

# Human Serum Albumin ELISA kit

**Easy-to-use ELISA kit for detecting the presence of Human Serum Albumin contamination during manufacturing of biologics**

The Human Serum Albumin ELISA kit is a quantitative sandwich enzyme-linked immunosorbent assay technology with results in 2.5 hours.

Human Serum Albumin is the serum albumin found in human blood. It constitutes about half of serum proteins, and is produced in the liver. Albumin transports hormones, fatty acids, and other compounds, buffers pH, and maintains oncotic pressure, among other functions. Human Serum Albumin (HSA) is a highly water-soluble globular monomeric plasma protein with a relative molecular weight of 67 KDa. Among nanoparticulate carriers, HSA nanoparticles have long been the center of attention in the pharmaceutical industry due to their ability to bind to various drug molecules, great stability during storage and in vivo usage, no toxicity and antigenicity, biodegradability, reproducibility, scale up of the production process and a better control over release properties. In addition, significant amounts of drug can be incorporated into the particle matrix because of the large number of drug binding sites on the albumin molecule.

## Ordering Information

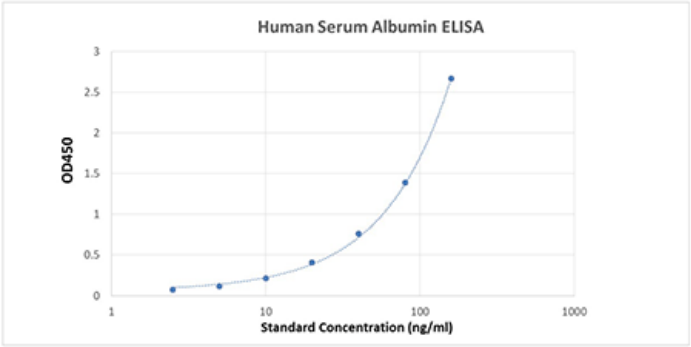
[Order Online »](#)

ENZ-KIT221-0001	96 wells
-----------------	----------

## Manuals, SDS & CofA

[View Online »](#)

- Specifically recognizes Human Serum Albumin
- Measures as little as 1ng/ml
- Results in 2.5 hours



# Handling & Storage

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## Product Details

Alternative Name	HSA
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of Human Serum Albumin in plasma, serum, cell culture supernatant and other samples.
Assay Time	2.5 hours
Compatibility	This product is compatible with the <a href="#">Absorbance 96 Plate Reader</a> .
Contents	Microtiter plate pre-coated with anti-Human Serum Albumin Antibody, HSA Standard, Standard Diluent buffer, Anti-Human Serum Albumin:HRP Conjugate, Sample Diluent buffer, 20 × Wash Buffer, TMB Substrate, Stop Solution
Crossreactivity	Does not detect Bovine Albumin or Mouse Serum
Sensitivity	1 ng/ml (range 1.56 – 500 ng/ml)
Species Reactivity	Human
Wavelength	450 nm

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-  
fr@enzolifesciences.com](mailto:info-fr@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)