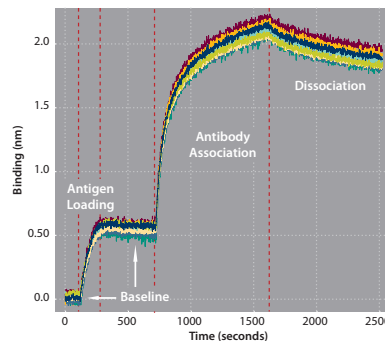


ForteBio's Octet System

Label-free, real-time detection of molecular interactions

KEY FEATURES OF BLI

- **Label-free detection**
BLI technology does not require the use of detection labels to measure the binding interaction at the biosensor surface.
- **Real-time analysis**
BLI continuously measures the protein binding at the biosensor surface throughout the interaction. Raw data is displayed in real time and the rapid analysis fits with process workflow.
- **Minimal interference**
BLI only detects binding at the sensor surface, so there is minimal interference from biological sample media. Proteins can be assayed in cell culture media or crude lysates without interference.
- **Automated**
Octet is automated to perform multi-step experimental protocols and complete data analysis. The system runs parallel, and up to 96 samples in unattended operation.
- **Easy to use**
Octet is fully automated and simple to operate. Flexible protocols using crude sample media are programmed in advance and do not require additional processing such as washing, centrifugation or long incubations to produce real-time results.



ForteBio's Octet systems incorporate a dedicated instrument for reading proprietary Bio-Layer Interferometry (BLI) plus a complete range of ready-to-use biosensors for real-time, label-free detection of protein:protein interactions.

By combining ready-to-use biosensors with sensitive optical detection, Octet can be broadly applied to a variety of research applications for protein:protein binding, affinity and kinetics. To meet your research and process development requirements, Octet is available in two configurations:

- Octet-Q: *quantitation only*
- Octet-QK: *quantitation and kinetics*

Contact ForteBio today to see what the Octet System can do for you!



fortéBIO™

ForteBio, Inc.
1360 Willow Road, Suite 205
Menlo Park, CA 94025

t: 888-OCTET-QK or 650-322-1360
f: 650-242.8706
web: www.fortebio.com

OCTET SYSTEM SPECIFICATIONS*

Technical Information and Specifications

| | |
|----------------------------|--|
| Detection technology | BioLayer Interferometry (BLI) based on fiber optic biosensors |
| Sensor type | Disposable, single-use sensors, application specific |
| Information provided | <ul style="list-style-type: none"> • Kinetic and affinity analysis (k_{obs}, k_a, k_d, K_D) • Kinetic screening for k_a or k_d • Binding specificity and binding cooperativity • Concentration monitoring of real-time binding (with no need for background subtraction) • Automated concentration determinations |
| Data presentation | <ul style="list-style-type: none"> • Plots displaying kinetic binding, equation fits, and residuals of fits • Tabulated kinetic data and data charts |
| Automation | 96 samples unattended operation |
| Sample types include: | Proteins, antibodies, peptides (>900 da), serum containing media (up to 10%), DMSO containing buffers, periplasmic fractions, untreated cell culture supernatants and crude cell lysates |
| Sample volume | 200 μ L at final dilution; non-destructive testing, easily recovered |
| Sample plate | Standard, 96-well, black, flat-bottom microplate |
| Orbital flow capacity | Static, or 100–1500 rpm |
| Analysis temperature range | Ambient + 2 $^{\circ}$ C to 40 $^{\circ}$ C, 1 $^{\circ}$ C increments |
| Sample refractive index | Independent of index changes in biological samples |

Instrument

| | |
|-------------------------|--|
| Dimensions (H x W x D) | 18.6" x 17" x 20.8" (47.5 cm x 43 cm x 53 cm) |
| Net weight | 50 lbs (23 kg) |
| Mains requirement/power | AC 100–240 V, 2.0–0.9 A, 50/60 Hz, single phase/120 w (240 peak) |

Data Handling and Storage

| | |
|---------------------|------------|
| PC operating system | Windows XP |
| Interfaces | RS232, USB |

Compliance

| | |
|------------------|---------|
| Safety standards | CE, CSA |
|------------------|---------|

Kinetics

| | |
|----------------------------------|---|
| Workflow | Up to 8 assays in parallel; 32 assays unattended operation |
| Analysis time per sample | Real-time kinetic binding experiments from 5 minutes to 3 hours |
| Association rate constant k_a | 10^3 to 10^7 $M^{-1} s^{-1}$ |
| Dissociation rate constant k_d | 10^{-5} to $10^{-1} s^{-1}$ |
| Sample concentration | 1 mM to 5 μ M |
| Molecular weight detection | >900 da |
| Baseline noise | 0.05 nm (RMS) |
| Baseline drift | < 0.1 nm/hour |

Kinetic Screening

| | |
|--------------------------|---|
| Workflow | 1–96 samples off-line batch immobilization for increased throughput |
| Analysis time per sample | Kinetic screening for 96 samples from 1 hour (depending on application) |
| Kinetic reporting | Rank ordering of clones for either k_a or k_d |

Human IgG Quantitation

| | |
|---------------------------|--|
| Workflow | Up to 8 assays in parallel; 96 assays unattended operation |
| Analysis time per sample | Quantitation in <2 minutes for 8 samples, <20 minutes for 96 samples |
| Quantitation range | 1 μ g/ml–200 μ g/ml |
| Precision range | 1 μ g/ml to 100 μ g/ml (CV <10%) application specific |
| Specificity (Fc specific) | Minimal cross reactivity with mouse, bovine, rat, goat and rabbit. Binds all IgG subclasses, minimal cross reactivity to IgM and IgE |

*All specifications are subject to change without notice.

