

# Bovine Serum Albumin ELISA kit

Easy-to-use BSA ELISA kit for optimizing purification processes during bioproduction, as well as testing the final product

The BSA ELISA kit is a colorimetric competitive enzyme immunoassay kit with results in 2 hours.

Bovine serum albumin (BSA) is a serum albumin protein derived from cows. The mature BSA protein contains 583 amino acids. BSA, like other serum albumins, is critical in providing oncotic pressure within capillaries, transporting fatty acids, bilirubin, minerals and hormones, and functioning as both an anticoagulant and an antioxidant. BSA has numerous biochemical applications including ELISAs (Enzyme-Linked Immunosorbent Assay), immunoblots, and immunohistochemistry. BSA is a major component of fetal bovine serum (FBS), which is commonly used in mammalian cell culture media for the production of recombinant biologics.

## Ordering Information

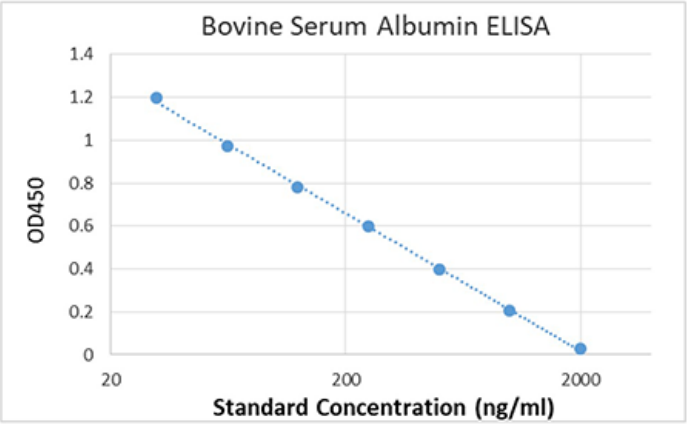
[Order Online »](#)

ENZ-KIT220-0001	96 wells
-----------------	----------

## Manuals, SDS & CofA

[View Online »](#)

- Competitive enzyme-linked immunosorbent assay for specific quantification of Bovine Serum Albumin
- Minimum detection limit - 80 ng/ml
- Results in less than 1h



## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name BSA

Application Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of Bovine Serum Albumin in plasma, serum, adherent and suspension cell culture

Assay Time 45 minutes

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

**Contents** Anti-BSA antibody coated Microtiter plate, BSA Standard, BSA:HRP Conjugate (concentrated), Sample Diluent, HRP Conjugate Dilution Buffer, Wash buffer concentrate (20x), TMB Substrate, Stop solution

**Crossreactivity** No significant cross-reactivity or interference between BSA and analogues was observed.

**Sensitivity** 94 ng/ml (range 156 – 5000 ng/ml)

**Species Reactivity** Bovine

**Wavelength** 450 nm



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)