

Recombinant Human Vitamin D-Binding Protein Conjugates

Vitamin D-Binding Protein (Vitamin D-BP, also known as GC Globulin) binds monomeric G-actin in fixed and permeabilized cells. Available conjugated to a selection of bright and photostable CF® Dyes for fluorescence microscopy.



Product Description

Recombinant Human Vitamin D-Binding Protein (Vitamin D-BP, also known as GC Globulin) binds monomeric G-actin (1-3). Fluorescent conjugates of Vitamin D-BP can be used to stain monomeric G-actin in fixed and permeabilized cells.

- Labels monomeric G-actin in fixed and permeabilized cells
- Can be co-stained with fluorescent phalloidin for comparing G-actin and F-actin distribution
- Developed and optimized for immunofluorescence
- Available with bright and photostable CF® Dyes, including near-infrared CF®740

Vitamin D-BP staining may also be used in combination with fluorescent phalloidin staining of F-actin to visualize the distribution of unpolymerized G-actin relative to actin filaments. Vitamin D-BP is more specific for staining G-actin compared to fluorescent conjugates of DNase-I, which binds to DNA in addition to G-actin. Biotium's Recombinant Human Vitamin D-BP Conjugates are labeled with a selection of our bright and photostable CF® Dyes for fluorescence microscopy.

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](#).

Recombinant Human Vitamin D-Binding Protein Conjugates

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	1 mL (100 ug)	70087	CF®488A Features
CF®594	593/615 nm	1 mL (100 ug)	70088	CF®594 Features
CF®640R	642/663 nm	1 mL (100 ug)	70089	CF®640R Features
CF®647	652/668 nm	1 mL (100 ug)	70090	CF®647 Features
CF®740	742/767 nm	1 mL (100 ug)	70091	CF®740 Features

For staining F-actin, Biotium recommends [ActinBrite™ High Affinity Phalloidin Conjugates](#) which were designed to preserve strong F-actin binding over conventional phalloidin conjugates. With ActinBrite™, samples can be imaged after for a month or more (depending on the conjugate and mounting method)—making delayed imaging easier and more dependable.

Browse Biotium's comprehensive catalog of [fluorescent bioconjugates](#) as well [novel and classic organelle stains](#).

References

1. J Biol Chem, 255(6), 2270(1980), [S0021-9258\(19\)85885-4](#)
2. J Cell Biol, 123(1), 1(1993), [10.1083/jcb.123.1.1](#)
3. Biochim Biophys Acta, 1452(1), 12(1999), [10.1016/s0167-4889\(99\)00119-6](#)

This datasheet was generated on August 31, 2025 at 02:16:49 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/recombinant-human-vitamin-d-binding-protein-conjugates/>

Product attributes

Probe cellular localization	G-Actin, Cytoskeleton
For live or fixed cells	For fixed cells
Assay type/options	Tissue staining
Detection method/readout	Fluorescence microscopy
Cell permeability	Membrane impermeant
Fixation options	Fix before staining (formaldehyde), Permeabilize before staining
Colors	Green, Red, Far-red, Near-infrared
Antibody/conjugate formulation	100 ug/mL in PBS/0.01% rBSA/0.05% sodium azide
Product origin	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant Vitamin D Binding Protein (His tag) produced in HEK293 cells
Shipping condition	Room temperature