

Biosensors & Reagents

Item	Part No.	UOM	Qty.	Description
Anti-human IgG Biosensors (FC Specific)	18-5001	Tray	1	Tray of 96 biosensors coated with anti-human IgG (Includes 1 bottle of diluent.)
Anti-human IgG Biosensors (FC Specific)	18-5004	Pack	1	Five trays of 96 biosensors coated with anti-human IgG (Includes 5 bottles of diluent.)
Anti-human IgG Biosensors (FC Specific)	18-5005	Case	1	Twenty trays of 96 biosensors coated with anti-human IgG (Includes 20 bottles of diluent.)
Protein A Biosensors	18-5010	Tray	1	Tray of 96 biosensors coated Protein A (Includes 1 bottle of diluent.)
Protein A Biosensors	18-5012	Pack	1	Five trays of 96 biosensors coated with Protein A (Includes 5 bottles of diluent.)
Protein A Biosensors	18-5013	Case	1	Twenty trays of 96 biosensors coated with Protein A (Includes 20 bottles of diluent.)
Streptavidin FA Biosensors	18-5003	Tray	1	One tray of 96 biosensors coated with streptavidin for kinetic screening. Off rate ≤ 15 minutes. (Includes 2 bottles of diluent.)
Streptavidin FA Biosensors	18-5008	Pack	1	Five trays of 96 biosensors coated with streptavidin for kinetic screening. Off rate ≤ 15 minutes. (Includes 10 bottles of diluent.)
Streptavidin FA Biosensors	18-5009	Case	1	Twenty trays of 96 biosensors coated with streptavidin for kinetic screening. Off rate ≤ 15 minutes. (Includes 40 bottles of diluent.)
Streptavidin SBC Biosensors	18-5002	Tray	1	One tray of 32 biosensors coated with streptavidin for standard kinetic analysis. (Includes 1 bottle of diluent.)
Streptavidin SBC Biosensors	18-5006	Pack	1	Five trays of 32 biosensors coated with streptavidin for standard kinetic analysis. (Includes 5 bottles of diluent.)
Streptavidin SBC Biosensors	18-5007	Case	1	Twenty trays of 32 biosensors coated with streptavidin for standard kinetic analysis. (Includes 20 bottles of diluent.)
Amine Reactive Biosensors, Standard	18-5011	Tray	1	One tray of 32 biosensors compatible with amine reactive coupling chemistry for kinetic analysis. (Reagents not included)
Amine Reactive Biosensors, Standard	18-5014	Pack	1	Five trays of 32 biosensors compatible with amine reactive coupling chemistry for kinetic analysis. (Reagents not included)
Amine Reactive	18-5015	Case	1	Twenty trays of 32 biosensors compatible with amine reactive coupling chemistry for kinetic

Biosensors, Standard				analysis. (Reagents not included)
----------------------	--	--	--	---

Biosensors & Reagents, continued

Item	Part No.	UOM	Qty.	Description
Sample Diluent	18-1000	Bottle	1	Diluent for sample dilution and biosensor pre-wetting, 50 ml
IgG Calibrators	18-1030	Set	1	Set of 8 IgG Calibrators filled at 10 mL/vial. (plus a negative control)

Instruments

Item	Part No.	Description
Octet Q System	30-5002	Octet, Quantitation only instrument for label-free, real-time analysis. Includes computer, monitor and accessory kit.
Octet QK System	30-5003	Octet, Quantitation and kinetics instrument for label-free, real-time analysis. Includes computer, monitor and accessory kit.

Software

Item	Part No.	Qty.	Description
Seat License - Q Analysis, Single	50-5001	1	Single Seat License for offline Q Analysis
Seat License - Q Analysis, Double	50-5002	2	2 Seat License for offline Q Analysis
Seat License - QK Analysis, Single	50-5003	1	Single Seat License for offline QK Analysis
Seat License - QK Analysis, Double	50-5004	2	2 Seat License for offline QK Analysis

Ordering

To order biosensor consumables, a purchase order referencing a quote number should be faxed or emailed to the following:

Email: ordering@fortebio.com

eFax: 866-327-9115

Outside of the US:

Fax: 650-322-1370

Please supply the following information on your purchase order:

- Full delivery address with contact name and telephone number
- Invoice address
- Purchase order number
- ForteBio quote #
- Fax number or e-mail address for receiving confirmation



Product Listing

Limitations on Use

The Octet System and its biosensor consumables are sold for Research Use Only.