



Sensitive and Reliable Measurement of Protein A

Protein A ELISA Kit (ADI-900-057)

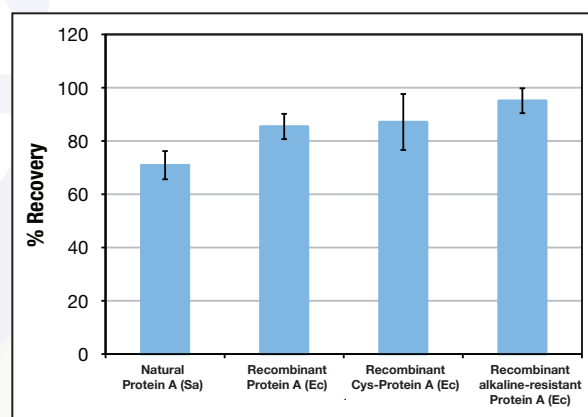
Sensitive, accurate, and reproducible assay to measure natural and recombinant Protein A

Staphylococcus aureus Protein A is a 42 kDa cell wall constituent that is characterized by its binding affinity to the Fc portion of some immunoglobulins, especially the IgG class. Protein A is typically used to purify antisera and in commonly used immunodetection and visualization techniques. Protein A participates in a number of protective biological functions including anti-tumor, toxic and carcinogenic activities, thus necessitating the removal of potential Protein A contaminants from antibody preparations for therapeutic use. The Protein A ELISA kit is a sensitive, accurate, and reproducible sandwich assay used to measure natural and recombinant Protein A residuals in monoclonal antibody preparations. The Protein A ELISA kit has been extensively validated in humanized IgG preparations to ensure accurate recovery and sensitivity.

Key Features

- Superior sensitivity
 - Detect 1ppm Protein A residuals in human IgG preparations
 - Lower standard curve range versus conventional methods
- Flexible detection
 - Detect 4 different Protein A constructs
 - Efficiently detect natural and recombinant Protein A variants
- Great recovery and reproducibility
 - Up to 100% recovery in the presence of human IgG
 - Reproducible results with 4-5% CV
- Time and money savings
 - Fast results in < 3 hours
 - Low cost per test

Assay Recognition of Different Protein A Constructs



RELATED PRODUCTS

PRODUCT NAME	PRODUCT #	PRODUCT NAME	PRODUCT #
BIOLIKE® Conserved mAb Sequence Conformational ELISA Kit	ENZ-KIT123	PROTEOSTAT® Protein Aggregation Assay	ENZ-51023
CHO HCP ELISA Kit	ENZ-KIT128	PROTEOSTAT® Protein Aggregation Standards	ENZ-51039
<i>E. coli</i> HCP ELISA Kit	ENZ-KIT127	PROTEOSTAT® Protein Refolding and Aggregation Sensing Kit	ENZ-51040
PEGylated Protein ELISA Kit	ADI-900-213	PROTEOSTAT® Thermal Shift Stability Assay Kit	ENZ-51027

Accelerate Your Research:
www.enzolifesciences.com

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Phone: 800.942.0430
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen, Switzerland
Phone: +41 61 926 8989
Fax: +41 61 926 8979
info-eu@enzolifesciences.com