

SARTORIUS

AAV Capsid Titer on the Octet® 80% Faster Compared to ELISA

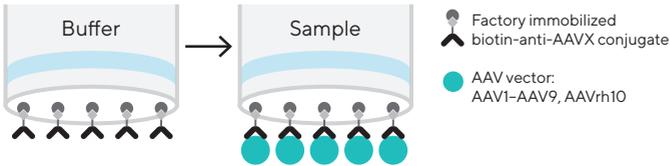


Bioprocess Five
Additional Batches
of AAV Samples
Annually

Your Key Benefits

- Reduce time-to-results by **> 80%** compared to ELISA for AAV capsid titer quantification
- Enable **near real-time** decisions on the bioprocess
- Reduce operator hands-on time by **50%**
- Reduce sample preparation time by **15–25%**

AAV Quantitation Workflow Using Octet® AAVX Biosensors



Octet® AAVX Biosensors can be used to quantitate 10 different AAV serotypes (AAV1 to AAV9 and AAVrh10) in both purified and crude cell culture samples with titer in the range of 8.5E8 – 1.0E13 vp/mL.

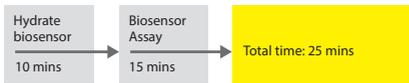
Further, Octet® AAVX Biosensors can be efficiently regenerated up to 10 – 20 times, which makes them an extremely useful and cost-effective solution for AAV capsid titer measurement.

Octet® AAVX Biosensors offer:

- An at-line, high-throughput one step AAV capsid titer quantification
- Rapid analysis of both crude and purified samples with minimal matrix interference
- Wide dynamic range: 8.5E8 – 1.0E13 vp/mL for most AAV serotypes
- Good correlation with ELISA

Flow chart of process steps and step time summary

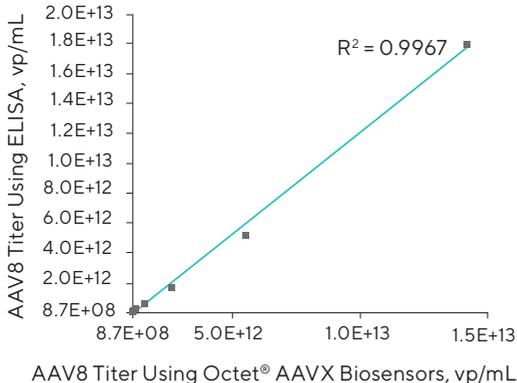
The Octet® BLI System



ELISA



Correlation of AAV8 Titer Measured Using the Octet® AAVX Biosensors vs. ELISA.



Octet® BLI systems are a complete solution for bioprocess analytics and are the industry standard for:

- Antibody titer determination
- Clone selection
- Protein and peptide quantitation
- Dynamic binding capacity determination
- Biotherapeutic drug activity assays
- Chromatography optimization
- Residual Protein A contaminant testing
- Host cell protein contaminant testing