brochures, you will receive a FREE MetaMorph Basic offline software package!

The free MetaMorph®  
Software Basics  
Training Course is  
April 20 & 21, 2010!

Email:  
[Mary.David@moldev.com](mailto:Mary.David@moldev.com)  
for more information on  
incentive programs.  
For training courses click here:  
[MetaMorph Software Training](#)

The MetaMorph®  
Software Advanced  
Training Course is  
April 22 & 23, 2010!

### FOCUS: Graphing Astrocytic Calcium Intensity with MetaMorph® Software

---

*Sarah E. Crowe and Graham Ellis-Davies  
Department of Pharmacology and Physiology, Dr. Graham Ellis-Davies Laboratory  
Drexel University College of Medicine*

Changes in intracellular calcium are one of the most widely used mechanisms of signaling in cells and are involved in processes such as cell movement, neurotransmitter release, cell division, and more. Many efforts have been made to monitor these calcium fluctuations both *in vivo* and *in vitro* with fluorescent calcium indicators. We have used a non-ratiometric indicator, X-Rhod-1, and 2-photon microscopy to study intrinsic *in vivo* astrocytic calcium signaling in a mouse model of Alzheimer's disease. Dysfunctional astrocytic Ca<sup>2+</sup> signaling has been found in several diseases or disorders associated with the brain, such as epilepsy, traumatic brain injury, and Alzheimer's disease. Due to the significance of Ca<sup>2+</sup> in astrocytic signaling, the observed dysfunctions are believed to play an important role in the pathology of such disorders.

We have used the "graph intensities" MetaMorph application to analyze the astrocytic calcium signaling observed in a mouse model of Alzheimer's disease. An advantage to using MetaMorph software is the ability to import T-Series sets as their native 12-bit format, thus maintaining all information collected. Our data are collected as .tiff files, which are easily handled by MetaMorph software, and opened as a tiff image sequence. Once opened in the software, we applied pseudocoloring to the series. (**con't pg.3**)

---

# FOCUS: Graphing Astrocytic Calcium Intensity with MetaMorph® Software

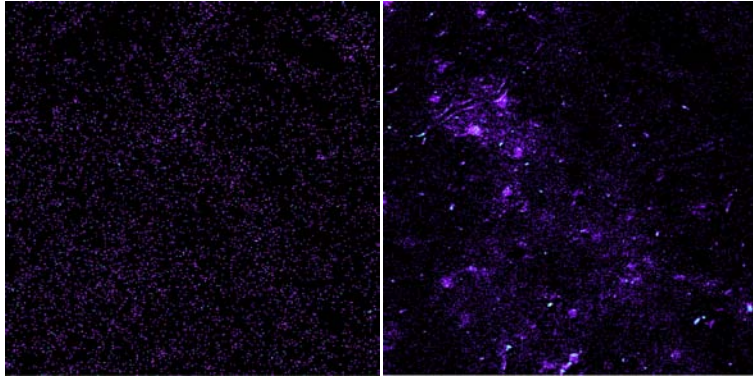


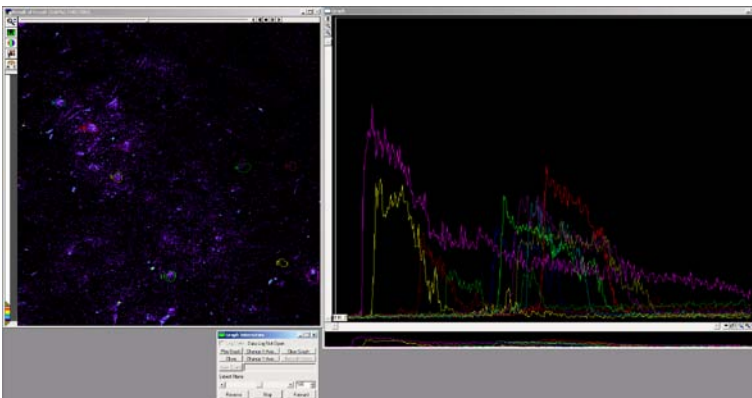
Figure 1- The  $\Delta F/F_0$  for a non-signaling frame (A), and a frame containing multiple astrocytic signals (B).

(continued from page 2)

This displays an increase in fluorescence, which indicates an increase in intracellular calcium, as warmer colors. An important step in analyzing any type of intensity over time data is the removal of background noise and the generation of the  $\Delta F/F_0$  signal. For our data background,  $F_0$ , we selected the first five frames of the T-Series and compiled an averaged stack image. Using the MetaMorph arithmetic tools, we then subtracted the averaged image from all frames of the T-Series ( $F-F_0$ ), producing  $\Delta F$  for each frame. Next, we divided all  $\Delta F$  frames by the averaged image to generate the  $\Delta F/F_0$  for each image frame (Fig 1). It is from the  $\Delta F/F_0$  T-Series that we analyze all calcium signal intensities.

The fluorescence intensity over time for each astrocyte must be individually examined, rather than the general change in intensity for the entire image frame. MetaMorph software allows for the selection of multiple astrocytes, or simply regions, within the same frame that can be analyzed simultaneously, which is significantly more efficient than single cell analysis. Individual astrocytes were manually selected with the "trace regions" tool. The region of the astrocyte chosen was the soma, with the trace following the edge of the soma as closely as possible (Fig 2A).

Figure 2, A & B



After the astrocytes of interest were selected, we opened the "graph intensities" module from the "apps" menu. This opened a dialogue box of options entitled "Configure Graph Intensities". For a T-Series such as ours, the options "stack", "plane number", and "average intensity" were selected. After clicking the "ok" button, a new box of parameters, named "Graph Intensities" appeared (Fig 2B). To start graphing the intensities of the selected astrocytes, we hit the "begin" button and the application simultaneously graphed each selected astrocyte into one window (Fig 2C). A right click of the mouse within the graph window will bring up a menu of options, including graph

settings. It is in here that the graph can be customized. Along with graph settings, the right click options menu also includes "show graph data", which will list the intensity values for each astrocyte for each image frame. These data can be logged into Excel for any further analysis.

The "graph intensities" application in MetaMorph software allows us to analyze calcium signals from multiple astrocytes with ease and efficiency. Simultaneous graphing of many cells provides information on dynamics of calcium in the astrocytic network. From this information, it is possible to examine propagation speed of intercellular signaling, distance of signal propagation, and basic kinetics of the intracellular signal. The importance of astrocytic calcium signaling is quickly emerging in a variety of different states of pathology and analysis of such may provide great insights into the mechanisms of diseases and disorders.

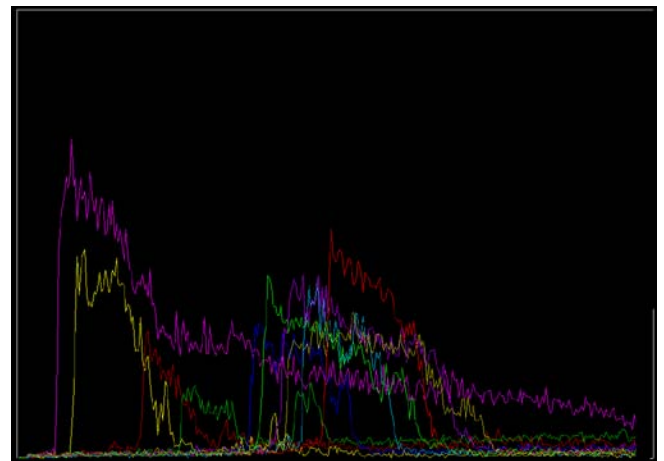


Figure 2C



402 Boot Road  
Downingtown, PA 19335

Phone: 800-635-5577  
Fax: 610-873-5499  
meta.admin@moldev.com  
support.dtn@moldev.com  
training.dtn@moldev.com

**MetaMorph® Software...  
the gold standard in  
research imaging**

**We're on the web!  
MetaMorph.com**

## → NEW 64-bit Camera Drivers

- Photometrics PVCAM
- Hamamatsu DCAM
- QImaging QCAM
- Andor

## → NEW 64-bit Hardware Drivers

- Generic serial port devices that do not require an SDK (such as Uniblitz, Lambda-10, etc)
- Nikon Eclipse Ti microscope
- Zeiss COM MTB microscopes (current models)
- Sutter USB controllers (Lambda 10-3, etc.)
- Prior USB controllers (ProScan II, ProScan III, etc.)
- TILL Polychrome USB

→ Note on supported hardware:  
Latest SDK from vendor must be downloaded and installed.

## Upcoming Training, Courses and Conferences

APRIL 20 – 21, 2010

Fundamentals of  
MetaMorph Software  
Downingtown, PA

APRIL 22 – 23, 2010

Advanced Topics of  
MetaMorph Software  
Downingtown, PA

SEPTEMBER 21 – 22, 2010

Fundamentals of  
MetaMorph Software  
Downingtown, PA

SEPTEMBER 23 – 24, 2010

Advanced Topics of  
MetaMorph Software  
Downingtown, PA

MAY, 2010

Analytical and Quantitative  
Light Microscopy  
Woods Hole, MA

JUNE 13 - 19, 2010

Quantitative Fluorescence  
Microscopy  
Bar Harbor, ME

OCTOBER 13 – 26, 2010

Immunocytochemistry, In Situ  
& Live Cell Imaging  
Cold Spring Harbor

NOVEMBER 13 – 17, 2010

Society for Neuroscience  
San Diego, CA

DECEMBER 11 – 15, 2010

American Society for Cell  
Biology  
Philadelphia, PA



MetaMorph® products @ QFM,  
2008