

Using Internet Explorer 8 to access MetaMorph, MetaView, MetaFluor FTP Installation Files

Overview

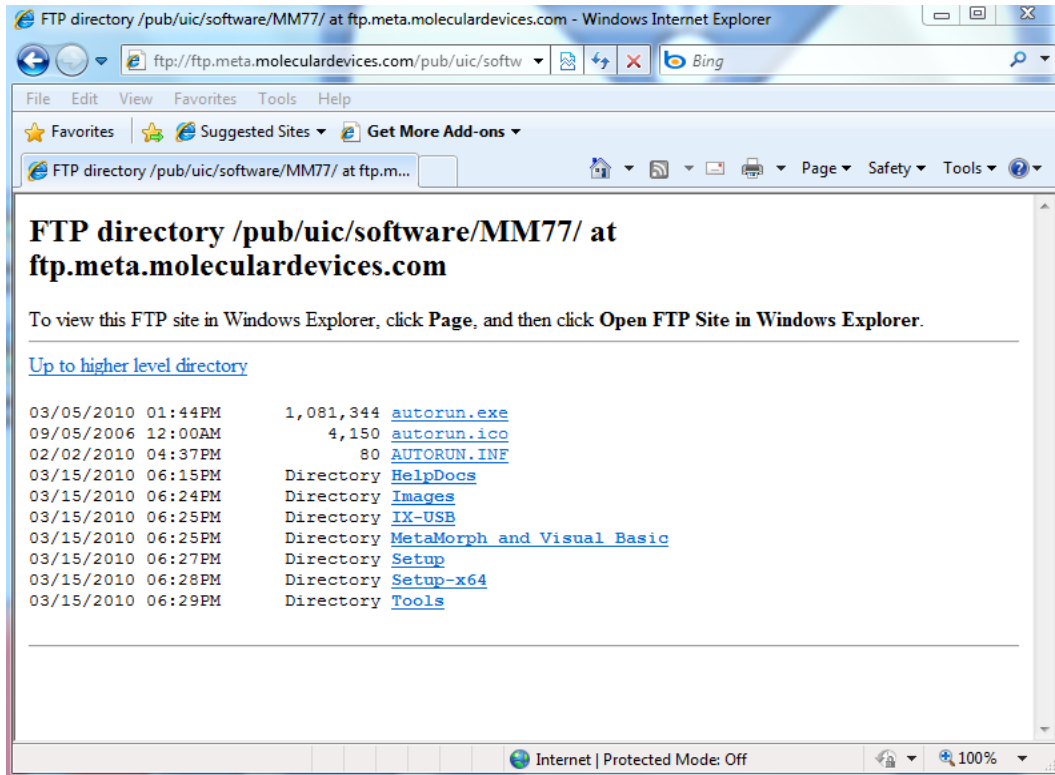
- Due to the changes that Microsoft has implemented in the newer versions of Internet Explorer the following extra steps are required to download the installation files for your software.
- Each step will be outlined below and all the current FTP CD links have been provided

Links

1. <ftp://ftp.meta.moleculardevices.com/pub/uic/software/MM45/>
2. <ftp://ftp.meta.moleculardevices.com/pub/uic/software/MM46/>
3. <ftp://ftp.meta.moleculardevices.com/pub/uic/software/MM63/>
4. <ftp://ftp.meta.moleculardevices.com/pub/uic/software/MM70/>
5. <ftp://ftp.meta.moleculardevices.com/pub/uic/software/MM71/>
6. <ftp://ftp.meta.moleculardevices.com/pub/uic/software/MM75/>
7. <ftp://ftp.meta.moleculardevices.com/pub/uic/software/MM76/>
8. <ftp://ftp.meta.moleculardevices.com/pub/uic/software/MM77/>

Instructions

1. Open Internet Explorer
2. Navigate to one of the FTP CD links listed above corresponding to the files you require for your installation. For this document we will be using the MM77 files.



As you can see the files and folders listed are not very accessible as links.

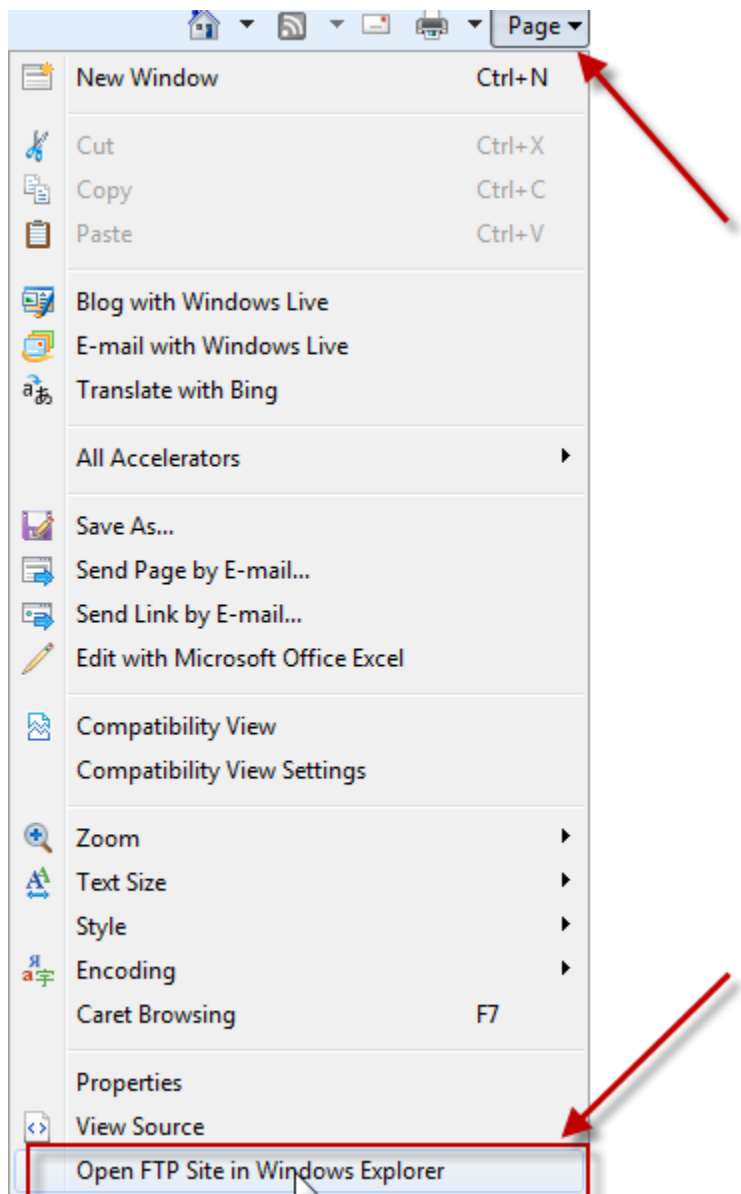
3. Navigate to the Page option in the upper left hand corner of your window and select (Open FTP Site in Windows Explorer)

ARTICLE #
T20141

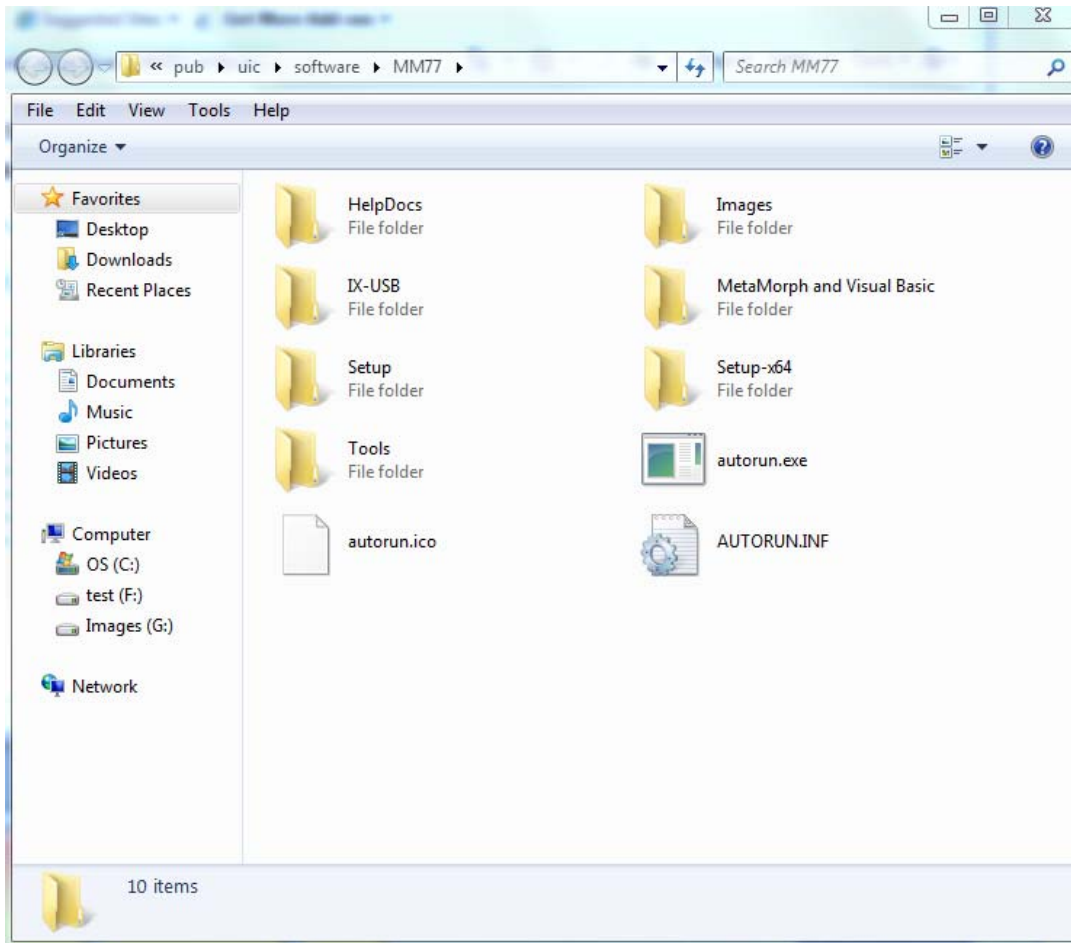
PRODUCTS
MetaMorph®
MetaFluor®
MetaVue®

CREATED
06-July-2010

LAST UPDATED
07-July-2010



4. Your newly opened window should look like this.



5. You may now copy all the files and folders listed by highlighting them and pasting them to a folder on your local computer. (Your desktop or a folder on your desktop will do just fine)
6. Once you have completed the file copy you can execute or double click the (autorun.exe) to start your MM installation.

If there are any further questions, please contact Technical support at 800-635-5577 x1820 or support.dtn@moldev.com.