Bovine Serum Albumin 30% Solution

Can be used as a stabilizing agent for proteins and enzymes, or added to blocking buffers to minimize non-specific antibody binding in immuno-detection procedures such as western blotting and cell staining.



Product attributes

Storage Conditions

Store at 2 to 8 °C

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

Bovine serum albumin (BSA) has uses in many applications such as stabilizing proteins in solution, blocking non-specific binding in immuno-assays, and as a protein standard in various protein assays.

- Stabilizer for proteins and enzymes in solution
- Blocking agent for immuno-assays (ie, western blot, IHC, IF)
- Minimizes non-specific antibody binding
- IgG- and protease-free
- 100 mL solution of 30% BSA

Bovine serum albumin (BSA) is commonly used as a stabilizing agent for proteins and enzymes, including dilute solutions of antibody. It is also used as a blocking agent to reduce non-specific antibody binding in immuno-detection procedures such as western blotting, immunofluorescence, and IHC. It is also suitable for use in protease sensitive assays such as RIA and EIA, as well as nucleic acid hybridization. This BSA is IgG- and protease-free.

Note: This product contains <0.1% sodium azide. When diluted to a typical working concentration of BSA, the final sodium azide concentration is not expected to inhibit HRP or affect most applications. However, if you are concerned about the presence of sodium azide affecting your experiment, we recommend preparing a BSA stock solution from powder using <u>Bovine Serum Albumin Fraction V</u>.

This item is only available for customers in the USA.

This datasheet was generated on November 7, 2025 at 01:50:54 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/bovine-serum-albumin-30-solution/