

## OCTET BIOSENSORS QUANTITATION GRADE

Grade	Item	Part No.	UOM	Qty.	Description
Quant	Anti-human IgG Biosensors (FC Specific)	18-5001	Tray	1	Tray of 96 biosensors coated with anti-human IgG (Includes 1 bottle of diluent.)
Quant	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.)
Quant	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.)
Quant	Protein A Biosensors	18-5010	Tray	1	Tray of 96 biosensors coated Protein A (Includes 1 bottle of diluent.)
Quant	Protein A Biosensors	18-5012	Pack	1	Five trays of 96 biosensors coated with Protein A (Includes 5 bottles of diluent.)
Quant	Protein A Biosensors	18-5013	Case	1	Twenty trays of 96 biosensors coated with Protein A (Includes 20 bottles of diluent.)
Quant	Anti-Murine IgG Biosensors	18-5022	Tray	1	Tray of 96 biosensors coated with goat anti-mouse IgG antibody (Includes 1 bottle of diluent.)
Quant	Anti-Murine IgG Biosensors	18-5023	Pack	1	Five trays of 96 biosensors coated with goat anti-mouse IgG antibody (Includes 5 bottles of diluent.)
Quant	Anti-Murine IgG Biosensors	18-5024	Case	1	Twenty trays of 96 biosensors coated with goat anti-mouse IgG antibody (Includes 20 bottles of diluent.)

## OCTET BIOSENSORS SCREENING GRADE

- screening applications with off-rates < 15 minutes

Grade	Item	Part No.	UOM	Qty.	Description
Screen	Streptavidin High Binding FA Biosensors	18-5019	Tray	1	One tray of 96 biosensors coated with streptavidin for custom quantitation (Includes 1 bottle of kinetics buffer.)
Screen	Streptavidin High Binding FA Biosensors	18-5020	Pack	1	Five trays of 96 biosensors coated with streptavidin for custom quantitation (Includes 1 bottle of kinetics buffer.)
Screen	Streptavidin High Binding FA Biosensors	18-5021	Case	1	Twenty trays of 96 biosensors coated with streptavidin for custom quantitation. (Includes 4 bottles of kinetics buffer.)
Screen	Streptavidin FA Biosensors	18-5003	Tray	1	One tray of 96 biosensors coated with streptavidin for kinetic screening. (Includes 2 bottles of diluent.)
Screen	Streptavidin FA Biosensors	18-5008	Pack	1	Five trays of 96 biosensors coated with streptavidin for kinetic screening. (Includes 10 bottles of diluent.)
Screen	Streptavidin FA Biosensors	18-5009	Case	1	Twenty trays of 96 biosensors coated with streptavidin for kinetic screening. (Includes 40 bottles of diluent.)

## OCTET BIOSENSORS – KINETICS GRADE

- baseline stable for 1 hour

Grade	Item	Part No.	UOM	Qty.	Description
<b>THESE BIOSENSORS ARE PACKAGED AS 32 BIOSENSORS/ TRAY.</b>					
Kinetics	Streptavidin SBC Biosensors	18-5002	Tray	1	One tray of 32 biosensors coated with streptavidin for standard kinetic analysis. <i>(Includes 1 bottle of diluent.)</i>
Kinetics	Streptavidin SBC Biosensors	18-5006	Pack	1	Five trays of 32 biosensors coated with streptavidin for standard kinetic analysis. <i>(Includes 5 bottles of diluent.)</i>
Kinetics	Streptavidin SBC Biosensors	18-5007	Case	1	Twenty trays of 32 biosensors coated with streptavidin for standard kinetic analysis. <i>(Includes 20 bottles of diluent.)</i>
Kinetics	Amine Reactive Biosensors, Standard	18-5011	Tray	1	One tray of 32 biosensors compatible with amine reactive coupling chemistry for kinetic analysis. <i>(Reagents not included)</i>
Kinetics	Amine Reactive Biosensors, Standard	18-5014	Pack	1	Five trays of 32 biosensors compatible with amine reactive coupling chemistry for kinetic analysis. <i>(Reagents not included)</i>
Kinetics	Amine Reactive Biosensors, Standard	18-5015	Case	1	Twenty trays of 32 biosensors compatible with amine reactive coupling chemistry for kinetic analysis. <i>(Reagents not included)</i>
<b>THESE BIOSENSORS ARE PACKAGED AS 96 BIOSENSORS/ TRAY.</b>					
Kinetics	Amine Reactive Biosensors, Standard	18-5029	Tray	1	One tray of 96 biosensors compatible with amine reactive coupling chemistry for kinetic analysis. <i>(Reagents not included)</i>
Kinetics	Amine Reactive Biosensors, Standard	18-5030	Pack	1	Five trays of 96 biosensors compatible with amine reactive coupling chemistry for kinetic analysis. <i>(Reagents not included)</i>
Kinetics	Amine Reactive Biosensors, Standard	18-5031	Case	1	Twenty trays of 96 biosensors compatible with amine reactive coupling chemistry for kinetic analysis. <i>(Reagents not included)</i>

### Reagents & companion products

Item	Part No.	UOM	Qty.	Description
Sample Diluent	18-5028	Bottle	1	Diluent for sample dilution and biosensor hydration, 50 ml, packaged <i>(not for use with Amine Reactive Biosensors)</i>
IgG Calibrators	18-1030	Set	1	Set of 8 IgG Calibrators filled at 10 mL/vial. (plus a negative control)
Amine Coupling Reagent Kit	18 5017	Kit	1	Amine Coupling Reagent Kit is sufficient for 96 immobilizations onto Amine Reactive Biosensors
MES Buffer, pH 4.0	18-5025	Bottle	1	MES, 100mM, pH 4.0, packaged
MES Buffer, pH 5.0	18-5026	Bottle	1	MES, 100mM, pH 5.0, packaged
MES Buffer, pH 6.0	18-5027	Bottle	1	MES, 100mM, pH 6.0, packaged
Biosensor Dispenser	18-5016	Ea	1	Octet biosensor transfer tool

## Instruments

Item	Part No.	Description						
Octet Q System	30-5002	<p>Octet Q, Quantitation only instrument for label-free, real-time analysis. System includes the following:</p> <table border="0"> <tr> <td>(1) Octet Q instrument</td> <td>(1) Accessory Kit</td> </tr> <tr> <td>(1) Desktop computer</td> <td>(1) Installation</td> </tr> <tr> <td>(1) 19" LCD Monitor</td> <td>(1) Basic Operator Training</td> </tr> </table> <p>Includes Warranty for Octet Q System, 1 year coverage</p>	(1) Octet Q instrument	(1) Accessory Kit	(1) Desktop computer	(1) Installation	(1) 19" LCD Monitor	(1) Basic Operator Training
(1) Octet Q instrument	(1) Accessory Kit							
(1) Desktop computer	(1) Installation							
(1) 19" LCD Monitor	(1) Basic Operator Training							
Octet QK System	30-5003	<p>Octet QK, Quantitation and kinetics instrument for label-free, real-time analysis. System includes the following:</p> <table border="0"> <tr> <td>(1) Octet QK instrument</td> <td>(1) Accessory Kit</td> </tr> <tr> <td>(1) Desktop computer</td> <td>(1) Installation</td> </tr> <tr> <td>(1) 19" LCD Monitor</td> <td>(1) Basic Operator Training</td> </tr> </table> <p>Includes Warranty for Octet QK System, 1 year coverage</p>	(1) Octet QK instrument	(1) Accessory Kit	(1) Desktop computer	(1) Installation	(1) 19" LCD Monitor	(1) Basic Operator Training
(1) Octet QK instrument	(1) Accessory Kit							
(1) Desktop computer	(1) Installation							
(1) 19" LCD Monitor	(1) Basic Operator Training							

## 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
Octet Software Upgrade	50-5005	1	Octet Software Upgrade

## Operator Trainings

Item	Part No.	UOM	Qty.	Description
Training, Basic - Octet Q	30-6008	Ea	1	One day training course to provide basic overview of the Octet Q System and its main applications
Training, Basic - Octet QK	30-6005	Ea	1	One day training course to provide basic overview of the Octet QK System and its main applications
Training, Advanced - Octet QK	30-6007	Ea	1	Two days training course to provide basic overview of the Octet QK System and its main applications

## Ordering

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

Email: [sales@fortebio.com](mailto:sales@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

## Limitations on Use

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