

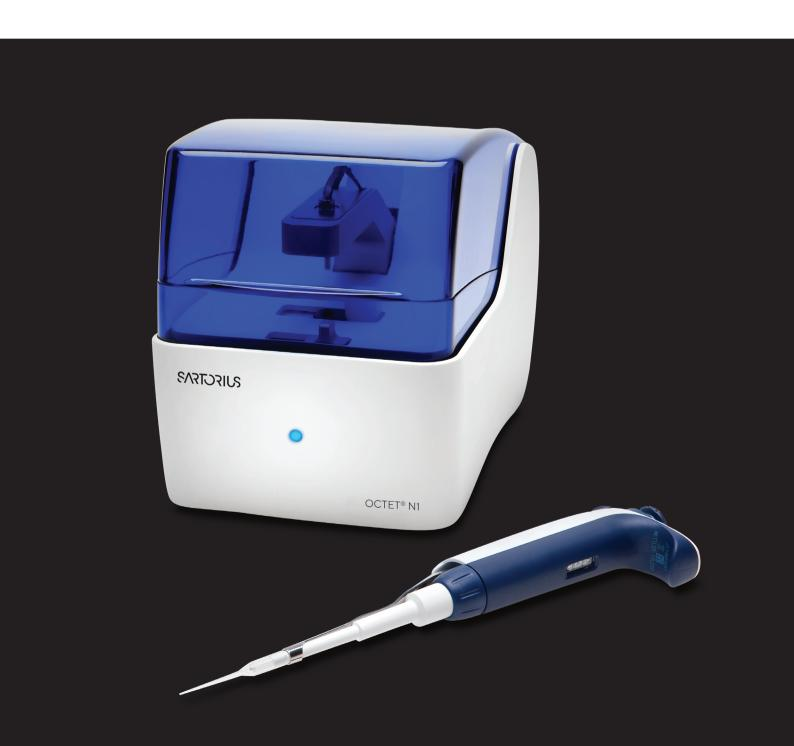
Label-Free Assays in a Drop

Protein Presence. Quantitation. Kinetics. Assay Development. Simplifying Progress



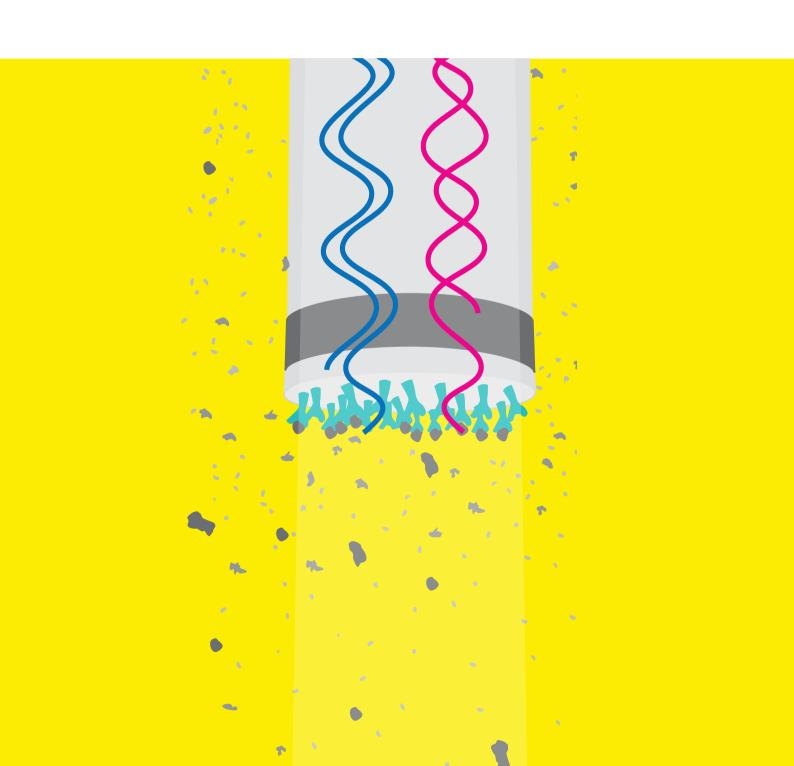
Drop. Read. Done.

The Octet® N1 System brings label-free analysis to individual researchers and is affordably priced so everyone can have one. With this personal assay system, you can run protein quantitation and kinetics experiments in only 4 μL of sample—right at your own lab bench.



It's Label-Free.

The Octet® N1 system detects molecules as they bind to the surface of its Dip and Read Octet® Biosensors using label-free BLI technology. This enables sensitive, realtime detection even in crude media. Octet® Biosensors are ready-to-use, disposable and available with a variety of surface chemistries to meet all your application needs.

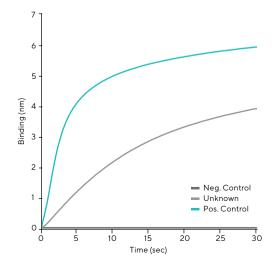


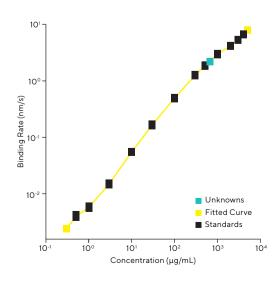
Protein Presence: Instantly.

Detect the presence or absence of active proteins in just a drop of sample, and easily identify specific proteins in complex solutions. Data from the Octet® N1 system complements existing methods like blots and gels, and its real-time binding curves give instant visual confirmation.

Protein Quantitation: Faster.

Measure protein concentration without labeling or secondary reagents—even in crude media. With a wide dynamic range and sensitivity down to a few ng/mL, the Octet® N1 system gives you accurate quantitation in seconds compared to traditional methods that can often take hours.



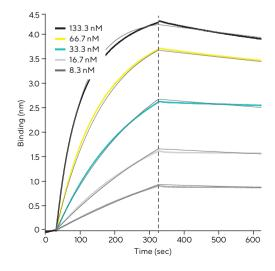


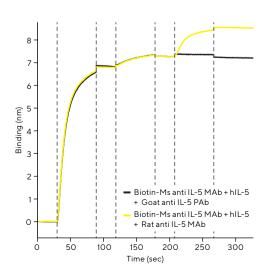
Kinetics: Effortless.

Get rate and affinity constants for binding interactions ($k_{\rm a}$, $k_{\rm d}$, $K_{\rm D}$) in minutes using 4 μ L of sample. With both a small price and a small size, the Octet® N1 system can be your own personal label-free kinetics system. Ready to go right out of the box, its simple workflows let you run assays right at your lab bench with ease.

Assay Development: Smarter.

Monitoring the effects of changing conditions on binding interactions in real-time, the Octet® N1 system lets you develop better immunoassays in only a few minutes. Find the best antibody pairs, identify the best assay parameters, and even build and optimize an ELISA step-by-step.





About the Octet® N1 System

Technology	Bio-Layer Interferometry (BLI)	
Sample volume	4 μL	
Dimensions	6.8" H × 6.0" W × 8.7" D (17.4 cm H × 15.3 cm W × 22.2 cm D)	
Weight	7.2 lb (3.3 kg)	
Power	100 – 240 V AC, 0.2 – 0.1 A, 50/60 Hz, 8 W (18 W peak)	

Biosensors

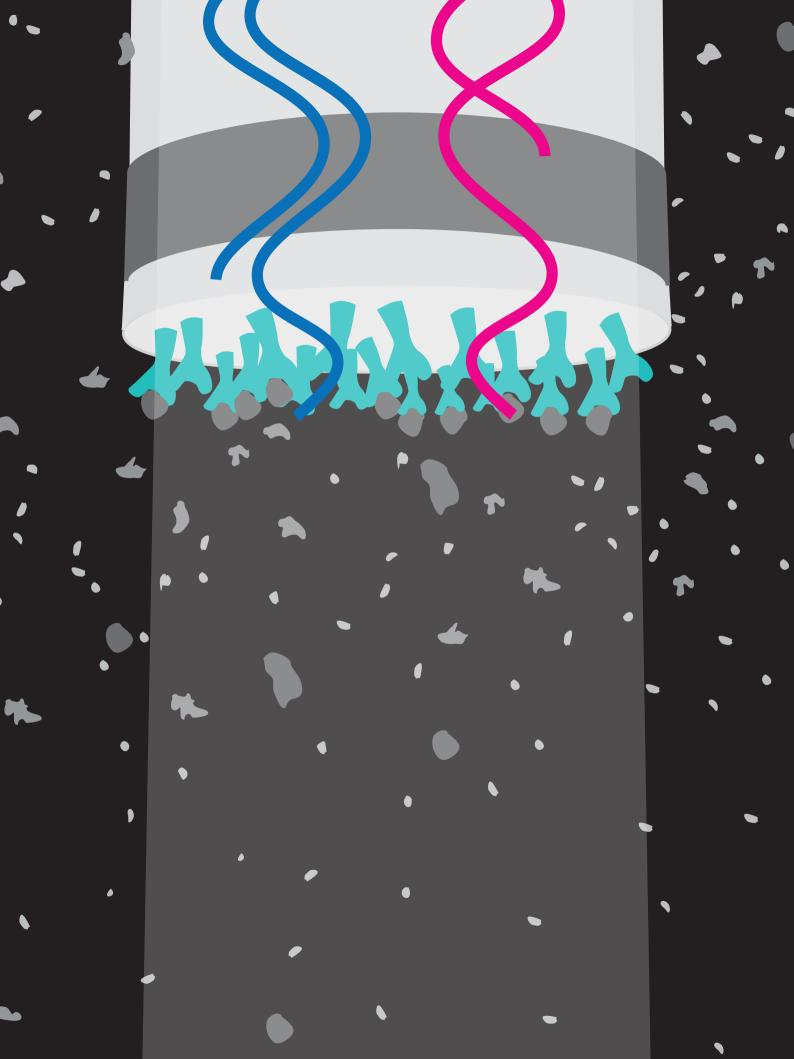
Octet® APS Biosensors	Octet® SAX Biosensors	Octet® ProA Biosensors
Octet® AR2G Biosensors	Octet® SAX2 Biosensors	Octet® ProG Biosensors
Octet® SSA Biosensors	Octet® AHQ Biosensors	Octet® ProL Biosensors
Octet® AHC Biosensors	Octet® AMQ Biosensors	Octet® FAB2G Biosensors
Octet® AMC Biosensors	Octet® HIS1K Biosensors	Octet® GST Biosensors
Octet® SA Biosensors	Octet® HIS2 Biosensors	Octet® NTA Biosensors

To Order

Octet® N1 system with Octet® N1 Software and starter kit: PN 45-5000 Visit www.sartorius.com/Octet-N1

Learn More

Visit www.sartorius.com/octet-support to get more information on the system, applications or BLI technology.



Germany

Sartorius Lab Instruments GmbH & Co. KG Otto-Brenner-Strasse 20 37079 Goettingen Phone +49 551 308 0

USA

Sartorius Corporation 565 Johnson Avenue Bohemia, NY 11716 Phone +1888 OCTET 75 Or +1 650 322 1360



For further contacts, visit www.sartorius.com/octet-support