

Antibody Wave 6 | Installation Guide

Introduction

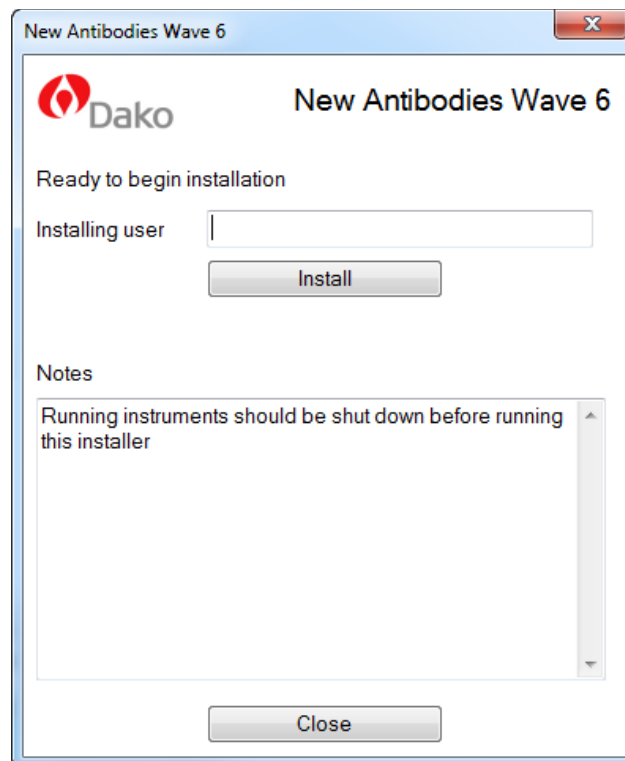
The Antibody Wave 6 installer will:

- Add the Antibody Wave 6 to the DakoLink™ database.
- Define the antibody rules for the added antibodies

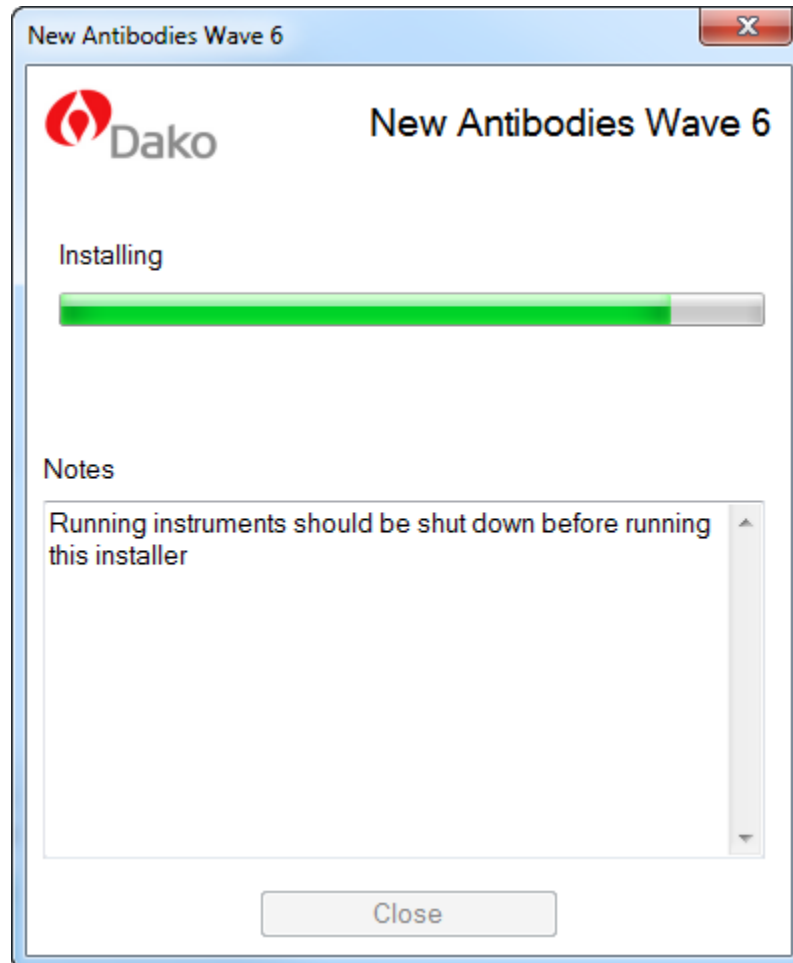
Section 1 | Installing the software

Follow the instructions below to add the Wave 6 Antibodies to DakoLink:

1. Download the installer file (**NewAntibodyWave6.exe**) from www.dako.com/installer to a portable device; e.g., a USB flash drive.
2. Copy the file to a DakoLink™ Instrument Workstation Computer **or** the DakoLink server computer.
3. Before starting the installation, close DakoLink™ and **ensure** that no DakoLink instruments have staining runs in progress. This includes all Autostainer Link, PT Link and Artisan™ Link/Link Pro instruments.
4. Run the **NewAntibodyWave6.exe** file. This will open the installer. The following dialog will be displayed:



5. Enter your name in the **Installing User** field and click the **Install** button. The note stating that “Running instruments should be shut down before running this installer” will be displayed if any ACS (Autostainer Control Software) or ArtiCS (Artisan Control Software) are not shut down.
6. Wait for the installation to complete:



7. The dialog below is shown when the installation is completed:

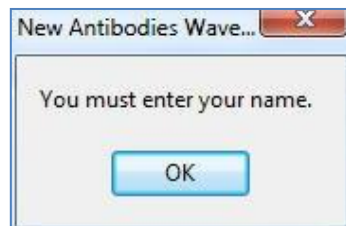


8. Click the **Close** button to exit.
9. LIS/TPID users should contact your local Dako representative for instructions on mapping the new protocols.

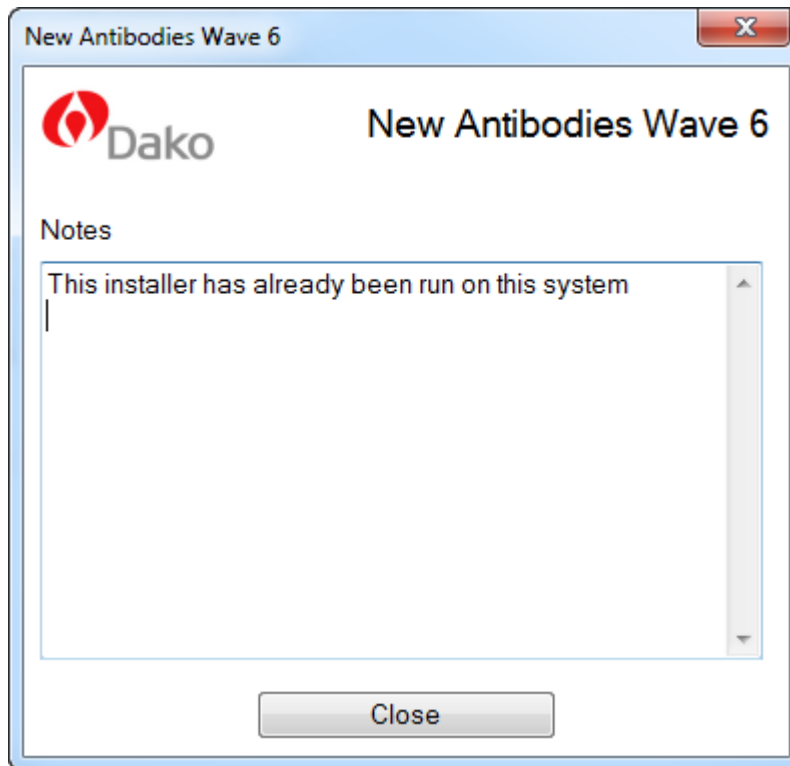
Section 2 | Troubleshooting

The installer may display dialogs with instructions which are not covered by the normal installation procedure. These dialogs are explained below. If you have questions or need assistance with the installation, please contact your local Dako representative.

1. If you do not enter a name in the Installing user field, the installation will not start and the error below is shown. You must enter a name to start the installation.



2. The dialog below is displayed if the protocols have already been installed on the system.



3. If the following dialog is displayed, please ensure that you are running the installer on a DakoLink™ instrument workstation computer or a DakoLink™ server computer. Click the **Close** button and attempt the installation on a different computer.



4. If the installation fails, the following dialog will be shown. Ensure all DakoLink software (client and instrument) is closed and click the **Close** button. Restart from Step 3 in [Section 1 | Installing the software](#).



The text displayed may also be:

Could not reconnect with Dako Link after backup.