

BLI Selection Guide

	BLitz System	Octet QK [®] System	Octet QK384 System	Octet RED96 System	Octet RED384 System	Octet HTX System
	Simple-to-use single-sample protein detection	Automated large molecule characterization	High throughput large molecule characterization	Automated small and large molecule characterization	High throughput small and large molecule characterization	Highest throughput small and large molecule characterization
Ideal Uses						
Small Molecule Assays				● ●	● ●	● ● ●
VLP Assays	●	● ●	● ●	● ● ●	● ● ●	● ● ●
Screening Applications	●	● ●	● ● ●	● ●	● ● ●	● ● ●
Measuring Weak Binding Affinities	●	● ●	● ●	● ● ●	● ● ●	● ● ●
Measuring Strong Binding Affinities	●	● ● ●	● ● ●	● ● ●	● ● ●	● ● ●
Multi-step Quantitation Assays	●	● ●	● ●	● ● ●	● ● ●	● ● ●
Low-volume Sample Measurements	● ● ●	●	● ●	●	● ●	● ● ●
At-line Bioprocess Monitoring	● ● ●	● ●	●	● ●	●	●
Quantitation in Media and Lysates	● ● ●	● ● ●	● ● ●	● ● ●	● ● ●	● ● ●
Determining Antibody Titer	● ● ●	● ● ●	● ● ●	● ● ●	● ● ●	● ● ●
Immunoassay Development and QC	● ● ●	● ● ●	● ● ●	● ● ●	● ● ●	● ● ●
Performance						
Affinity Range	1 μM – 0.1 nM	0.1 mM – 0.1 nM	0.1 mM – 0.1 nM	1 mM – 10 pm	1 mM – 10 pM	1 mM – 10 pM
Molecular Weight	> 10,000 Da	>5,000 Da	>5,000 Da	>150 Da	>150 Da	>150 Da
Minimum Sample Volume	4 μL	100 μL†	40 μL*	200 μL	40 μL*	40 μL*
Acquisition Rate	5 Hz	0.3, 0.6 Hz	0.3, 0.6 Hz	2, 5, 10 Hz	2, 5, 10 Hz	2, 5, 10 Hz
Simultaneous Reads	1	8	16	8	16	96
Specifications						
Number of Spectrometers	1	1	2	8	16	16
Temperature Control	No	Ambient +4°C – 40°C	Ambient +4°C – 40°C	Ambient +4°C – 40°C	Ambient +4°C – 40°C	Ambient +4°C – 40°C
Microplate Positions	NA	1 (96-well)	2 (96- or 384-well)	1 (96-well)	2 (96- or 384-well)	2 (96- or 384-well)
Integration with Automation	No	No	Yes	No	Yes	Yes
Dimensions						
Size H x W x D (cm)	17.4 x 15.3 x 22	47 x 43 x 53	77 x 80 x 80	47 x 43 x 53	77 x 80 x 80	77 x 80 x 80
Weight (Kg)	3.3	23	68.2	24.5	68.2	90

* in a 384-well tilted-bottom microplate (ForteBio, part no.18-5080)

† in a 96-well half-area microplate (VWR, part no. 82050-044)