

Dako Omnis Import of Protocols

To import protocols to the Dako Omnis follow the following five steps:

1. Check whether the new protocol name, the one you want to import, already exists in the **Protocol Library**.
2. If the new protocol name exists in the **Protocol Library**, you will need to rename the current custom protocol (see the **Rename protocol** instruction below).
3. Import protocols (see the **Import protocol** instructions below).
4. Create test as needed (see the **Create Tests** and **Adding New Tests** instructions below).
5. Map protocol to test as needed (see the **Map protocol** instruction below).



Rename Protocol

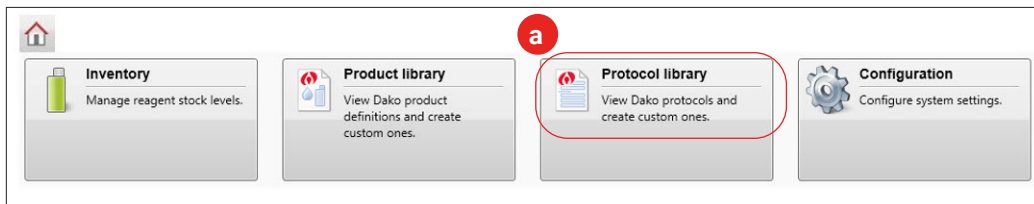
- a. Click the **Protocol library** panel on the **Home** screen to access the **Protocol library** screen.
- b. Highlight the custom protocol you want to rename, CD23 in the example below:

Protocol na...	Test category	Protocol category	Origin	Version	Approved for
CD20cy	IHC	EnVision FLEX prot...	Dako	27-08-2013 16:57	Diagnostic use
CD23	IHC	EnVision FLEX prot...	Custom	01-01-2013 14:58	Diagnostic use
CD3	IHC	EnVision FLEX prot...	Dako	29-04-2013 07:52	Diagnostic use

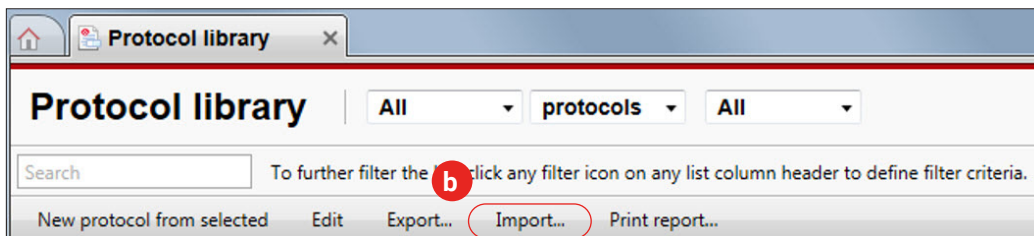
- c. Click the **New protocol from selected** button.
- d. Rename the protocol and click **Save**.
- e. If the custom protocol was mapped to a test, this must be updated through **Configuration** to the new protocol name to continue using the custom protocol.

Import Protocol

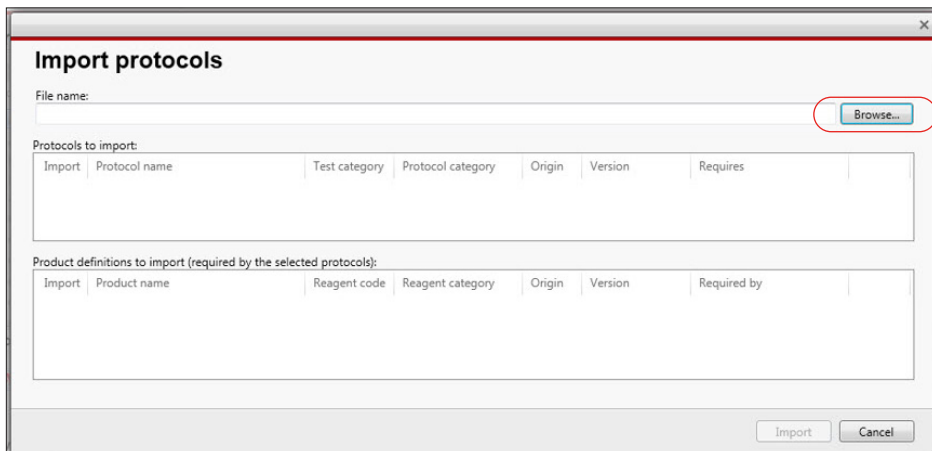
- a. Click the **Protocol library** panel on the Home screen to access the **Protocol library** screen



- b. Click the **Import...** button in the **Toolbar**



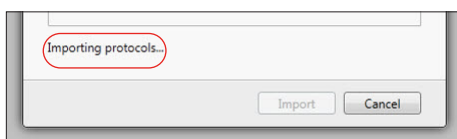
This opens the **Import protocols** dialog:



It allows you to browse for new protocols in the .ddx (Dako Data Exchange) file format to import into your **Protocol library**, e.g., when Dako releases new protocols.

- c. Click Browse... and select the data file(s) of interest
- d. Select the protocols you wish to import
- e. Click the Import button

During import, the following message appears:



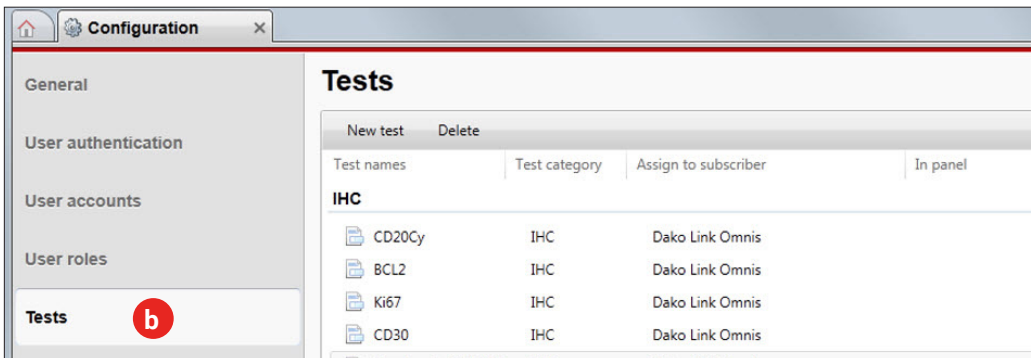
Note: Related products are also imported as part of the process.

Create Tests

- a. Click the **Configuration** panel on the **Home** screen to access the **Configuration** screen



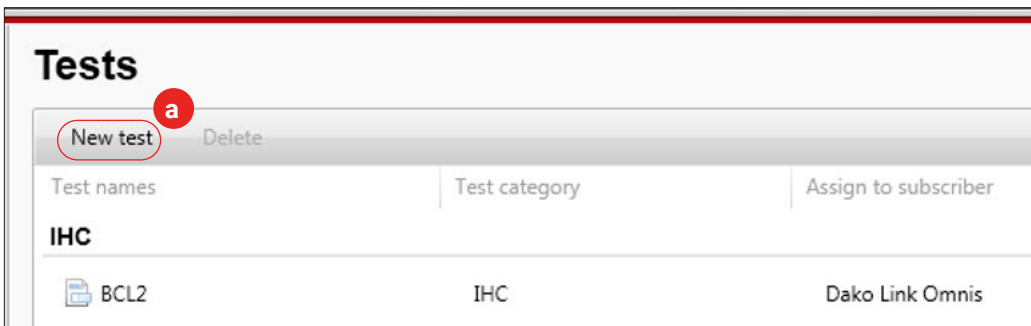
- b. Click the **Tests** button to open the **Tests** screen



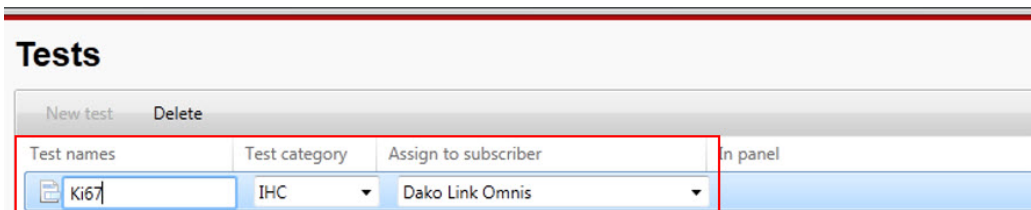
Adding New Tests

If your system is set up to LIS, tests from LIS will automatically be transferred to the Dako Omnis Workstation. To manually add tests, follow the steps below:

- a. Click the **New test** button



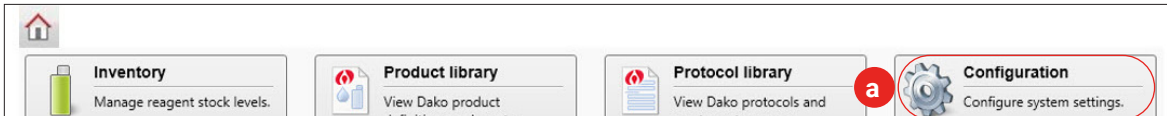
A blank field appears at the top of the Tests list:



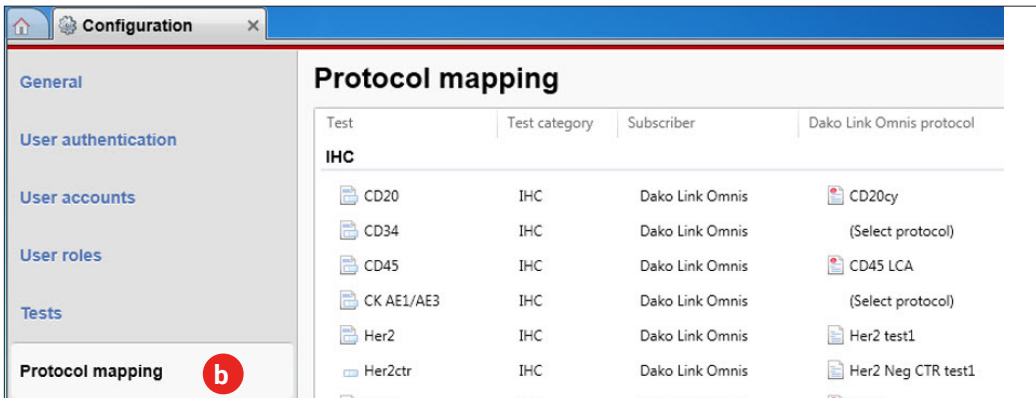
- b. Enter the specific test name in the **Test names** field
- c. Select the category from the **Test category** drop-down list:
HE / SS / IHC / CISH / FISH
- d. Select the instrument for distribution from the **Assign to subscriber** drop-down list
- e. Repeat until you have entered all tests
- f. Click **Save**

Map Protocol

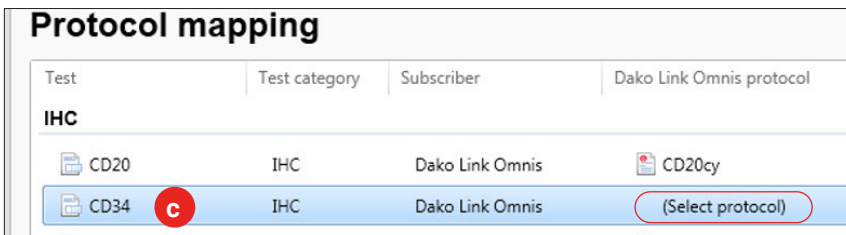
- a. Click the **Configuration** panel on the **Home** screen to access the **Configuration** screen.



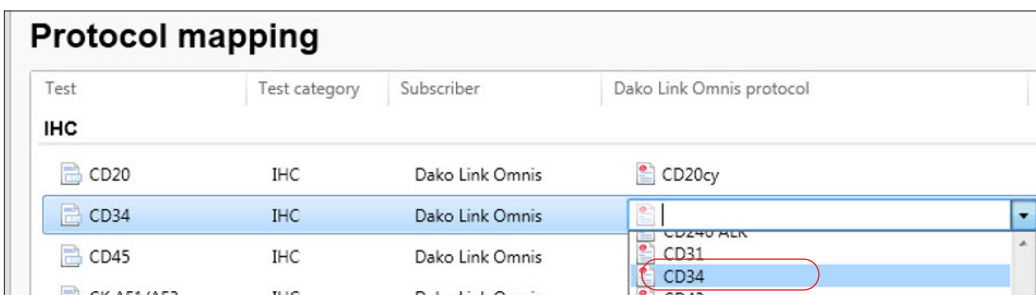
- b. Click the **Protocol mapping** button to open the **Protocol mapping** screen.



- c. Select the new test without protocol ('Select protocol'). Through the **Tests** screen, the new test has been named and subscribed to Dako Omnis.



- d. Select the relevant protocol from the **Dako Link Omnis protocol** drop-down list. This list contains all protocols.



In this manner, you point to the specific CD34 protocol that is approved for diagnostic use and do not by accident use a non-approved protocol.

- e. Repeat steps 3 and 4 until all new tests are mapped to protocols.
- f. Click **Save**

www.agilent.com

This information is subject to change without notice.

© Agilent Technologies, Inc. 2020
D59736 2020APR16