

Octet Software version 4.0

Product Overview

Software Upgrade

The latest software upgrade from FortéBio provides minor bug fixes in both the assay setup and data analysis portions of the FortéBio Octet Software. The bug fixes were implemented to provide the user with increased stability and productivity while using the Octet System.

Octet User Manual

There are no changes to the Octet System User Manual, *part no. 41-0006 Rev. C* in support of Octet Software version 4.0.

Bug Fixes from Version 3.2 to 4.0

The following Octet Software bug fixes were implemented in Octet Software version 4.0:

GENERAL SETUP, RESULTS AND DEFAULT OPTIONS

- Function to disable call-out boxes in data analysis
- Units are displayed in results tables

KINETICS ASSAY SET-UP AND DATA ANALYSIS

- Global Fitting
 - Improved association fitting to eliminate sporadic negative k_a
 - Improved dissociation fitting
- Import sample information button re-implemented

QUANTITATION ASSAY SET-UP AND DATA ANALYSIS

- Anti-Murine IgG Assay included in assay selection
- Programmed units selection is preserved in the method file
- Default options of well data for unknowns or standards reverts to "by value"
- De-selection is functional when analyzing both standard curve data and imported standard curves
- De-selection of single points is functional when analyzing an imported standard curve
- Improved error messages when samples are out of range instead of "NaN" (not a number)
- Inconsistencies between saved method file concentration units and data analysis are eliminated
- Improved creation of new assays under the Assay Development Mode

Known Issues in the Octet 4.0 Software

There are known issues in Octet Software 4.0 that we have provided a work-around solution.

REFERENCE SUBTRACTION FOR 2 DIFFERENT DATA SETS DURING DATA ANALYSIS IN THE KINETICS MODE

During data analysis, reference subtraction for 2 unmatched data sets will result in an intended error message. However upon closing the error message, the user can still proceed with data analysis with improperly subtracted data.

Workaround: It is recommended when comparing multiple data sets that require reference subtraction that they match in timing for the steps being analyzed.

STANDARD CURVE IN QUANTITATION GETS DRAWN INCORRECTLY AFTER MANIPULATIONS OF THE DATA SET

A standard curve that contains data files that are removed and then re-applied will not get refreshed properly while calculating unknowns. The user will see the standard curve with all data files during the calculation of the unknowns. *Although the screen does not get refreshed during the calculation of unknowns, the calculations will not be affected and are of high confidence.*

Workaround: If the standard curve is not displayed the same way as in the standard curve tab when the calculation of the unknowns are finished, then click on the recalculate button at the top of the data analysis screen to properly refresh the visualization of the data.

LOW CONFIDENCE DATA IN QUANTITATION

During data analysis of quantitation data, the user may see an asterisk (*) for some of the calculated binding rates. This will occur when the binding rate is below the programmed threshold to be reliably calculated.

CREATING NEW METHODS WHILE AN EXPERIMENT IS RUNNING

Refrain from setting up a new method with a different units selection while an experiment is running. This will alter the unit selection for the ongoing experiment.

continued on next page

NEW Octet System: Knowledge Base is now live
In ForteBio's continuing effort to provide a high level of service and support, we introduce the launch of the online Knowledge Base for the Octet System and its consumables located at <http://www.fortebio.com/pages/support.html>

Content

- Searchable user manuals
- Website material
- Downloadable documents only available to Octet Users
 - Product Inserts
 - Technical Notes
 - Product Bulletins
 - Product Listings
 - MSDSs
- FAQs
- Data analysis troubleshooting

Access to the Octet System Knowledge Base and support library requires a one time product registration.

Technical and Product Support

Contact support for more information and support of any technical issues.

Email: support@fortebio.com
Phone: 888.OCTET-QK
650-322-1360

NEW PRODUCTS

Part No.	UOM	Item/ Description
18-5042	Tray	Streptavidin High Binding Biosensors <i>Streptavidin coated biosensors with approx. 3x more binding capacity – Kinetics grade for baselines up to 1 hour</i>
18-5043	Pack	Streptavidin High Binding Biosensors <i>Streptavidin coated biosensors with approx. 3x more binding capacity – Kinetics grade for baselines up to 1 hour</i>
18-5044	Case	Streptavidin High Binding FA Biosensors <i>Streptavidin coated biosensors with approx. 3x more binding capacity – Kinetics grade for baselines up to 1 hour</i>
18-5045	Bottle	Kinetics grade buffer for sample dilution and biosensor hydration, 50 ml, packaged (10X).
18-5016	Each	Biosensor Dispenser <i>Octet biosensor transfer tool</i>

For more information about FortéBio's Octet System for label-free, real-time detection, applications and services, visit our website at www.fortebio.com or contact us directly.