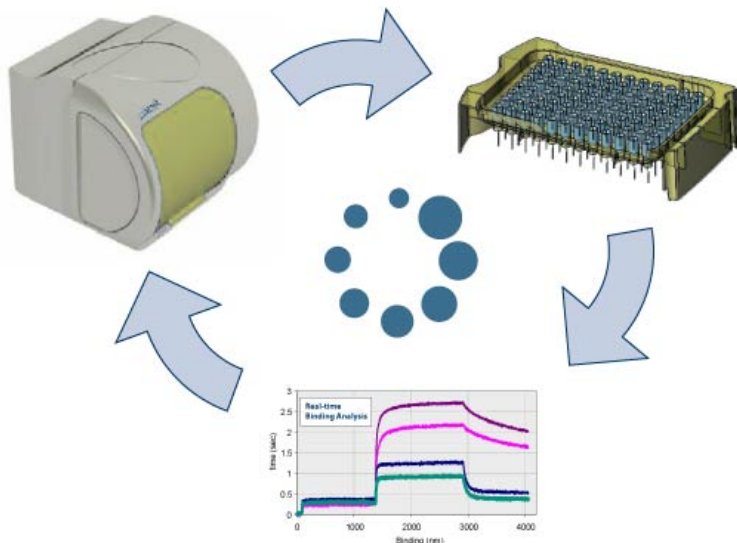




## ForteBio Octet System

Label-free, Real-time detection of molecular interactions



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

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

- Octet-Q                      Quantitation Only
- Octet-QK                    Quantitation and Kinetics

### 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**  
Because BLI only detects binding at the sensor surface, 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's system 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.

### For more information, contact:

ForteBio, Inc.  
1360 Willow Road, Suite 101  
Menlo Park, CA 94025

(650) 322-1360  
info@fortebio.com  
www.fortebio.com

**forté**BIO™