



MACSQuantify™ App

Frequently asked questions

Lab instrument

(How) Can I connect my lab instrument to a network?

Each MACSQuant® Analyzer can be connected to a network. In order to use the cloud solution together with the MACSQuantify™ App, an internet connection is required. The initial software configuration to be able to gain cloud access needs to be done by your local admin or a service technician.

Local systems / local IT

Can I integrate my own/local data within the cloud databases?

It is possible to upload experiment settings and sample annotations into the cloud database by using our Microsoft® Excel Add-In.

Can I connect to the cloud via my proxy?

Yes, this is possible. Please specify the correct proxy/port during the MACSQuantify installation process on your MACSQuant Analyzer and analysis computer.

Cloud

Where is the data physically stored/where is the location of the server?

The cloud server we are using is located in Ireland. Therefore, the European data protection law applies.

Which data is stored?

Data that is entered by the user is stored as "user data". In order to enable tracking, each action and data entry is connected with the user account and a time stamp. In detail, the data that is stored includes: sample information, reagent information, panels, rack configuration with samples, reagents, and settings. In order to enable future services, additional data (instrument-specific data, instrument logs) will also be collected once the services are available. The results (raw data, or analyzed data) are not synchronized to the cloud.

For how long is the data stored?

There is no limitation in terms of time for data storage. Data that is not used anymore can be moved to an archive section, from which it can be retrieved if required.

Is my data used for any other purposes than planning experiments?

When using the cloud, your data is solely used for planning experiments, which means that your experiment configurations are stored in order to transfer them to your MACSQuant Analyzer. Miltenyi Biotec is not using your user data for analysis or re-distribution.

MACSQuantify App

How can I install the app on my device?

To register for a MACSQuantify App test license, please visit www.miltenyibiotec.com/MACSQuantifyApp and fill in the contact form. Our development team will contact you with detailed information on how to download and install the app. Upon official release, the app will be made available in public app stores (Apple® App Store, Google Play™, Windows® Store). After registering with our cloud backend, your user account will be activated. You will receive a message as soon as you can connect to the cloud and use the app.

How can I update the app on my device?

During the test-phase, the app can be updated via the test store from which the app was downloaded. After the official release, updates are provided via the common app stores. The app can be updated following the standard update routines.

Are there any hardware prerequisites for using the app?

The app will only be available on tablet devices (minimum 10 inch display) with operating systems iOS 9, Android 4.1, Windows® 8.1, or higher.

Is a special WLAN or data transmission rate required?

A slow internet connection may result in delayed performance speed when it comes to data synchronization to the cloud.

Microsoft® Excel Add-In

How can I install the add-in on my computer?

To download the add-in, please visit:

www.miltenyibiotec.com/MACSQuantifyApp.

You will also find an installation guide there. In order to use the add-in together with the app, you have to log-in with your (cloud/app) user account.

Are there any software prerequisites for using the add-in?

The add-in requires Excel 2013 (Service Pack1) or higher.

Are there any hardware prerequisites for using the add-in?

No special hardware prerequisites apply.

Is a special WLAN or data transmission rate required?

A slow internet connection may result in delayed performance speed when it comes to data synchronization to the cloud.



Miltenyi Biotec

Miltenyi Biotec GmbH | Phone +49 2204 8306-0 | Fax +49 2204 85197 | macs@miltenyibiotec.de | **www.miltenyibiotec.com**

Miltenyi Biotec provides products and services worldwide. Visit **www.miltenyibiotec.com/local** to find your nearest Miltenyi Biotec contact.

Unless otherwise specifically indicated, Miltenyi Biotec products and services are for research use only and not for therapeutic or diagnostic use. MACS, MACSQuantify and the MACS logo are registered trademarks or trademarks of Miltenyi Biotec GmbH and/or its affiliates in countries worldwide. All other trademarks mentioned in this document are the property of their respective owners and are used for identification purposes only. Copyright © 2017 Miltenyi Biotec GmbH and/or its affiliates. All rights reserved.