

Pall ForteBio Octet Data Acquisition and Data Analysis software is rapidly evolving to better address the growing applications and needs of our customers. The table below provides a general guide to the features included in various versions of the software. The latest release, Version 9, is available for use on Octet QK, QK^e, RED, RED96, QK384, RED384, HTX and the new Octet K2 system. For more information or for upgrade options, please contact your local sales representative or fill out our [software download request form](#).

Feature	Version 7.x	Version 8.x	Version 9.x
Data Acquisition			
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			✓
Latest offering of all biosensor types			✓
Visualization of all assay steps in advanced quantitation module		✓	✓
Regeneration steps in kinetics module, similar to quantitation		✓	✓
Multiple assay steps can be added at one time in "Step Data List" in kinetics for quick assay setup		✓	✓
All assay steps in multiple assays within an experiment can be displayed in a grid for better visualization		✓	✓
User selectable read-head (Octet HTX system only)		✓	✓
Automation compatible (Octet QK384, RED384 and HTX instruments only)	✓	✓	✓
Remote monitoring during experimental run	✓	✓	✓
Quantitation Data Analysis			
Edit cells in table such as known concentration (standard wells), dilution factor, replicate group and sample ID			✓
Mask all proprietary information (sample ID, biosensor info, etc.) in data			✓
View and analyze all assay steps in advanced quantitation		✓	✓
View all graphs in either logarithmic or linear scale		✓	✓
Multiple read-head configurations for Load and Detect assay steps (Octet HTX system only)		✓	✓
Interactive data visualization of plate map, results table and trace data		✓	✓
Spacebar shortcut key includes/excludes data in results table	✓		✓
Quantitation of association step from kinetic experiments	✓	✓	✓
Import/export multiple external standard curves to quantitate new datasets	✓	✓	✓
Kinetics Data Analysis			
New Octet Data Analysis HT Software for Epitope Binning Analysis			
Append data from multiple experiments, plates and biosensor trays (for example: Antibody 1 loading is performed separately from the rest of the epitope binning assay antigen and Antibody 2 associations)			✓
Overlay data from multiple experiments, plates and biosensor trays for analysis of large binning datasets (64x64, 96x96 and more)			✓
Automatically find binning cycle pattern, assign Ab 1 and Ab 2 and create 2-D traffic light matrix			✓
Highlight unidirectional binning pairs			✓
Flag Antibody 1 traces with low response or no binding in the 2-D matrix			✓
Traces with low response or high off-rates are highlighted with flag icons			✓
Octet System Data Analysis Software			
Mask all proprietary information (sample ID, biosensor info, etc.) in data			✓
Assign % K_{D1} and K_{D2} contribution to 2:1 model interaction			✓
Transfer reference sensor(s) and reference well(s) info from Data Acquisition to Data Analysis software			✓
Expanded color palette: up to 144 colors for biosensor traces			✓
Edit cells, customize and sort results table			✓
Cursors to add report points to kinetics data		✓	✓
Two new cursors to add multiple report points to processed data		✓	✓
Multiple read-head configurations within an experiment (Octet HTX system only)		✓	✓
View all graphs in either logarithmic or linear scale		✓	✓
Spacebar shortcut key includes/excludes data in results table	✓		✓
Order Columns option enables reordering of columns from the default order in saved report	✓		✓