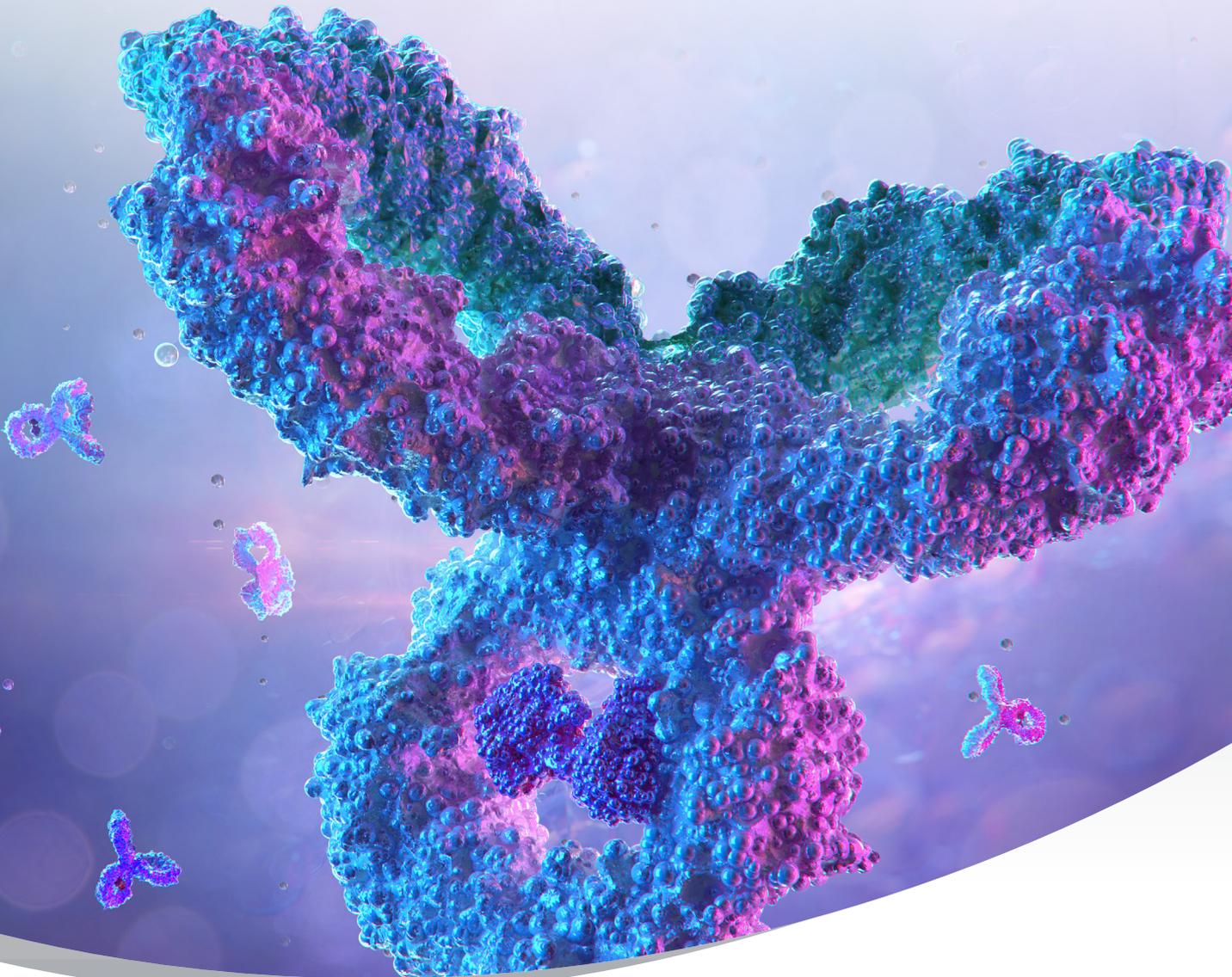




# Valita Aggregation Pure

## PROTEIN AGGREGATION DETECTION AND QUANTIFICATION ASSAY

SPEED • PERFORMANCE • SIMPLICITY  
AUTOMATION • 96-WELL FORMAT



ACCELERATING  
*answers*

 **BECKMAN  
COULTER**  
*Life Sciences*

# Rapid Determination of Protein Aggregation in Pure Solution

## Results in minutes not hours

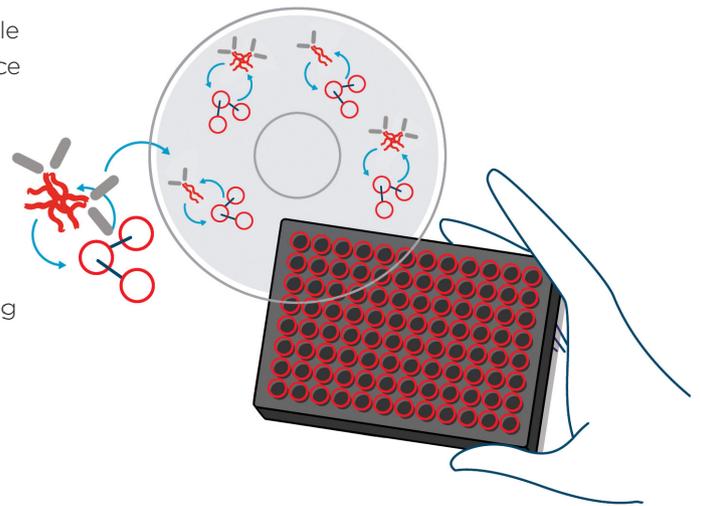
The Valita Aggregation Pure assay is a plate-based, 96-well screening tool that offers rapid, high throughput protein aggregation detection and quantification. The Valita Aggregation Pure assay is among the fastest and simplest aggregation detection tools on the market, and is compatible with high-performance microplate readers with Fluorescence Polarization functionality.

## Workflow benefits

- No additional reagents (just the plate!)
- No wash steps required
- Reduce project timelines with rapid in-lab/in-house testing

## Assay Benefits

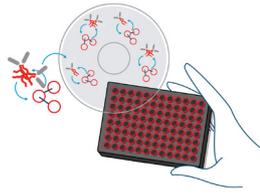
- **Simple:** Add, mix, read
- **Rapid:** Results in less than 15 minutes
- **Automation friendly:** 3 steps in standard plates
- **Cost effective:** All that's needed is a plate reader



## VALITA AGGREGATION PURE ASSAY TECHNICAL FEATURES

- **Limit of detection:**  $\geq 0.5\%$  at 1 g/L
- **Plate format:** 96-well format
- **Accuracy:**  $> 95\%$  for human IgG1 samples with levels of aggregation  $> 0.5\%$  at 1 g/L
- **Assay time:** Add, mix, read  $< 15$  minutes
- **Fluorescent binding peptide:** Proprietary molecule (Excitation max = 560 nm, Emission max = 590 nm)
- **Validated protein constructs:** Human IgG1, human IgG2, human IgG4, Fab, F(ab')<sub>2</sub>, IgG hybrid, IgG Fc fusion, murine IgG, rabbit IgG
- **Recommended sample concentration:** 200 mg/L to 2 g/L
- **Validated assay matrix:** Purified solution

# PROCESS DEVELOPMENT WORKFLOW



Quantify aggregation in minutes, not hours with the Valita Aggregation Pure assay

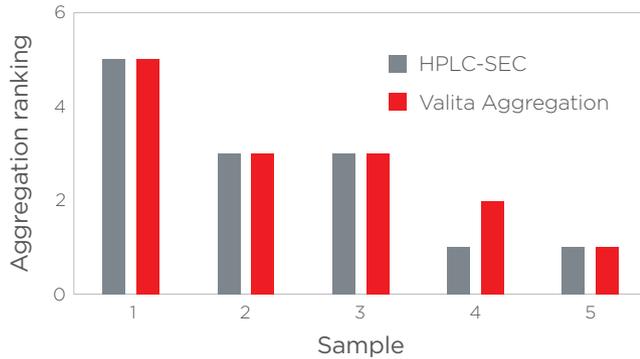


Other possible applications include: late cell line development, stability studies, formulation screening

|                     | Valita Aggregation Pure Assay | DLS/SLS          | HPLC - SEC |
|---------------------|-------------------------------|------------------|------------|
| Assay Time          | 15 mins                       | 1-2 hours        | >6 hours   |
| Throughput          | ++++                          | ++               | +          |
| Automation Friendly | Yes                           | Semi-automatable | No         |
| Number of Reagents  | 1                             | >1               | >2         |
| CAPEX               | \$                            | \$\$             | \$\$\$     |

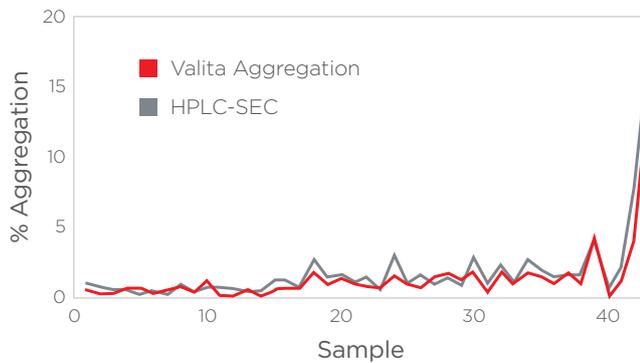
# BENCHMARK DATA

Ranking: Valita Aggregation Pure assay vs HPLC-SEC



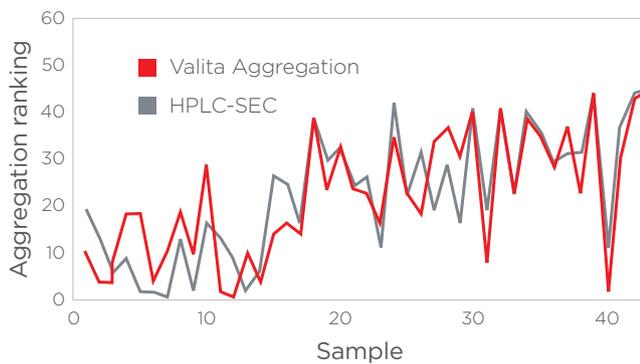
Protein Aggregation ranking, comparing 5 samples with < 1% aggregation measured on both Valita Aggregation Pure and HPLC-SEC

% Aggregation: Valita Aggregation Pure assay vs HPLC-SEC



Percentage Protein Aggregation, comparing 43 samples measured on both Valita Aggregation Pure and HPLC-SEC.

Ranking: Valita Aggregation Pure assay vs HPLC-SEC



Protein Aggregation ranking, comparing 43 samples measured on both Valita Aggregation Pure and HPLC-SEC.