

Microsoft Windows® File Privileges Required for MetaMorph® Administrators and Users

OVERVIEW

This document describes which Microsoft Windows directories must be accessible to MetaMorph administrators and users. A MetaMorph administrator's tasks, such as creating users and groups, and defining hardware profiles, typically require greater file privileges than the tasks of a normal MetaMorph user.

NOTES:

- This document assumes that MetaMorph has been installed under **C:\MM** (the default location).
- You can install MetaMorph on any drive or location on your computer. For example, there is nothing that prevents you from installing the software to the path: **D:\Program Files\MM**.

READ/WRITE FILE PRIVILEGES REQUIRED FOR A METAMORPH SOFTWARE ADMINISTRATOR

MetaMorph software administrators need read/write access for everything under the **C:\MM** directory.

READ-ONLY FILE PRIVILEGES REQUIRED FOR METAMORPH USERS

MetaMorph users require a minimum of read-only access to the following directories:

- **C:\MM**
- **C:\MM\app\mmproc**
- **C:\MM\app\mmproc\Dropins**
- **C:\MM\Help**
- **C:\MM\Groups**

If the computer is an acquisition computer, users will also need read-only access to **C:\MM\Hardware** and all subdirectories.

READ/WRITE FILE PRIVILEGES REQUIRED FOR METAMORPH USERS

If MetaMorph is configured to run in single-user mode, the user will require read/write access to the following directory:

C:\MM\Groups\MetaMorph

If MetaMorph is configured to run in multi-user mode, the user will require read/write access to the following directory:

C:\MM\Groups\MetaMorph\Users\USERNAME

NOTE: If you are storing images, journals, or other MetaMorph files in directories outside the MetaMorph root directory (**C:\MM** by default), users must have Read/Write access to those directories in addition to the directories listed in this document.

ARTICLE #
T20053

PRODUCTS
MetaMorph® 6.0 and
above

CREATED
16-March-2006

LAST UPDATED
22-March-2006