



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Bovine Serum Albumin CF® Dye Conjugates



Product attributes

Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Green, Red, Far-red, Near-infrared

Bright and photostable fluorescent BSA conjugates for applications in which well-defined molecular weight fluid phase tracers are needed. These applications include studying the rate of endocytosis and exocytosis and following vesicle trafficking.

Product Description

CF® dye bovine serum albumin (BSA) are bright and photostable fluorescent BSA conjugates for applications in which well-defined molecular weight fluid phase tracers are needed.

- Choice of 4 CF® dyes from green to near-infrared
- Superior CF® dyes are bright, photostable, and water-soluble
- For endocytosis and exocytosis studies

These applications include studying the rate of endocytosis and exocytosis and following vesicle trafficking.

Superior CF® Dyes

Biotium's next-generation CF® dyes were designed to be highly water-soluble with advantages in brightness and photostability compared to Alexa Fluor®, DyLight®, and other fluorescent dyes. Learn more about [CF® Dyes](#).

Bovine Serum Albumin CF® Dye Conjugates

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@488A	490/515 nm	5 x 1 mg	20289	CF@488A Features
CF@594	593/614 nm	5 x 1 mg	20290	CF@594 Features
CF@640R	642/662 nm	5 x 1 mg	20291	CF@640R Features
CF@680	681/698 nm	5 x 1 mg	20292	CF@680 Features

CF is a registered trademark of Biotium, Inc. Alexa Fluor, Texas Red, and DyLight are registered trademarks of Thermo Fisher Scientific.

References

J Biomed Mater Res A. (2019), [doi: 10.1002/jbm.a.36726](https://doi.org/10.1002/jbm.a.36726).

Download a list of [CF® dye references](#).

This datasheet was generated on June 10, 2026 at 03:24:15 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/bovine-serum-albumin-conjugate-cf-dye-conjugates/>