



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 12, is available for use on all current Octet systems. 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	V10.x	V11.x	V12.x
Increased flexibility in Kinetic method generation when entering Dilution Factors and designating Controls			✓
Increased flexibility in Advanced Quantitation method generation – more step types and reuse step types		✓	✓
Customize default settings - plate type, plate definition (row, columns or quadrants) and data input		✓	✓
Support for new Sialic Acid glycoanalysis biosensor (GlyS)		✓	✓
Shortcuts to minimize number of clicks		✓	✓
Support for the Biosensor Mount Cleaning Tool for Octet 384 and HTX systems	✓	✓	✓
Support for Octet RED96e system and its microplate evaporation cover consumable	✓	✓	✓
Experiment Wizard includes built-in protocol templates for kinetics, quantitation and epitope binning	✓	✓	✓

Octet Data Analysis HT software version feature comparison

Feature	V10.x	V11.x	V12.x
Automated analysis of datasets			
Available as a 64-bit application			✓
Ability to sort multiple fields in data tables			✓
Pre-defined analysis settings import to maximize speed		✓	✓
Automated kinetic analysis of multiple plates	✓	✓	✓
Automated epitope binning analysis of multiple plates	✓	✓	✓
Automated quantitation analysis of multiple plates	✓	✓	✓
Data preprocessing*			
Cross-plate referencing			✓
Epitope Binning: ability to exclude full sensor trace			✓
Epitope Binning: ability to export data to a csv file			✓
New buttons for reference assignment for better usability		✓	✓

Feature	V10.x	V11.x	V12.x
More editable columns for flexibility - loading sample ID, loading concentration, dilution factor		✓	✓
Perform data processing for a combined data from multiple experiments, plates and/or biosensor trays	✓	✓	✓
More flexible referencing options – by column, row, selected wells, pairs, pattern and between biosensor trays	✓	✓	✓
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	✓	✓	✓
Epitope binning analysis			
NEW BinChart for automated binning of antibodies			✓
NEW color schemes for Matrix display			✓
Ability to export BinChart, original and corrected matrix, table with clustering info			✓
Enable sample ID and information correction, if necessary		✓	✓
Copy epitope binning matrix image		✓	✓
Subtract self-binding in the matrix	✓	✓	✓
Subtract the matrix against a select row or column	✓	✓	✓
Append or overlay data from multiple experiments, plates and biosensor trays	✓	✓	✓
Automatically create 2D traffic light matrix	✓	✓	✓
Kinetic analysis			
Ability to use multiple parameters in global fitting groups			✓
Analyze kinetic cycles with different association/disassociation times in the same experiment file			✓
Batch export of group graphs to image files			✓
Pause auto-fitting of kinetic datasets option		✓	✓
Flip data option		✓	✓
Enable kinetic analysis of multiple experiments that can be visualized together in All Assays tab	✓	✓	✓
More graphical options – thicken line, tooltip, legends, custom y-scaling	✓	✓	✓
Quantitation analysis			
Overlay fitted traces on top of raw binding traces			✓
Exportable Fitted Graph that shows binding signal and curve fit for each sample well			✓
NEW Advanced option to choose the model classifier for binding rate determination			✓
Pause auto-fitting of quantitation datasets option		✓	✓
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	✓	✓	✓
More advanced sample alerts for data quality	✓	✓	✓

Feature	V10.x	V11.x	V12.x
Custom report generation			
Multi-page table support for large experiments			✓
Formatting and sorting capability for data tables			✓
Add residual, steady-state, iso-affinity, XY scatter and grouped graphs to report		✓	✓
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)	✓	✓	✓
Export the report in PDF format or print a hard copy	✓	✓	✓
Save report template and load to generate similar reports for other datasets	✓	✓	✓

21 CFR Part 11 support

Feature	V10.x	V11.x	V12.x
Ability to authorize users to remove signatures from an analysis file			✓
Ability to inactivate the default Administrator account			✓
More detailed audit log of changes made within ForteBio Server Monitor			✓
Ability to choose time zone and date/time format			✓
New Security tab in Octet Data Analysis HT software		✓	✓
Controlled access with multiple user levels and expanded privileges		✓	✓
Electronic signatures to lock dataset and prevent changes after analysis is complete		✓	✓
Digitally signed data files that are invalidated if any modification is made outside the software		✓	✓
Log all critical user actions in Data Acquisition and Data Analysis HT software Audit Trail with old versus new values		✓	✓
Audit Trail can be filtered by project, machine, user, experiment and date interval		✓	✓
Exporting Audit Trail and user information to PDF reports		✓	✓
Option to add customized comments to each analysis event in Audit Trail		✓	✓
Full control over running and analyzing routine assays in regulated labs		✓	✓
Lock all Microsoft® Excel® export reports		✓	✓
Software Validation Package product support		✓	✓

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



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