

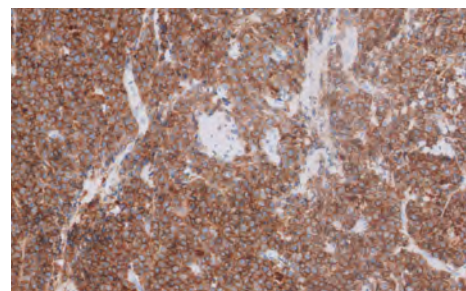
CD10 DAK-CD10 (IR786)

Protocol Installation Guide for DakoLink

Purpose

This document provides instructions for the self-installation of the CD10 DAK-CD10 protocol to support the launch of the new CD10 FLEX RTU (IR786). The protocol must be installed for DakoLink Software to recognize the new product FLEX RTU IR786.

The protocol is compatible with Dakolink 4.1 and versions thereafter.



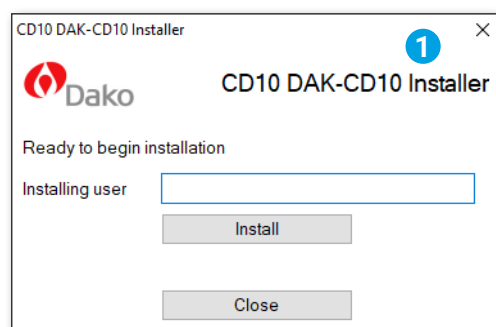
Scope

This procedure covers the following:

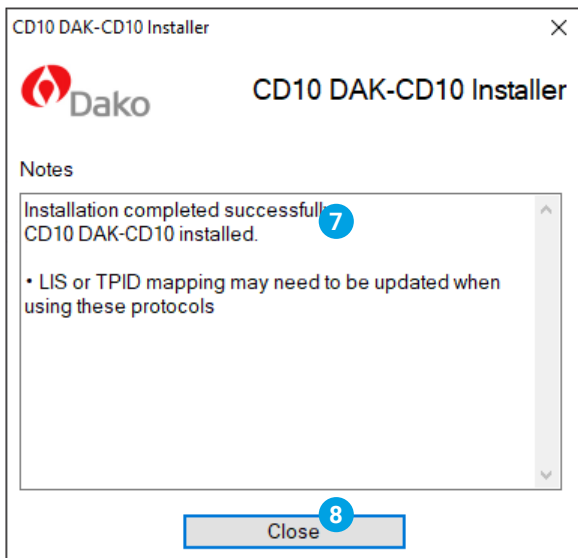
- Installing the Protocol
- Protocol Name for LIS Connected Systems
- Troubleshooting

Installing the Protocol

1. Download the installer file (**CD10 DAK-CD10 Installer.exe**) from <http://www.agilent.com/pathology/installer> to a portable device; e.g., a USB flash drive
2. Before inserting the USB drive into your DakoLink workstation or Server ensure the USB is free from viruses or malware
3. Copy the file to a DakoLink Instrument Workstation Computer or the DakoLink server computer if the system is connected to a network
4. 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
5. Run the **CD10 DAK-CD10 Installer.exe** file. This will open the installer. The following dialog will be displayed:



- 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.
- Wait for the installation to complete
- Click the **Close** button to exit.



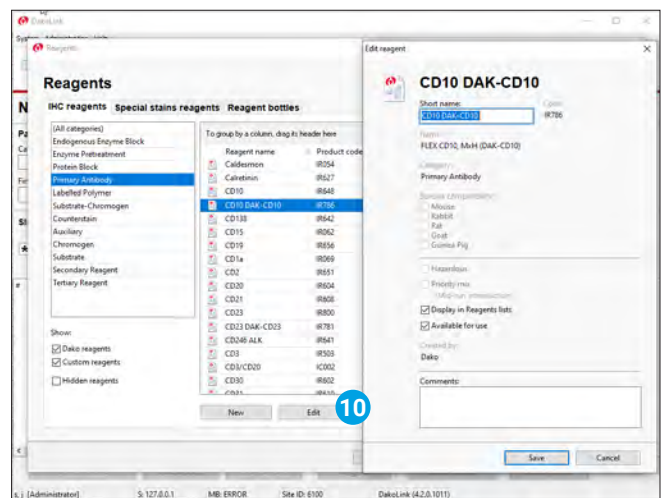
- Log on to DakoLink software and from the **New Slides** tab confirm that the new protocol appears on the **Protocol** list



Protocol Name for LIS Connected Systems

The protocol short name "CD10 DAK-CD10" will likely need to be updated to match the name ordered from the LIS as it likely exceeds the name length that many LIS systems can handle. Determine what name will be sent from the LIS and modify the short name accordingly. Note that if another CD10 product is currently used by the LIS there cannot be two protocols with the same short name. Rename the other CD10 or alter the new CD10 name as necessary.

- From **Administration** → **Reagents** → **Primary Antibody** scroll down to the new CD10 DAK-CD10 antibody and select **Edit**



- Change the **Short name** to match the name that will be ordered from the LIS system and select **Save**

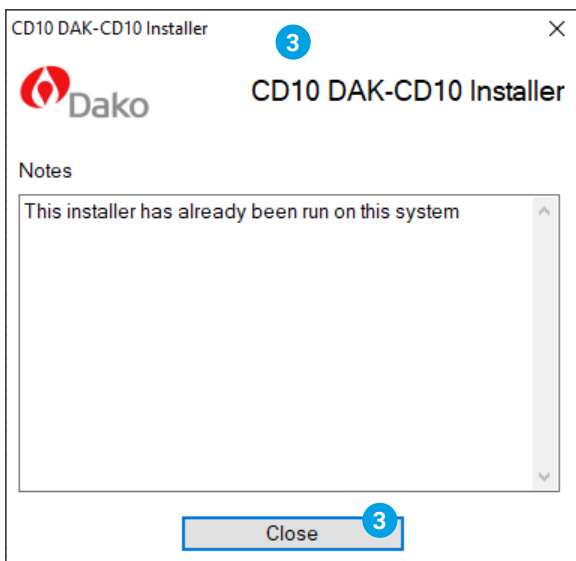
Troubleshooting

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.

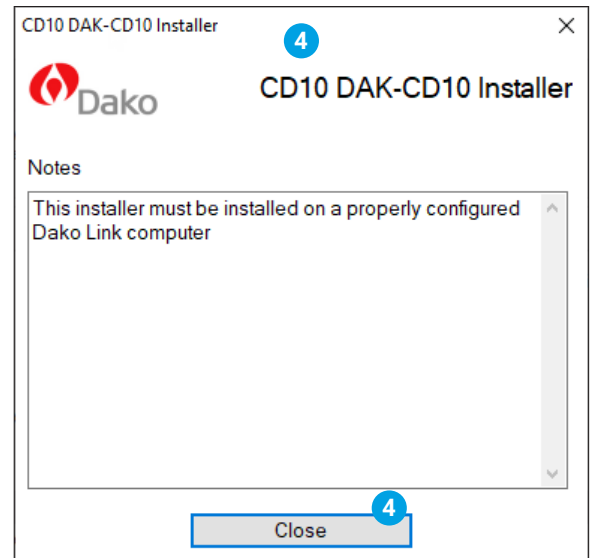
1. If the USB device you have used to download the installer does not appear in Windows Explorer on your Autostainer workstation contact your local Agilent Technical Support as the ability to recognize USB storage devices may be temporarily disabled on your system
2. If you do not enter a name in the Installing user field, the installation will not start and the error below is shown. You must enter a name to start the installation



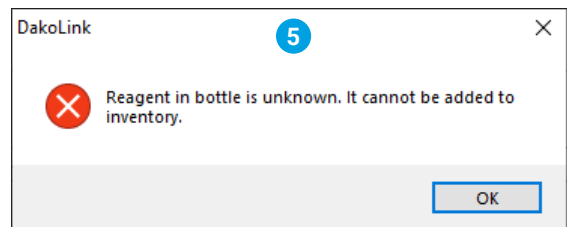
3. The dialog below is displayed if the protocols have already been installed on the system



4. 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



5. If the new reagent is scanned using the hand-held Barcode Scanner before the protocol has been installed, the following message will appear. Confirm that the protocol installer has been run and that the protocol is shown in the Protocols list.



6. If the new reagent is scanned by placing the reagent on the stainer before the protocol has been installed, the following message will appear. Confirm that the protocol installer has been run and that the protocol is shown in the Protocols list.

211170 **6** The reagent in rack <reagent rack number> position <reagent number in rack> cannot be found in DakoLink's database.

For any issues or questions which may arise during this process please contact your local Agilent Technical Support representative.

www.agilent.com

This information is subject to change without notice.

© Agilent Technologies, Inc. 2021
Printed in the USA, September 24, 2021
D68396_01 2021SEP23

