

What is your policy on installing MetaMorph® / MetaFluor® / MetaVue™ software programs on multiple computers?

Question:

We have multiple users who wish to use MetaMorph®, MetaFluor® or MetaVue™ software programs on their own computers or laptops. What do we need to do? What are your policies?

Answer:

The current policy is that you can install the software on as many computers as you want. However, the number of licenses you have purchased limits the number of open copies. We offer two types of Meta Imaging Series® security keys. If using the standard security key the software may only be run on the computers that have the security key (dongle) installed. You can move the security key to any computer (e.g., between your desktop and your laptop) but beware that you do not damage or lose it. We also offer a network security key. The network security key can be purchased with a specific number of software licenses, installed on one computer, and copies of the software can be opened and run on as many computers as the number of software licenses you have purchased. You can also purchase additional security keys (standard or network) to run the software on multiple computers. Please contact your distributor to receive a quote for this. You can find your distributor on our website at: <http://www.moleculardevices.com/pages/distributors.html>.

ARTICLE #T20107

PRODUCTS

MetaMorph®
MetaFluor®
MetaVue®

CREATED

17-March-2009

LAST UPDATED

17-March-2009