

SARTORIUS



Powered and Prepared

Octet® SF3 SPR

Maintenance Kits
and Sensor Chips

**Octet® SPR maintenance solutions
for maximizing instrument run-time
and optimizing data generation**

- Octet® SPR Maintenance kit for instrument cleaning, desorb and decontaminate protocols
- Octet® SPR Maintenance Sensor Chip for use during all cleaning protocols

Octet® SPR Maintenance Solutions and Sensor Chips

Regular maintenance of the Octet® SF3 is critical in ensuring consistent and reliable results. Day to day use can result in the adsorption of material to the microfluidic tubing and potential microbial growth within the system, which may impact instrument performance.

The Octet® SPR Maintenance Kit has been developed to be used during each weekly instrument desorb protocol, as well as during each monthly desorb and decontaminate program, to remove any absorbed proteins or other materials from the system and ensure long-term instrument performance.

In addition, a new inert Octet® SPR Maintenance Chip has been developed for use during all predefined Octet® SF3 cleaning protocols within the Octet® SPR Discovery Software, to ensure that your system is maintained in excellent working order.

The combined use of both the Octet® SPR Maintenance Kit and Sensor Chip will help maximize instrument run time, provide total confidence in all results generated, and prolong the overall longevity of the instrument.

Octet® SPR Maintenance Chip

Pack of 1
Product Number: **19-0136**

Fully inert glass sensor chip with no associated surface chemistry. For use in all Octet® SF3 SPR maintenance and cleaning protocols.



Octet® SPR Maintenance Kit

Sufficient for 6 months standard instrument use
Product Number: **19-0137**

Contains

- 250 mL Octet® SPR 0.5% SDS
- 250 mL Octet® SPR 50 mM Glycine pH 9.5
- 1 x Octet® SPR Maintenance Chip

For use with the Octet® SF3 platform. Packaged in polythene bags.

