

iCycler iQ™ Real-Time PCR Detection System Software Version 3.1 Installation Instructions

This installation disk is for the following operating systems: Windows 98, ME, XP, NT and 2000.

1. Insert the iCycler installation CD in a CD-ROM drive.
2. If the installation program does not begin automatically, select **Run...** from the Start menu and then type **X:\Setup**, where X is the drive letter of the CD-ROM drive. For example, if the CD-ROM is the E drive, type **E:\Setup**.
3. You must type a number in the serial number field in order for the installation program to continue. It does not matter what number you type in this field.
4. Follow all screen prompts to finalize the installation.
5. If you have a previous version of iCycler software, the report templates will be copied to a new folder called Old Report Templates. New report templates will be installed in a folder called Report Templates. We recommend that you do not use the old report templates with version 3.1 of the iCycler software.

NOTES for Windows NT, 2000 and XP users

Administrator-level privileges are required to install the software. If this is a first-time installation of the software, the administrator must also start the application initially.

Certain configurations of Windows NT, 2000 and XP initialize new folders by assigning Read and Execute permission for the members of the 'Users' group. If you have this type of operating system, and this is a first-time installation, the administrator must change the protection on the Program Files/Bio-Rad folder or the Program Files/Bio-Rad/iCycler folder so that you can save protocol, plate setup and data files. If, after changing the protection on either of these folders, it is still not possible to write to the folders beneath Program Files/Bio-Rad/iCycler, check the Properties of each individual folder to be sure that under the Securities tab, the box is checked that allows inheritable permissions to propagate to that folder.

Allow inheritable permissions from parent to propagate to this object