

Biacore X100 Plus Package Software v2.1

Release Notes

Introduction

Scope of the document

This document describes added and changed functionality of Biacore™ X100 Plus Package Software v2.1 compared to Biacore X100 Plus Package Software v2.0.3. Known remaining software issue is also described.

No user functionality has been added or changed for Biacore X100 Plus Package Software v2.1 compared to v2.0.3. For changes to Biacore X100 Software v2.1, see the *Biacore X100 Software v2.1 Release Notes*.

The *Release Notes* is also accessible on the web using the product key. See [cytiva.com](https://www.cytiva.com) for the latest information.

Software included

Biacore X100 Plus Package Software v2.1.

System requirements

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

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

Upgrading from previous software versions

Before Biacore X100 Plus Package Software v2.1 is installed, Biacore X100 Control Software and Evaluation Software version 2.1 needs to be installed. Read the *Biacore X100 Software v2.1 Release Notes* for further details.

Functionality

The optional Biacore X100 Plus Package is a combined hardware and software package that provides enhanced functionality and flexibility.

There is no new functionality, and known remaining issues are listed in the section below.

Biacore X100 Plus Package Software v2.1

Functionality

See the table below for a brief description of functionalities. For detailed information refer to *Biacore X100 Handbook (BR100810)*.

Functionality	Description
Solvent correction	Analysis of small molecules requiring presence of dimethyl sulfoxide (DMSO) for aqueous solubility is supported through a number of pre-defined wizard templates that are used with the Custom Assay Wizard (see below).
Concentration analysis	Analyte concentration related to specific binding in samples can be determined either with the help of a calibration curve using known concentrations, or using the method Calibration-free concentration analysis (CFCA), which determines concentration without the need for calibration samples.
Variable analysis temperature control	The analysis temperature can be set between 4°C to 40°C.
Custom assay wizard	Offers flexibility in instrument control and provides support for applications that cannot be handled conveniently with the other wizards. Templates from application-specific wizards can be opened in the Custom assay wizard to provide a starting point for development of customized applications.
Custom immobilization	Supports custom immobilization methods, which can be used if none of the predefined immobilization methods is suitable for the application.
User-defined fitting models	New kinetics/affinity models can be created.

New functionality

There is no new functionality in this version.

Changed functionality

No functionality has changed compared to v2.0.3.

Remaining issue

No	Description
1	Chi2 is undefined for four-parameter fits of concentration calibration curves with four data points. Despite this, the Evaluation software reports Chi2 to be 0 in this case instead of N/A which would be the correct value.

**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.

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

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

28930054 AH V:8 04/2023