

Product Information

Bovine Serum Albumin, 30% Solution

Catalog Number: 22014

Unit Size: 100 mL

Color and Form: Amber solution

Storage and Handling

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

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.

Cat.#	Product Name	Unit Size
22010	Fish Gelatin Blocking Buffer, 10X	100 mL
22011	Gelatin, cold water fish	100 g
22012	Non-fat dry milk powder	4 x 25 g
22002	Tween® 20	50 mL
23001	EverBrite™ antifade mounting medium	10 mL
23002	EverBrite™ antifade mounting medium with DAPI	10 mL

TWEEN is a registered trademark of Uniqema Americas LLC.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.