

SARTORIUS



Powered and Prepared

Octet® SF3 SPR

Powering Opportunity

A robust, high-throughput, low maintenance SPR solution

- State-of-the-art microfluidics and optimized electronic designs
- Complete kinetics and affinity for up to 768 samples in a single unattended assay
- Eliminate the need to prepare multiple dilution series using OneStep® Injection Technology
- Determine full kinetics and affinity in the presence of multiple competitors

Octet® SF3 Ordering

Description	Product Number
Octet® SF3 SPR System	Octet-SF3
Octet® SF3 SPR System Installation	84IAFF3

Octet® SF3 Specifications

Baseline Properties	
Refractive Index Range	1.33 - 1.40
Baseline Noise	Typically < 0.025 RU (RMS)
Baseline Drift	Typically < 0.3 RU/min
Molecular Weight Cut-Off	No lower limit for organic molecules
Variable Data Rate	1, 2, 5, 10, 20 and 40 Hz
Affinity Range	fM to mM
Working Ranges	
Association Rate Constant (k_a)	$10^2 - 10^9 \text{ M}^{-1} \text{ s}^{-1}$
Dissociation Rate Constant (k_d)	$10^{-6} - 2.5 \text{ s}^{-1}$
Global Affinity Constant (K_D)	$10^{-3} - 10^{-12} \text{ M}$
Concentration	Sample concentration >1 pM
Sample Acquisition	
Unattended Run Time	>72 hours with no built in runtime limit
Sample Capacity	Any 2 sample racks plus 2 reagent racks
Sample Rack Options	96 vial, deep well and PCR formats, 384-well microplates, custom high volume
System and Sample Temperature Control	4 to 40°C (max 15° below ambient)
Buffer Line Selection	Automatic switching between 3 independent lines
Inline Buffer Degasser	Yes
Flow Cell and Injections	
Number of Flow Channels	3
Flow Path	1, 1-2, 1-2-3, 3, 3-2, 3-2-1
Flow Channel Volume	<90 nL
Channel-Channel Dead Volume	<20 nL
Injection Volume	2-700 µL
Flow Rate	0.1-200 µL/min
Injection Rise and Fall Time	<0.75 second @ 25 µL/min
Gradient Injections	OneStep® (including in High-Throughput Mode), NeXtStep™
Simultaneous Injections	Yes
Software	
Octet® SPR Analysis Hit Selection	Yes
Dimensions	
Width x Height x Depth	24" x 24" x 20" (61 cm x 61 cm x 51 cm)
Weight	66 kg