

spectroquant® prove connect

to Dashboard



1	Introduction	3
2	Quickstart	3
	System Requirements	3
3	Create Account	3
4	Login.....	4
5	Register a Spectroquant® Prove Instrument.....	5
6	Connect a Spectroquant® Prove Instrument	7
7	Add New User	9
8	Register as User	10
9	Add New Groups (Admin)	11
10	Manage the Measurement Tasks	12
11	Perform a Measurement	13
12	Plot the Data	15
13	Get Audit Trails.....	16

1 Introduction

Thank you for choosing Prove Connect. Prove Connect is a cloud application to connect your Spectroquant® Prove instruments to the cloud and share results with your team. You can access your data from everywhere at any time, using any device. Just open a web browser, go to www.proveconnect.com and log in. The software receives and saves measurement data from Spectroquant® Prove instruments over the internet.

2 Quickstart

System Requirements

System requirements:

- Spectroquant® Prove 100
- Spectroquant® Prove 300
- Spectroquant® Prove 600

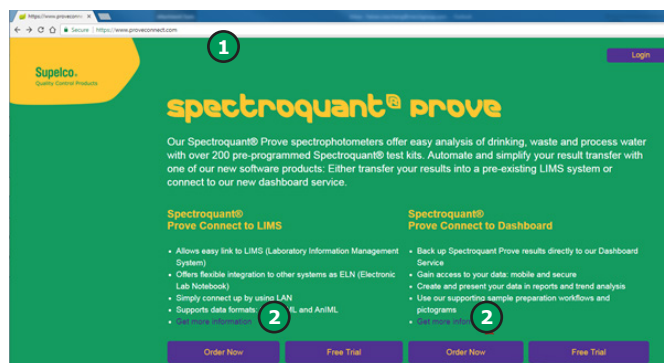
The software requires a computer with the following specification:
Windows Version 7, 8 or 10 or Mac OS X, 32bit or 64bit
Google Chrome Version 45+,
Mozilla Firefox 45.9+, or
Safari 10+
Network connection

NOTE

Other configurations may work, but are not supported.

3 Create Account

1. Open <https://www.proveconnect.com> in a web browser ①.
2. Click on Order now ②. You will be redirected to our webshop where you can purchase your Prove Connect to Dashboard account



3. After you followed the order process you receive a confirmation email containing your login details ③. It is recommended, to change the password immediately, refer to section 8.



Welcome

Hi Tim Test,

An account has been created for you with the following credentials:

Login: Tim.Test@Tester.com
Password: [Tim@Test](#)

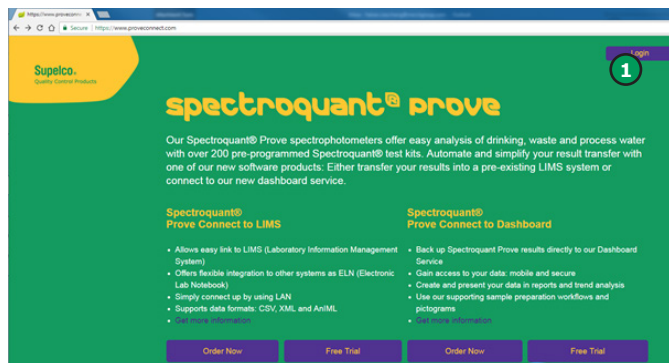
③

[Change password](#)

4 Login

4 Login

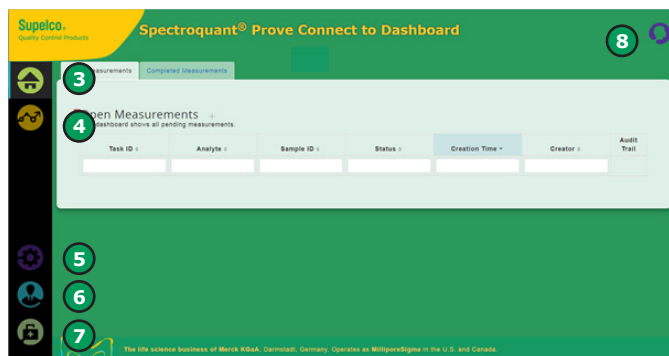
1. Log in to your Prove Connect to Dashboard account on www.proveconnect.com using the Login button **1**.



2. After successful login the Prove connect to dashboard start page **2** appears.

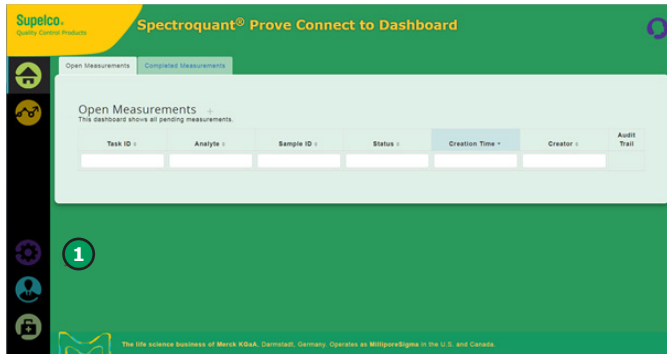
On the left hand side you find the menu bar containing all the icons for the available functions:

- 3**: Home
- 4**: Trending
- 5**: Administration
- 6**: Contact us
- 7**: Help?
- 8**: Profile

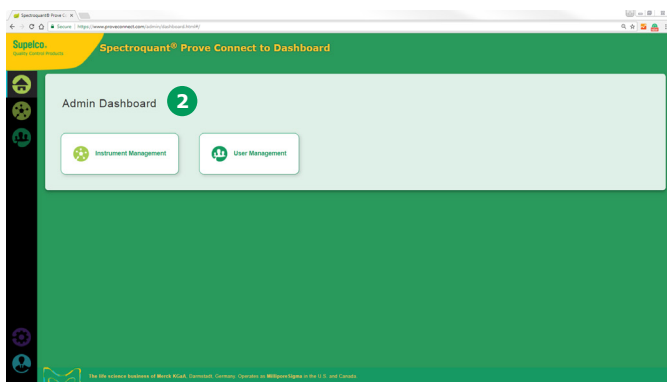


5 Register a Spectroquant® Prove Instrument

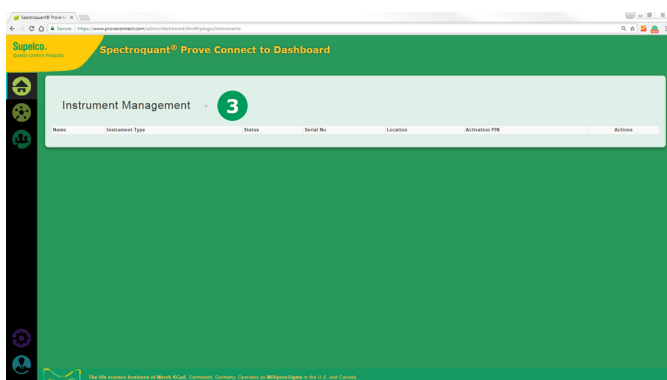
1. Log in to Prove Connect.
2. Select Administration **1**.



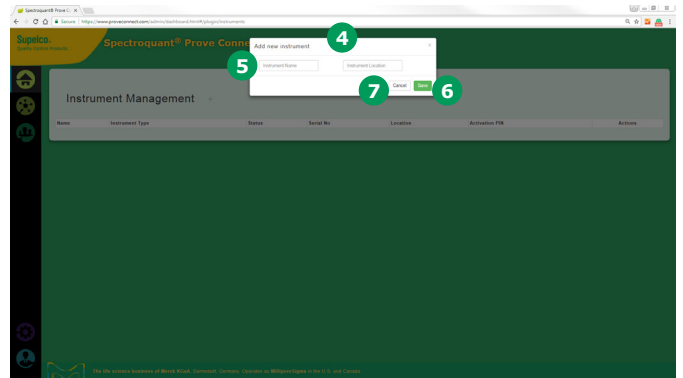
3. On the Admin Dashboard select Instrument Management **2**.



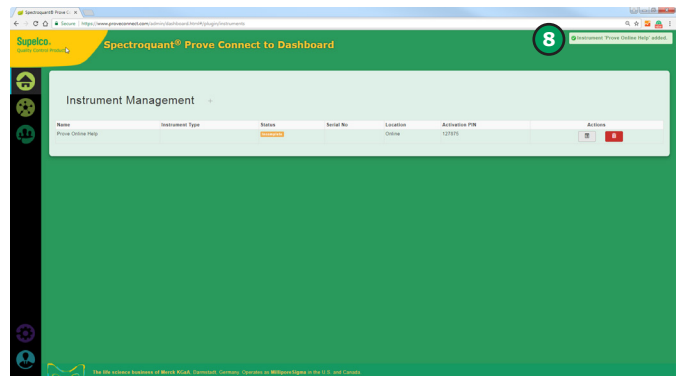
4. In the Instrument Management click on the + symbol **3** next to Instrument Management.



5. A pop up "Add a new instrument" **4** appears.
6. To register the instrument insert the instrument name and the instrument location **5**.
7. Click the Save button **6** to apply the entries. To abort click the Cancel button **7**.

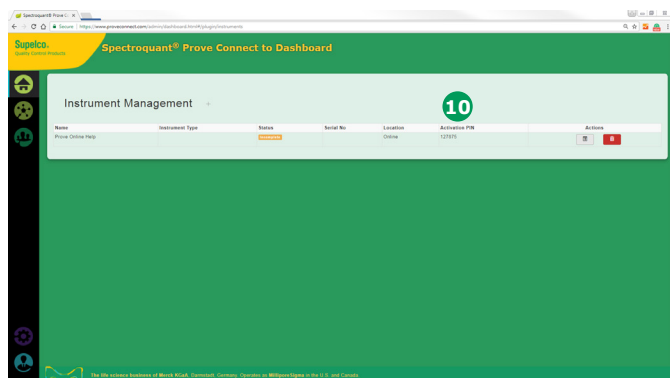


8. A notification informs the user that the instrument has been successfully registered **8**.



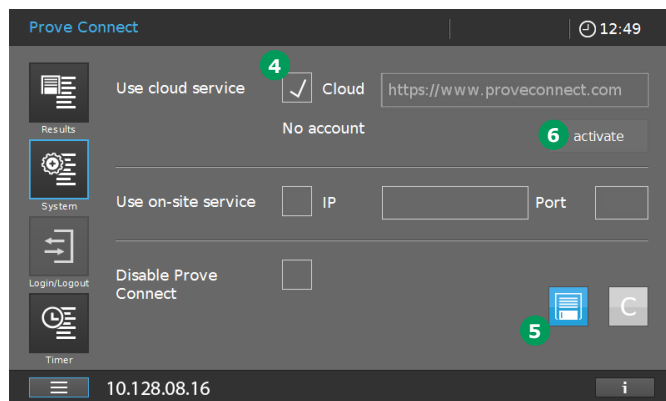
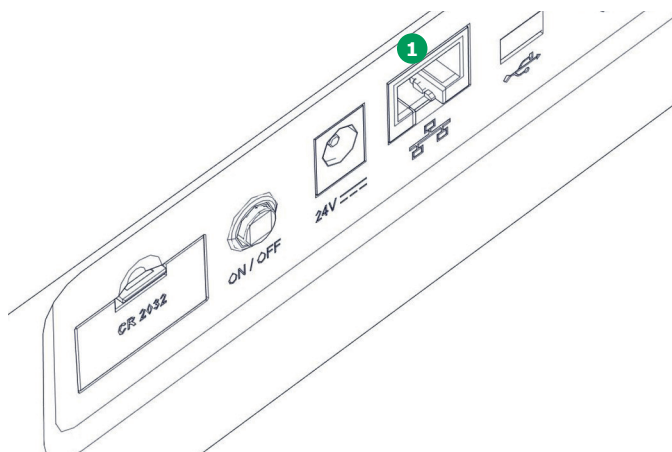
5 Register a Spectroquant® Prove Instrument

9. The instrument is now listed in the updated instrument management list.
10. Prove Connect generates an Activation PIN **10**, which needs to be entered into the Prove instrument (see section 6: Connecting a Spectroquant® Prove Instrument).

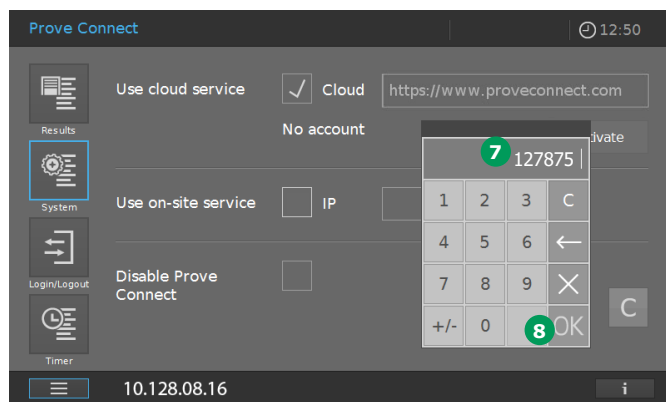


6 Connecting a Spectroquant® Prove Instrument

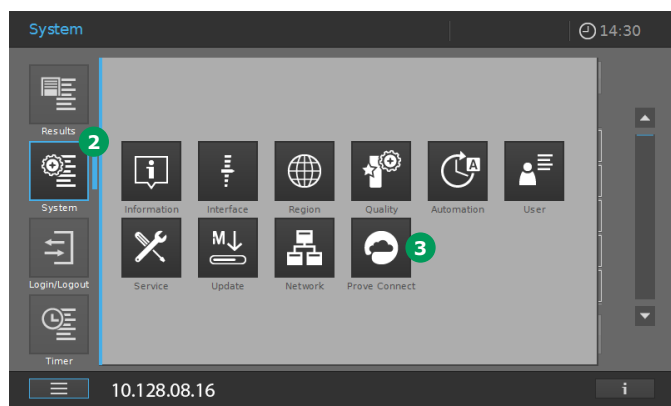
1. Use the Ethernet port **1** on the back side of the instrument to connect the Spectroquant® Prove to the internet and turn on the instrument. A network cable (RJ45) is required (not included).
2. To configure the Prove instrument open the setup tab in the instrument operation system **2**.
3. Select Prove Connect **3**.
4. Select Enable Use Cloud Service **4** in the Prove Connect menu.
5. Click on the Save button **5**.
6. Click on Activate **6**.



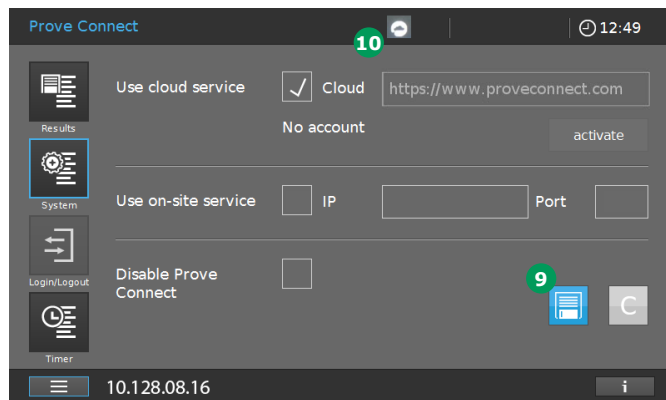
7. Enter PIN **7** and confirm with OK **8**.



2. To configure the Prove instrument open the setup tab in the instrument operation system **2**.
3. Select Prove Connect **3**.

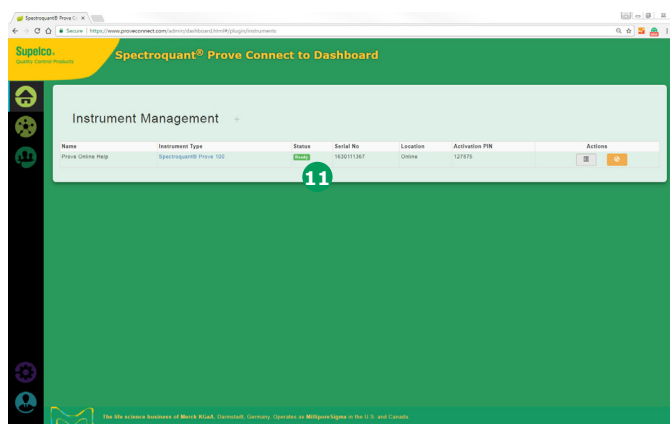


8. Save the settings by clicking on the save button **9**.
9. A cloud icon **10** appears in the upper status bar. The instrument is now connected.



5 Connecting a Spectroquant® Prove Instrument

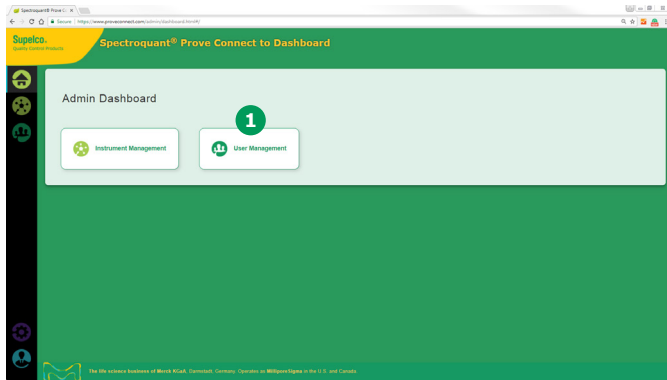
10. The status in the instrument list changes to "Ready" **11** in the Prove Connect to Dashboard Instrument Management tab.



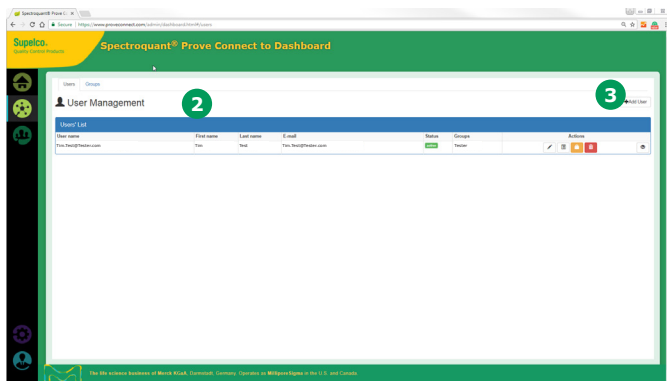
7 Add New User

To add users to your group go to the "User Management" menu in the Admin Dashboard.

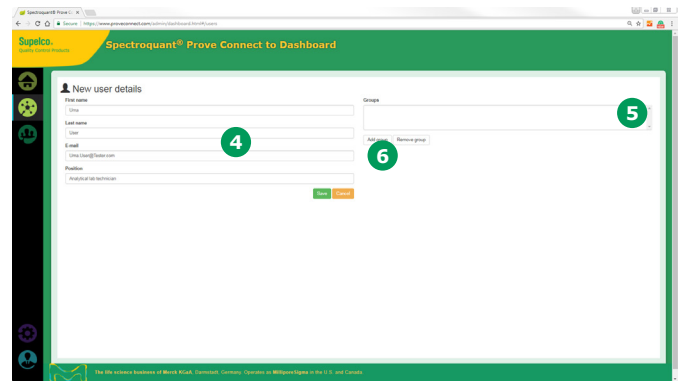
1. Click on "User Management" **1**.



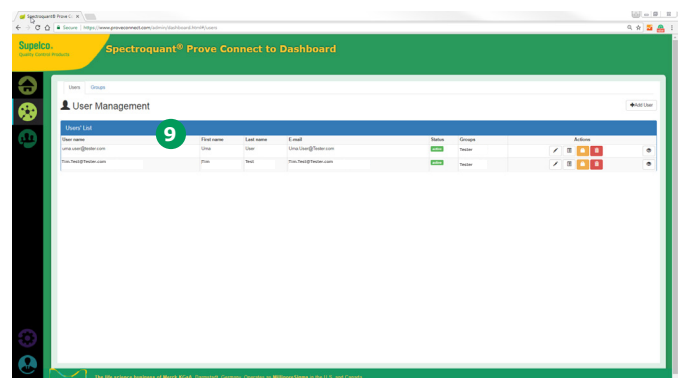
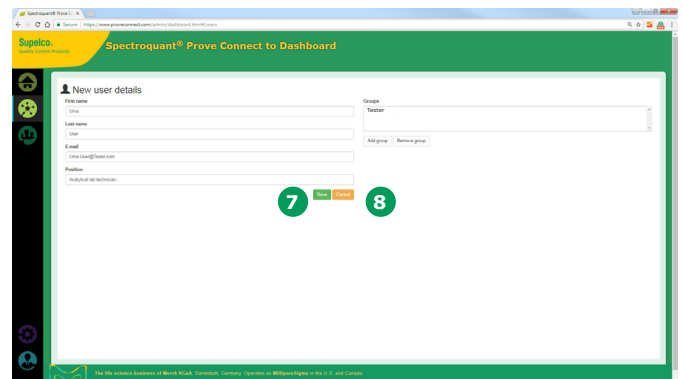
2. An overview of all current users appears **2**.
3. To add a new user click on the "Add User" button **3**.



4. Fill in the user information into the registration form on the left side **4**.
5. To assign the new user to a user group click on "Add group" and select the correct group from the drop down menu **5**. Click "Add to groups" **6** to confirm the selection and to assign the user to the group. You can assign a user to multiple groups.



6. Click on the Save button **7** to create the new user account. To discard all entries and abort click the Cancel button **8**.
7. The new user **9** has been created and receives an email to register.



8 Register as User

8 Register as User

As soon as the Admin has added you as new user in his organization, you will receive a registration email:

1. In your registration email click on the "Change Password" button **1**. Ensure you have turned off your pop up blocker in the standard web browser.



Welcome

Hi Tim Test,

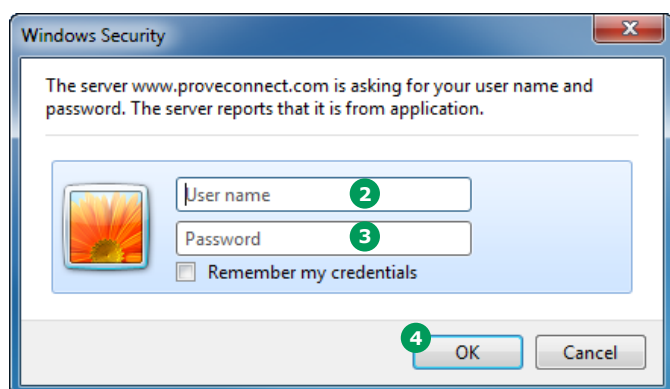
An account has been created for you with the following credentials:

Login: Uma.User@Tester.com

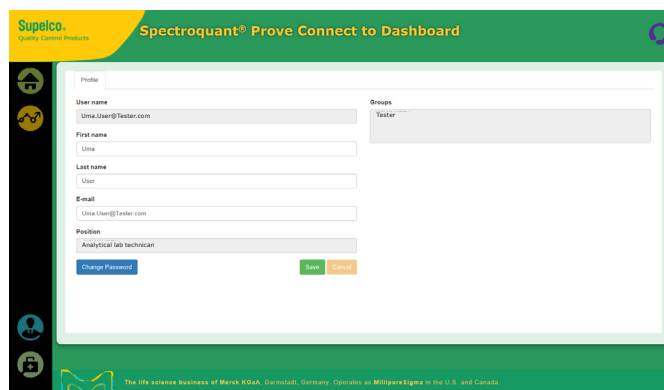
Password: XVEUnu80

Change password **1**

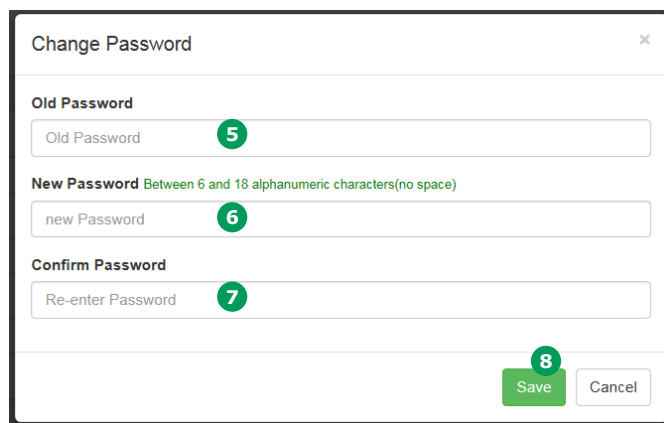
2. Enter your login details (user name/login **2** and password **3**) into the dialog pop up and click OK **4**. The Prove Connect user settings web page opens.



3. The form on the user profile page is prepopulated with the user data provide during the account creation. Click on change password to set a new password.



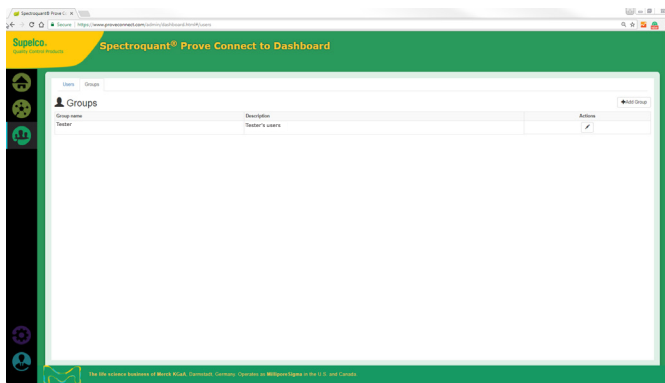
4. Enter you initial password **5** and set a new password **6**. Confirm the new password **7**. Press Save **8** and the new password is effective immediately.



5. If you have used any browser other than the ones given in section 2.1, please close the browser and all its instances to log off and start i. e. "Google Chrome" to work with Prove Connect.

9 Add New Groups (Admin)

If you have more than one team in your organization, you can generate separate groups for each team. Users can only access the data and instruments the admin has assigned to the group the user belongs to. Click on the "groups" tab in the user administration tab to add new groups, delete them or change the name and description of an existing group. All groups that you generate belong also to a basic group that is automatically created with the account.

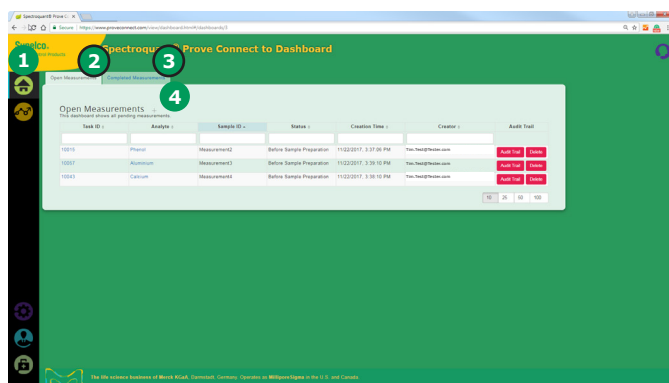


10 Manage the Measurement Tasks

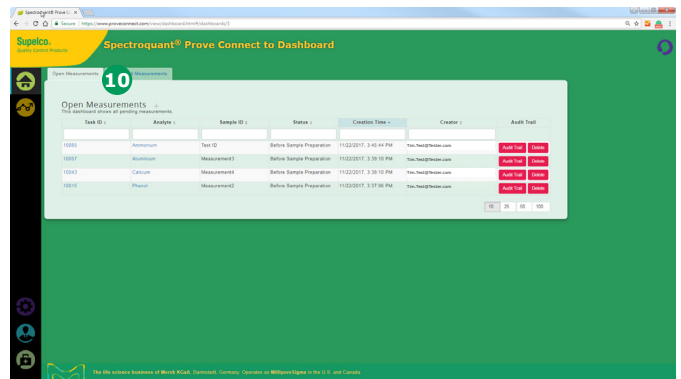
10 Manage the Measurement Tasks

Prove Connect to Dashboard offers you an easy to use measurement & sample administration system to reduce your documentation work.

Click the Home button **1** in the menu bar. The dashboard appears. It contains two lists in two individual tabs. One for open measurement tasks **2**, and a second one for completed measurements **3**.



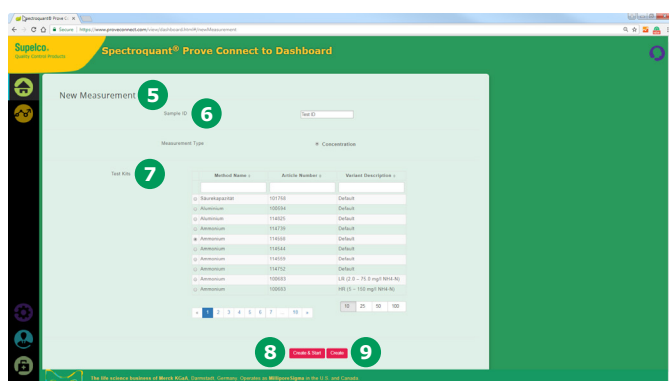
- The newly created measurement is listed under the Open Measurement tab **10**.



NOTE

You can use the open measurements tab like a sample receipt book.

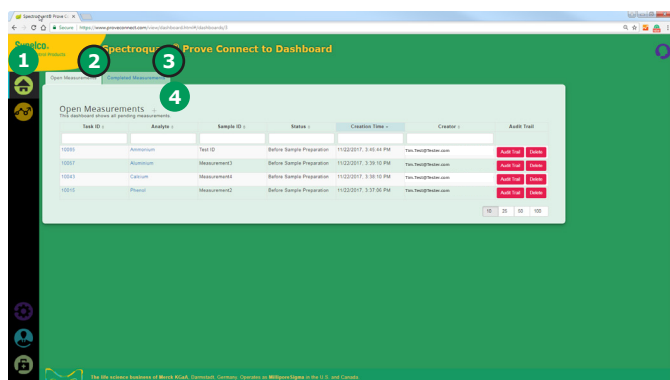
- Click on "+" next to **4** "Open measurements" to add new samples.
- A New Measurements form opens **5**.
- Enter the Sample ID **6** and choose the Test Kit **7**.
- Click on "Create" **8**, to create a new entry in the "Open Measurements" list or click on "Create & Start" **9** to create a new entry in the "Open Measurements" list and start the workflow.



11 Perform a Measurement

Prove Connect to Dashboard offers you an easy to use measurement & sample administration system to reduce your documentation work.

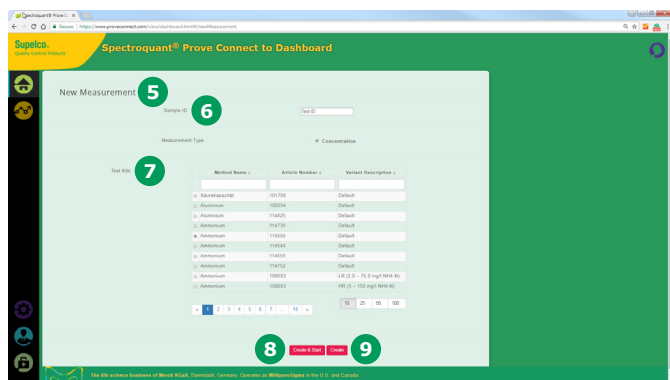
Click the Home button **1** in the menu bar. The dashboard appears. It contains two lists in two individual tabs. One for open measurements tasks **2**, and a second one for completed measurements **3**.



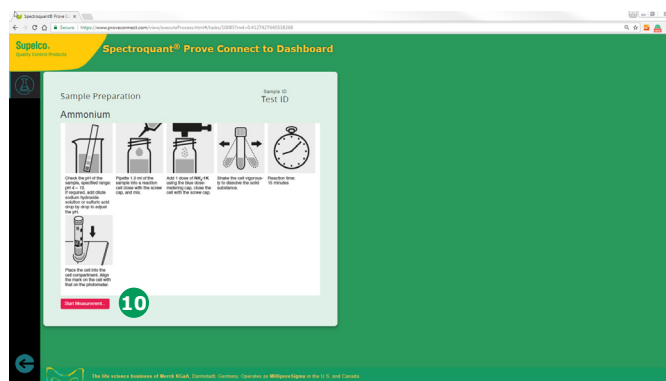
NOTE

You can use the open measurements tab like a sample receipt book.

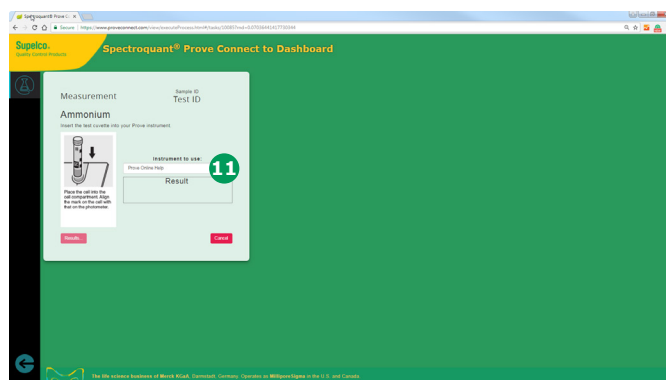
1. Click on "+" next to **4** "Open measurements" to add new samples.
2. A New Measurements form opens **5**.
3. Enter the Sample ID **6** and choose the Test Kit **7**.
4. Click on "Create" **8**, to create a new entry in the "Open Measurements" list or click on "Create & Start" **9** to create a new entry in the "Open Measurements" list and start the workflow.



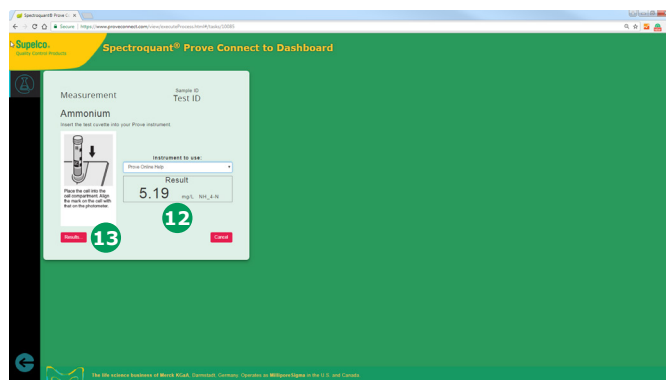
5. When you start your measurement Prove Connect will guide you through the sample preparation with explanatory pictograms.
6. Click "Start measurement" **10**.



7. Select an instrument from the drop down menu **11**.
8. Start the measurement on the instrument.

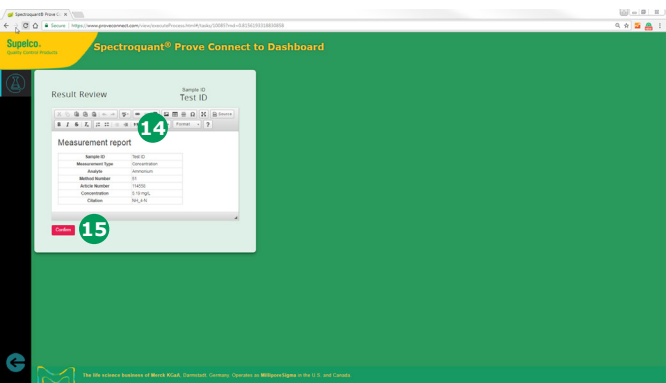


9. The result is automatically transferred from the instrument and presented on the screen **12**.

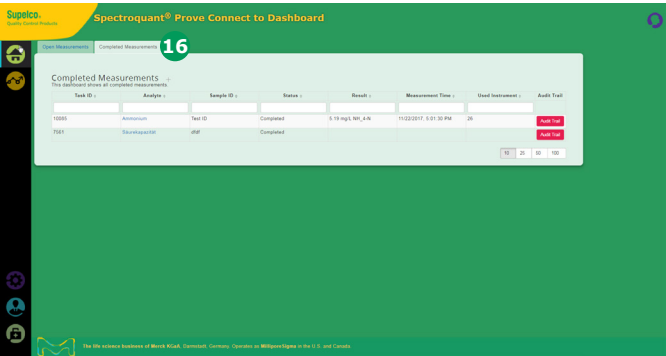


10. Click on "Results" **13** to review and edit the measurement report **14**.

11 Perform a Measurement



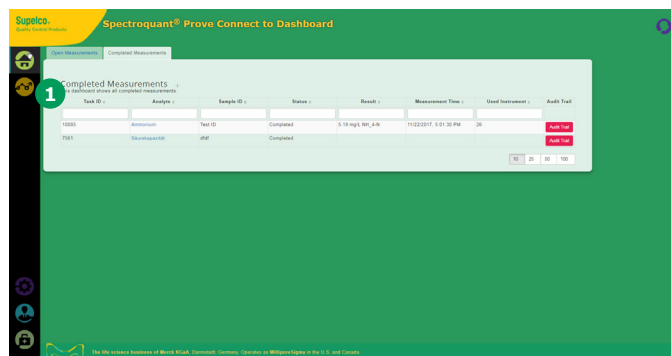
- 11. Click "Confirm" 15 to end the measurement.
- 12. The completed measurements are listed under in Completed Measurements tab 16.



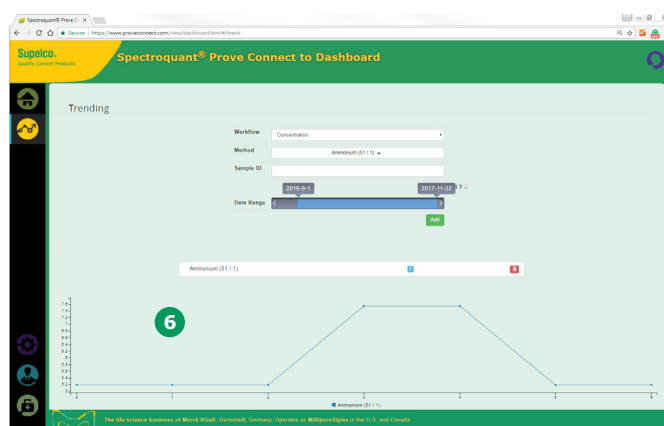
12 Plot the Data

Prove Connect to Dashboard offers the option to create trend charts based on your measurement results.

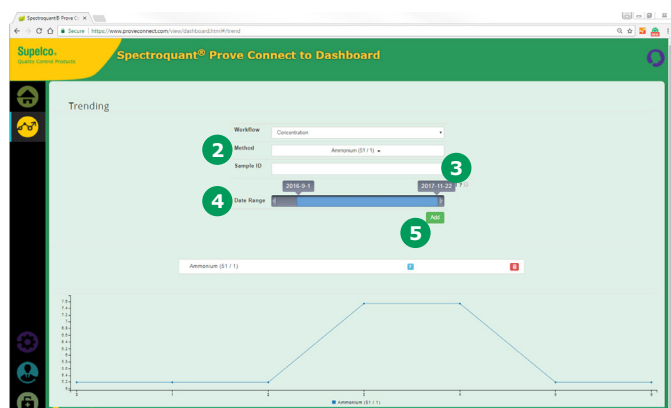
1. Click on the "Trend" icon **1** in the menu bar.



6. Add further data series by choosing another method or sample ID. The data series are overlaid **6**.



2. The trending dashboard opens.
3. Select the method **2**.
4. You can narrow the result included into the trend analysis by filtering for sample ID **3** and/or by selecting a time period **4**.

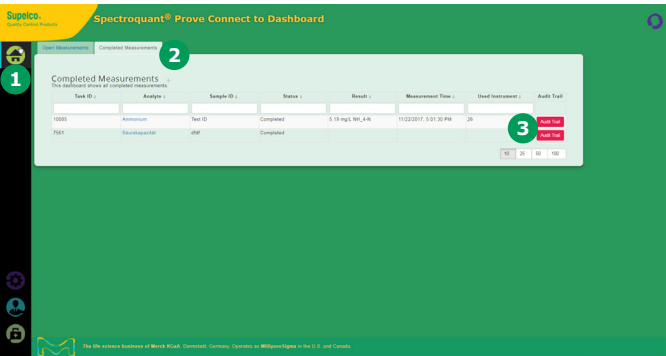


5. Click on "Add" **5**.

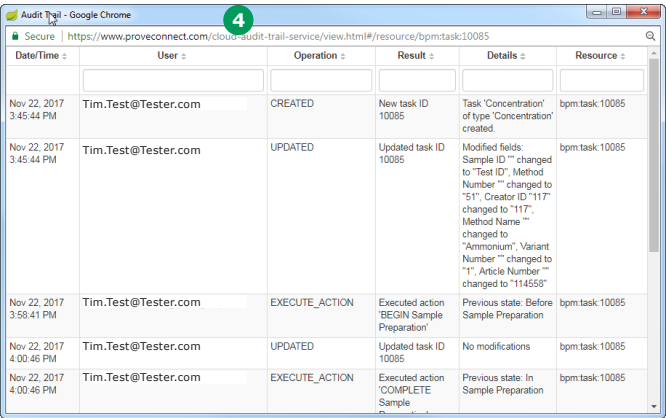
13 Get Audit Trails

To view audit trails for all your instruments, users and tasks, click on the button "Audit Trail" at the end of each row in the instrument, user or task list. You get an overview of all activities that were performed.

- 1. Click on the Home Button 1.
- 2. Switch Tab to "Completed Measurements" 2.
- 3. Click on "Audit Trail" 3 .



- 4. A list with all available information is presented 4.



We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.

Manufactured by Merck KGaA, 64271 Darmstadt, Germany, Tel. +49(0)6151 72-2440
www.analytical-test-kits.com

Distributed by EMD Millipore Corporation, 290 Concord Road, Billerica,
MA 01821, USA, Tel. +1-978-715-4321

