



Octet Software v7.0 for Data Acquisition and Data Analysis

Ease of Use, Fast Assay Setup and Powerful Data Analysis

Key Features

- Acquisition “Wizards” for easy setup
- Pre-defined label-free assay protocols
- Multiple binding models for curve fitting
- Affinity, kinetics and concentration analysis
- FDA 21 CFR Part 11 tools
- Screening capabilities for proteins and small molecules
- Robotic integration (acquisition and analysis)
- Customizable data reports

FortéBio’s Octet software provides an intuitive and easy to use interface for data acquisition and analysis on all Octet instruments, enabling label-free kinetic, affinity, concentration and screening capabilities. An “Experiment Wizard” guides the operator through a four-step data acquisition process that facilitates both assay setup and efficient experimental design. Sample and plate maps are represented by interactive graphics that allow rapid plate definition through user-friendly click and drag mouse movements, while sample wells, biosensors and regeneration steps are descriptively color-coded.

During acquisition, data are displayed in real time with a detailed graphical interface that depicts experimental progress. Interactive zoom tools and advanced legends enable rapid association of important elements of a dataset moments after the data is acquired.

Post-acquisition, a clutter-free environment enables a focused and efficient analysis of the data set through data grouping, auto-alignment and a multi-level search tool. Multiple analyses can be applied to raw data, including single, double and parallel reference subtraction. Curve fitting can be highly customized by both assay step and binding model, and, additionally, any two numerical acquisition or processing parameters can be plotted in an interactive X-Y map.

Through an interface that makes no assumptions about communication medium or language, Octet instruments can be fully integrated with third-party robotic software for automat-

ed data acquisition and analysis. Optional FDA 21 CFR Part 11 GxP tools are also available.

Octet software continues to rapidly evolve to address the needs of FortéBio customers. Version 7.0 enhances workflow with improved data viewing, simplified navigation of large data sets and the ability to remotely monitor experiment on the Octet platform. To upgrade, contact your local sales representative or e-mail sales@fortebio.com.



OCTET SOFTWARE FEATURE COMPARISON — DATA ACQUISITION

Feature	v6.1–6.4	v7
21 CFR Part 11 tools available	✓*	✓
Pre-defined templates	✓	✓
Automation compatible (Octet 384 instruments)	✓	✓
Real-time reference subtraction during data acquisition	✓	✓
Threshold option to control loading density	✓	✓
Individual control of sample and reagent plates (Octet 384 instruments)	✓	✓
Review Experiment tab available to “dry run” experiments	✓	✓
Multi-analyte data acquisition		✓
Remote monitoring		✓
Replicate groups		✓
“Flip” quantitation data		✓

*Available for software versions 6.1,6.2 and 6.3 only.

NEW DATA ACQUISITION FEATURES IN VERSION 7.0

Replicate Groups

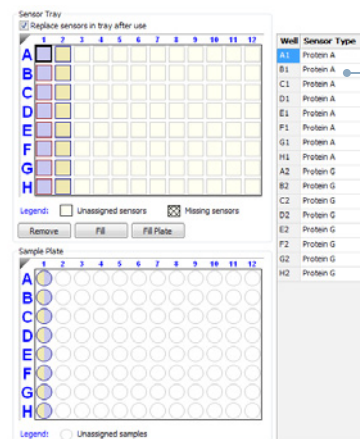
Organize replicate samples into groups for convenient calculation of average concentration, average binding rate, standard deviation and CV%. Statistics are automatically calculated during data analysis and can be plotted in the new X-Y graphing tool for custom viewing.

Enter replicate group names during acquisition for automated calculation of statistics during analysis

Well	Sample ID	Replicate Group	Type	Conc (µg/ml)	Dilution Factor	Information
A4	Ab1	Ab1	Unknown	n/a	2	Sample Diluent
A5	Ab1	Ab1	Unknown	n/a	2	Sample Diluent
A6	Ab1	Ab1	Unknown	n/a	2	Sample Diluent
B4	Ab2	Ab2	Unknown	n/a	2	Sample Diluent
B5	Ab2	Ab2	Unknown	n/a	2	Sample Diluent
B6	Ab2	Ab2	Unknown	n/a	2	Sample Diluent
C4	Ab3	Ab3	Unknown	n/a	2	Sample Diluent
C5	Ab3	Ab3	Unknown	n/a	2	Sample Diluent
C6	Ab3	Ab3	Unknown	n/a	2	Sample Diluent
D4	Ab4	Ab4	Unknown	n/a	2	Sample Diluent
D5	Ab4	Ab4	Unknown	n/a	2	Sample Diluent
D6	Ab4	Ab4	Unknown	n/a	2	Sample Diluent
E4	Ab5	Ab5	Unknown	n/a	2	Sample Diluent
E5	Ab5	Ab5	Unknown	n/a	2	Sample Diluent
E6	Ab5	Ab5	Unknown	n/a	2	Sample Diluent

Multi-analyte Data Acquisition

Assign multiple biosensor types during a quantitation experiment for a new level of versatility and increase the walk-away time of experiments. Screen multiple biosensor chemistries against a single analyte for biosensor selection (i.e. Protein A, Protein G and Protein L biosensors against one antibody). Alternatively, use different biosensor chemistries to serially multiplex the quantitation of an array of analytes.

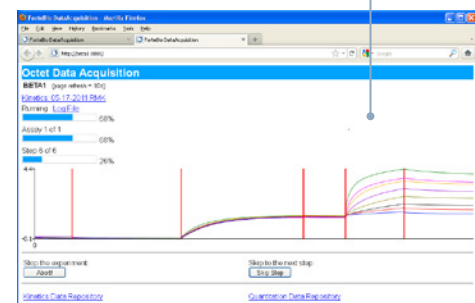


Use different biosensor chemistries (Protein A and Protein G shown) to multiplex quantitation experiments

Remotely monitor real-time data collection through a standard web browser — even on a smartphone

Remote Monitoring

Want to observe your data in real time, but unable to stay in front of the Octet instrument? Monitor experiments remotely from any networked computer, smartphone or mobile device using a standard web browser using remote monitoring.



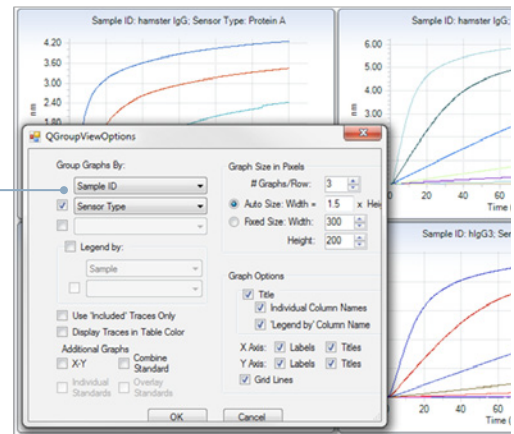
OCTET SOFTWARE FEATURE COMPARISON — QUANTITATION DATA ANALYSIS

Feature	v6.1–6.4	v7
Import/export standard curves	✓	✓
Import kinetic experiments for quantitation analysis	✓	✓
Thresholds for R ² and residual data flagging	✓	✓
Save analysis settings	✓	✓
Automation and batch processing of multiple data sets	✓	✓
Save processing and analysis settings	✓	✓
Advanced legend display	✓	✓
Grouped view		✓
X-Y graph		✓
Accelerated binding calculation		✓
“Flip” quantitation data		✓
Sample alerts		✓

NEW QUANTITATION DATA ANALYSIS FEATURES IN VERSION 7.0

Grouped View (Quantitation)

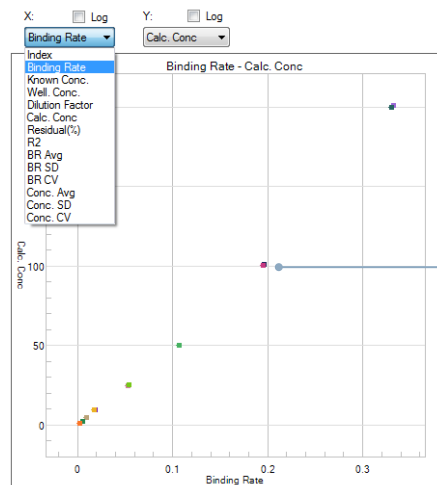
Simplify the visual analysis of datasets with highly tailored displays organized by groups. Groups can be multi-level with customizable dimensions and layouts, including options for coloring, legends and the display of standard curves.



Organize datasets into customizable groups for rapid visual analysis and navigation

X-Y Graph

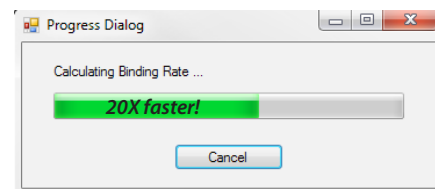
Analyze for trends and correlations between parameters by plotting any two numerical values from the analysis table in an X-Y Cartesian map. Toggle between linear and logarithmic scales, customize the dimensions with click-and-drag and copy to the clipboard for easy export.



Plot any two numerical values to analyze for data trends and correlations

Accelerated Binding Calculation

Process datasets in a fraction of the time and view final data sooner. An improved and validated algorithm provides the same quality quantitation data — just faster.



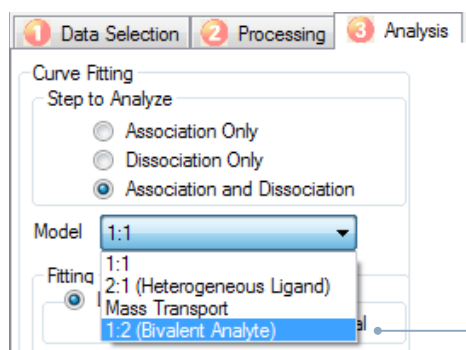
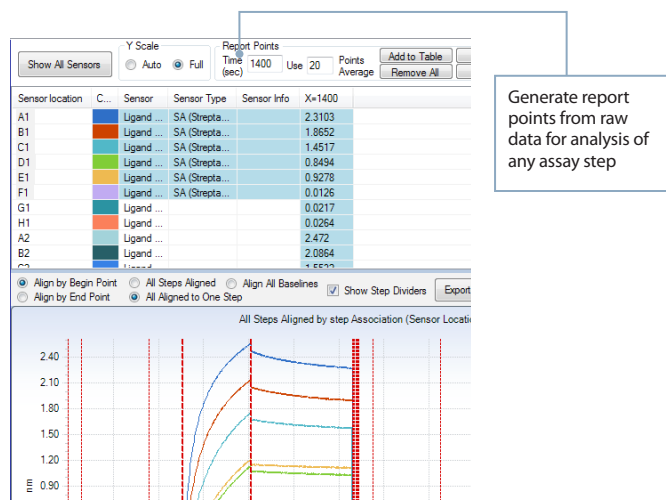
OCTET SOFTWARE FEATURE COMPARISON — KINETICS DATA ANALYSIS

Feature	v6.1–6.4	v7
Global and local fitting with 2:1 (HL) model	✓	✓
Grouped view visualization	✓	✓
Fit data with an unlinked R_{max}	✓	✓
Mass transport model	✓	✓
Customizable data reports	✓	✓
Advanced search function	✓	✓
Batch processing of multiple data sets	✓	✓
1:2 Bivalent analyte model		✓
Report points with raw data		✓
Save and load report points		✓

NEW KINETICS DATA ANALYSIS FEATURES IN VERSION 7.0

Report Points for Raw Data

Generate report points from raw data and analyze any assay step, such as baseline, loading or association within a concatenated experiment. Save the report point settings to a file and recall those settings during future experiments to streamline analysis.



1:2 Bivalent Analyte Model

Perform a full kinetic analysis of data for bivalent analytes interacting with an immobilized ligand. The bivalent model provides terms for two association constants and two dissociation constants.

Perform full kinetic analyses of bivalent analytes with the 1:2 binding model