

Biacore™ X100 Software v2.1

Release Notes

Introduction

Scope of the document

This document describes new and changed functionality of Biacore™ X100 Software v2.1 compared to Biacore X100 Software v2.0.3. Known software issues are also listed.

The difference in v2.1 compared to v2.0.3 is that v2.1 has improved security and employs a SQL Server database. The functionality for database administration and backup procedures has changed from version 2.0.3.

The release notes is also accessible on the web using the product key. See [cytiva.com](https://www.cytiva.com) for the latest available version.

Included software

This document includes information about Biacore X100 Software v2.1 that consists of the following parts:

- Biacore X100 Control Software v2.1
- Biacore X100 Evaluation Software v2.1
- Biacore X100 Database v2.1

System requirements

Biacore X100 Software v2.1 requires Windows 10 or 11 Professional or Enterprise operating system, 64-bit US English version.

The Biacore X100 Software v2.1 does not support Windows 32-bit operating system.

The Biacore X100 Software includes installation of a local SQL Server database.

Upgrading to v2.1

It is recommended to upgrade to v2.1 because of its improved security. An upgrade to v2.1 requires installation of SQL Server Express 2022 database and a migration of data from the previous Oracle database to the version 2.1 SQL Server database. Refer to the *Biacore X100 Software v2.1 Installation Instructions (29288193)* for more information. It is critical that any data migration is done after the installation of v2.1 and before any other use of the new Biacore X100 system. Any run data acquired by the Biacore X100 v2.1 system will otherwise be deleted by the data migration.

Biacore X100 Software v2.1 can not be installed on the same computer as a Biacore X100 Software v2.0.1 or earlier version due to difference in supported Windows operating system (64-bit instead of 32-bit).

Because of the new type of user accounts, an upgrade to v2.1 may lead to duplicates of user names in the folder structure. Refer to the *Biacore X100 Software v2.1 Installation Instructions* for recommended strategies to handle this.

Note: *It is recommended to backup the current database before installing the v2.1 database.*

Functionality

Biacore X100 instrument is run by Biacore X100 Control Software. It is mainly used on a system controller connected to an instrument, but can also be used with limited functionality on a computer not connected to an instrument. The software provides workflows to guide the user through the steps needed to develop and run an assay. The wizards used in the workflows can also be accessed directly outside a workflow. It is also possible to operate the system via manual run, which provides interactive control of the instrument operation.

Biacore X100 Evaluation Software offers general functions for presentation of results as sensorgrams, report point plots and/or bar charts and application specific evaluation functions for common applications. Evaluation of assay development steps is included in Biacore X100 Control Software.

Both software components come with integrated support including methodology recommendations, tips and troubleshooting. The support also includes direct links to more in-depth information from Biacore's extensive knowledge base (these links require access to Internet).

Biacore X100 Control Software v2.1

The functionality is described in more detail in *Biacore X100 Handbook (BR100810)*.

New functionality

No	Description
1	<p>Biacore X100 Control Software v2.1 is accessed with Windows or Active Directory accounts. It is no longer possible to log in with the Biacore X100 user accounts from previous software versions.</p> <p>Note: <i>It is not possible to log in with user accounts that have no password.</i></p> <p>Note: <i>If you migrate the data from the old Oracle database to the new SQL Server database any saved data produced with the old Biacore X100 user accounts remain in the database in the same folder structure as before.</i></p>
2	<p>Web links in Support Navigator does not open through Internet explorer as the browser is no longer supported by Microsoft. Instead, from v2.1 all links are opened in default web browser.</p>
3	<p>Updates in Support Navigator are in line with updates in the respective Instructions for use:</p> <ul style="list-style-type: none">• The recommendation to use hydrazine for immobilization with aldehyde coupling has been replaced with recommendation to use carbonylhydrazide that is less toxic.• A warning about the toxicity of cyanoborohydride has been included.

Corrected issues

No	Description
1	<p>The Biotin Capture Reagent was only injected over the active flow cell even though the Binding Wizard for CAP chips stated that it should be injected over both. From v2.1, the reagent is injected over both flow cells.</p>
2	<p>New models based on the Two State Reaction model created in model editor were not correct. This now fixed.</p> <p>Note: <i>Limited to the model editor. The Two State Reaction model used for curve fitting never contained this error.</i></p>

Known remaining issues

There are no known issues in this version.

Biacore X100 Evaluation Software v2.1

The functionality is described in more detail in *Biacore X100 Handbook (BR100810)*.

New functionality

No	Description
1	<p>Biacore X100 Evaluation Software v2.1 is accessed with Windows or Active Directory accounts. It is no longer possible to log in with the Biacore X100 user accounts from previous software versions.</p> <p>Note: <i>It is not possible to log in with user accounts that have no password.</i></p> <p>Note: <i>If you migrate the data from the old Oracle database to the new SQL Server database any saved data produced with the old Biacore X100 user accounts remains in the database in the same folder structure as before.</i></p>
2	<p>Web links in Support Navigator does not open through Internet explorer as the browser is no longer supported by Microsoft. Instead, from v2.1 all links are opened in default web browser.</p>

Corrected issues

No issues are corrected in this version.

Known remaining issues

There are no known issues in this version.

Biacore X100 Database version 2.1

The Biacore X100 Software v2.1 includes a SQL Server 2022 Express database.

New functionality

No.	Description
1	<p>The Preferences tool has been removed from Biacore X100 Control Software and Biacore X100 Evaluation Software. The following database solutions are used instead:</p> <ul style="list-style-type: none">• users are deleted from the Quick Filter tab• user roles are defined through Windows Computer Management• backup routines are managed through the Biacore X100 Backup and Restore application• database restoration is done from the Biacore X100 Backup and Restore application

Identical functionality with v2.0.3

No	Description
1	Manual backup and restore cannot be started while a scheduled backup is ongoing.
2	If the installation is done by user other than a Windows administrator, a dialog appears, prompting the user to start the installation as a Windows administrator.

Corrected issues

There are no issues corrected in this version.

Important information

No	Description
1	<p>Shutting down the computer: The database files may be corrupted if the Microsoft database software is not shut down properly, e.g. if the computer is turned off by switching off the power. If the files become corrupt, it will not be possible to run the software.</p> <p>To avoid corrupt files, always make sure to turn off the computer by clicking on the Shut down button in the operating system. The database files can be repaired. If the problem occurs, contact your local service representative for more information.</p>

No	Description
2	<p>The backup files are saved locally. It is the responsibility of the users to follow their local backup policies and save the files off-site if required.</p> <p>Note: <i>The folder where the backup files are saved is hidden by default. Hidden items can be made visible from settings in Windows Explorer.</i></p>

Page intentionally left blank



Give feedback on this document

Visit cytiva.com/techdocfeedback or scan the QR code.



cytiva.com

Cytiva and the Drop logo are trademarks of Life Sciences IP Holdings Corp. or an affiliate doing business as Cytiva.

Biacore is a trademark of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

Active Directory, Microsoft, SQL Server and Windows are registered trademarks of Microsoft group of companies.

Oracle is a registered trademark of Oracle or its affiliates..

Any other third-party trademarks are the property of their respective owners.

© 2020–2023 Cytiva

Any use of software may be subject to one or more end user license agreements, a copy of, or notice of which, are available on request.

For local office contact information, visit cytiva.com/contact

28930042 AK V:7 04/2023