SARTURIUS



Powered and Prepared

Octet® SPR Sensor Chips and Consumables

Accurate, Robust and Reproducible

SPR Sensor Chips and Consumables for every application

- Highly accurate and robust SPR sensor chips for all applications, ranging from small molecules to whole cells
- Extensive range of SPR consumables to optimize assay reproducibility and throughput

Octet® SPR Consumables

Octet® SF3 SPR Consumables	Product Number
Octet® SPR Glassware	
Octet® SPR Vial, 2 mL (100 vials)	19-0140
Octet® SPR Vial, 0.9 mL (100 vials)	19-0141
Octet® SPR Vial, 7.5 mL (3 vials)	19-0142
Octet® SPR Vial, 7.5 mL (1 vial)	19-0143
Octet® SPR Sample Rack and Sample Rack Plate Adapters	
Octet® SPR Sample Rack, 96 well, 0.9 mL (pack of 1)	19-0145
Octet® SPR Sample Rack Plate Adaptor, Axygen 96 deep well, 1 mL (pack of 1)	27-0448
Octet® SPR Sample Rack Plate Adaptor, Axygen 96 well, 450 μL (pack of 1)	27-0449
Octet® SPR Sample Rack Plate Adaptor, Axygen 96 well PCR (pack of 1)	27-0450
Octet® SPR Sample Rack Plate Adaptor, Axygen 384 well, 240 μL (pack of 1)	27-0451
Octet® SPR Sample Rack Plate Adaptor, Greiner 384 well, 120 μL (pack of 1)	27-0453
Octet® SPR Sample Rack Plate Adaptor, Greiner 384 deep well, 225 μL (pack of 1)	27-0454
Octet® SPR Sample Rack Plate Adaptor, Greiner 96 well, 280 μL (pack of 1)	27-0455
Octet® SPR Sample Rack Plate Adaptor, Greiner 96 well, 300 µL (pack of 1)	27-0456
Octet® SPR Microfoils	
Octet® SPR Plate Seal, 96 well PCR (pack of 5)	19-0102
Octet® SPR Sample Rack Seal, 96 well (pack of 5)	19-0103
Octet® SPR Plate Seal, 384 well (pack of 5)	19-0104
Octet® SPR Plate Seal, Varied (pack of 5)	19-0105
Octet® SPR Rear Reagent Rack Seal (pack of 5)	19-0107
Octet® SPR Front Reagent Rack Seal (pack of 5)	19-0108
Octet® SPR Plate Seal, 384 square well (pack of 20)	19-0110
Octet® SPR Plate Seal, 96 well (pack of 20)	19-0111
Octet® SPR Maintenance and Performance	
Octet® SPR Clean Station Cover (pack of 1)	19-0106
Octet® SPR Buffer Filter Replacement Cartridges, 10 µm (pack of 2)	19-0109
Octet® SPR Sensor Cleaning Cards (pack of 6)	19-0114
Octet® SPR Maintenance Chip (pack of 1)	19-0136
Octet® SPR Maintenance Kit (6 months)	19-0137

Octet® SPR Sensor Chips

Octet® SF3 SPR Sensor Chips	Product Number
Octet® SPR COOH1 Sensor Chip (pack of 3)	19-0053
Octet® SPR CDL Sensor Chip (pack of 3)	19-0127
Octet® SPR CDH Sensor Chip (pack of 3)	19-0128
Octet® SPR PCH Sensor Chip (pack of 3)	19-0129
Octet® SPR HisCap Sensor Chip (pack of 3)	19-0058
Octet® SPR SADH Sensor Chip (pack of 3)	19-0130