

Product Information

Bovine Serum Albumin, CF™ dye Conjugates

Catalog No.	Product Description	Unit Size
20289	Bovine Serum Albumin, CF488A Conjugate	5x 1 mg
20290	Bovine Serum Albumin, CF594 Conjugate	5x 1 mg
20291	Bovine Serum Albumin, CF640R Conjugate	5x 1 mg
20292	Bovine Serum Albumin, CF680 Conjugate	5x 1 mg

Concentration: 2 mg/mL upon addition of 0.5 mL PBS.

Form: lyophilized powder.

Spectral Property

$\lambda_{abs}/\lambda_{em}$ (in pH 7.4 PBS buffer)

Product Description	Abs _{max} nm	Em _{max} nm
Bovine Serum Albumin, CF488A Conjugate	490	515
Bovine Serum Albumin, CF594 Conjugate	593	614
Bovine Serum Albumin, CF640R Conjugate	642	662
Bovine Serum Albumin, CF680 Conjugate	681	698

Storage and Handling

Product is stable for 2 years at -20°C with dessicant. Upon reconstitution of the lyophilized powder in 0.5 mL PBS, store at 4°C and protect from light.

Product Description

Biotium offers Bovine serum albumin (BSA) conjugates that labeled with our superior CF dyes. Bovine serum albumin (BSA) conjugates are commonly used as tracers in applications where physical dimensions are important. The conjugates can also be used for quantitative studies of electroporation and measurement of plasma volume in rats. Other applications include studying the rate of receptor-mediated endocytosis and exocytosis and following intracellular protein processing.

BSA conjugates are supplied as lyophilized powder in a 1 mg size. Solution can be made by dissolving the powder in 1X PBS buffer. It is good practice to centrifuge the protein conjugate solutions briefly in a microcentrifuge before use; only the supernatants should then be added to the experiment. This step will eliminate any protein aggregates that may have formed during storage, thereby reducing nonspecific background staining.

A full selection of CF™ dye labeled products including secondary antibodies, antibody labeling kits, and other bioconjugates such as phalloidins, annexin V and a-bungarotoxin are also available. Please visit the Biotium website at www.biotium.com for details.

Other Related Products

You may also be interested in the following related products:

CF Dye Product	Application	Unit Size
NHS esters and maleimide	Labeling antibodies and other biomolecules	1 μmole
CF™ dye aminoxy and hydrazide	Labeling oligonucleotides and other biomolecules	1 mg
CF™ dye protein labeling kits	Labeling antibodies and other biomolecules	3 labelings
Anti-Mouse, Rabbit, Goat, Rat, Sheep, Guinea Pkg, Human or Swine IgG (H+L), 2 mg/mL, whole antibody	Microscopy, flow cytometry and Western blotting	0.25 or 0.5 mL
Anti-Chicken IgY* (H+L), 2 mg/mL, whole antibody [* IgY is the original designation for the IgG like protein found in serum and egg yolk]	Microscopy, flow cytometry and Western blotting	0.25 or 0.5 mL
F(ab') ₂ fragment, Anti-Mouse or Anti-Rabbit, 2 mg/mL	Microscopy, flow cytometry and Western blotting	0.25 mL
EverBrite™ Mounting Medium (with or without DAPI)	Microscopy	10 mL
Annexin V Conjugates	Labeling apoptotic cells for microscopy or flow cytometry	0.5 mL (50 μg/mL)
Phalloidin Conjugates	Microscopy	300 U
α-Bungarotoxin	Labeling Ach-R for microscopy	0.5 mg