

# PROTEOSTAT<sup>®</sup>

## Protein aggregation standards

### The only *commercially available* protein aggregation standards assay!

The PROTEOSTAT<sup>®</sup> Protein Aggregation standards are ideal for reliable and accurate quantification of protein aggregation in solution. The standards are used to determine low levels of aggregated protein in a sample by comparing the assay response of a test sample to that of the standard curve; comprised of standards with known concentrations of aggregated IgG. Protein test samples and protein aggregation standards should be analyzed in the same manner, by mixing them with detection reagent and using a fluorescence microplate reader to measure the fluorescence intensity values. As with any protein assay, different protein aggregates will elicit slightly greater or lesser fluorescence intensity response based upon their inherent amino acid composition and sequence. It is recommended that a standard curve be prepared each time the PROTEOSTAT<sup>®</sup> Protein Aggregation assay (Prod. No. ENZ-51023) is performed.

The PROTEOSTAT<sup>®</sup> Protein Aggregation standards contain stabilized, high-quality reference samples for generating an accurate standard curve. The standard curve is formulated from different amounts of aggregated IgG, combined with monomeric IgG. Once reconstituted, the samples contain 12.5%, 6.25%, 3.13%, 1.56%, 0.78%, 0.39%, 0.20% and 0% aggregated protein, while the total protein concentration of each standard is maintained at 1mg/ml. The total protein concentration of the IgG standard has been calibrated by direct comparison to an internal protein standard, in order to ensure lot-to-lot consistency and accuracy. Orthogonal characterization methods, including laser-based nanoparticle tracking and micro-flow imaging, have been employed to validate monomer integrity during manufacturing.

Citations: 2

[View Online »](#)

### Ordering Information

[Order Online »](#)

ENZ-51039-KP002

2x96 wells

Manuals, SDS & CofA

[View Online »](#)

- Stabilized, high-quality reference IgG in known concentrations
- Developed for use with the PROTEOSTAT<sup>®</sup> Protein Aggregation Assay
- Easy to use - *Simply add water!*

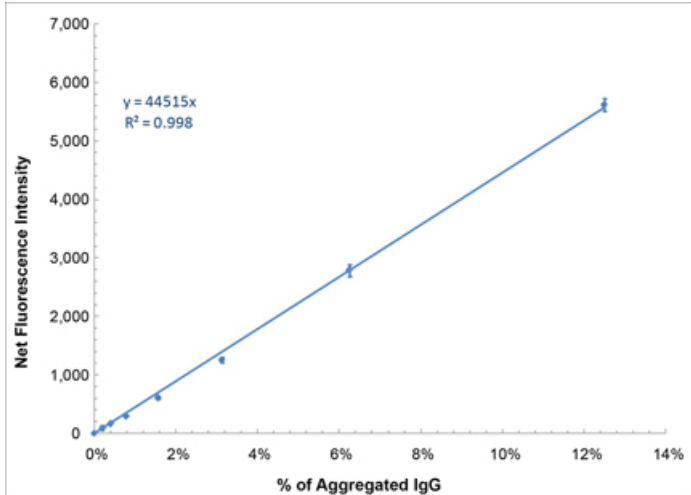


Figure 1: Typical standard curve produced using the PROTEOSTAT<sup>®</sup> Protein Aggregation standards.



[Particle analysis of therapeutic protein formulations with ImageStreamX<sup>®</sup> Imaging Flow Cytometry and the PROTEOSTAT<sup>®</sup> Protein Aggregation Assay](#)

[Prediction of Aggregation Propensity and Monitoring of Aggregation of Antibody-Drug Conjugates \(ADC\) using ProteoStat<sup>®</sup> Reagents](#)

This product is manufactured and sold by ENZO LIFE SCIENCES, INC. for research use only by the end-user in the research market and is not intended for diagnostic or therapeutic use. Purchase does not include any right or license to use, develop or otherwise exploit this product commercially. Any commercial use, development or exploitation of this product or development using this product without the express prior written authorization of ENZO LIFE SCIENCES, INC. is strictly prohibited.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)