

Product Information

Bovine Serum Albumin, Fraction V

Catalog Number: 22013

Unit Size: 50 g

Color and Form: Off-white solid

Storage and Handling

Store at 4°C. Stable for at least 12 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. BSA is typically used at 1-5% final concentration for blocking. IgG- and protease-free.

Related Products

Cat.#	Product Name	Unit Size
40061-T	RedDot™2 Far Red Nuclear Counterstain, 200X in DMSO, Trial Size (15-20 tests)	25 uL
23001	EverBrite™ Mounting Medium	10 mL
23002	EverBrite™ Mounting Medium with DAPI	10 mL
23003	EverBrite™ Hardset Mounting Medium	10 mL
23004	EverBrite™ Hardset Mounting Medium with DAPI	10 mL
23005	CoverGrip™ Coverslip Sealant	15 mL
22005	Mini Super ^{HT} Pap Pen 2.5 mm tip, ~400 uses	1 pen
22006	Super ^{HT} Pap Pen 4 mm tip, ~800 uses	1 pen
22015	Fixation Buffer	100 mL
22016	Permeabilization Buffer	100 mL
22017	Permeabilization and Blocking Buffer	100 mL
22010	10% Fish Gelatin Blocking Buffer	100 mL
22011	Fish Gelatin Powder	2 x 50 g
22014	30% Bovine Serum Albumin Solution	100 mL
22002	Tween®-20	50 mL

Please visit www.biotium.com to view our full selection of products featuring bright and photostable fluorescent CF™ dyes, including secondary antibodies and Mix-n-Stain™ antibody labeling kits, and R-PE dye conjugates. Biotium also offers a variety of apoptosis and cell viability assays for flow cytometry analysis, including mitochondrial membrane potential dyes, fluorescent Annexin V conjugates, and NucView™488 Caspase-3 Substrate for live cells.

TWEEN is a registered trademark of Uniqema Americas LLC.

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