

# Product Information

## Bovine Serum Albumin, 30% Solution

**Catalog Number:** 22014

**Unit Size:** 100 mL

### Storage and Handling

Store at 4°C. Product is stable for at least 6 months from date of receipt when stored as recommended.

**Color and Form:** Amber solution

### Product Description

Bovine serum albumin (BSA) is commonly used as a stabilizing agent for proteins such as antibodies and enzymes, and as a blocking agent to reduce non-specific protein binding in immunoassays such as ELISA, immunofluorescence staining, and western blotting. Supplied as a 30% solution of IgG- and protease-free BSA in water. BSA is typically used at 1-5% final concentration for blocking.

**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 (Cat. no. 22013).

### Related Products

Catalog number	Product
22013	Bovine Serum Albumin Fraction V
23013	TrueBlack® WB Blocking Buffer Kit
23007	TrueBlack® Lipofuscin Autofluorescence Quencher
23014	TrueBlack® Plus Lipofuscin Autofluorescence Quencher
23012	TrueBlack® IF Background Suppressor System (Permeabilizing)
22010	Fish Gelatin Blocking Agent, 10X
22011	Fish Gelatin Powder
22012	Dry Milk Powder
22002	Tween® 20

Please visit our website at [www.biotium.com](http://www.biotium.com) for information on our life science research products, including our line of TrueBlack® background reducers, fluorescent CF® Dye antibody conjugates, cellular stains, and more.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.