

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

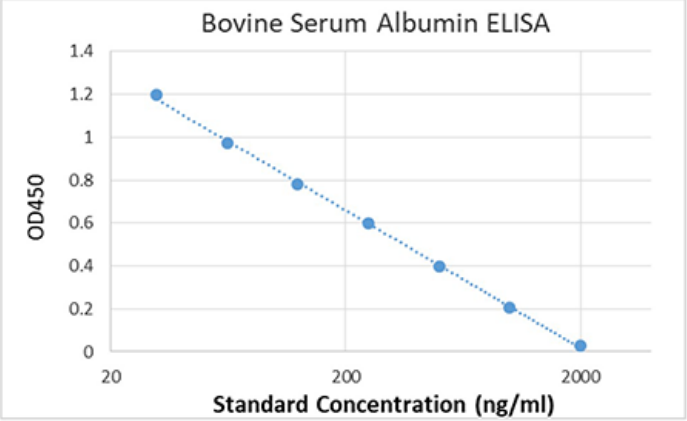
[Order Online »](#)

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

Last modified: June 4, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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