
Bovine Serum Albumin Polyclonal Antibody

(Catalog # A55766)

Background

Serum albumin, the main protein of plasma, has a good binding capacity for water, Ca²⁺, Na⁺, K⁺, fatty acids, hormones, bilirubin and drugs. Its main function is the regulation of the colloidal osmotic pressure of blood. Major zinc transporter in plasma, typically binds about 80% of all plasma zinc.

Description

Bovine Serum Albumin Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300. 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Bovine

Isotype

IgG

Uniprot ID

P02769

Purification

Protein G purified

Immunogen

Serum Albumin purified from Bovine serum (Native Protein)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

BSA

Application

ELISA
