

New Octet® R Series of Systems

The aggressive pace of scientific advancements and breakthroughs demands that labs around the globe produce high quality, reproducible data with limited budgets and short timelines. The Octet® platform utilizing Bio-Layer Interferometry (BLI) has served such needs for over a decade, by providing microfluidics-free, low maintenance and sensitive systems. These systems increase lab productivity, reduce cost and shorten experimental timelines. This platform is now expanded by introducing the new Octet® R series with upgradeable throughput to retain the initial system's investment value and futureproof your lab and research.

Biomolecular Interaction Analysis: From Basic Research to Drug Discovery to Manufacturing



Label-Free Interaction
Analysis



One Platform,
Many Applications



Fast Time to Results



Easy-to-Use



Low Cost of Ownership



5000+ Publications

New! Modular Design for
Upgradeable Throughput



Introducing the New Octet® R Series: The Octet® That Grows With Your Needs

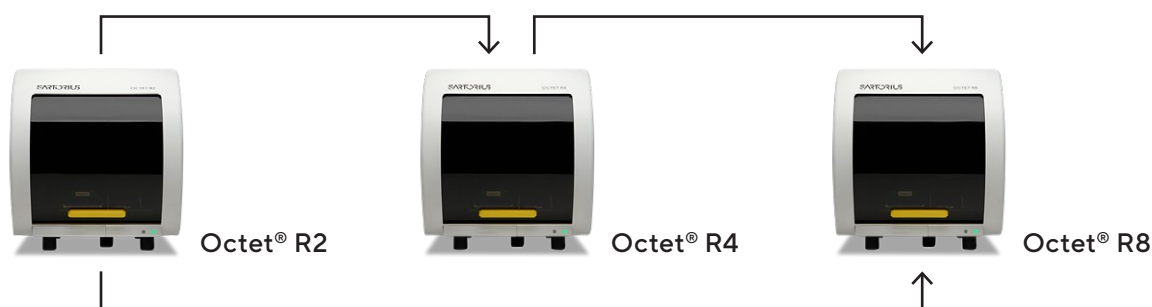
Find out more: www.sartorius.com/octet

Contact us for more information:
octet@sartorius.com

Upgrade Your Throughput With Minimal Downtime

The Octet® R series is ideal for academic start-up and low volume labs that are looking for high quality, label-free capabilities at varying price points, and the flexibility to upgrade upon increasing throughput needs.

Upgrade from a 2-channel to a 4-channel or 8-channel system, and a 4-channel to an 8-channel system within a single day by a field service engineer, without the hassle of trading in the existing model, thereby retaining the value of your current system.

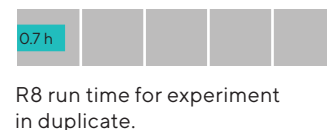
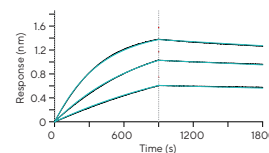
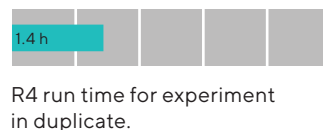
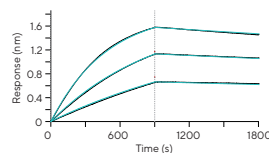
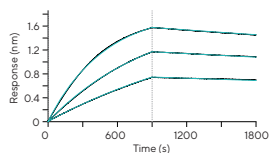


Octet® R Series Instrument Highlights

	Octet® R2	Octet® R4	Octet® R8
No. of channels	2	4	8
Analysis temperature range	15–40° C	15–40° C	15–40° C
Evaporation control	No	No	Yes
Sample plate format	96-well	96-well	96-well
Field upgradability	To R4, R8	To R8	–
GxP compatibility	No	No	Yes

Kinetics and affinity characterization of IgG1_k mAb binding to hCD64 across Octet R Series systems

$K_D = 0.15 \pm 0.02$ nM
 $k_a = (5.4 \pm 0.8) \times 10^5$ M⁻¹s⁻¹
 $k_d = (8.1 \pm 0.2) \times 10^{-5}$ s⁻¹



“Octet® provides the convenience of delivering a K_D in just 30 minutes. We cannot get that with ITC or SPR. We also love the ability to quickly reverse assay orientations to confirm binding or when one assay direction does not work well. Hard to do this in ITC as sample concentrations are required to be 10–200 fold above the K_D and are often not feasible in terms of required reagent amounts or reagent stabilities that will not withstand the duration of the experiment.”

–Dr. Arabela Grigorescu, Managing Director of Keck Biophysics Core Facility, Northwestern University, IL, USA