



Octet software comparison

ForteBio Octet Data Acquisition and Data Analysis software is rapidly evolving to better address the growing applications and needs of our customers. These comparison tables provide a general guide to the features included in various versions of the software. The latest release, Version 10, is available for use on all Octet systems. **Note:** Octet Data Analysis HT software is not available in the Octet CFR software suite installer. For more information or for upgrade options, please contact your local sales representative or fill out our [software download request form](#).

Octet Data Acquisition software version feature comparison

Feature	V8.x	V9.x	V10.x
Experiment Wizard includes built-in protocol templates for kinetics, quantitation and epitope binning		✓	✓
Capture Antibody step type in Advanced Quantitation for easier conversion of ELISA to Octet assays		✓	✓
Visualization of all assay steps in advanced quantitation module	✓	✓	✓
Add multiple assay steps at one time in "Step Data List" in kinetics for quick assay setup	✓	✓	✓
In kinetics, all assay steps in multiple assays within an experiment can be displayed in a grid for better visualization	✓	✓	✓
Remote monitoring during experimental run	✓	✓	✓
Support for the Octet RED96e system			✓
Support for the Biosensor Mount Cleaning Tool for Octet 384 and HTX systems			✓
Support for microplate evaporation cover consumable on Octet RED96e system only			✓

Octet Data Analysis software version feature comparison

Feature	V8.x	V9.x	V10.x
Quantitation			
Mask all proprietary information (sample ID, sensor info, etc.) from data		✓	✓
Editable cells, customize and sort Results table		✓	✓
Import standard curve of experiments run at the same temperature (Octet RED96e system only)			✓
Show temperature trace for the experiment			✓
Kinetics			
View all graphs in either logarithmic or linear scale	✓	✓	✓
Mask all proprietary information (sample ID, sensor info, etc.) from data		✓	✓
Editable cells, customize and sort Results table		✓	✓
Assign % KD1 and KD2 contribution to 2:1 model interaction		✓	✓
Show temperature trace for the experiment			✓

Octet Data Analysis HT software version feature comparison

Feature	V8.x	V9.x	V10.x
Epitope binning analysis			
Append or overlay data from multiple experiments, plates and biosensor trays		✓	✓
Automatically create 2D traffic light matrix		✓	✓
Highlight unidirectional binning pairs		✓	✓
Flag Ab 1 traces with low response or no binding in the 2D matrix		✓	✓
Subtract self-binding in the matrix			✓
Subtract the matrix against a select row or column			✓
Data preprocessing for kinetics and quantitation assays*			
Perform data processing for a combined mega dataset consisting of data from multiple experiments, plates and/or biosensor trays			✓
More flexible referencing options – by column, row, selected wells, pairs, pattern and between biosensor trays			✓
Additional referencing for quantitation assays – reference well and reference load well when no or irrelevant capture molecule is immobilized			✓
Advanced inter-step correction to remove bulk shifts or assay artifacts			✓
Report point feature for whole sensor traces – can be added in any step in the trace			✓
More graphical options – thicken line, change trace colors, customize x- and y-axis scales			✓
Processed data exported per sensor in a single Microsoft® Excel® file			✓
Kinetic analysis			
Enable kinetic analysis of multiple experiments and assays that can be visualized together in All Assays tab			✓
Custom y-axis scaling available for all graphs – stacked, individual and grouped			✓
More intuitive group by dropdown menu options under global fitting			✓
Faster and more optimized 1:1 and 2:1 algorithms			✓
More graphical options – thicken line, tooltip, legends, etc.			✓
Quantitation analysis			
One step or one step type (when regeneration is performed) can be quantitated from a kinetic assay			✓
Enables quantitation analysis of multiple experiments and plates by forming a single, combined mega dataset			✓
Standard curve sorted automatically by sensor type if multiple types of sensors are used in an experiment			✓
Import standard curve for experiments run at the same temperature (Octet RED96e system only)			✓
More advanced sample alerts for data quality			✓
Custom report generation			
Add experiment summary, sensor tray layout, sample plate layout, graphs and tables from any analysis module – quantitation, kinetics, preprocessing or epitope binning			✓
Add text boxes and any image(s)			✓
Format report by adding sections, headers, footers, gridlines, border, page size, etc.			✓
Export the report in PDF format or print a hard copy			✓
Save report template and load to generate similar reports for other datasets			✓

*Most features from Octet Data Analysis software plus additional ones shown.

Feature	V8.x	V9.x	V10.x
Automated analysis of datasets			
Automated quantitation analysis of multiple plates		✓	✓
Automated kinetic analysis of multiple plates			✓
Automated epitope binning analysis of multiple plates			✓



ForteBio
47661 Fremont Boulevard
Fremont, CA 94538
888.OCTET-75 or 650.322.1360
fortebio.info@moldev.com

ForteBio Analytics (Shanghai) Co., Ltd.
No. 88 Shang Ke Road
Zhangjiang Hi-tech Park
Shanghai, China 201210
salesops.china@moldev.com

Molecular Devices (UK) Ltd.
660-665 Eskdale
Winnersh Triangle
Wokingham, Berkshire
RG41 5TS, United Kingdom
+44 118 944 8000
uk@moldev.com

Molecular Devices (Germany) GmbH
Bismarckring 39
88400 Biberach an der Riss
Germany
+ 00800 665 32860